Breastmilk-promoted bifidobacteria produce aromatic lactic acids in the infant gut

3

4 Martin F. Laursen^{1#}, Mikiyasu Sakanaka^{1,2}, Nicole von Burg³, Daniel Andersen⁴, Urs Mörbe³,

5 Aymeric Rivollier³, Ceyda T. Pekmez⁵, Janne Marie Moll⁴, Kim F. Michaelsen⁵, Christian

6 Mølgaard⁵, Mads Vendelbo Lind⁵, Lars O. Dragsted⁵, Takane Katayama^{2,6}, Henrik L. Frandsen¹,

7 Anne Marie Vinggaard¹, Martin I. Bahl¹, Susanne Brix⁴, William Agace³, Tine R. Licht^{1*} and

- 8 Henrik M. Roager^{1,5#*}
- 9 [#] These authors contributed equally

10 Affiliations

- ¹ National Food Institute, Technical University of Denmark, Kgs. Lyngby, Denmark
- ² Faculty of Bioresources and Environmental Sciences, Ishikawa Prefectural University, Ishikawa
 921-8836, Japan
- ³ Department of Health Technology, Technical University of Denmark, Kgs. Lyngby, Denmark
- ⁴ Department of Biotechnology and Biomedicine, Technical University of Denmark, Kgs. Lyngby,
 Denmark
- ⁵ Department of Nutrition, Exercise and Sports, University of Copenhagen, Frederiksberg C,
- 18 Denmark
- ⁶ Graduate school of Biostudies, Kyoto University, Kyoto 606-8502, Japan

20 Correspondence

- * Tine R. Licht, Tel. + 45 35887186, <u>trli@food.dtu.dk</u>, Kemitorvet, 2800 Kongens Lyngby,
- 22 Denmark
- * Henrik M. Roager, Tel. + 45 35324928, <u>hero@nexs.ku.dk</u>, Rolighedsvej 30, 1958 Frederiksberg
 C, Denmark
- 25

ABSTRACT 26

Breastfeeding profoundly shapes the infant gut microbiota, which is critical for early life immune 27 development. However, few breastmilk-dependent microbial metabolites mediating host-microbiota 28 interactions are currently known. We here demonstrate that breastmilk-promoted Bifidobacterium 29 30 species convert aromatic amino acids (tryptophan, phenylalanine and tyrosine) into their respective aromatic lactic acids (indolelactate, phenyllactate and 4-hydroxyphenyllactate) via a previously 31 unrecognised aromatic lactate dehydrogenase. By longitudinal profiling of the gut microbiota 32 33 composition and metabolome of stool samples of infants obtained from birth until 6 months of age, we show that stool concentrations of aromatic lactic acids is determined by the abundance of human 34 milk oligosaccharide degrading Bifidobacterium species containing the aromatic lactate 35 dehydrogenase. Finally, we demonstrate that stool concentrations of Bifidobacterium-derived 36 indolelactate are associated with the capacity of infant stool samples to activate the aryl 37 hydrocarbon receptor, a receptor important for maintenance of intestinal homeostasis and immune 38 system development. These findings open up new directions towards understanding the role of 39 breastmilk-promoted *Bifidobacterium* in mediating host-microbiota interactions in early life. 40

INTRODUCTION 41

Human breastmilk is a perfectly adapted nutritional supply for the infant¹. Breastfeeding provides 42

children with important short-term protection against infections, and may also provide long-term 43

- metabolic benefits^{1,2}. These benefits may partly be mediated through the gut microbiota, since 44
- breastfeeding is the strongest determinant of gut microbiota composition and function during 45

infancy^{3–5}. Human breastmilk contains human milk oligosaccharides (HMOs), which are complex, 46

- highly abundant sugars serving as substrates for specific microbes including certain species of 47
- *Bifidobacterium*⁶. This co-evolution between bifidobacteria and the host, mediated by HMOs, to a 48
- large extent directs the colonization of the gut in early life, which has critical impact on the immune 49
- system⁷. Depletion of specific microbes, including *Bifidobacterium*, in early life has been associated 50
- with increased risk of allergy and asthma development in childhood^{8,9}, and is suggested to 51
- compromise immune function and lead to increased susceptibility to infectious disease^{10,11}. Despite 52
- Bifidobacterium dominating the gut of breastfed infants and being widely acknowledged as 53
- beneficial, mechanistic insights on the contribution of these bacteria and their metabolites to 54 immune development during infancy remain limited. Recent studies show that microbial aromatic
- 55 amino acid metabolites including tryptophan-derived indoles¹², via activation of the aryl 56
- hydrocarbon receptor (AhR), can fortify the intestinal barrier^{13,14}, protect against pathogenic
- 57
- infections^{15,16} and influence host metabolism^{13,17,18}, which makes this group of microbial 58
- metabolites of particular interest in the context of early life. 59
- 60 Here, we show that breastmilk-promoted Bifidobacterium species, via a previously unrecognised
- aromatic lactate dehydrogenase, produce aromatic lactic acids including indolelactate (ILA) in 61
- substantial amounts in the infant gut, and that stool concentrations of this metabolite are associated 62
- with the capacity of infant stool to activate AhR, which is known to impact immune development in 63
- 64 early life.

65 **RESULTS**

66 Bifidobacterium species associate with aromatic amino acid catabolites during late infancy

To explore interactions between breastfeeding status, gut microbial composition and metabolism of 67 aromatic amino acid in early life, we employed 16S rRNA amplicon sequencing to infer gut 68 microbiota composition and a targeted ultra-performance liquid chromatography mass spectrometry 69 (UPLC-MS) metabolomics approach to quantify 19 aromatic amino acids and derivatives thereof 70 71 (Supplementary Table 1) in stool samples from 59 healthy Danish infants from the SKOT I cohort¹⁹. The SKOT I infants included were born full term, 9.1 ± 0.3 (mean \pm SD) months of age at 72 sampling, and 40.7% were still partially breastfed (Supplementary Data 1a,b). After stratification 73 of the 9 months old infants based on breastfeeding status (partially breastfed versus weaned), 74 75 Principal Coordinates Analysis (PCoA) of weighted UniFrac distances showed a significant separation across the first PC-axis ($r^2 = 0.093$, p < 0.001, Adonis test; Fig. 1a), which was due to an 76 increasing gradient in relative abundance of *Bifidobacterium* in breastfed infants ($r^2 = 0.397$, p < 77 0.001, Adonis test; Fig. 1b). Other metadata (age, gender, mode of delivery, current formula intake 78 and age of introduction to solid foods) did not explain gut microbiota variation to the same degree 79 as breastfeeding status (Supplementary Data 1c.d). Principal Component Analysis (PCA) of faecal 80 81 aromatic amino acid metabolite concentrations also revealed separation according to breastfeeding status, which was largely driven by three aromatic lactic acids, 4-hydroxyphenyllactic acid (4-OH-82 83 PLA), phenyllactic acid (PLA) and indolelactic acid (ILA) (Fig. 1c). Correlation analysis revealed that Bifidobacterium, but no other bacterial genera, were significantly associated with faecal 84 concentrations of all three aromatic lactic acids (4-OH-PLA, PLA and ILA), in addition to 85 indolealdehyde (IAld) (Fig. 1d and Supplementary Data 1e). Bifidobacterium species 86 (Supplementary Fig. 1a and Supplementary Data 1f) significantly enriched in the breastfed 87 88 infants; B. longum, B. bifidum, and B. breve, were positively associated with the faecal concentrations of aromatic lactic acids (4-OH-PLA, PLA and ILA) and IAld (cluster 1 in Fig. 1e), 89 but negatively associated with the faecal concentrations of aromatic propionic acids, aromatic 90 amino acids and to a lesser degree with aromatic acetic acids (cluster 2 in Fig. 1e). In contrast, post 91 92 weaning type Bifidobacterium species, including B. adolescentis, B. animalis/pseudolongum and B. *catenulatum* group^{20,21}, were not significantly associated with aromatic lactic acids nor 93 breastfeeding status (Fig. 1e). These associations were in agreement with the observation that the 94 concentrations of the three aromatic lactic acids were higher in the faeces of breastfed than in 95 96 weaned infants (Supplementary Fig. 1b). Furthermore, the abundances of the three aromatic lactic acids in infant urine (Supplementary Fig. 2-4) showed similar positive associations between 97 relative abundances of breastmilk-promoted Bifidobacterium species (Supplementary Fig. 1c). In 98 addition, faecal and urinary levels of ILA were positively correlated (rho = 0.68, p < 0.0001), 99 showing that faecal levels of this metabolite are reflected systemically. Consistently, urine 100 abundance of ILA, but not of PLA and 4-OH-PLA, were significantly higher in breastfed compared 101 to weaned infants (Supplementary Fig. 1b). Together, this suggests that specific breastmilk-102 promoted Bifidobacterium species in the gut of infants convert aromatic amino acids to the 103 corresponding aromatic lactic acids (Fig. 1f). 104

105



107

108 Figure 1. Breastfeeding associates with gut microbiota composition and aromatic amino acid catabolism in 9 109 months old infants

- a-b, Principal coordinate analysis plots of weighted UniFrac distances based on OTUs from faecal samples of 9 months 110
- 111 old infants participating in the SKOT cohort (n = 59). Samples are coloured according to (a) breastfeeding status with
- 112 ellipses indicating 80%CI of data points for partially breastfed (red, n = 24) and weaned (blue, n = 35) infants or (b) 113 relative abundance of the genus Bifidobacterium.
- 114 c, Principal component analysis plot of concentrations (µmol/mg faeces) of aromatic amino acids and their catabolites
- 115 in SKOT faecal samples, coloured according to breastfeeding status with ellipses indicating 80%CI of data points for
- 116 breastfed (red, n = 24) and weaned (blue, n = 35) infants. Loadings are shown with arrows. See abbreviations in (d).
- 117 d, Heatmap illustrating Spearman's Rank correlation coefficients between the relative abundance of Bifidobacterium
- 118 and concentrations of aromatic amino acids and their catabolites in SKOT faecal samples (n=59). The concentration of
- 119 each metabolite (µmol/mg faeces) is presented as the median with 95%CI.
- 120 e, Heatmap illustrating hierarchical clustering of Spearman's Rank correlation coefficients between the relative
- 121 abundance of the different Bifidobacterium species and selected microbial-derived aromatic amino acid catabolites. Bar
- 122 plots are showing relative abundance (mean+SD) of the Bifidobacterium spp., stratified according to breastfeeding 123 status, with statistical significance evaluated by Mann-Whitney U test.
- 124 f, The pathway of aromatic amino acid catabolism by gut microbes (modified from Smith and Macfarlane 1996⁹⁵, Smith
- and Macfarlane 1997%, and Zelante et al. 2013¹⁶). Asterisks indicate statistical significance: * p<0.05, ** p<0.01, *** 125
- 126 p<0.001, **** p<0.0001.
- 127 See also Supplementary Figure 1-4.

128 Bifidobacterium species produce aromatic lactic acids in vitro

129 To confirm the ability of *Bifidobacterium* species detected in infants to produce aromatic lactic

- acids, *Bifidobacterium* type strains were grown anaerobically in a medium containing all three
- aromatic amino acids with either glucose or HMOs as sole carbohydrate sources. Analyses of
- 132 culture supernatants revealed that ILA, PLA and 4-OH-PLA were produced mainly by *B. bifidum*,
- 133 *B. breve, B. longum* subsp. *longum, B. longum* subsp. *infantis* and *B. scardovii* (Fig. 2a), in
- accordance with the associations observed in the 9 months old infants (Fig. 1e). Other
- 135 Bifidobacterium species, namely B. adolescentis, B. animalis subsp. lactis, B. animalis subsp.
- 136 animalis, B. dentium, B. catenulatum, B. pseudocatenulatum, B. pseudolongum subsp.
- 137 *pseudolongum* produced only low amounts of these metabolites (Fig. 2a). The ability of
- 138 *Bifidobacterium* species to produce high levels of the aromatic lactic acids was generally
- 139 convergent with the ability to utilize HMOs as carbohydrate source (**Fig. 2a**), supporting the link
- 140 between breastmilk-promoted bifidobacteria and production of aromatic lactic acids. None of the
- 141 downstream products of the aromatic lactic acids (**Fig. 1f**) were detected in any of the culture
- supernatants.

143 An aromatic lactate dehydrogenase is responsible for the aromatic lactic acid production

- 144 Since it has been reported that an L-lactate dehydrogenase (LDH) in *Lactobacillus* spp. can convert
- 145 phenylpyruvic acid to PLA^{22} , we hypothesized that a corresponding enzyme was present in
- phenylpyruvic acid to PLA²², we hypothesized that a corresponding enzyme was present in *Bifidobacterium* species. Alignment and phylogenetic analysis of all genes annotated as *ldh* in the *Bifidobacterium* type strains included in this study, revealed four clusters (Fig. 2b). Whereas all *Bifidobacterium* genomes contain an *ldh* responsible for conversion of pyruvate to lactate (here
- 149 designated as type 1 *ldh*) in the bifidobacterial fructose-6-phosphate shunt^{23,24}, some species have
- an extra *ldh*, here designated as type 2, type 3 and type 4, respectively. In agreement with the *in*
- *vitro* fermentations (**Fig. 2a**), all prominent aromatic lactic acid-producing *Bifidobacterium* species contain the type 4 *ldh*. Interestingly, genomic analysis of the *Bifidobacterium* type strains revealed
- that the type 4 *ldh* gene is part of a genetic element containing an amino acid transaminase gene
- 154 (suspected to be responsible for converting the aromatic amino acids into aromatic pyruvic acids)
- and a haloacid dehydrogenase gene (of unknown importance) (**Supplementary Fig. 5**), which has been indicated to constitute an operon in *B. breve*²⁵. Cloning of the type 4 *ldh* gene from *B. longum*
- subsp. *infantis*^T into a vector transformed into *E. coli* revealed that the expression of the type 4 ldh
- gene indeed resulted in the appearance of PLA, 4-OH-PLA and ILA in the culture supernatant (Fig.
- **2c**). To verify the type 4 *ldh* dependent production of aromatic lactic acids in *Bifidobacterium*
- species, we generated a type 4 *ldh* insertional mutant strain by homologous recombination in *B*.
- 161 *longum* subsp. *longum* 105-A (**Supplementary Fig. 6**), a genetically tractable strain containing the 162 type 4 LDH (**Supplementary Fig. 7**)^{26,27}. Cultivation of the wild-type (WT), the type 4 *ldh* mutant
- type 4 LDH (**Supplementary Fig. 7**)^{26,27}. Cultivation of the wild-type (WT), the type 4 *ldh* mutar strain and a complemented type 4 *ldh* mutant strain in a medium containing the three aromatic
- amino acids (**Supplementary Fig. 6**) confirmed that type 4 *ldh* disruption did not impair growth
- (Fig. 2d). ILA, PLA and 4-OH-PLA accumulated in the supernatant of the WT and of the
- 166 complemented type 4 *ldh* mutant strains, but not in the type 4 *ldh* mutant (**Fig. 2e**). Importantly, the
- type 4 *ldh* mutant was not significantly compromised in its ability to convert pyruvate to lactate
- 168 (Fig. 2e), supporting the distinct role of type 4 *ldh* in converting aromatic pyruvic acids.

Purification and characterization of the recombinant type 4 LDH enzyme revealed that it had a mass 169 of 33.9 kDa (Supplementary Fig. 8a), while the native molecular mass was estimated to be 72.9 170 kDa by size exclusion chromatography, indicating dimer formation in solution (Supplementary 171 172 Fig. 8b). No metal requirement was observed, the optimal pH was 8.0-8.5 and the enzyme was most stable at 37°C (Supplementary Fig. 8c-e). Heterotrophic effects were neither observed for 173 fructose-1.6-bisphosphate (an allosteric effector for type 1 LDH) nor for several intermediates for 174 aromatic amino acid synthesis^{23,24} (Supplementary Fig. 9). However, we found that phosphate 175 served as a positive effector suggesting that type 4 LDH is an intracellular enzyme (Supplementary 176 Fig. 10a-c). Assay at the different phosphate concentrations revealed the type 4 LDH is a K-type 177 allosteric enzyme (Supplementary Fig. 10b,c). The catalytic rate (k_{cat}) was moderate to high for 178 the aromatic pyruvic acid substrates, but very low for pyruvate (Fig. 2f,g), in accordance with the 179 non-impaired lactate production observed for the type 4 *ldh* mutant (Fig. 2e). Production of ILA, 180 PLA and 4-OH-PLA from the respective aromatic pyruvic acid substrates was verified by high-181 performance liquid chromatography (HPLC) (Supplementary Fig. 10d). The enzyme showed 182 highest affinity (lowest $K_{0.5}$) for indole pyruvic acid, but highest catalytic rate for 4-OH-phenyl 183 pyruvic acid in the presence of 100 mM phosphate (Fig. 2f,g). However, maximal catalytic 184 efficiency ($k_{cat}/K_{0.5}$) was observed for indole pyruvic acid (190 s⁻¹ mM⁻¹), followed by 4-OH-phenyl 185 pyruvic acid (16 s⁻¹ mM⁻¹) and phenyl pyruvic acid (11 s⁻¹ mM⁻¹), suggesting preference for indole 186 pyruvic acid. The observed Hill coefficient ($n_{\rm H} = 1-1.4$) for all substrates indicate weak positive 187 cooperativity under the conditions tested. Collectively, these results show that the type 4 *ldh* gene 188 encodes an aromatic lactate dehydrogenase responsible for the production of ILA, PLA and 4-OH-189 PLA in Bifidobacterium species associated with breastfeeding. We therefore suggest that the type 4 190 *ldh* gene should be re-classified as an aromatic lactate dehydrogenase gene (*aldh*). 191

192





193

a, *In vitro* production of indolelactic acid (ILA), phenyllactic acid (PLA) and 4-hydroxyphenyllactic acid (4-OH-PLA)

- by *Bifidobacterium* spp. type strains in modified MRS medium (MRSc) with 2% (w/v) glucose or a mix human milk
- 197 oligosaccharides as sole carbohydrate source. For the type strains of *B. adolescentis*, *B. animalis* subsp. animalis, *B.*
- animalis subsp. *lactis, B. dentium* and *B. catenulatum* no or very poor growth ($OD_{600nm} < 0.4$) was observed with HMOs
- as carbohydrate source. Mean of three biological replicates is shown.
- **b**, Neighbor-Joining phylogenetic tree of all genes in the *Bifidobacterium* spp. type strains annotated as L-lactate
- 201 dehydrogenases (*ldh*). The four clusters are designated type 1-4.
- 202 c, Production of ILA, PLA and 4-OH-PLA by E. coli LMG194 cells transformed with an inducible vector lacking
- 203 (empty vector) or containing the type 4 *ldh* (Type4_*ldh*⁺) from *B. longum* subsp. *infantis* DSM 20088^T in LB-medium 5
- h post-induction of gene expression by addition of L-arabinose and supplementation with the aromatic pyruvic acids
- 205 (indolepyruvate, phenylpyruvate and 4-OH-phenylpyruvate). Bars show mean ± SD of three biological replicates.
- **d**, Growth curves of *Bifidobacterium longum* subsp. *longum* 105-A (wild type), its isogenic insertional type 4 *ldh*
- 207 mutant (Type4_*ldh* mutant) and the type 4 *ldh* mutant strain complemented with the type 4 *ldh* gene (Complemented). 208 Curves show mean \pm SD of three biological replicates and doubling times reported as mean \pm SD.
- e, Production of ILA, PLA, 4-OH-PLA and lactate by wild type, Type4 *ldh* mutant and the complemented strain in
- early stationary phase cultures. Bars show mean \pm SD of three biological replicates. Statistical significance was
- 211 evaluated by one-way ANOVA, with * p<0.05, ** p<0.01, *** p<0.001, **** p<0.0001.
- 212 f-g, Enzyme kinetics of the type 4 lactate dehydrogenase. f, substrate saturation curves of type 4 LDH with indole
- 213 pyruvate, phenylpyruvate, 4-OH-phenylpyruvate or pyruvate as substrates. Data is one representative of two
- independent assays. g, kinetic parameters of type 4 LDH with the different substrates. Data are mean \pm SD of two
- 215 independent assays.
- ND, not detected
- 217 See also Supplementary Figure 5-10.

218 Bifidobacterium species govern aromatic lactic acid profiles during early infancy

To study the dynamics of *Bifidobacterium* species establishment and aromatic lactic acids in 219 infants, we established the Copenhagen Infant Gut (CIG) cohort including 25 healthy breast- or 220 mixed fed infants, which were sampled every 2-4 weeks from birth until the age of 6 months 221 (Supplementary Data 2a) for microbiome profiling and targeted metabolite quantification 222 including aromatic lactic acids (Supplementary Fig. 11 and Supplementary Table 2). A total of 223 145 operational taxonomic units (OTUs) were detected by 16S rRNA amplicon sequencing, 224 however, collapsing of OTUs with identical taxonomic classifications and using a cut-off of average 225 226 relative abundance of 0.1%, resulted in the identification of 39 bacterial species/taxa, representing 97.5% of the total community (Supplementary Fig. 11a and Supplementary Data 2b). As 227 expected, the gut microbiota was highly dominated by *Bifidobacterium* (average of 64.2 %) and 228 among the top 10 dominating taxa, *B. longum* (38.5 %), *B. breve* (9.1 %), *B. bifidum* (7.9 %), *B.* 229 catenulatum group (6.4 %) and B. dentium (1.7 %) were found (Supplementary Fig. 11a), with the 230 remaining *Bifidobacterium* spp. being assigned to *B. scardovii* (0.24 %), *B. adolescentis* (0.15 %) 231 and B. animalis/pseudolongum (0.10%) (Supplementary Data 2b,c). Although the relative 232 abundance of Bifidobacterium increased with time, on average the microbial composition and 233 Shannon diversity did not change dramatically during the six months (Supplementary Fig. 11b). 234 235 However, the subject specific gut microbiota profiles revealed a highly individual species composition (Supplementary Fig. 11c) and 48% of the variation in community structure was 236 explained by subject (weighted UniFrac, $r^2 = 0.48$, p < 0.001, Adonis, Supplementary Fig. 12a,b). 237 Indeed, the difference in microbial community structure between faecal samples were primarily 238 239 driven by the abundance of *Bifidobacterium* as assessed by PCoA of weighted UniFrac distances (r²

240 = 0.48, p < 0.001, Adonis, Fig. 3a). The non-phylogenetic Bray-Curtis dissimilarity analysis revealed a separation of the communities based on abundance of the five dominating 241 Bifidobacterium species (B. longum, B. bifidum, B. breve, B. catenulatum group and B. dentium) 242 (Fig. 3b and Supplementary Fig. 12c-h). Community abundance of *B. longum*, *B. bifidum* and *B.* 243 breve, but not B. catenulatum group and B. dentium (Fig. 3b) matched the measured faecal 244 concentrations of aromatic lactic acids (Fig. 3c). Consistently, faecal concentrations of ILA, PLA 245 and 4-OH-PLA increased concurrently with an increase in absolute abundance of infant type 246 247 Bifidobacterium species (defined as the summarized abundance of B. longum, B. bifidum, B. breve and *B. scardovii*) from birth to around 6 months of age (Fig. 3d, upper panels). The gut microbiota 248 of individuals dominated by infant type *Bifidobacterium* spp. was more stable over time than in 249 individuals with a gut microbiota not dominated by infant type *Bifidobacterium* spp. (p < 0.0001, 250 Mann-Whitney U test) (Supplementary Fig. 12i-j). Using solely samples from breastfed infants, 251 faecal abundances of HMO residuals showed a progressive decline with age, concurrent with the 252 progressive increase in infant type *Bifidobacterium* species (Fig. 3d, lower panels). By repeated 253 measure correlation analyses²⁸ (Supplementary Fig. 13) and partial Spearman's Rank correlation 254 analyses²⁹ adjusted for age, we confirmed that the abundance of infant type *Bifidobacterium* species 255 were positively associated with faecal levels of ILA, PLA and 4-OH-PLA and negatively associated 256 with abundances of HMOs in faeces (Fig. 3e). Both subspecies of *B. longum* were associated with 257 the aromatic lactic acids, but mainly B. longum subsp. infantis was associated with the HMO 258 residuals in faeces (Fig. 3e). We have thus established a link between breastfeeding, degradation of 259 260 HMOs, abundance of specific infant type Bifidobacterium spp. and concentrations of aromatic lactic acids in early infancy. 261

Examination of the Bifidobacterium and aromatic lactic acid dynamics in each of the 25 infants 262 during the first 6 months of life (Fig. 3f-h and Supplementary Fig. 14-15) revealed that breastfed 263 infants early colonised by infant type Bifidobacterium species consistently showed high 264 concentrations of aromatic lactic acids in faeces (Fig. 3f and Supplementary Fig. 15a,b). In 265 contrast, infants with delayed infant type Bifidobacterium species colonization showed considerably 266 lower concentrations of the aromatic lactic acids, in particular of ILA, despite breastfeeding (Fig. 267 3g and Supplementary Fig. 15c). Among the latter, CIG08 and CIG09 were twins, born late 268 preterm, and dominated by an OTU assigned to *Clostridium neonatale* (Supplementary Fig. 11c 269 and Supplementary Data 2b) in accordance with previous reports on C. neonatale overgrowth³⁰ 270 and delayed *Bifidobacterium* colonization³¹⁻³⁴ in preterm infants. CIG07 who also showed delayed 271 colonization with infant type Bifidobacterium, was mixed fed throughout the whole period and 272 predominantly colonised with E. coli and Clostridium spp. (Supplementary Fig. 11c). CIG18 had 273 274 relatively low faecal concentrations of aromatic lactic acids until age 172 days, when B. breve replaced *B. dentium* (Fig. 3g), consistent with the fact that *B. dentium* lacks the *aldh* gene while *B.* 275 *breve* contains it (**Fig. 2a,b**). Finally, in the three infants treated with antibiotics during our study, 276 Bifidobacterium species abundance were temporarily decreased simultaneously with reduced 277 278 concentrations of the aromatic lactic acids (Fig. 3h). This indicates that bifidobacterial aromatic lactic acid production is compromised by pre-term delivery, exposure to antibiotics and formula 279 supplementation. 280



284 Fig. 3. Infant type *Bifidobacterium* species determine aromatic lactic acid concentrations during early infancy

- **a-c**, Principal coordinate analysis plots of weighted UniFrac (a) or Bray-Curtis (b-c) distances/dissimilarities (n=234[#]),
- coloured according to relative abundance of *Bifidobacterium* (a-b) or log₁₀ transformed absolute concentration
- 287 [µmol/mg faeces] of aromatic lactic acids (sum of indolelactic acid (ILA), phenyllactic acid (PLA) and 4-
- 288 hydroxyphenyllactic acid (4-OH-PLA)) in the Copenhagen Infant Gut (CIG) cohort. Dashed lined circles indicate
- communities dominated (relative abundance > 50%) either by *B. longum*, *B. bifidum*, *B. breve*, *B. catenulatum* group or
- 290 *B. dentium (B. adolescentis, B. scardovii and B. animalis/pseudolongum* never dominated any of the communities, see
- **Supplementary Fig 11-12**). #6 samples were omitted from the analyses due to low read counts (<8000).
- **292 d**, Temporal development in absolute abundance of infant type *Bifidobacterium* spp. (defined as the sum of absolute
- abundances of *B. longum*, *B. breve*, *B. bifidum* and *B. scardovii*), faecal concentrations of aromatic lactic acids (ILA,
- 294 PLA and 4-OH-PLA, n=240) and relative faecal abundance of HMOs (2'FL/3FL, 2'/3-O-fucosyllactose; 3'SL/6'SL,
- 295 3'/6'-O-sialyllactose; LNT/LNnT, lacto-N-tetraose/ lacto-N-*neo*tetraose, n=228) during the first 6 months of life in the
- 296 CIG cohort. A local polynomial regression (LOESS) fit is shown with 95% CI shaded in grey. Statistical significance is
- $\label{eq:297} evaluated by repeated measures correlations (r_{rm}).$
- e, Heatmap illustrating partial Spearman's Rank correlation coefficients (adjusted for age) between the absolute
 abundance of *Bifidobacterium* species/subspecies and absolute faecal concentrations of aromatic lactic acids (n=240) or
 relative abundances of HMOs (n=228) in the CIG cohort. Infant type *Bifidobacterium* spp. is the sum of absolute
 abundances of *B. longum*, *B. breve*, *B. bifidum* and *B. scardovii*. Statistical significance is indicated by asterisks with *
 p<0.05, **p<0.01, ***p<0.001 and ****p<0.0001.
- 303 f-h, Absolute abundance of *Bifidobacterium* spp. (>1% of total community) and concentrations of ILA, PLA and 4-OH-
- 304 PLA in selected individuals from the CIG cohort. **f**, Fully breastfed infants and early colonised with high abundances of
- 305 one or more of the infant *Bifidobacterium* spp. (*B. longum*, *B. breve* or *B. bifidum*) and concurrent high absolute
 306 concentrations of ILA, PLA and 4-OH-PLA through the first 6 months of life. g, Infants with delayed colonization of
- 307 infant type *Bifidobacterium* spp. and concurrent low concentrations of ILA, PLA and 4-OH-PLA, **h**. Infants with
- 308 recorded oral antibiotics intake during the first 6 months of life. Similar dynamics of the remaining infants can be seen
- 309 in Supplementary Fig. 15.
- 310 See Supplementary Fig 11-15

311 Indolelactic acid is a relevant early life AhR agonist

The tryptophan-derived metabolite ILA was the most abundant aromatic lactic acid, as well as the 312 most abundant tryptophan catabolite measured in the faeces of infants at 0-6 months 313 (Supplementary Table 2) and 9 months of age (Fig. 1d). Microbial tryptophan catabolites have 314 been found to contribute to intestinal and systemic homeostasis, in particular by their ability to bind 315 the aryl hydrocarbon receptor $(AhR)^{12}$. In accordance with previous reports^{16,35}, we observed 316 modest but significant dose-dependent increases in agonistic activity of ILA in both rat and human 317 AhR reporter gene cell lines (Supplementary Fig. 16). Notably, ILA was more potent in the human 318 AhR compared to the rat AhR assay (Supplementary Fig. 16). To investigate the relationship 319 between gut microbiota, aromatic amino acid metabolites and AhR signalling, the AhR activity 320 induced by sterile-filtered faecal water from selected CIG infants (Fig. 3f-h) was associated to the 321 most abundant bacterial taxa and all quantified aromatic amino acid metabolites (n=20) in the same 322 samples (Fig. 4a). This revealed that particularly the infant type *Bifidobacterium* spp. were 323 positively associated with AhR activity (Fig. 4b & Supplementary Fig. 17a). B. dentium also 324 correlated positively with AhR activity despite the absence of an *aldh* gene in this species (**Fig. 4a**), 325 which might be due to its association to the AhR agonist IAld (Fig. 1e), showing similar AhR 326 reporter activity as ILA (Supplementary Fig. 16a). Of all the aromatic amino acid metabolites 327 measured, only faecal concentrations of three known AhR agonists indoleacetic acid (IAA), IAld 328 329 and in particular ILA were significantly associated with AhR activity (Fig. 4a,c & Supplementary Fig. 17b). Temporal development in AhR activity was assessed in the CIG samples, and 330 stratification of infants into those colonised early (Fig. 3f) and late (Fig. 3g) with infant type 331 *Bifidobacterium* spp. revealed a higher AhR activity of the faecal water of early colonised infants 332 (Fig. 4d-e). These data show that ILA, produced by Bifidobacterium spp. of the infant gut, is a 333

highly relevant AhR agonist in early life.

335





338 Fig. 4. Faecal concentrations of infant type *Bifidobacterium* spp. and indolelactic acid associate with AhR activity

a, Scatterplot of age adjusted Spearman's Rank correlation coefficients (versus associated FDR-adjusted p-values)
 between AhR activity (Luminescence Units, LU) of faecal water (n=119) from selected Copenhagen Infant Gut (CIG)
 infants (Fig. 3f-h, n=11) in a reporter cell line assay and absolute abundance of gut bacterial taxa (relative
 abundance>0.1%, n=40, circles) or quantities of aromatic amino acid catabolites (n=20, triangles) measured in the same

samples. Coloured circles/triangles mark taxa/metabolites measures that are significantly positively (red) or negatively
(blue) associated with AhR activity, within an FDR-adjusted p-value of 0.1. Labels are coloured brown for metabolites
and black for bacterial taxa. Asterisks indicate known AhR agonists¹⁶.

b-c, Scatterplots showing the associations from (a) between infant type *Bifidobacterium* spp. (b) or ILA (c) and AhR
 activity in faecal water from selected CIG infants, assessed by age adjusted Spearman's Rank correlation analysis.

d, Temporal development (11 samplings points) of *in vitro* AhR activity in CIG infants early (green; CIG04, CIG06,

CIG11 and CIG14) or late (orange; CIG07, CIG08, CIG09, CIG18) colonised with infant type *Bifidobacterium* spp (See
 Fig. 3f-g). Locally weighted regression scatterplot smoothing (LOWESS) curves were fitted to the data points.

e, Scatterplots (mean \pm SD) of AhR activity for each sampling point, stratifying samples into individuals early and late colonised with infant type *Bifidobacterium*. Statistical significance was evaluated by unpaired *t*-test.

353 See Supplementary Fig 16-17

354

355 DISCUSSION

The importance of intestinal commensal bacteria in regulation of the intestinal barrier function and 356 immune development during infancy is well established^{36,37}. Yet, specifically the symbiotic role of 357 the breastmilk-promoted *Bifidobacterium* species, which are highly abundant in breastfed infants, 358 remains largely unknown. Here we identified an aromatic lactate dehydrogenase, which catalyses 359 the last step of the conversion of aromatic amino acids into their respective aromatic lactic acids in 360 the infant gut. We show that only the infant type Bifidobacterium species contain the aromatic 361 lactate dehydrogenase, explaining why Bifidobacterium species commonly isolated from the infant 362 gut *in vitro* produce relative higher levels of ILA compared with other *Bifidobacterium* species³⁸. 363 The infant type *Bifidobacterium* species are adapted to breastfeeding by their HMO-transport and 364 degradation genes, which provide them with a colonization advantage in infant gut^{39–43}. Unlike 365 HMO-utilization genes, the *aldh* gene appeared not to be important for bifidobacterial proliferation 366 since disruption of the *aldh* gene did not compromise growth. However, acquisition of *aldh* genes in 367 breastmilk-promoted *Bifidobacterium* species may provide an evolutionary advantage to the host. 368 Here, our data suggest that the production of the AhR agonist ILA by breastmilk-promoted 369 *Bifidobacterium* is a key determinant of AhR-dependent signalling in the gut during infancy. 370 Indeed, our enzymatic assays show strong adaptation of ALDH towards indolepyruvate, resulting 371 372 preferentially in the formation of ILA, the most abundant of the aromatic lactic acids and microbial tryptophan catabolites detected in our two cohorts. This is potentially of fundamental importance, 373 since AhR signalling is essential for gut barrier function^{44,45} and immune development^{15,46–48}. 374 Microbiota-produced tryptophan-derived AhR ligands (e.g. indolealdehyde) have been shown to 375 confer colonization resistance against gastrointestinal pathogens¹⁶, improve intestinal barrier 376 function⁴⁴, attenuate induced colitis⁴⁹ and improve features of metabolic syndrome⁴⁴. Further, 377 although AhR is conserved across invertebrates and vertebrates, microbiota-derived tryptophan 378 agonists have shown human-AhR selectivity⁵⁰. In support of this, we found that ILA is a more 379 potent ligand in the human compared to the rat AhR assay. Considering the strong dominance of 380 ILA producing Bifidobacterium spp. in the gut of breastfed infants, we speculate that ILA provides 381 an advantage to the infant host by contributing to AhR-mediated early life intestinal and metabolic 382 homeostasis. A putative mechanism relates to the development of innate lymphoid cells (ILCs), 383 which are important mediators of innate immunity to bacterial, viral and fungal infections⁵¹. A 384 specific subset of ILCs, group 3 ILCs, shows AhR agonist dependency for survival and function 385 and are reported to be crucial for resistance against gastrointestinal pathogens through the 386 production of IL-22^{16,48}. This effect has been shown to be mediated by microbiota-produced AhR-387 agonistic indoles¹⁶. Thus, we find it plausible that high levels of ILA produced by *Bifidobacterium* 388 389 spp. in the infant gut contribute to innate protection against gastrointestinal pathogens. Further, AhR ligands (including ILA) have been shown to modulate the adaptive immune system by promoting 390 IL-10+ T_{reg} induction and suppressing IL-17 and IFN γ production by T-cells, potentially reducing 391 the risk of autoimmunity⁵²⁻⁵⁴. Non-AhR mediated mechanisms may add to the protective effect, as 392 ILA and PLA have been shown *in vitro* to have direct anti-bacterial^{55,56} and anti-fungal 393 properties^{57,58}. Furthermore, ILA and PLA were recently identified as potent agonists of human 394 hydroxycarboxylic acid receptor 3 (HCA3)⁵⁹, a receptor expressed in adipocytes, immune cells and 395 in the intestinal epithelium⁶⁰, involved in the regulation of immune functions and energy 396

homeostasis 61,62 . Thus, the production of aromatic lactic acids by human breastmilk-promoted

398 *Bifidobacterium* during infancy may represent an evolutionary explanation as to why HCA3 is only

expressed in humans and hominids 59 . In summary, the production of aromatic lactic acids via the

400 newly identified aromatic lactate dehydrogenase found specifically in breastmilk-promoted species

401 of *Bifidobacterium* might be an unprecedented feature contributing to intestinal homeostasis and

402 immune development during infancy. These findings represent a major progress in the

- 403 understanding of the beneficial effects of breastmilk, *Bifidobacterium* species and their metabolites
- 404 for human health in early life.

405 METHODS

406 Human study populations and metadata

407 SKOT cohort

The discovery cohort consisted of a random subset of 59 healthy infants of the observational SKOT 408 I cohort¹⁹. These infants were originally recruited from Copenhagen and Frederiksberg by random 409 selection from the National Danish Civil Registry⁶³. Inclusion criteria were single birth and full 410 term delivery, absence of chronic illness and age of 9 months \pm 2 weeks at inclusion. Mode of 411 delivery, gender, age at sampling, use of medication, breast- and formula feeding prevalence as well 412 as exclusive and total breastfeeding duration and age of introduction to solid foods was recorded by 413 parental questionnaires (Supplementary Data 1a,b). Anthropometrics, full dietary assessment and 414 other relevant metadata have been published previously^{3,64}. Faecal samples were obtained at 9 415 months ± 2 weeks of age and were stored at -80°C until DNA extraction, as described previously³. 416 Urine samples were collected by the use of cotton balls placed in the infants' disposable nappies 417 418 from which the urine was squeezed into a sterile tube and stored at -80°C. In case of faeces in the nappy, the urine sample was discarded. The study protocol was approved by the Committees on 419 Biomedical Research Ethics for the Capital Region of Denmark (H-KF-2007-0003) and The Data 420 Protection Agency (2002-54-0938, 2007-54-026) approved the study. Informed consent was 421

422 obtained from all parents of infants participating in the SKOT I study.

423 CIG cohort

424 The validation cohort, CIG, consisted of 25 healthy infants, vaginally born (23/25) and full-term

425 (23/25) delivered. Infants in CIG were recruited through social media and limited to the

426 Copenhagen region. Parents collected faecal samples approximately every second week, starting

- from the first week of life until 6 months of age (i.e. within week 0, 2, 4, 6, 8, 10, 12, 16, 20 and
- 428 24). Parents were instructed to collect faecal samples from nappies into sterile faeces collection

tubes (Sarstedt, Nümbrecht, Germany) and immediately store them at -18°C in a home freezer until

- transportation to the Technical University of Denmark where the samples were stored at -80°C until
- 431 sample preparation. Gender, pre-term vs full-term birth, mode of delivery, infant/maternal
- antibiotics, feeding patterns (breastmilk vs formula), introduction to solid foods and consumption of
- 433 probiotics was recorded (**Supplementary Data 2a**). The Data Protection Agency (18/02459)
- 434 approved the study. The office of the Committees on Biomedical Research Ethics for the Capital
- 435 Region of Denmark confirmed that the CIG study was not notifiable according to the Act on

Research Ethics Review of Health Research Projects (Journal nr.: 16049041), as the study only
concerned the faecal microbial composition and activity and not the health of the children. Informed
consent was obtained from all parents of infants participating in the CIG study.

439 Gut microbiota analysis

440 16S rRNA gene amplicon sequencing

Sample preparation and sequencing was performed as previously described³ using a subset of 59 441 faecal samples originating from infants participating in the SKOT I cohort and 241 faecal samples 442 from 25 infants participating in the CIG cohort (data from a total of 28 samples were missing due to 443 444 insufficient DNA extraction, lack of PCR product, very low number of sequencing reads or resemblance of community to negative controls). Briefly, DNA was extracted from 250 mg faeces 445 (PowerLyzer® PowerSoil® DNA isolation kit, MoBio 12855-100) and the V3 region of the 16S 446 rRNA gene was amplified (30s at 98°C, 24-30 cycles of 15s at 98°C and 30s at 72°C, followed by 5 447 min at 72°C) using non-degenerate universal barcoded primers⁶⁵ and then sequenced with the Ion 448 OneTouchTM and Ion PGM platform with a 318-Chip v2. Sequences from SKOT and CIG were 449 analyzed separately. Briefly, they were de-multiplexed according to barcode and trimmed as 450 previously described^{65,66} in CLC Genomic Workbench (v8.5. CLCbio, Qiagen, Aarhus, DK). 451 Ouality filtering (-fastq filter, MAX $EE_{ISKOTI} = 2.0$, MAX $EE_{ICIGI} = 1.0$), dereplication, OTU 452 clustering (-cluster_otus, minsize 4), chimera filtering (-uchime_ref, RDP_gold database), mapping 453 of reads to OTUs (-usearch_global, id 97%) and generation of OTU tables (python, uc2otutab.py) 454 was done according to the UPARSE pipeline⁶⁷. In QIIME⁶⁸, OTU tables ($n_{OTUs[SKOT]} = 545$, 455 $n_{OTUs[CIG]} = 478$) was filtered to include only OTUs with abundance across all samples above 456 0.005% of the total OTU counts ($n_{OTUs[SKOT]}$ = 258, $n_{OTUs[CIG]}$ = 145). OTU relative abundances 457 within samples were estimated by total sum scaling. Taxonomy was assigned to the OTUs using the 458 rdp classifier with confidence threshold 0.5^{69} and the GreenGenes database v13.8⁷⁰. Estimating 459 species composition in the CIG cohort, the OTUs detected with identical taxonomy were collapsed 460 and using a cutoff of average relative abundance of 0.1%, only 39 bacterial species/taxa remained. 461 representing 97.5% of total community (Supplementary Data 2b & Supplementary Fig. 11). 462 Based on PvNAST alignment of representative OTU sequences from each cohort separately, a 463 phylogenetic tree was created with FastTree, as described previously⁶⁶. Alpha diversity (Shannon 464 index, Observed OTUs, Pielou's evenness index) and beta diversity (weighted and unweighted 465 UniFrac distances, abundance weighted and binary Bray-Curtis and abundance weighted Jaccard 466 dissimilarities) measures were calculated in QIIME, with the sequencing depth rarefied to 2,000 467 (SKOT) - 8,000 (CIG) sequences per sample. Jaccard similarity index (1- abundance weighted 468 Jaccard distance) was computed by calculating the median of all Jaccard similarity index values 469 between adjacent time points within each individual of the CIG cohort. In order to investigate 470 Bifidobacterium species composition OTUs sequences classified as Bifidobacterium according to 471 the GreenGenes database v13.8 were filtered to remove low abundant OTUs (cutoff 0.1% of total 472 *Bifidobacterium*) and the taxonomy of these resulting OTUs ($n_{OTUs[SKOT]} = 23$, $n_{OTUs[CIG]} = 8$) was 473 confirmed by BLAST⁷¹ search against the 16S rRNA gene sequence database at NCBI. The top 474 BLAST hit indicated species annotation (Supplementary Data 1f and 2c). OTUs were collapsed 475

476 into *Bifidobacterium* species (*B. longum*, *B. bifidum*, *B. breve*, *B. catenulatum* group, *B.*

477 adolescentis, B. scardovii, B. dentium, B. animalis/pseudolongum) based on the top BLAST hit

478 (Supplementary Data 1f and 2c).

479 Quantitative PCR

Total bacterial load (universal primers) and absolute abundances of *B. longum* subsp. *longum* and *B.* 480 longum subsp. infantis (subspecies specific primers) were estimated by quantitative PCR (qPCR), 481 using the primers listed in **Supplementary Table 3**. Each reaction was performed (in triplicates) 482 with 5 µl PCR-grade water, 1.5 µl forward and reverse primer, 10 µl SYBR Green I Master 2X 483 (LightCycler[®] 480 SYBR Green I Master, Roche) and 2 µl template DNA, in a total volume of 20 484 ul. Standard curves were generated from 10-fold serial dilutions of linearized plasmid (containing 485 10^8 - 10^0 gene copies/µl), constructed by cloning a PCR amplified 199bp fragment of the 16S rRNA 486 gene (V3-region) of *E. coli* (ATCC 25922) or a 307bp fragment of the Blon0915 gene⁷² of *B.* 487 longum subsp. infantis (DSM 20088) or a 301bp fragment of the BL0274 gene⁷³ of B. longum 488 subsp. longum (DSM 20219) into a pCR4-Blunt-TOPO (Invitrogen) or pCRII-Blunt-TOPO vector 489 (Invitrogen). Plates were run on the LightCycler[®] 480 Instrument II (Roche) with the program 490 including 5 min pre-incubation at 95°C, followed by 45 cycles with 10 sec at 95°C, 15 sec at 50-491 492 60°C and 15 sec at 72°C and a subsequent melting curve analysis including 5 min at 95°C, 1 min at 65°C and continuous temperature increase (ramp rate 0.11 °C/s) until 98°C. Data were analyzed 493 with the LightCycler[®] 480 Software (v1.5) (Roche). Bacterial load data (using the universal 494 primers) were used to estimate absolute abundances of each microbial taxa by multiplying with 495 relative abundances (derived from 16S rRNA gene amplicon sequencing). 496

497 Bifidobacterium strains and growth experiments

498 Aromatic lactic acid production by Bifidobacterium type strains

Bifidobacterium type strains (Supplementary Table 4) were cultivated on MRSc (MRS containing 499 2% (w/v) glucose and supplemented with 0.05% (w/v) L-cysteine) agar plates for 48h at 37°C 500 anaerobically. Single colonies were dissolved in 5.0 mL pre-reduced MRSc broth and incubated for 501 502 24h at 37°C anaerobically with shake. The overnight cultures were washed (10000xg, room temperature, 5 min) and resuspended in sterile 0.9% NaCl water, diluted 1:20 (in triplicates) in pre-503 504 reduced MRSc or MRSc+HMOs (MRS broth without glucose, but supplemented with 2.0% (w/v) HMO mixture and 0.05% (w/v) L-cysteine) and re-incubated at 37°C anaerobically for 72h, after 505 which OD_{600nm} was measured and the culture supernatants (16000xg, 5 min, 4°C) were analysed by 506 UPLC-MS. The individual HMOs were kindly donated by Glycom A/S (Hørsholm, Denmark); 2'-507 O-fucosyllactose (2'FL), 3-O-fucosyllactose (3FL), lacto-N-tetraose (LNT), lacto-N-neotetraose 508 (LNnT), 6'-O-sialyllactose (6'SL), 3'-O-sialyllactose (3'SL), together representing the three 509 structures found in human breastmilk (fucosylated, sialylated, and neutral core). Based on the HMO 510 composition in breastmilk^{74,75}, these were mixed in a ratio of 53% 2'FL, 18% 3FL, 13% LNT, 5% 511 LNnT, 7% 6'SL and 4% 3'SL in sterile water to obtain a representative HMO mix used in the in 512 513 vitro experiments at 2% (w/v).

Growth experiment with B. longum subsp. longum 105-A strains 514

- B. longum subsp. longum 105-A (JCM 31944) was obtained from Japan Collection of 515
- Microorganisms (RIKEN BioResource Research Center, Tsukuba, Japan). B. longum subsp. 516
- longum 105-A strains (wild type [WT], insertional mutant [type4 ldh::pMSK127] and 517
- complemented insertional mutant [type 4 *ldh*::pMSK127 / pMSK128 (Pxfp-type4 *ldh*)]; 518
- Supplementary Table 4) were cultivated on MRSc or MRSc-Chl (MRSc supplemented with 2.5 519
- µg/mL chloramphenicol) agar plates for 48h at 37°C anaerobically. Single colonies were dissolved 520
- 521 in 5.4 mL MRSc or MRSc-Chl broth, 10-fold serially diluted and incubated for 15h at 37°C
- 522 anaerobically with shake. The most diluted culture (exponential phase) was washed in same
- medium (10000xg, room temperature, 5 min) and resuspended in MRSc or MRSc-Chl broth to 523
- yield $OD_{600nm} = 1$ and subsequently diluted 1:40 in prewarmed and reduced MRSc or MRSc-Chl 524
- broth (in triplicates), before incubation at 37°C, anaerobically with shake. The cultures were 525
- sampled (500 μ L) every hour for OD_{600nm} measurements and the culture supernatants (16000xg, 526 4°C, 5 min) from early (13h) stationary phase was analyzed by UPLC-MS for aromatic amino
- 527
- metabolites and by GC-MS for lactate. 528

Alignments and construction of phylogenetic trees 529

- 530 From the full genome sequences (available at https://www.ncbi.nlm.nih.gov/genome/) of
- Bifidobacterium type strains included in this study (Supplementary Table 4) all genes annotated as 531
- L-lactate dehydrogenases were aligned (gap cost 10, gap extension cost 1) and subsequently a 532
- phylogenetic tree (Algorithm = Neighbor-Joining, Distance measure = Jukes-Cantor, 100 bootstrap 533
- replications) was constructed in CLC Main Workbench (v7.6.3, CLCbio, Qiagen, Aarhus, DK). The 534
- tree was visualized by use of the FigTree software v1.4.3 (http://tree.bio.ed.ac.uk/software/figtree/). 535
- In addition, the type 4 LDH amino acid sequence (translated from the type 4 ldh nucleotide 536
- sequence) of *B. longum* subsp. *longum* 105-A was aligned (gap cost 10, gap extension cost 1) with 537
- the type 4 LDH amino acid sequences of the B. longum subsp. longum, B. longum subsp. infantis, 538
- B. bifidum, B. breve and B. scardovii type strains. Comparison of type 4 ldh gene cluster/operon in 539
- 12 Bifidobacterium type strains (Supplementary Fig. 5) was basically conducted by pairwise 540
- alignments in MBGD (Microbial Genome Database for Comparative Analysis; 541
- http://mbgd.genome.ad.jp/). The amino acid sequences of the gene cluster from *B. pseudolongum* 542
- subsp. pseudolongum type strain was collected from NCBI database 543
- (https://www.ncbi.nlm.nih.gov/genome/) and was used for comparison with that from B. animalis 544
- 545 subsp. animalis type strain.

Recombinant expression of type 4 *ldh* 546

547 Chemically competent cells for recombinant expression

- E. coli LMG 194 ON culture (200µL) was inoculated into 5 ml Luria-Bertani (LB) medium and 548
- incubated at 37°C, 250 rpm until $OD_{600nm} = 0.5$, at which the culture was centrifuged 5 min at 549
- 10000xg at 4°C and supernatant discarded. Cell pellet was resuspended in ice-cold 1.8 ml 10mM 550
- MgSO₄ (Sigma, M2643) and centrifuged for 2 min at 5000xg at 0°C. Supernatant was discarded 551

and cell pellet resuspended in 1.8 ml ice-cold 50 mM CaCL₂ (Merck, 1.02083.0250), incubated on
ice for 20 min and centrifuged for 2 min at 5000xg at 0°C. Cell pellet was resuspended in 0.2 ml
ice-cold 100mM CaCl₂, 10mM MgSO₄ and placed on ice until transformation.

555 Cloning and recombinant expression

Genomic DNA was extracted (PowerLyzer® PowerSoil® DNA isolation kit, MoBio 12855-100) 556 from colony material of *B. longum* subsp. *infantis* DSM 20088^T. In order to amplify the type 4 *ldh* 557 gene, 50 ng template DNA was mixed with 5 µL 10X PCR buffer, 0.5 µL (50 mM) dNTP mix, 1 µL 558 (10µM) forward primer (ldh4_F, 5'-ACCATGGTCACTATGAACCG-3'), 1 µL (10µM) reverse 559 primer (ldh4 R, 5'-AATCACAGCAGCCCCTTG-3') and 1 µL (1 U/µL) Platinum Taq DNA 560 polymerase (Invitrogen, 10966-018) in a 50 µL total reaction volume. The PCR program included 2 561 min at 94°C, 35 cycles of 30 sec at 94°C, 30 sec at 55°C, 60 sec at 72°C, followed by a final 562 extension 10 min at 72°C. The PCR product was purified (MinElute PCR purification kit, Qiagen, 563 28004) and 4 µL was mixed with 1 µL Salt solution (1.2M NaCl, 0.06M MgCl₂) and 1 µL pBAD-564 TOPO® plasmid (Invitrogen, K4300-01) and incubated for 5 minutes at room temperature. 2 µL of 565 the cloning mixture was transformed into 50 µL One Shot® TOP10 Competent Cells (Invitrogen, 566 K4300-01) by gentle mix, incubation 15 min on ice and heat-shock for 30 sec at 42°C. 250µL 567 568 S.O.C medium (Invitrogen, K4300-01) was added and incubated at 37°C for 1 h at 200 rpm and subsequently spread on LB-AMP (LB supplemented with 20 µg/ml Ampicillin (Sigma, A9518)) 569 570 agar plates and incubated at 37°C ON. Transformants were picked and clean streaked on LB-AMP agar plates, incubated 37°C, ON and afterwards single colonies of each transformant was inoculated 571 into 5 ml LB-AMP broth and incubated at 37°C for 15 h at 250 rpm. Plasmid DNA was isolated 572 (QIAprep Spin Miniprep Kit, Qiagen, 27104) from each transformant and subsequently 5 µL 573 plasmid DNA (80-100 ng/µL) was mixed with 5µL (5pmol/µL) pBAD forward (5'-574 575 ATGCCATAGCATTTTTATCC-3') or reverse (5'-GATTTAATCTGTATCAGG-3') sequencing primers (5pmol/µL) and shipped for sequencing at GATC (GATC-biotech, Koln, Germany). In 576 order to remove the leader peptide in pBAD-TOPO, 10 µL plasmid (0.1 µg) with correct insert was 577 cut with FastDigest Ncol (Thermo Scientific, FD0563) for 10 min at 37°C and the enzyme 578 579 inactivated 15 min at 65°C. Plasmid was ligated using 1 µL (1U/µL) T4 DNA Ligase (Invitrogen, 15224-017) for 5 min at room temperature and subsequently 2µl plasmid was transformed into 580 100µl chemically competent E. coli LMG194 cells by incubation on ice for 30 min, followed by 581 heat-shock at 43°C for 3 min and incubation on ice for 2 min. 900µL LB medium was added and 582 cells were incubated at 37°C for 1h at 250 rpm, before plating on LB-AMP agar plates and 583 incubation at 37°C ON. Transformants were picked, clean streaked and plasmid DNA isolated and 584 sequenced as described above. A transformant with correct insert was selected for recombinant 585 expression of the type 4 ldh gene; 2 ml LB-AMP broth was inoculated with a single recombinant 586 colony or the non-transformed E. coli LMG194 (negative control) and grown at 37°C ON at 587 250rpm. In 3x triplicates, 100 µL of the ON cultures (2x3x 100 µL transformant culture + 1x3x 100 588 µL non-transformed E. coli LMG194 culture) were diluted 100-fold into 9.9 mL prewarmed LB-589 590 AMP/LB broth and grown at 37°C, 250 rpm until $OD_{600nm} \approx 0.5$, at which 9 mL culture was added 1 mL mix of indolepyruvic acid, phenylpyruvic acid and 4-hydroxyphenylpyruvic acid (1mg/mL 591 each). The cultures were sampled (time zero) and subsequently 100 µL 20% L-arabinose (or 100 uL 592

- sterile water; control for induction) was added to induce gene expression and the cultures were re-
- incubated at 37°C, 250 rpm, before sampling at 1 h and 5 h post-induction for OD_{600nm}
- 595 measurements and assessment of production of aromatic lactic acids. For the latter, samples were
- centrifuged at 16000xg, 5 min, 4° C and supernatants were stored at -20° C for UPLC-MS analyses.

597 **Construction of type 4** *ldh* **insertional mutant**

598 Transformation of B. longum subsp. longum 105-A

B. longum subsp. *longum* 105-A cells were grown to exponential phase at 37°C in Gifu anaerobic liquid medium (Nissui Pharmaceutical Co., Ltd., Tokyo, catalog no. 05422), harvested by centrifugation, and washed twice with ice-cold 1 mM ammonium citrate buffer containing 50 mM sucrose (pH = 6.0). The cells were concentrated 200 times with the same buffer and used for electroporation with settings of 10 kV/cm, 25 μ F, and 200 Ω . After recovery culturing in Gifu anaerobic liquid medium at 37°C for 3 h, the cells were spread onto Gifu anaerobic agar containing antibiotics (i.e. 30 μ g/mL spectinomycin and/or 2.5 μ g/mL chloramphenicol) for selection.

606 Insertional mutant construction and plasmid complementation

The type 4 ldh gene (BL105A 0985) of B. longum subsp. longum 105-A was disrupted by a 607 plasmid-mediated single crossover event as described previously⁷⁶. The plasmid used for disruption 608 was constructed using the In-Fusion cloning kit (Clontech Laboratories, Inc., Mountain View, CA, 609 USA, catalog no. 639649). Escherichia coli DH5a was used as a host. In brief, the internal region of 610 the *ldh* gene (position 142-638 of the nucleotide sequence of BL105A_0985, see Supplementary 611 Fig. 6) was amplified by PCR using a primer pair Pr-580/581 (Supplementary Table 5) and 612 ligated with the BamHI-digested pBS423 fragment carrying pUC ori and a spectinomycin 613 resistance gene²⁶. The resulting plasmid pMSK127 was introduced into *B. longum* subsp. *longum* 614 105-A by electroporation to be integrated into type 4 *ldh* locus by single crossover recombination 615 (type 4 *ldh*::pMSK127). Type 4 *ldh* disruption was confirmed by PCR with a primer pair designed 616 to anneal outside of the gene (Supplementary Fig. 6 and Supplementary Table 5). The amplified 617 fragment was also sequenced to ensure the correct recombination event. Complementation plasmid 618 pMSK128 was constructed by ligating PCR-amplified xfp (xylulose 5-phosphate/fructose 6-619 phosphate phosphoketolase) promoter region (Pxfp) and the type 4 ldh coding region with PstI- and 620 SalI-digested pBFS38⁷⁷ using the In-Fusion cloning kit, by which type 4 *ldh* was placed under the 621 control of Pxfp. Primer pairs of Pr-598/Pr-599 and Pr-600/Pr-601 were used for amplifying Pxfp 622 from pBFS48⁷⁷ and the type 4 *ldh* gene from the *B. longum* subsp. *longum* 105-A genome, 623 respectively (Supplementary Table 5). The resulting plasmid was electroporated into type 4 624 *ldh*::pMSK127 to give type 4 *ldh*::pMSK127 / pMSK128 (Pxfp-type4_*ldh*) (Supplementary Fig. 625 6). 626

627

628 Biochemical characterization of type 4 LDH

629 Recombinant expression and purification

630 Type 4 LDH (BL105A_0985) was recombinantly expressed as a non-tagged form. The gene was amplified by PCR using the genomic DNA of *B. longum* subsp. *longum* 105-A as a template and a 631 632 635 (5'-AAGGAGATATACATATGGTCACTATGAACCGC-3'). Underlined bases indicate 15-633 bp for In-Fusion cloning (Clontech). The amplified DNA fragment was inserted into the NdeI and 634 XhoI site of pET23b(+) (Novagen) using an In-Fusion HD cloning kit (Clontech). The resulting 635 plasmid was introduced into E. coli BL21 (DE3) $\Delta lacZ$ carrying pRARE2⁷⁶, and the transformant 636 was cultured in LB medium supplemented with ampicillin (100 μ g ml⁻¹) and chloramphenicol (7.5 637 μ g ml⁻¹). When OD_{600nm} reached 0.5, isopropyl β -D-thiogalactopyranoside was added at a final 638 639 concentration of 0.02 mM to induce the protein expression. The culture was incubated for four days at 18°C, harvested by centrifugation, and resuspended in 50 mM potassium phosphate buffer (KPB; 640 pH 7.0) supplemented with 1 mM 2-mercaptoethanol (2-ME) and 200 µM phenylmethane sulfonyl 641 fluoride. Following cell disruption by sonication, the cleared lysate was saturated with ammonium 642 sulfate (40–60%). The resulting precipitate was dissolved, dialyzed against 20 mM KPB (pH 7.0) 643 containing 1 mM 2-ME, and concentrated by Amicon Ultra 10K centrifugal device (Merck 644 Millipore). The sample was then loaded onto an Affigel blue column (Bio-Rad) preequilibrated with 645 20 mM KPB (pH 7.0) containing 1 mM 2-ME, and eluted by the same buffer containing 1 M NaCl. 646 The protein was further purified by a Mono Q 5/50 (GE Healthcare; a linear gradient of 0-1 M 647 648 NaCl in 20 mM Tris-HCl (pH 8.0) containing 1 mM 2-ME) and Superdex 200 Increase 10/300 GL column (GE Healthcare; 10 mM KPB [pH 7.0] containing 50 mM NaCl and 1 mM 2-ME). Protein 649 concentration was determined by measuring the absorbance at 280 nm based on a theoretical 650 extinction coefficient of 26,470 M⁻¹ cm⁻¹. 651

652 Enzyme assay

The standard reaction mixture contained 100 mM KPB (pH 8.0), 1 mM 2-ME, 0.1 mM β-NADH, 653 and the substrate. The reaction was initiated by adding the enzyme, and the mixture was incubated 654 at 37°C for an appropriate time, in which the linearity of the reaction rate was observed. The 655 substrate concentrations were varied between 0.01 and 0.25 mM for IPA, 1.5 and 12.75 mM for 656 PPA, 2 and 24 mM for 4-OH-PPA, and 2.5 and 40 mM for pyruvic acid. The enzyme was used at 657 the concentrations of 0.22 nM for IPA, 1.47 nM for PPA, 0.12 nM for 4-OH-PPA, and 88.50 nM 658 for pyruvic acid. The reducing reactions of PPA and pyruvic acid was continuously monitored by 659 measuring the decrease of the absorbance at 340 nm (NADH consumption). When 4-OH-PPA and 660 IPA were used as the substrates, the reaction products 4-OH-PLA and ILA were quantified by 661 HPLC after the termination of the reactions by adding 5 % (w/v) trichloroacetic acid. HPLC 662 analysis was performed using a Waters e2695 separation module (Waters) equipped with a 663 LiChrospher 100 RP-18 column (250×4 mm, $\varphi = 5 \mu m$; Merck Millipore) at 50°C. Following 664 equilibration with a mixture of 10% solvent A (50% methanol, 0.05% trifluoroacetic acid) and 90% 665 666 solvent B (0.05% trifluoroacetic acid) at a flow rate of 1 mL/min, the concentration of solvent A was linearly increased to 100 % for 25 min and maintained at 100 % for additional 15 min. 4-OH-667

668 PLA and ILA were detected by a Waters 2475 Fluorescence Detector with λ_{ex} 277 nm and λ_{em} 301 nm and λ_{ex} 282 nm and λ_{em} 349 nm, respectively. The standard curves were created using the 669 known concentrations of both compounds. The kinetic parameters (k_{cat} , $K_{0.5}$, and Hill coefficient $n_{\rm H}$) 670 671 were calculated by curve-fitting the experimental data to the Hill equation, using KaleidaGraph version 4.1 (Synergy Software). Experiments were performed at least in duplicate. Physicochemical 672 property of the enzyme was examined by using 1 mM PPA as a substrate. The effects of metal ions 673 (0.1 mM each) on the enzyme activity was examined using 50 mM MES (2-(N-674 675 morpholino)ethanesulfonic acid) buffer (pH 7.0). EDTA (ethylenediaminetetraacetic acid) was added at the final concentration of 0.1, 0.5, or 1 mM. The optimal pH was determined using 50 mM 676 KPB (pH 6.0-8.5) and TAPS (N-Tris(hydroxymethyl)methyl-3-aminopropanesulfonic acid) buffer 677 (pH 8.0–9.0). The thermostability was evaluated by the residual activities after incubating the 678 enzyme (1.0 mg/ml in 10 mM KPB [pH 7.0] containing 50 mM NaCl and 1 mM 2-ME) at the 679 indicated temperatures for 30 min prior to the assay. Fructose-1,6-bisphosphate, shikimate-3-680 phosphate, D-erythrose-4-phosphate, and phosphoenolpyruvate were added to the reaction mixtures 681 at the concentrations of 0.1 and 1 mM to examine their heterotropic effects. KPB, TAPS buffer, or 682 683 HEPES (4-(2-hydroxyethyl)-1-piperazineethanesulfonic acid) buffer (pH 8.0 each) containing 1 and 4 mM PPA as a substrate was used. The effect of phosphate ion was analyzed by adding various 684 concentration of KPB (pH 8.0) into 10 mM HEPES buffer (pH 8.0). All experiments were 685 conducted at least in duplicate. In the subsequent kinetic analysis, we used phosphate ion at the 686 concentration of 100 mM because (i) no saturation was obtained for phosphate under the tested 687 688 conditions (Supplementary Fig. 10a), (ii) the intracellular phosphate concentration in Grampositive bacteria is known to be 130 mM at maximum⁷⁸, and (iii) the strong homotrophic effect of 689 the substrate PPA was observed only in the presence of 10 mM phosphate ion. 690

691 Metabolomics

692 Chemicals

693 Authentic standards of the aromatic amino acids and derivatives (**Supplementary Table 1**) were 694 obtained from Sigma Aldrich (Germany), whereas isotope-labelled aromatic amino acids used as

- 695 internal standards (L-Phenylalanine (ring-d5, 98%), L-Tyrosine (ring-d4, 98%), L-Tryptophan
- (indole-d5, 98%) and indoleacetic acid (2,2-d2, 96%)) of the highest purity grade available were
- 697 obtained from Cambridge Isotope Laboratories Inc. (Andover, MA).

698 Extraction of metabolites from faecal samples

Faecal samples (100-500 mg) were diluted 1:2 with sterile MQ water, vortexed for 10 seconds and 699 centrifuged at 16.000xg, 4°C for 5 minutes. Subsequently, the supernatant liquor was transferred to 700 a new tube and centrifuged again at 16.000xg, 4°C for 10 minutes. Finally, an aliquot of 150-300 701 µL was stored at -20°C. All samples were later thawed at 4°C, centrifuged at 16.000xg, 4°C for 5 702 minutes, and diluted in a total volume of 80 µL water corresponding to a 1:5 dilution of the faecal 703 sample. To each sample, 20 µL internal standard mix (4 µg/mL) and 240 µL of acetonitrile were 704 added. The tubes were vortexed for 10 seconds and left at -20°C for 10 minutes to precipitate the 705 proteins. The tubes were then centrifuged at 16.000xg, 4°C for 10 minutes and each supernatant 706

($320 \ \mu$ L) was transferred to a new tube, which was dried with nitrogen gas. Subsequently, the residues were reconstituted in 80 μ L water (equalling a 1:5 dilution of the faecal sample with internal standards having a concentration of 1 μ g/mL), vortexed for 10 seconds, centrifuged at 16.000xg, 4°C for 5 minutes, and transferred to LC vial, which was stored at -20°C until analysis.

- 711 Extraction of metabolites from urine samples
- 712 Urine samples (n=49) from the SKOT cohort were thawed in a refrigerator and all procedures
- during the sample preparation were carried out at 0-4°C using an ice bath. The subjects were
- randomized between analytical batches by placing all the samples from the each subject in the same
- 715 96 well-plate. The run order of the samples was randomized within the analytical batch. Urine
- samples were centrifuged at 3000xg for 2 minutes at 4°C. 150 μ L of each urine sample were added
- 717 to separate wells and diluted with 150 μL of diluent (MQ water: Formic acid (99.9:0.1, v/v) /
- 718 Internal standard mixture (100 μ g/mL) (90:10, v/v). A blank sample (diluent), standard mixture of
- external standard containing 44 biologically relevant metabolites (metabolomics standard)⁷⁹ and
- pooled sample containing equal amount of each sample $(20 \,\mu L)$ were added to spare wells as
- quality control samples. The plates were stored at -80°C until the analysis. Immediately prior to
- analysis, the plates were thawed and mixed by vortex stirring for 10 minutes.

723 Extraction of metabolites from in vitro fermentation samples

- Supernatants from *in vitro* fermentations were thawed at 4°C, centrifuged at 16.000xg, 4°C for 10
- minutes, before 80 μ L was transferred to a new tube. To each sample, 20 μ L internal standard (40
- $\mu g/mL$) and 300 μL of acetonitrile were added. The tubes were vortexed for 10 seconds and left at -
- 20°C for 10 minutes to precipitate the proteins. Following, the tubes were centrifuged at 16.000xg,
- 4°C for 10 minutes before 50 μ L of each sample was diluted with 50 μ L of sterile water and
- transferred to a LC vial (equalling a 1:10 dilution of the sample with internal standards having a
- 730 concentration of $1 \mu g/mL$).

731 Metabolic profiling of faecal and in vitro samples using UPLC-MS

- Aromatic amino acids and derivatives (**Supplementary Table 1**) of faecal and *in vitro* samples were quantified by ultra performance liquid chromatography mass spectrometry (UPLC-MS) as
- 734 previously published⁸⁰.
- In brief, samples were analysed in random order. For the analysis of the CIG faecal samples, a
 pooled quality control (OC) sample was injected for every 10 sample. In all cases, five standard mix
- solutions (0.1 μ g/mL, 0.5 μ g/mL, 1 μ g/mL, 2 μ g/mL and 4 μ g/mL) were analysed once for every 10
- solutions (0.1 μ g/mL, 0.5 μ g/mL, 1 μ g/mL, 2 μ g/mL and 1 μ g/mL) were unarysed once for every 1 samples to obtain a standard curve for every 10 samples. For each sample, a volume of 2 μ L was
- 739 injected into a ultra-performance liquid chromatography quadrupole time-of-flight mass
- 740 spectrometry (UPLC-QTOF-MS) system consisting of Dionex Ultimate 3000 RS liquid
- 741 chromatograph (Thermo Scientific, CA, USA) coupled to a Bruker maXis time of flight mass
- spectrometer equipped with an electrospray interphase (Bruker Daltonics, Bremen, Germany)
- operating in positive mode. The analytes were separated on a Poroshell 120 SB-C18 column with a
- dimension of 2.1x100 mm and 2.7 μ m particle size (Agilent Technologies, CA, USA) as previously

published⁸⁰. Aromatic amino acids and derivatives were detected by selected ions and quantified by

- isotopic internal standards with similar molecular structures as listed in **Supplementary Table 1**.
- 747 Data were processed using QuantAnalysis version 2.2 (Bruker Daltonics, Bremen, Germany) and
- bracket calibration curves for every 10 lumen samples were obtained for each metabolite. The
- calibration curves were established by plotting the peak area ratios of all of the analytes with respect
- to the internal standard against the concentrations of the calibration standards. The calibration
- 751 curves were fitted to a quadratic regression.
- For untargeted metabolomics, the raw UPLC-MS data, obtained by analysis of the CIG faecal
- samples in positive ionization mode, were converted to mzXML files using Bruker Compass
- 754 DataAnalysis 4.2 software (Bruker Daltonics) and pre-processed as previously reported⁸¹ using the
- 755 R packpage XCMS $(v.1.38.0)^{82}$. Noise filtering settings included that features should be detected in
- minimum 50% of the samples. A data table was generated comprising mass-to charge (m/z),
- retention time and intensity (peak area) for each feature in the every sample. The data were
- normalized to the total intensity. Subsequently, features with a coefficient of variation above 0.3 in
- the QC samples and features with a retention time below 0.5 min were excluded from the data.
- Parent ion masses of compounds of interest (2'FL/3FL, LNT/LN*n*T, 3'SL/6'SL) were searched in
- the cleaned dataset with 0.02 Da m/z and 0.02 min retention time tolerance. Subsequently, the
- identities of the features of interest were confirmed at level 1^{83} by tandem mass spectrometry and
- 763 comparison to authentic standards (**Supplementary Table 6**). Of notice, HMO isomers could not
- be distinguished with the method applied due to identical retention times.

765 Metabolic profiling of urine samples using UPLC-MS

The samples were analysed by UPLC coupled with a quadrupole-Time of Flight Mass Spectrometer 766 (q-TOF-MS) equipped with an electrospray ionization (ESI) (Waters Corporation, Manchester, UK). 767 Reverse phase HSS T3 C₁₈ column (2.1x100 mm, 1.8 µm) coupled with a pre-column (VanGuard 768 769 HSS T3 C18 column (2.1x5 mm, 1.8 µm)) were used for chromatographic separation. Five µl of each 770 well was injected into the mobile phase A (0.1 % formic acid in MQ water), mobile phase B (10% 1M ammonium acetate in methanol), mobile phase C (methanol) and mobile phase D (isopropanol). 771 Mobile phase gradient during the run time of 10 min was as follows: start condition (100 % A), 0.75 772 min (100 % A), 6 min (100 % C), 6.5 min (70 % B, 30 % D), 8 min (70 % B, 30 % D), 8.1 min (70 773 % B, 30 % D), 9 min (100 % A), 10 min (100 % A). The flow rate gradient was as follows: start 774 condition (0.4 mL/min), 0.75 min (0.4 mL/min), 6 min (0.5 mL/min), 6.5 min (0.5 mL/min), 8 min 775 (0.6 mL/min), 8.10 min (0.4 mL/min), 9 min (0.4 mL/min), 10 min (0.4 mL/min). ESI was operated 776 in negative mode with 3.0 kV capillary probe voltage. The cone voltage and the collision energy were 777 778 set at 30 kV and 5 eV, respectively. Ion source and desolvation gas (nitrogen) temperature were 120 and 400 °C while sampling cone and desolvation gas flow rates were 50 and 1000 l/hr. Scan time set 779 as 0.08 s with 0.02 sec interscan time for both modes. Data were acquired in centroid mode with mass 780 range between 50 to 1500 Da. Leucine-enkephalin (500 ng/ml) was infused as the lock-spray agent 781 to calibrate the mass accuracy every 5 sec with 1 sec scan time. Quality control samples were used to 782 evaluate possible contamination, monitoring the changes in mass accuracy, retention time and 783 instrumental sensitivity drifts^{79,84}. 784

The raw data were converted to netCDF format using DataBridge Software (Waters, Manchester, UK) and imported into MZmine version 2.28^{85} . A subset of samples was used to optimize the preprocessing parameters for the positive and negative mode data separately. Optimized pre-processing parameters are listed in **Supplementary Table 7**. Data pre-processing was employed with the following steps: mass detection, chromatogram builder, chromatogram deconvolution, deisotoping, peak alignment and gap filling. After the pre-processing, each detected peak was represented by a feature defined with a retention time, m/z and peak area.

The data matrix was imported into MATLAB R2015b (The MathWorksInc., Natick, MA). Features that were present in the blanks, were very early and late eluting (rt<0.30 and rt>9.46 min), potential isotopes, duplicates as well as features with masses indicating multiple charges were removed from the dataset using an in-house algorithm. The data were normalized using unit length normalization to correct the variation in urine concentration. Parent ion masses of the aromatic lactic acids (ILA, PLA

and 4-OH-PLA) were searched in the cleaned dataset with 0.02 Da m/z and 0.02 s retention time tolerance. A linear regression model were employed feature wise to correct for batch differences and instrumental sensitivity drifts⁸⁶. The aromatic lactic acids were confirmed at level 1⁸³ by comparison to authentic standards and by tandem mass spectrometry using the same experimental conditions (**Supplementary Fig. 2-4**).

802 Lactate production by B. longum subsp. longum 105-A strains using GC-MS

The lactate production of the *B. longum* subsp. *longum* 105-A wild-type, type 4 *ldh* mutant and type 803 804 4 *ldh* complemented strains were assessed in supernatants obtained after 13h of growth (early stationary phase) by gas-chromatography mass spectrometry (GC-MS) upon methyl chloroformate 805 (MCF) derivatisation using a slightly modified version of the protocol previously described⁸⁷. All 806 samples were analysed in a randomized order. Analysis was performed using GC (7890B, Agilent 807 808 Technologies, Inc., Santa Clara, CA) coupled with a quadrupole detector (59977B, Agilent Technologies, Inc., Santa Clara, CA). The system was controlled by ChemStation (Agilent 809 Technologies, Inc., Santa Clara, CA). Raw data was converted to netCDF format using 810 Chemstation, before the data was imported and processed in Matlab R2014b (Mathworks, Inc.) 811 using the PARADISe software⁸⁸. 812

813 Rat Aryl hydrocarbon receptor (AhR) reporter gene assay

814 Stably transfected rat hepatoma (H4IIE-CALUX) cells provided by Dr. Michael Denison

(University of California, USA) were used. The assay was conducted as previously described⁸⁹,

816 where cells were incubated for ~ 22 h in Minimum Essential Medium (MEM) α with 1% foetal

bovine serum (FBS) and 1% penicillin/streptomycin/fungizone. Chemical exposure was performed

818 for 24 h, and successively luminescence was measured. Cell viability was analyzed by measuring

819 ATP levels with the CellTiter-Glo® Luminescense Assay according to the manufacturer's

820 instruction (Promega, Denmark). 2,3,7,8-Tetrachlorodibenzo-p-dioxin (TCDD) was used as a

positive control. Three experiments in triplicates were conducted with five 2-fold dilutions of ILA

- and IAld ranging from 12.5 to 200 μ M with a constant vehicle concentration in all wells. Further,
- sterile filtered fecal water (10 mg faeces/ml MQ water) obtained from all samples (n=119) of 11

selected CIG infants (Fig. 3f-h) were run in technical triplicates in the assay. Only mild toxicity that
 did not correlate with AhR-luminesence signal was observed for some fecal water samples.

826 Human AhR reporter gene assay

827 ILA and IAld (positive control)¹⁶ were tested for activation of the human AhR. AhR Reporter Cells

from Indigo Biosciences (PA, USA) that include the luciferase reporter gene functionally linked to

an AhR-responsive promoter were used. The assay was run according to the instructions of the

830 manufacturer (technical manual version 6.0) with the reference agonist MeBIO as the positive

control. Three experiments in triplicates were conducted with five 2-fold dilutions of ILA and IAld

ranging from 12.5 to $200 \,\mu\text{M}$ with a constant vehicle concentration in all wells. No cytotoxicity was

833 observed for any of the tests as determined by a resazurin toxicity assay.

834 Statistical analyses

Statistical analyses were performed using QIIME v1.9⁶⁸, R v3.1⁹⁰ and GraphPad Prism v8.1

836 (GraphPad Software, Inc. CA). PCoA, ADONIS and PERMDISP tests (permutations = 999) of

837 OTU distance/dissimilarity matrices were performed in QIIME, and PCoA plots were illustrated in

838 R using the ggplot2 package⁹¹. PCA was performed in R using the ggbiplot package⁹². Spearman's

839 Rank correlations were performed in GraphPad Prism, whereas partial Spearman's Rank correlation

analyses with adjustments for age and repeated measures correlation analyses were performed in R using the $ppcor^{29}$ and rmcorr packages²⁸, respectively. Heatmaps and hierarchically clustering of

using the $ppcor^{29}$ and rmcorr packages²⁸, respectively. Heatmaps and hierarchically clustering of correlation coefficient were generated in R using the *gplots* package⁹³ and visualized in GraphPad

Prism. Longitudinal metabolite and taxonomic abundance were modelled using LOESS regression,

and implemented and plotted with 95% confidence intervals in R using the *ggplot2* package⁹¹.

Longitudinal AhR activity was modelled used coarse LOWESS curve fits within the GraphPad

Prism software. Two-tailed Student's *t* test (if normally distributed, evaluated by D'Agostino-

847 Pearson test) or two-tailed non-parametric Mann-Whitney U test (if not normally distributed) were

848 performed when comparing two groups. For comparison of more than two groups, statistical

significance was evaluated by one-way ANOVA (if normally distributed) or the non-parametric

850 Kruskal-Wallis test (if not normally distributed). P-values < 0.05 were considered statistically

significant. When applicable p-values were corrected for multiple testing by the Benjamini–

Hochberg false discovery rate $(FDR)^{94}$ using a cutoff of 0.1.

853 ACKNOWLEDGMENTS

We thank the children and families participating in the SKOT I study, which was supported by the

B55 Danish Directorate for Food, Fisheries and Agribusiness (Grant no. 3304-FSE-06-0503), as well as

the children and parents participating in the CIG cohort. Furthermore, we thank Aarstiderne A/S for

857 providing a small gift for the CIG participants. We thank Marlene Danner Dalgaard at the Technical

858 University of Denmark in-house facility (DTU Multi-Assay Core, DMAC) for performing the 16S

rRNA gene sequencing, MS-Omics (Hørsholm, Denmark) for performing the lactate analyses,

860 Glycom A/S (Hørsholm, Denmark) for kindly donating the human milk oligosaccharides, and

861 Anette Schnipper for her efforts in supporting the work. We also thank Satoru Fukiya and Atsushi

862 Yokota for providing *Bifidobacterium* gene manipulation tools and for technical suggestions, Yuta

- 863 Sugiyama for helpful discussions on HPLC analysis, and Shingo Maeda for technical support on
- enzyme purification. This work was supported by Augustinus Fonden (grant no. 17-2003 to
- H.M.R.), Hørslev Fonden (grant no. 203866 to H.M.R.), Beckett Fonden (grant no. 17-2-0551 to
- H.M.R.), Aase og Ejnar Danielsens Fond (grant no. 10-002019 to H.M.R.), the Innovation Fund
- B67 Denmark (grant no. 11-116163/0603-00487B; Center for Gut, Grain and Greens to T.R.L.), JSPS-
- KAKENHI (18K14379 to M.S., 19K22277 to T.K.), JSPS Overseas Research Fellowships
- 869 (201860637 to M.S.) and the Institute for Fermentation, Osaka (to M.S. and T.K.) and "Diet-
- 870 induced Arrangement of the gut Microbiome for improvement of Cardiometabolic health"
- 871 (DINAMIC) under the Joint Programming Initiative, "A Healthy Diet for a Healthy Life" (JPI-
- HDHL), supported by the Innovation Fund Denmark (grant no. 5195-00001B to L.O.D.).

873 AUTHOR CONTRIBUTIONS

H.M.R. and M.F.L. conceived and designed the experiments. M.F.L. prepared the samples for

- sequencing/qPCR and analyzed the sequencing/qPCR data. H.M.R. prepared the samples for faecal
- and *in vitro* metabolome analyses and performed together with H.L.F. the targeted and untargeted
- metabolomics experiments. C.T.P. and L.O.D. performed the urine metabolomics. M.F.L. and M.S.
- performed the *in vitro* growth and mutant construction experiments. M.S. and T.K. performed
- enzyme kinetics. K.F.M. and C.M. designed the SKOT I study and M.V.L. managed the data.
- 880 H.M.R. and M.F.L. designed the CIG cohort, recruited the study participants and managed the data.
- A.M.V. performed the AhR assays. N.B., D.A., U.M., A.R., and J.M.M. contributed to the
- interpretation of the results. S.B., W.A., T.K., M.I.B. and T.R.L. contributed with expert
- supervision. H.M.R. and M.F.L. led the work, undertook the integrative data analyses and drafted
- the manuscript. All authors contributed to and approved the final manuscript.

885 Data statement

Partial (V3 region) 16S rRNA gene amplicon sequencing data is deposited in the Sequence Read

- 887 Archive (SRA) under the BioProjects PRJNA273694 (SKOT) and PRJNA554596 (CIG).
- 888 Competing Financial Interests statement
- 889 The authors declare no competing financial interests.

890 Corresponding authors

- 891 # Henrik M. Roager
- 892 # Tine R. Licht

893 **REFERENCES**

- Victora, C. G. *et al.* Breastfeeding in the 21st century: epidemiology, mechanisms, and
 lifelong effect. *Lancet (London, England)* 387, 475–90 (2016).
- 896 2. Horta, B. L., Loret de Mola, C. & Victora, C. G. Long-term consequences of breastfeeding
 897 on cholesterol, obesity, systolic blood pressure and type 2 diabetes: a systematic review and
 898 meta-analysis. *Acta Paediatr.* 104, 30–7 (2015).
- Baursen, M. F. *et al.* Infant Gut Microbiota Development Is Driven by Transition to Family
 Foods Independent of Maternal Obesity. *mSphere* 1, e00069-15 (2016).
- 901 4. Bäckhed, F. *et al.* Dynamics and stabilization of the human gut microbiome during the first
 902 year of life. *Cell Host Microbe* 17, 690–703 (2015).
- 5. Stewart, C. J. *et al.* Temporal development of the gut microbiome in early childhood from
 the TEDDY study. *Nature* 562, 583–588 (2018).
- Bode, L. The functional biology of human milk oligosaccharides. *Early Hum. Dev.* 91, 619–622 (2015).
- 907 7. Gensollen, T., Iyer, S. S., Kasper, D. L. & Blumberg, R. S. How colonization by microbiota in early life shapes the immune system. *Science (80-.).* 352, 539–544 (2016).
- 8. Stokholm, J. *et al.* Maturation of the gut microbiome and risk of asthma in childhood. *Nat. Commun.* 9, 141 (2018).
- 9. Fujimura, K. E. *et al.* Neonatal gut microbiota associates with childhood multisensitized atopy and T cell differentiation. *Nat. Med.* 22, 1187–1191 (2016).
- 913 10. Khosravi, A. & Mazmanian, S. K. Disruption of the gut microbiome as a risk factor for
 914 microbial infections. *Curr. Opin. Microbiol.* 16, 221–227 (2013).
- 915 11. Fukuda, S. *et al.* Bifidobacteria can protect from enteropathogenic infection through
 916 production of acetate. *Nature* 469, 543–547 (2011).
- P17 12. Roager, H. M. & Licht, T. R. Microbial tryptophan catabolites in health and disease. *Nat.*P18 *Commun.* 9, 3294 (2018).
- 919 13. Natividad, J. M. *et al.* Impaired Aryl Hydrocarbon Receptor Ligand Production by the Gut
 920 Microbiota Is a Key Factor in Metabolic Syndrome. *Cell Metab.* (2018).
 921 doi:10.1016/j.cmet.2018.07.001
- 922 14. Dodd, D. *et al.* A gut bacterial pathway metabolizes aromatic amino acids into nine
 923 circulating metabolites. *Nature* 551, 648–652 (2017).
- 924 15. Guo, X. *et al.* Innate Lymphoid Cells Control Early Colonization Resistance against
 925 Intestinal Pathogens through ID2-Dependent Regulation of the Microbiota. *Immunity* 42,
 926 731–743 (2015).
- 927 16. Zelante, T. *et al.* Tryptophan catabolites from microbiota engage aryl hydrocarbon receptor

928		and balance mucosal reactivity via interleukin-22. Immunity 39, 372–385 (2013).		
929 930	17.	Hoyles, L. <i>et al.</i> Molecular phenomics and metagenomics of hepatic steatosis in non-diabet obese women. <i>Nat. Med.</i> 24 , 1070–1080 (2018).		
931 932	18.	Krishnan, S. <i>et al.</i> Gut Microbiota-Derived Tryptophan Metabolites Modulate Inflammato Response in Hepatocytes and Macrophages. <i>Cell Rep.</i> 23 , 1099–1111 (2018).		
933 934 935	19.	Madsen, A. L., Schack-Nielsen, L., Larnkjaer, A., Mølgaard, C. & Michaelsen, K. F. Determinants of blood glucose and insulin in healthy 9-month-old term Danish infants; the SKOT cohort. <i>Diabet. Med.</i> 27 , 1350–7 (2010).		
936 937	20.	Kato, K. <i>et al.</i> Age-Related Changes in the Composition of Gut Bifidobacterium Species. <i>Curr. Microbiol.</i> 74 , 987–995 (2017).		
938 939	21.	Koga, Y. <i>et al.</i> Age-associated effect of kestose on Faecalibacterium prausnitzii and symptoms in the atopic dermatitis infants. <i>Pediatr. Res.</i> 80 , 844–851 (2016).		
940 941 942	22.	Li, X., Jiang, B., Pan, B., Mu, W. & Zhang, T. Purification and partial characterization of Lactobacillus species SK007 lactate dehydrogenase (LDH) catalyzing phenylpyruvic aci (PPA) conversion into phenyllactic acid (PLA). <i>J. Agric. Food Chem.</i> 56 , 2392–2399 (2)		
943 944 945	23.	Koide, S., Iwata, S., Matsuzawa, H. & Ohta, T. Crystallization of allosteric L-lactate dehydrogenase from Thermus caldophilus and preliminary crystallographic data. <i>J. Biochem.</i> 109 , 6–7 (1991).		
946 947 948	24.	Takashi, M., So, I., Hiroshi, S., Haruhiko, M. & Takahisa, O. Sequence and characteristics of the Bifidobacterium longum gene encoding l-lactate dehydrogenase and the primary structure of the enzyme: a new feature of the allosteric site. <i>Gene</i> 85 , 161–168 (1989).		
949 950	25.	Bottacini, F. <i>et al.</i> Global transcriptional landscape and promoter mapping of the gut commensal Bifidobacterium breve UCC2003. <i>BMC Genomics</i> 18 , 991 (2017).		
951 952 953	26.	Hirayama, Y. <i>et al.</i> Development of a double-crossover markerless gene deletion system in Bifidobacterium longum: Functional analysis of the ??-galactosidase gene for raffinose assimilation. <i>Appl. Environ. Microbiol.</i> 78 , 4984–4994 (2012).		
954 955 956	27.	Matsumura, H., Takeuchi, A. & Kano, Y. Construction of Escherichia coli-Bifidobacterium longum shuttle vector transforming B. longum 105-A and 108-A. <i>Biosci. Biotechnol. Biochem.</i> 61 , 1211–2 (1997).		
957 958	28.	Bakdash, J. Z. & Marusich, L. R. Repeated Measures Correlation. <i>Front. Psychol.</i> 8 , 456 (2017).		
959 960	29.	Kim, S. ppcor: An R Package for a Fast Calculation to Semi-partial Correlation Coefficients. <i>Commun. Stat. Appl. methods</i> 22 , 665–674 (2015).		
961 962 963	30.	Alfa, M. J. <i>et al.</i> An Outbreak of Necrotizing Enterocolitis Associated with a Novel Clostridium Species in a Neonatal Intensive Care Unit. <i>Clin. Infect. Dis.</i> 35 , S101–S105 (2002).		

964	31.	Butel, MJ. et al. Conditions of Bifidobacterial Colonization in Preterm Infants: A
965		Prospective Analysis. J. Pediatr. Gastroenterol. Nutr. 44, (2007).

- 32. Moles, L. *et al.* Bacterial Diversity in Meconium of Preterm Neonates and Evolution of Their
 Fecal Microbiota during the First Month of Life. *PLoS One* 8, e66986 (2013).
- 33. Arboleya, S. *et al.* Establishment and development of intestinal microbiota in preterm
 neonates. *FEMS Microbiol. Ecol.* **79**, 763–772 (2012).
- 34. Korpela, K. *et al.* Intestinal microbiota development and gestational age in preterm neonates. *Sci. Rep.* 8, 2453 (2018).
- 972 35. Cervantes-Barragan, L. *et al.* Lactobacillus reuteri induces gut intraepithelial
 973 CD4(+)CD8αα(+) T cells. *Science* eaah5825 (2017). doi:10.1126/science.aah5825
- 36. Tamburini, S., Shen, N., Wu, H. C. & Clemente, J. C. The microbiome in early life:
 implications for health outcomes. *Nat. Med.* 22, 713–722 (2016).
- 37. Dominguez-Bello, M. G., Godoy-Vitorino, F., Knight, R. & Blaser, M. J. Role of the
 microbiome in human development. *Gut* 68, 1108–1114 (2019).
- 38. Sakurai, T., Odamaki, T. & Xiao, J.-Z. Production of Indole-3-Lactic Acid by
 Bifidobacterium Strains Isolated fromHuman Infants. *Microorganisms* 7, 340 (2019).
- 39. Yamada, C. *et al.* Molecular Insight into Evolution of Symbiosis between Breast-Fed Infants
 and a Member of the Human Gut Microbiome Bifidobacterium longum. *Cell Chem. Biol.* 24,
 515-524.e5 (2017).
- 40. Katayama, T. Host-derived glycans serve as selected nutrients for the gut microbe: human
 milk oligosaccharides and bifidobacteria. *Biosci. Biotechnol. Biochem.* 80, 621–632 (2016).
- 41. Thomson, P., Medina, D. A. & Garrido, D. Human milk oligosaccharides and infant gut
 bifidobacteria: Molecular strategies for their utilization. *Food Microbiol.* 75, 37–46 (2018).
- 987 42. Sakanaka, M. *et al.* Evolutionary adaptation in fucosyllactose uptake systems supports
 988 bifidobacteria-infant symbiosis. *Sci. Adv.* 5, eaaw7696 (2019).
- Vatanen, T. *et al.* The human gut microbiome in early-onset type 1 diabetes from the
 TEDDY study. *Nature* 562, 589–594 (2018).
- 44. Natividad, J. M. *et al.* Impaired Aryl Hydrocarbon Receptor Ligand Production by the Gut
 Microbiota Is a Key Factor in Metabolic Syndrome. *Cell Metab.* 28, 737-749.e4 (2018).
- Metidji, A. *et al.* The Environmental Sensor AHR Protects from Inflammatory Damage by
 Maintaining Intestinal Stem Cell Homeostasis and Barrier Integrity. *Immunity* 49, 353-362.e5
 (2018).
- 46. Gomez de Aguero, M. *et al.* The maternal microbiota drives early postnatal innate immune
 development. *Science (80-.).* 351, 1296–1302 (2016).
- 998 47. Kiss, E. A. et al. Natural Aryl Hydrocarbon Receptor Ligands Control Organogenesis of

999 Intestinal Lymphoid Follicles. Science (80-.). 334, 1561–1565 (2011). 1000 48. Qiu, J. et al. The Aryl Hydrocarbon Receptor Regulates Gut Immunity through Modulation of Innate Lymphoid Cells. Immunity 36, 92-104 (2012). 1001 49. Lamas, B. et al. CARD9 impacts colitis by altering gut microbiota metabolism of tryptophan 1002 1003 into aryl hydrocarbon receptor ligands. Nat. Med. (2016). doi:10.1038/nm.4102 50. Hubbard, T. D. et al. Adaptation of the human aryl hydrocarbon receptor to sense 1004 microbiota-derived indoles. Sci. Rep. 5, 12689 (2015). 1005 Artis, D. & Spits, H. The biology of innate lymphoid cells. *Nature* 517, 293–301 (2015). 1006 51. Dant, T. A. et al. T-cell expression of AhR inhibits the maintenance of pT reg cells in the 1007 52. gastrointestinal tract in acute GVHD. Blood 130, 348-359 (2017). 1008 1009 53. Wilck, N. et al. Salt-responsive gut commensal modulates TH17 axis and disease. Nature 551, 585–589 (2017). 1010 Quintana, F. J. et al. Control of T(reg) and T(H)17 cell differentiation by the aryl 1011 54. hydrocarbon receptor. Nature 453, 65-71 (2008). 1012 55. 1013 Ning, Y. et al. Antibacterial activity of phenyllactic acid against Listeria monocytogenes and Escherichia coli by dual mechanisms. Food Chem. 228, 533-540 (2017). 1014 1015 56. Rodríguez, N., Salgado, J. M., Cortés, S. & Domínguez, J. M. Antimicrobial activity of d-3phenyllactic acid produced by fed-batch process against Salmonella enterica. Food Control 1016 1017 25, 274–284 (2012). 1018 57. Narayanan, T. K. & Rao, G. R. Beta-indoleethanol and beta-indolelactic acid production by Candida species: their antibacterial and autoantibiotic action. Antimicrob. Agents Chemother. 1019 9, 375-80 (1976). 1020 Honoré, A. H. et al. Metabolic footprinting for investigation of antifungal properties of 1021 58. 1022 Lactobacillus paracasei. Anal. Bioanal. Chem. 408, 83-96 (2016). 59. Peters, A. et al. Metabolites of lactic acid bacteria present in fermented foods are highly 1023 potent agonists of human hydroxycarboxylic acid receptor 3. PLOS Genet. 15, e1008145 1024 1025 (2019). 1026 60. Offermanns, S. Hydroxy-Carboxylic Acid Receptor Actions in Metabolism. Trends Endocrinol. Metab. 28, 227–236 (2017). 1027 1028 61. Ahmed, K. et al. Deorphanization of GPR109B as a Receptor for the β-Oxidation Intermediate 3-OH-octanoic Acid and Its Role in the Regulation of Lipolysis. J. Biol. Chem. 1029 1030 **284**, 21928–21933 (2009). 1031 62. Irukayama-Tomobe, Y. et al. Aromatic D-amino acids act as chemoattractant factors for human leukocytes through a G protein-coupled receptor, GPR109B. Proc. Natl. Acad. Sci. 1032 106, 3930–3934 (2009). 1033

1034 1035 1036	63.	Madsen, A. L., Larnkjær, A., Mølgaard, C. & Michaelsen, K. F. IGF-I and IGFBP-3 in healthy 9 month old infants from the SKOT cohort: Breastfeeding, diet, and later obesity. <i>Growth Horm. IGF Res.</i> 21 , 199–204 (2011).		
1037 1038	64.	Laursen, M. F. <i>et al.</i> Having older siblings is associated with gut microbiota development during early childhood. <i>BMC Microbiol.</i> 15 , 154 (2015).		
1039 1040 1041	65.	Laursen, M. F., Dalgaard, M. D. & Bahl, M. I. Genomic GC-Content Affects the Accuracy 16S rRNA Gene Sequencing Based Microbial Profiling due to PCR Bias. <i>Front. Microbiol.</i> 8 , 1934 (2017).		
1042 1043 1044	66.	Laursen, M. F. <i>et al.</i> Administration of two probiotic strains during early childhood does no affect the endogenous gut microbiota composition despite probiotic proliferation. <i>BMC Microbiol.</i> 17 , 175 (2017).		
1045 1046	67.	Edgar, R. C. UPARSE: highly accurate OTU sequences from microbial amplicon reads. <i>Na Methods</i> 10 , 996–8 (2013).		
1047 1048	68.	Caporaso, J. G. <i>et al.</i> QIIME allows analysis of high-throughput community sequencing data. <i>Nat. Methods</i> 7 , 335–6 (2010).		
1049 1050 1051	69.	Wang, Q., Garrity, G. M., Tiedje, J. M. & Cole, J. R. Naive Bayesian classifier for rapid assignment of rRNA sequences into the new bacterial taxonomy. <i>Appl. Environ. Microbiol.</i> 73 , 5261–7 (2007).		
1052 1053	70.	DeSantis, T. Z. <i>et al.</i> Greengenes, a chimera-checked 16S rRNA gene database and workbench compatible with ARB. <i>Appl. Environ. Microbiol.</i> 72 , 5069–72 (2006).		
1054 1055	71.	Altschul, S. F., Gish, W., Miller, W., Myers, E. W. & Lipman, D. J. Basic local alignment search tool. <i>J. Mol. Biol.</i> 215 , 403–410 (1990).		
1056 1057	72.	Frese, S. A. <i>et al.</i> Persistence of Supplemented Bifidobacterium longum subsp. infantis EVC001 in Breastfed Infants. (2017). doi:10.1128/mSphere		
1058 1059	73.	Lawley, B. <i>et al.</i> Differentiation of <i>Bifidobacterium longum</i> subspecies <i>longum</i> and <i>infantis</i> by quantitative PCR using functional gene targets. <i>PeerJ</i> 5 , e3375 (2017).		
1060 1061 1062	74.	Sprenger, N., Lee, L. Y., De Castro, C. A., Steenhout, P. & Thakkar, S. K. Longitudinal change of selected human milk oligosaccharides and association to infants' growth, an observatory, single center, longitudinal cohort study. <i>PLoS One</i> 12 , e0171814 (2017).		
1063 1064	75.	Xu, G. <i>et al.</i> Absolute Quantitation of Human Milk Oligosaccharides Reveals Phenotypic Variations during Lactation. <i>J. Nutr.</i> 147 , 117–124 (2017).		
1065 1066 1067	76.	Sakurama, H. <i>et al.</i> Lacto-N-biosidase encoded by a novel gene of bifidobacterium longum subspecies longum shows unique substrate specificity and requires a designated chaperone for its active expression. <i>J. Biol. Chem.</i> 288 , 25194–25206 (2013).		
1068 1069	77.	Sakanaka, M. <i>et al.</i> Functional analysis of bifidobacterial promoters in Bifidobacterium longum and Escherichia coli using the α -galactosidase gene as a reporter. <i>J. Biosci. Bioeng.</i>		

1070 118, 489–495 (2014).

1071 78. Poolman, B., Nijssen, R. M. J. & Konings, W. N. Dependence of Streptococcus lactis
1072 phosphate transport on internal phosphate concentration and internal pH. *J. Bacteriol.* 169, 5373–5378 (1987).

- 1074 79. Barri, T., Holmer-Jensen, J., Hermansen, K. & Dragsted, L. O. Metabolic fingerprinting of high-fat plasma samples processed by centrifugation- and filtration-based protein
 1076 precipitation delineates significant differences in metabolite information coverage. *Anal.*1077 *Chim. Acta* **718**, 47–57 (2012).
- 1078 80. Nielsen, L. N. *et al.* Glyphosate has limited short-term effects on commensal bacterial
 1079 community composition in the gut environment due to sufficient aromatic amino acid levels.
 1080 *Environ. Pollut.* 233, (2018).
- 1081 81. Roager, H. M. *et al.* Colonic transit time is related to bacterial metabolism and mucosal
 1082 turnover in the gut. *Nat. Microbiol.* 1, 16093 (2016).

Smith, C. A., Want, E. J., O'Maille, G., Abagyan, R. & Siuzdak, G. XCMS: Processing Mass
Spectrometry Data for Metabolite Profiling Using Nonlinear Peak Alignment, Matching, and
Identification. *Anal. Chem.* 78, 779–787 (2006).

- 1086 83. Sumner, L. W. *et al.* Proposed minimum reporting standards for chemical analysis Chemical
 1087 Analysis Working Group (CAWG) Metabolomics Standards Initiative (MSI). *Metabolomics*1088 3, 211–221 (2007).
- 1089 84. Gürdeniz, G., Kristensen, M., Skov, T. & Dragsted, L. O. The Effect of LC-MS Data
 1090 Preprocessing Methods on the Selection of Plasma Biomarkers in Fed vs. Fasted Rats.
 1091 *Metabolites* 2, 77–99 (2012).
- Pluskal, T., Castillo, S., Villar-Briones, A. & Oresic, M. MZmine 2: modular framework for
 processing, visualizing, and analyzing mass spectrometry-based molecular profile data. *BMC Bioinformatics* 11, 395 (2010).
- 1095 86. van der Kloet, F. M., Bobeldijk, I., Verheij, E. R. & Jellema, R. H. Analytical error reduction
 1096 using single point calibration for accurate and precise metabolomic phenotyping. *J. Proteome*1097 *Res.* 8, 5132–41 (2009).
- 1098 87. Smart, K. F., Aggio, R. B. M., Houtte, J. V. R. Van & Villas-Bôas, S. G. Analytical platform
 1099 for metabolome analysis of microbial cells using methyl chloroformate derivatization
 1100 followed by gas chromatography-mass spectrometry. *Nat. Protoc.* 5, 1709–1729 (2010).
- 1101 88. Johnsen, L. G., Skou, P. B., Khakimov, B. & Bro, R. Gas chromatography mass
 1102 spectrometry data processing made easy. *J. Chromatogr. A* 1503, 57–64 (2017).
- 1103 89. Rosenmai, A. K. *et al.* Are structural analogues to bisphenol a safe alternatives? *Toxicol. Sci.*1104 139, 35–47 (2014).
- 1105 90. R Core Team. R: A language and environment for statistical computing. (2013).

1106 91. Wickhamm, H. ggplot2 Elegant Graphics for Data Analysis. (Springer, New York, 2009).

- 1107 92. Vu, V. ggbiplot: a ggplot2 based biplot. 2011
- 1108 93. Warnes, G. R. *et al.* gplots: Various R Programming Tools for Plotting Data. R package
 1109 version 2.16.0. *CRAN* (2015).
- 94. Benjamini, Y. & Hochberg, Y. Controlling the False Discovery Rate a Practical and
 Powerful Approach to Multiple Testing. *J. R. Stat. Soc. Ser. B-Methodological* 57, 289–300 (1995).
- Smith, E. A. & Macfarlane, G. T. Enumeration of human colonic bacteria producing phenolic and indolic compounds: effects of pH, carbohydrate availability and retention time on dissimilatory aromatic amino acid metabolism. *J. Appl. Bacteriol.* 81, 288–302 (1996).
- Smith, E. A. & Macfarlane, G. T. Formation of Phenolic and Indolic Compounds by
 Anaerobic Bacteria in the Human Large Intestine. *Microb. Ecol.* 33, 180–188 (1997).



С

b



Kinetic parameters (Hill equation) of type 4 LDH

Substrate	K _{0.5} (mM)	<i>k</i> _{cat} (s ⁻¹)	<i>k</i> _{cat} / <i>K</i> _{0.5} (s ⁻¹ mM ⁻¹)	n _H (Hill coefficient)
IPA	0.08 ± 0.01	15.21 ± 1.01	193.73	1.05 ± 0.02
PPA	5.96 ± 0.06	64.22 ± 2.38	10.77	1.20 ± 0.01
4-OH-PPA	6.61 ± 0.93	103.89 ± 4.32	15.83	1.39 ± 0.01
Pyruvate	14.88 ± 0.48	0.47 ± 0.03	0.03	1.28 ± 0.05

g



Age (days) 





