

Oxysterols are agonist ligands of ROR γ t and drive Th17 cell differentiation

Pejman Soroosh^{a,1}, Jiejun Wu^{a,1}, Xiaohua Xue^{a,1}, Jiao Song^a, Steven W. Sutton^a, Marciano Sablad^a, Jingxue Yu^a, Marina I. Nelen^b, Xuejun Liu^a, Glenda Castro^a, Rosa Luna^a, Shelby Crawford^{a,2}, Homayon Banie^a, Rose A. Dandridge^b, Xiaohu Deng^a, Anton Bittner^a, Chester Kuei^a, Mandana Tootoonchi^a, Natasha Rozenkrants^a, Krystal Herman^a, Jingjin Gao^a, Xia V. Yang^{a,3}, Kacey Sachin^{a,4}, Karen Ngo^a, Wai-Ping Fung-Leung^a, Steven Nguyen^a, Aimee de Leon-Tabaldo^a, Jonathan Blevitt^a, Yan Zhang^a, Maxwell D. Cummings^b, Tadimeti Rao^a, Neelakandha S. Mani^a, Changlu Liu^a, Murray McKinnon^b, Marcos E. Milla^a, Anne M. Fourie^a, and Siqian Sun^{a,5}

^aJanssen Research and Development, LLC, San Diego, CA 92121; and ^bJanssen Research and Development, Spring House, PA 19002

Edited by Peter S. Kim, Stanford University School of Medicine, Stanford, CA, and approved July 10, 2014 (received for review December 10, 2013)

The RAR-related orphan receptor gamma t (ROR γ t) is a nuclear receptor required for generating IL-17-producing CD4⁺ Th17 T cells, which are essential in host defense and may play key pathogenic roles in autoimmune diseases. Oxysterols elicit profound effects on immune and inflammatory responses as well as on cholesterol and lipid metabolism. Here, we describe the identification of several naturally occurring oxysterols as ROR γ t agonists. The most potent and selective activator for ROR γ t is 7 β , 27-dihydroxycholesterol (7 β , 27-OHC). We show that these oxysterols reverse the inhibitory effect of an ROR γ t antagonist, ursolic acid, in ROR γ - or ROR γ t-dependent cell-based reporter assays. These ligands bind directly to recombinant ROR γ ligand binding domain (LBD), promote recruitment of a coactivator peptide, and reduce binding of a corepressor peptide to ROR γ LBD. In primary cells, 7 β , 27-OHC and 7 α , 27-OHC enhance the differentiation of murine and human IL-17-producing Th17 cells in an ROR γ t-dependent manner. Importantly, we showed that Th17, but not Th1 cells, preferentially produce these two oxysterols. In vivo, administration of 7 β , 27-OHC in mice enhanced IL-17 production. Mice deficient in CYP27A1, a key enzyme in generating these oxysterols, showed significant reduction of IL-17-producing cells, including CD4⁺ and $\gamma\delta$ ⁺ T cells, similar to the deficiency observed in ROR γ t knockout mice. Our results reveal a previously unknown mechanism for selected oxysterols as immune modulators and a direct role for CYP27A1 in generating these ROR γ t agonist ligands, which we propose as ROR γ t endogenous ligands, driving both innate and adaptive IL-17-dependent immune responses.

IL-17-producing CD4⁺ Th17 cells are essential in protective immunity against extracellular bacterial and fungal infections but also, play key pathogenic roles in autoimmune diseases, such as psoriasis and multiple sclerosis. RAR-related orphan receptor gamma t (ROR γ t) is an orphan nuclear receptor expressed in several immune cell types, including CD4⁺ Th17 cells. Because of its essential role in driving IL-17 production, ROR γ t represents a potential target for therapeutic intervention, and a number of antagonists have been published that inhibited Th17 cell differentiation (1–3). However, the nature of ROR γ t endogenous ligands remains unknown, although accumulating evidence suggested the existence of such ligands. For example, crystal structures of the ligand binding domain of ROR γ clearly showed a well-defined pocket that could accommodate a cholesterol or 25-hydroxycholesterol (OHC) molecule (4). Purified naïve CD4⁺ T cells can be activated in culture to differentiate into IL-17-producing cells, suggesting that these cells have the capacity to produce endogenous ROR γ t ligand(s). In addition, IL-17-producing mouse ROR γ t⁺ innate lymphoid cells, believed to be essential in orchestrating immunity at mucosal sites, including intestine and lung, were shown to be generated in vivo under germfree conditions (5). Identification of the endogenous agonist ligand(s) and understanding of how the ligand engages

and activates ROR γ t may, therefore, offer critical insights into the rational design of ROR γ t modulators to block pathogenic IL-17-producing cells in disease.

Oxysterols are oxygenated derivatives of cholesterol or by-products of cholesterol synthesis (6–9). Some oxysterols serve as key intermediates for bile acid and steroid synthesis and function as signaling molecules or receptor ligands modulating cell proliferation and apoptosis, lipid and cholesterol synthesis, transportation and metabolism, and immune responses. For example, certain oxysterols have been shown to serve as endogenous ligands of EBI2 receptor (10, 11), activate the liver X receptor (12), function as a selective estrogen receptor modulator (13), or affect Hedgehog signaling by binding to the Smoothed molecule (14). In addition, several oxysterols were previously shown to function as ROR γ or ROR α inverse agonists (such as 7-hydroxylated oxysterols) or 24(S)-OHC agonists (such as 25-OHC) (4, 15, 16). However, their role as endogenous ROR γ , ROR γ t, or ROR α ligands in vivo was not clearly shown.

We have identified a number of naturally occurring oxysterols as ROR γ t agonists. Although we cannot exclude at present the existence of additional endogenous agonists of ROR γ t, here we show that two 7, 27-dihydroxycholesterols (i.e., 7 β , 27-OHC and 7 α , 27-OHC) are strong candidates as endogenous ROR γ t

Significance

Because of its essential role in driving IL-17 production, the orphan nuclear receptor RAR-related orphan receptor gamma t (ROR γ t) represents a potential therapeutic target for autoimmune diseases. Here, we present evidence that 7 β , 27-dihydroxycholesterol (7 β , 27-OHC) and 7 α , 27-OHC are ROR γ t agonist ligands and may serve as potential endogenous ROR γ t ligands in promoting the differentiation of mouse and human CD4⁺ Th17 cells. These findings may not only facilitate rational design of ROR γ t modulators but also, provide insights into additional targets for inhibiting IL-17 production.

Author contributions: P.S., J.W., X.X., J.S., T.R., N.S.M., M.M., A.M.F., and S.S. designed research; P.S., J.W., X.X., J.S., S.W.S., M.S., J.Y., M.I.N., G.C., R.L., S.C., H.B., R.A.D., X.D., A.B., C.K., M.T., N.R., K.H., J.G., X.V.Y., K.S., K.N., W.-P.F.-L., S.N., A.d.L.-T., J.B., Y.Z., and C.L. performed research; X.D., N.S.M., and C.L. contributed new reagents/analytic tools; P.S., J.W., X.X., J.S., S.W.S., J.Y., M.I.N., X.L., R.L., S.C., H.B., M.D.C., and S.S. analyzed data; and P.S., J.W., X.X., M.I.N., X.D., C.K., M.E.M., A.M.F., and S.S. wrote the paper.

The authors declare no conflict of interest.

This article is a PNAS Direct Submission.

Freely available online through the PNAS open access option.

¹P.S., J.W., and X.X. contributed equally to this work.

²Present address: Private address, San Diego, CA 92130.

³Present address: Regulus Therapeutics, San Diego, CA 92121.

⁴Present address: Takeda San Diego Inc., San Diego, CA 92121.

⁵To whom correspondence should be addressed. Email: ssun1@its.jnj.com.

This article contains supporting information online at www.pnas.org/lookup/suppl/doi:10.1073/pnas.1322807111/-DCSupplemental.

agonists driving IL-17 production in CD4⁺ Th17 cells as well as other IL-17-producing innate cells, such as $\gamma\delta^+$ T cells.

Results

ROR γ t Agonist Activity of Selected Oxysterols. To identify potential ROR γ t agonists, we screened a panel of naturally occurring oxysterols and a few other compounds (SI Appendix, Tables S1 and S2) for their agonist activity in reversing the inhibitory effect of ursolic acid (UA), an ROR γ t antagonist (3). We first used a cell-based ROR γ -dependent reporter assay, in which the DNA-binding domain of GAL4 was fused to the ligand binding domain (LBD) of human ROR γ . This construct was transiently expressed in HEK293T cells and showed constitutive activity in driving luciferase reporter gene expression. Because ROR γ t and the related isotype, ROR γ , share the same LBD, such an assay potentially allowed detection of agonists for both receptors. A number of oxysterols showed significant agonist activity, with 27-OHC, 7 β , 27-OHC, and 7-keto, 27-OHC being the most potent and efficacious (Fig. 1A and SI Appendix, Tables S1 and S2). In comparison, 7 α , 25-OHC, an endogenous ligand for EBI2 receptor, was inactive. Several previously reported ROR γ or ROR α inverse agonists or agonists including 7 α -OHC, 7 β -OHC,

7-KC, 25-OHC, 24(S), 25-EC, and 24(S)-OHC (15, 16) showed weak to moderate agonist activity. Cholestenic acid derivatives of the 27-hydroxylated sterols displayed minimal agonist activity (SI Appendix, Table S1), indicating that the hydroxyl group at carbon 27 is required for ROR γ t agonism. We next tested selected oxysterols, including all four 27-OHCs, in cell-based full-length human ROR γ - or ROR γ t-dependent reporter assays. One clear exception was 7-keto, 27-OHC, which was largely inactive in those two assays, whereas other 27-OHCs were active, and there was no apparent difference between ROR γ - and ROR γ t-dependent assays (SI Appendix, Fig. S1).

To establish the selectivity of the ROR γ t-active 27-OHCs, we tested them for potential agonist activity in several nuclear receptor (NR) reporter assays (SI Appendix, Fig. S2). While 27-OHC showed weak activity for LXR α and 27-OHC and 7-keto, 27-OHC showed moderate activity for LXR β , 7 β , 27-OHC and 7 α , 27-OHC were inactive up to 30 μ M. 27-OHC also showed weak activity for ER β , but 7 β , 27-OHC, 7 α , 27-OHC, and 7-keto, 27-OHC were inactive against ER α or ER β . In addition, none of the 27-OHCs showed agonist activity up to 30 μ M against other NRs, including PPAR α , PPAR β , PPAR γ , CAR, GR, and bile acid-sensing FXR. When tested in antagonist mode (SI Appendix, Fig. S2), 7-keto, 27-OHC showed inhibition of all NRs tested, except LXR α and LXR β . Neither 7 β , 27-OHC nor 7 α , 27-OHC showed any antagonist activity against all NRs tested except for ER β , against which all 27-OHCs showed some activity, particularly 7-keto, 27-OHC and 27-OHC. Taken together, 7 β , 27-OHC and 7 α , 27-OHC seemed to be the most selective ROR γ t agonists among the four 27-OHCs in cell-based reporter assays.

Direct Binding of Oxysterols to ROR γ LBD. We next measured direct binding of oxysterols to recombinant ROR γ LBD protein in the absence or presence of a peptide from the steroid receptor coactivator-1 (SRC1) using the thermal shift readout ThermoFluor (17–19) (SI Appendix, Fig. S3A). Direct binding of oxysterols to recombinant LBD of ROR α and ROR β in the presence of SRC1 peptide was also examined (SI Appendix, Fig. S3A). Consistent with their activities in the cell-based chimeric ROR γ -dependent reporter assay, 27-OHCs were among the most potent binders to ROR γ LBD, with a rank order of 7 β , 27-OHC > 7-keto, 27-OHC > 27-OHC > 7 α , 27-OHC. The presence of SRC1 peptide did not affect binding to ROR γ LBD by 27-OHCs except 27-OHC, which showed reduced affinity, suggesting that this most abundant oxysterol in the body is unlikely to be an endogenous ROR γ agonist. When tested against the LBDs of ROR α and ROR β in the presence of SRC1, some selectivity was observed: 7 α , 27-OHC was inactive in binding to ROR α and ROR β , and 7 β , 27-OHC was inactive in binding to ROR α . In contrast, 7-keto, 27-OHC bound to all three ROR LBD proteins and showed the highest affinity for ROR α in the presence of SRC1 peptide. Taken together, these findings suggest that 7 β , 27-OHC and 7 α , 27-OHC are more selective agonist ligands for ROR γ t than 27-OHC or 7-keto, 27-OHC.

25-OHC was previously shown to form a complex with ROR γ LBD, and ³H-labeled 25-OHC exhibited specific binding to recombinant ROR γ LBD in vitro (4), which showed a K_d of 10 nM in a saturation binding assay (SI Appendix, Fig. S3B). Consistent with their activities in the cell-based ROR γ reporter assay, the ROR γ -active oxysterols effectively competed with ³H-25-OHC for binding to recombinant ROR γ LBD (Fig. 1A). The same rank of potency for the four 27-OHCs was observed as that determined using the ThermoFluor readout, with 7 β , 27-OHC being most potent in competing with ³H-25-OHC for binding to ROR γ LBD. Moreover, we found that ROR γ -active oxysterols (i.e., 7 β , 27-OHC and 7 α , 27-OHC but much less for 7 α , 25-OHC) enhanced binding of ROR γ LBD to NCOA3 while decreasing binding to NCORNR (Fig. 1B) using surface plasmon resonance as readout.

7 β , 27-OHC and 7 α , 27-OHC Induce IL-17 Production by Mouse and Human Th17 Cells. Addition of 7 β , 27-OHC and 7 α , 27-OHC but not 7 α , 25-OHC (Fig. 2A and SI Appendix, Figs. S4A and S5) to

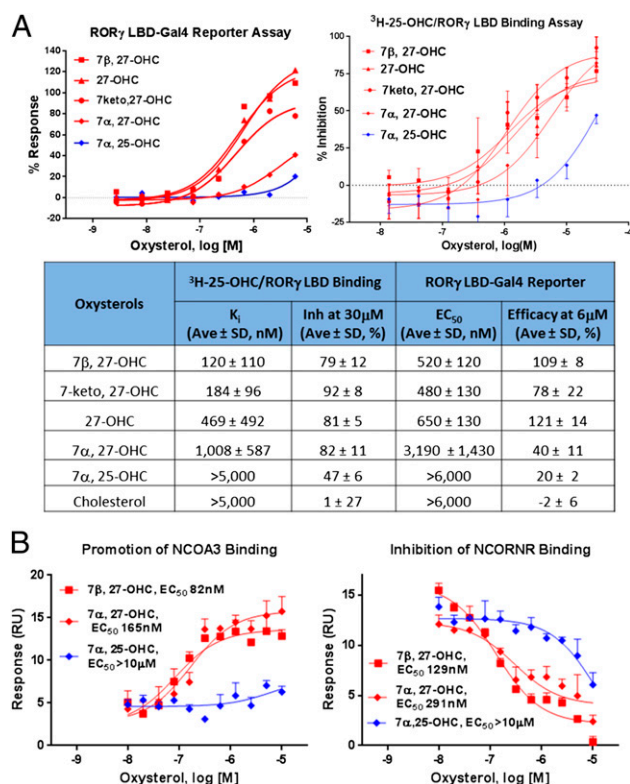


Fig. 1. Agonist activity of oxysterols in a cell-based ROR γ reporter assay and direct binding to ROR γ LBD. (A) Effect of 27-OHCs in reversing the inhibitory effect of UA in a cell-based chimeric ROR γ reporter assay and inhibition of binding of ³H-25-OHC to the ROR γ ligand binding domain. Shown in Upper are dose titration curves from representative experiments. For the reporter assay, oxysterols (starting at 6 μ M; 1:3 serial dilutions) were tested in duplicates in the presence (0% of response) or absence (100% of response) of 1 μ M UA. For the ³H-25-OHC/ROR γ LBD competition binding assay, oxysterols (starting at 30 μ M; 1:3 serial dilutions) were also tested in duplicates. Percent inhibition values were calculated based on without ROR γ LBD protein as 100% and DMSO only as 0%. EC_{50} or IC_{50} values were calculated using GraphPad Prism 5. K_i values were calculated based on $K_i = IC_{50}/([^3H-25-OHC]/K_d + 1)$. Average \pm SD ($n = 3$) of EC_{50} or K_i values and efficacy values at the highest tested concentration are shown in the table. (B) Effect of 7 β /27-OHC vs. 7 α , 25-OHC on binding of coactivator NCOA3 and corepressor NCORNR peptides by ROR γ LBD in a surface plasmon resonance binding assay.

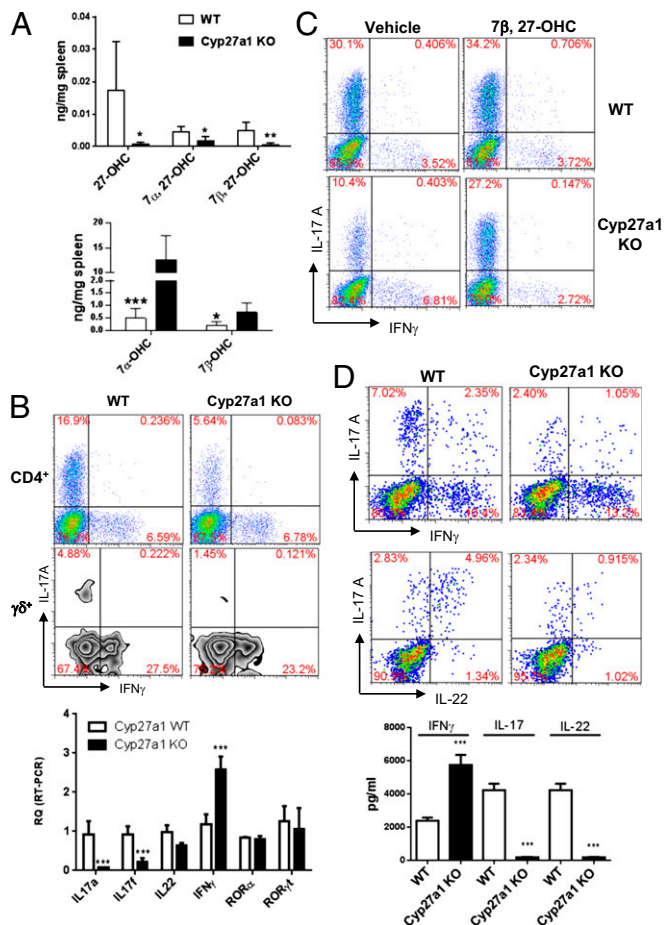


Fig. 3. Defective production of 27-OHCs and IL-17-producing cells in *Cyp27a1* KO mice. (A) Spleen levels of 27-OHCs and 7-OHCs of *Cyp27a1* KO vs. WT mice. Oxysterols in spleen samples ($n = 5$) were measured using deuterated oxysterols as internal controls as described in [SI Appendix](#). (B) Flow cytometry intracellular staining analysis for IL-17A and IFN- γ in spleen CD4 $^{+}$ and $\gamma\delta^{+}$ T cells from *Cyp27a1* KO vs. WT mice. Total splenocytes were activated under Th17-polarizing condition for 3 d. RT-PCR analyses were performed on selected IL-17 pathway genes from mRNA samples ($n = 2-3$). RQ, relative quantification. (C) Defective in vitro Th17 differentiation of naive CD4 $^{+}$ T cells from *Cyp27a1* KO mice. Purified naive CD4 $^{+}$ T cells of *Cyp27a1* KO and WT controls were activated under Th17-polarizing condition for 6 d. DMSO (vehicle) or 7β , 27-OHC (6 μ M) was added at the beginning of the culture. (D) Defective in vivo Th17 differentiation in *Cyp27a1* KO mice. *Cyp27a1* KO and WT mice were immunized with ovalbumin/complete Freund's adjuvant (OVA/CFA); 7 d later, draining lymph node (LN) cells were harvested and restimulated with OVA in vitro for 2 d. IL-17A, IFN- γ , and IL-22 production from cultures of draining LN cells were determined by flow cytometry intracellular staining analysis and ELISA. Statistics by two-tailed, unpaired Student t test. * $P < 0.05$; ** $P < 0.01$; *** $P < 0.001$.

were, in general, much weaker agonists. This observation may be because of conversion of 7-OHCs to 7, 27-OHCs by T cells in culture, because neither γ -OHC nor 7α -OHC enhanced IL-17 production in CYP27A1-deficient CD4⁺ T cells (*SI Appendix, Fig. S12B*). To directly test this hypothesis, we investigated whether mouse Th17 cells could convert exogenously added ^3H - 7α -OHC or γ -OHC into 7, 27-OHCs in culture. Mouse Th17 cells but not Th1 cells significantly converted ^3H - γ -OHC into more hydrophilic ones (Fig. 5A). Similarly, ^3H - 7α -OHC was converted by Th17 cells to a lower extent but not by Th1 cells. HPLC analysis of the product from ^3H - γ -OHC resulted in its separation into peaks A and B (Fig. 5B). Peak B is likely γ , 27-OHC, because it was absent from the converted product of cultured *Cyp27a1* KO CD4⁺ T cells; also, its identity was

confirmed by MS analysis. Peak A was complex as suggested by MS studies. Its generation seems to be partially dependent on CYP27A1: this peak was reduced but not absent from the converted product in *Cyp27a1* KO CD4⁺ T cells. Similar products were observed when ³H-7β-OHC was added to cultures of COS cells transiently expressing human CYP27A1 (Fig. 5C). These findings suggest that mouse Th17 but not Th1 cells preferentially produce 7β, 27-OHC from exogenously added 7β-OHC. To a lesser extent, they likely produced 7α, 27-OHC from exogenously added 7α-OHC in a CYP27A1-dependent manner. Interestingly, the expression of *Cyp27a1* was enriched in mouse Th17 cells, whereas the expression of *Ch25h* was higher in mouse Th1 cells (*SI Appendix, Fig. S13A*). Expected from this observation, higher levels of endogenous 25-OHCs were observed in vitro differentiated Th1 cells (~50% IFN-γ⁺ at harvest) (*SI Appendix, Fig. S13B*). Among 27-OHCs, in vitro differentiated Th17 cells (~30% IL-17⁺ at harvest) produced significantly higher levels of 7β, 27-OHC relative to Th1 cells. No apparent difference in the levels of 7α, 27-OHC and higher levels of 27-OHC were seen in Th1 cells. Whether the lower level of 27-OHC in Th17 was caused by conversion to 7β, 27-OHC is unclear, although neither Th17 nor Th1 cells showed much capacity for conversion of ³H-27-OHC in vitro (*SI Appendix, Fig. S14*). Taken together, our findings suggest that Th17 cells have higher capacity to produce 7β, 27-OHC, although it is less clear whether the same applies for 7α, 27-OHC.

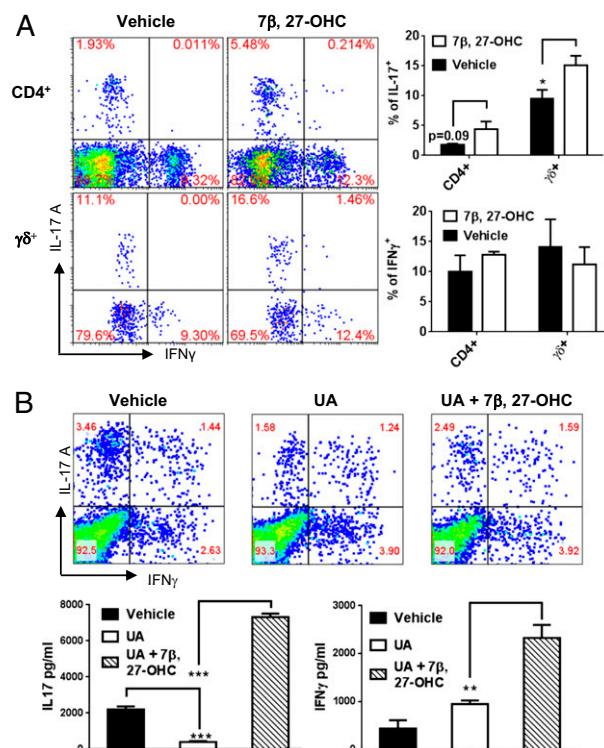


Fig. 4. Dosing of $\gamma\beta$, 27-OHC in mice enhanced IL-17 production in vivo. (A) Effect on LN IL-17⁺ CD4⁺ and $\gamma\delta$ ⁺ T cells. B6 mice ($n = 3$) were dosed with vehicle or $\gamma\beta$, 27-OHC [60 mg/kg s.c. bis in die (BID)] for 3 d. Inguinal lymph node (LN) cells were harvested and stimulated with phorbol myristate acetate/ionomycin for 4 h before intracellular staining. Data in the bar graph are averages \pm SEMs of three mice per group. (B) Effect on IL-17 production in the OVA/CFA priming model. Three groups of B6 WT mice ($n = 3$) were immunized with OVA/CFA and dosed with vehicle, UA, (150 mg/kg i.p. every other day), or UA + $\gamma\beta$, 27-OHC (60 mg/kg s.c. BID); 7 d later, draining LN cells were harvested and restimulated with OVA in vitro for 2 d. IL-17A and IFN- γ production from cultures of draining LN cells were determined by flow cytometry intracellular staining analysis and ELISA. Data in bar graph are averages \pm SEMs. Statistics by two-tailed, unpaired Student *t* test. * $P < 0.05$; ** $P < 0.01$; *** $P < 0.001$.

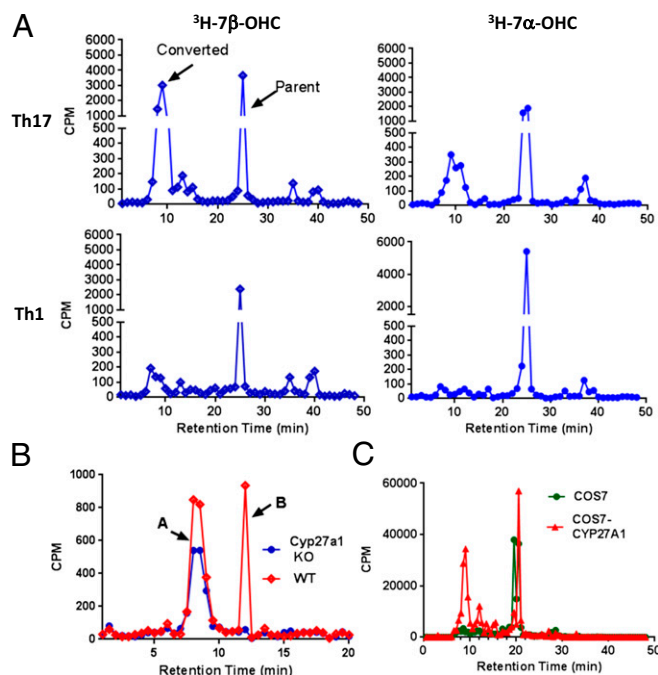


Fig. 5. ^3H - $7\beta/\alpha$ -OHC were preferentially converted into ^3H - $7\beta/\alpha$, 27-OHC in mouse Th17 but minimally in Th1 T cells. (A) Conversion of ^3H - 7β -OHC or ^3H - 7α -OHC by mouse Th17 and Th1 cells in culture (purified total CD4^+ T cells activated under Th17 or Th1 conditions for 3 d). Conditioned media were fractionated using HPLC as described in *SI Appendix*. Radio activities of 50 μL 1-min fractions were determined. Data shown were from a representative experiment of three independent experiments. (B) Conversion of ^3H - 7β -OHC by cultured Th17 cells differentiated from WT and *Cyp27a1* KO naive CD4^+ T cells. Conditioned media were purified as above. Fractions containing the converted products were pooled and further purified by another HPLC; 30-s fractions were collected to resolve converted species, which resulted in peaks A and B. MS analysis revealed that peak B was consistent with being 7β , 27-OHC. (C) HPLC fractionation of conditioned medium from ^3H - 7β -OHC-spiked culture of untransfected COS7 cells or COS7 cells transiently expressing human CYP27A1 protein. CPM, counts per minute.

Discussion

We have shown that 7β , 27-OHC and 7α , 27-OHC are ROR γ t agonist ligands and provided evidence suggesting their role as endogenous drivers of Th17 differentiation. Both bound to recombinant ROR γ LBD protein directly and were selective against other nuclear receptors tested. They were active in enhancing ROR γ t-dependent reporter expression and IL-17 production in primary mouse and human Th17 cells in vitro. In vivo dosing of 7β , 27-OHC in mice increased the number of IL-17 $^+$ cells. Production of these two oxysterols in vivo requires CYP27A1, a key bile acid synthesis enzyme. Mice deficient in CYP27A1 showed reduced levels of 7β , 27-OHC and 7α , 27-OHC and importantly, decrease in IL-17-producing cells. Thus, naive *Cyp27a1* KO CD4^+ T cells are defective in their ability to differentiate into Th17 cells in vitro, which could be rescued by addition of 7β , 27-OHC. In addition, in vivo antigen-specific generation of Th17 cells was also defective in CYP27A1-deficient mice in a priming model. The overall profile of 7α , 27-OHC seemed to be less favorable than that of 7β , 27-OHC; however, we cannot exclude it as a potential candidate endogenous ROR γ t agonist ligand.

Although a deficiency in IL-17-producing cells was consistently observed in younger *Cyp27a1* KO mice (<8 wk), one intriguing observation was that, as these KO mice age, the deficiency became less obvious. Some animals even seemed to develop a phenotype resembling autoinflammation (e.g., splenomegaly was observed on necropsy, and production of cytokines, including IL-17A and IFN- γ , increased). When tested in immune disease models of longer duration, such as collagen-induced arthritis, our

preliminary data indicated no difference between WT and *Cyp27a1* KO mice or even some exacerbation in KO mice. Although these observations seem to contradict our finding that these mice were deficient in IL-17 production, it is worth noting that certain IL-17- or IL-22-producing innate lymphocytes might play a regulatory role, and dysregulation of those cells could account, in part, for the apparent autoinflammatory phenotype in aged KO animals. Nonetheless, the observed IL-17 production in aged CYP27A1-deficient mice suggests the existence of endogenous ROR γ t agonists other than 27-OHCs, exogenous ROR γ t ligands, or ROR γ t-independent mechanisms. These alternative hypotheses require additional investigation.

Endogenous levels of 7β , 27-OHC and maybe 7α , 27-OHC seem to be in the same ranges of binding affinity values determined using the ^3H -25-OHC/ROR γ LBD competitive binding assay. Based on our LC/MS/MS studies, the levels of endogenous 7β , 27-OHC and 7α , 27-OHC in total spleen or in vitro-generated Th17 cells are in single to double digit nanomolar ranges (~ 3 –12 nM for 7β , 27-OHC and 1–35 nM for 7α , 27-OHC). In the ^3H -25-OHC/ROR γ LBD competitive binding assay, 7β , 27-OHC showed higher affinity ($K_i = 120$ nM) and 7α , 27-OHC showed lower affinity ($K_i = 1,008$ nM) than unlabeled 25-OHC ($K_i = 280$ nM) (Fig. 1A). Based on the K_d value of 10 nM for ^3H -25-OHC in this assay (*SI Appendix*, Fig. S3B), we estimate that K_d values of 7β , 27-OHC and 7α , 27-OHC are ~ 4 and ~ 36 nM, respectively. Additional studies will be needed to confirm this estimation.

We have further showed that cultured mouse Th17 cells but not Th1 cells showed enhanced capability to convert exogenously added 7β -OHC or 7α -OHC to 7, 27-OHCs in culture. This finding explained why 7-OHCs were also active in reversing the inhibitory effect of UA in mouse T cells and indicates that T cells themselves can synthesize these endogenous ROR γ t agonists. Addition of ^3H -labeled 27-OHC to Th17 or Th1 cells did not support generation of 7β , 27-OHC or 7α , 27-OHC from exogenously added 27-OHC (*SI Appendix*, Fig. S13). It may be due to low expression of CYP7B1 in mouse T cells, which is the enzyme required for converting 27-OHC into 7α , 27-OHC.

Because of its high cytotoxicity in primary mouse and human T cells in culture, we could not show whether 27-OHC had any agonist activity in enhancing IL-17 production. However, there are a number of findings suggesting that 27-OHC might not be an endogenous ROR γ t agonist. In the presence of SRC1, 27-OHC bound to ROR γ LBD with reduced affinity. In addition, 27-OHC also showed partial agonist activity against LXR α , LXR β , and ER β in cell-based reporter assays. We cannot rule out that 27-OHC may still have a role in modulating ROR γ t function in cells because of binding ROR γ LBD in the absence of SRC1 with relatively high affinity and because of its abundance at steady state.

Among other naturally occurring oxysterols, we showed that 25-OHCs are unlikely endogenous agonists for ROR γ t. First, they were less potent or efficacious than their 27-hydroxylated derivatives in the ROR γ /t reporter assays and the ROR γ LBD binding assay. Second, none of the 25-OHCs were active in enhancing IL-17 production in T cells. Furthermore, in *Cyp27a1* KO mice, which had decreased IL-17-producing cells, the levels of all 25-OHCs were increased. Finally, in *Ch25h* KO mice, in which 25-OHCs production was defective, normal or slightly increased numbers of IL-17-producing cells were observed.

After 27-OHC, 7α -OHC and 24(S)-OHC are the next two most abundant oxysterols in vivo. In the ROR γ reporter assay, both seemed to show partial efficacy at highest nontoxic concentrations compared with the efficacy observed for 27-OHCs. 7α -OHC failed to show significant agonist activity in enhancing IL-17 production in *Cyp27a1* KO CD4^+ T cells, and its level was highly increased in *Cyp27a1* KO mice, suggesting that it is not an ROR γ t agonist in vivo. Both 7α -OHC and 24(S)-OHC were previously reported by others (15, 16) to function as inverse agonists for ROR γ . It is possible that they may play a role in regulating the effect of full endogenous agonists, such as 7β , 27-OHC

and 7α , 27-OHC, by competing for binding to ROR γ or ROR γ t. We hypothesize that the balance between the abundant oxysterols, such as 27-OHC, 7α -OHC, and 24(S)-OHC, and the low levels of 7β , 27-OHC and 7α , 27-OHC at steady state maintains the homeostatic state of IL-17-producing innate cells. However, T-cell activation under Th17 conditions may favor generation of high levels of 7β , 27-OHC and 7α , 27-OHC, which serve as ROR γ t agonists to program Th17 differentiation and drive IL-17 production.

Nonetheless, we must point out that, although our data support that 7β , 27-OHC and 7α , 27-OHC are strong candidates as endogenous ROR γ t ligands, additional investigations are needed in many aspects. Because of the observations associated with aged *Cyp27a1* KO mice and the labile nature of oxysterols, such as 7β , 27-OHC, it has been difficult to explore these tools in more relevant in vivo IL-17-dependent models, such as experimental allergic encephalomyelitis or susceptibility to *Citrobacter* infection. Moreover, we cannot exclude the possibility that endogenous molecules other than the oxysterols that we studied or pathogen-derived factors may also act as exogenous ROR γ t ligands in activating or modulating ROR γ t in IL-17-producing cells. In addition, although we have focused on examining effects of oxysterols in ROR γ t-dependent functions in IL-17-producing cells, we also need to point out that the same oxysterols may serve as endogenous ROR γ agonists. To this end, it will be of interest to investigate whether some of the abnormalities observed in CYP27A1-deficient humans or mice might be partially caused by ROR γ functional defects.

Materials and Methods

Reagents. Oxysterols, antibodies, cytokines, and other chemicals are described in [SI Appendix](#).

Recombinant LBD of ROR γ , ROR α , and ROR β . The generation of the LBD proteins of human RORs are described in [SI Appendix](#).

ROR γ - or ROR γ t-Dependent Luciferase Reporter Assays. Cell-based reporter assays were performed by transiently transfecting HEK293T cells as described in [SI Appendix](#).

Nuclear Receptor Reporter Assays for Selectivity. Cell-based reporter assay kits were purchased from Indigo Biosciences, Inc. or Invitrogen and performed according to the manufacturer's instructions as described in [SI Appendix](#).

3 H-25-OH/ROR γ LBD Scintillation Proximity Assay. 3 H-25-OHC was purchased from Perkin-Elmer. The homogenous scintillation proximity assay was run as described in [SI Appendix](#).

ROR γ LBD Surface Plasmon Resonance Assay. Biacore studies were performed by Biosensor Tools as described in [SI Appendix](#).

ThermoFluor Binding Assay. ThermoFluor experiments were carried out as described in [SI Appendix](#).

Synthesis of $7\alpha/\beta$ /keto, 25-OHC, and 7β , 27-OHC. Detailed synthesis methods are described in [SI Appendix](#).

Extraction and Quantification of Oxysterols with LC/MS/MS. Detailed methods are described in [SI Appendix](#).

Mice. *Ch25h* KO mice, *Cyp27a1* KO mice, ROR γ t KO mice, and C57BL/6J control mice, as described in [SI Appendix](#), were purchased from The Jackson Laboratory.

Mouse Th17 Cell Activation and Differentiation in Vitro. Isolation and in vitro Th17 differentiation of total or naïve mouse CD4 $^+$ T cells were carried as described in [SI Appendix](#).

Gene Expression Studies in Mouse Splenocytes or Th17 Cells. RNA extraction using the RNeasy Plus Mini Kit (Qiagen) and real-time RT-PCR using the commercial Taqman RNA-to-Ct 1-Step Kit (Life Technologies) were carried out as described in [SI Appendix](#).

Conversion of 3 H- 7α -OHC, 3 H- 7β -OHC, or 3 H-27-OHC by Mouse Th1 or Th17 Cells in Culture. 3 H- 7α -OHC or 7β -OHC (American Radiolabeled Chemicals) or 3 H-27-OHC (synthesized by Quotient BioResearch) was added to in vitro differentiated mouse Th1 or Th17 cells, and conditioned media were collected and analyzed by HPLC as described in [SI Appendix](#).

Conversion of 3 H- 7α -OHC or 3 H- 7β -OHC to Dihydroxycholesterols by COS7 Cells Transiently Expressing Recombinant Human CYP27A1 Protein. COS7 cells transiently transfected with human CYP27A1 cDNA in pcDNA3.1 $^+$ (Invitrogen) or vector alone were spiked with 5 μ M 3 H- 7α -OHC or 3 H- 7β -OHC. Conditioned media of 24-h culture were collected and analyzed by HPLC described in [SI Appendix](#).

Human Th17 Cell Activation and Differentiation in Vitro. Total or naïve CD4 $^+$ T cells were isolated from peripheral blood mononuclear cells of healthy donor using the CD4 T-Cell Isolation Kit II or the Naïve CD4 T Cells Isolation Kit II, respectively (Miltenyi Biotec). Th17 cell differentiation in culture was carried out as described in [SI Appendix](#).

In Vivo Generation of Antigen-Specific Th17 Cells. Induction of Th17 cells in vivo and restimulation of draining lymph nodes cells in vitro are described in [SI Appendix](#).

ACKNOWLEDGMENTS. We thank Drs. Jonathan Sprent, David A. Fox, Timothy W. Lovenberg, Michael Jackson, and Anish Suri for their critical reading of this manuscript and valuable suggestions and James McDuffie, Antonio Guy, Jennifer Vegas, Brian Scott, and Leslie Nguyen for running the pharmacokinetics study and analysis of 7β , 27-OHC in mice.

- Solt LA, et al. (2011) Suppression of TH17 differentiation and autoimmunity by a synthetic ROR ligand. *Nature* 472(7344):491–494.
- Huh JR, et al. (2011) Digoxin and its derivatives suppress TH17 cell differentiation by antagonizing ROR γ t activity. *Nature* 472(7344):486–490.
- Xu T, et al. (2011) Ursolic acid suppresses interleukin-17 (IL-17) production by selectively antagonizing the function of ROR γ t protein. *J Biol Chem* 286(26):22707–22710.
- Jin L, et al. (2010) Structural basis for hydroxycholesterols as natural ligands of orphan nuclear receptor ROR γ . *Mol Endocrinol* 24(5):923–929.
- Sawa S, et al. (2010) Lineage relationship analysis of ROR γ t $^+$ innate lymphoid cells. *Science* 330(6004):665–669.
- Schroepfer GJ, Jr (2000) Oxysterols: Modulators of cholesterol metabolism and other processes. *Physiol Rev* 80(1):361–554.
- Olkonen VM, Béaslas O, Nissilä E (2012) Oxysterols and their cellular effectors. *Bio-molecules* 2(1):76–103.
- Sottero B, Gamba P, Gargiulo S, Leonarduzzi G, Poli G (2009) Cholesterol oxidation products and disease: An emerging topic of interest in medicinal chemistry. *Curr Med Chem* 16(6):685–705.
- Poli G, Biasi F, Leonarduzzi G (2013) Oxysterols in the pathogenesis of major chronic diseases. *Redox Biol* 1(1):125–130.
- Liu C, et al. (2011) Oxysterols direct B-cell migration through EBI2. *Nature* 475(7357):519–523.
- Hannedouché S, et al. (2011) Oxysterols direct immune cell migration via EBI2. *Nature* 475(7357):524–527.
- Janowski BA, et al. (1999) Structural requirements of ligands for the oxysterol liver X receptors LXRA and LXR β . *Proc Natl Acad Sci USA* 96(1):266–271.
- Umetani M, Shaul PW (2011) 27-Hydroxycholesterol: The first identified endogenous SERM. *Trends Endocrinol Metab* 22(4):130–135.
- Nachtergaele S, et al. (2012) Oxysterols are allosteric activators of the oncoprotein Smoothened. *Nat Chem Biol* 8(2):211–220.
- Wang Y, et al. (2010) Modulation of retinoic acid receptor-related orphan receptor alpha and gamma activity by 7-oxygenated sterol ligands. *J Biol Chem* 285(7):5013–5025.
- Wang Y, Kumar N, Crumbley C, Griffin PR, Burris TP (2010) A second class of nuclear receptors for oxysterols: Regulation of ROR α and ROR γ activity by 24S-hydroxycholesterol (cerebrosterol). *Biochim Biophys Acta* 1801(8):917–923.
- Pantoliano MW, et al. (2001) High-density miniaturized thermal shift assays as a general strategy for drug discovery. *J Biomol Screen* 6(6):429–440.
- Matulis D, Kranz JK, Salemme FR, Todd MJ (2005) Thermodynamic stability of carbonic anhydrase: Measurements of binding affinity and stoichiometry using ThermoFluor. *Biochemistry* 44(13):5258–5266.
- Clemente JC, et al. (2012) Screening and characterization of human monoglyceride lipase active site inhibitors using orthogonal binding and functional assays. *J Biomol Screen* 17(5):629–640.
- Cali JJ, Russell DW (1991) Characterization of human sterol 27-hydroxylase. A mitochondrial cytochrome P-450 that catalyzes multiple oxidation reaction in bile acid biosynthesis. *J Biol Chem* 266(12):7774–7778.
- Brown AJ, Watts GF, Burnett JR, Dean RT, Jessup W (2000) Sterol 27-hydroxylase acts on 7-ketocholesterol in human atherosclerotic lesions and macrophages in culture. *J Biol Chem* 275(36):27627–27633.
- Honda A, et al. (2001) Differences in hepatic levels of intermediates in bile acid biosynthesis between *Cyp27* $^{-/-}$ mice and CTX. *J Lipid Res* 42(2):291–300.