Bioorganic & Medicinal Chemistry Letters xxx (2014) xxx-xxx

Contents lists available at ScienceDirect



Bioorganic & Medicinal Chemistry Letters

journal homepage: www.elsevier.com/locate/bmcl

Identification of the first potent, selective and bioavailable PPAR α antagonist

Yalda Bravo^{*}, Christopher S. Baccei, Alex Broadhead, Richard Bundey, Austin Chen, Ryan Clark, Lucia Correa, Jason D. Jacintho, Daniel S. Lorrain, Davorka Messmer, Karin Stebbins, Peppi Prasit, Nicholas Stock

Inception Sciences, 5871 Oberlin Drive Suite 100, San Diego, CA 92121, USA

ARTICLE INFO

Article history: Received 20 January 2014 Revised 25 March 2014 Accepted 26 March 2014 Available online xxxx

Keywords: Antagonist PPAR alpha Nuclear hormone receptor Fatty acid oxidation Cancer

ABSTRACT

The discovery and SAR of a novel series of potent and selective PPARα antagonists are herein described. Exploration of replacements for the labile acyl sulfonamide linker led to a biaryl sulfonamide series of which compound **33** proved to be suitable for further profiling in vivo. Compound **33** demonstrated excellent potency, selectivity against other nuclear hormone receptors, and good pharmacokinetics in mouse. © 2014 Elsevier Ltd. All rights reserved.

Cancer cells are known to exhibit distinct characteristics from their 'normal' counterparts with regards to their energy requirements and metabolism. This divergence in turn, presents researchers with a variety of approaches to selectively apply metabolic stress on cancerous cells without adversely affecting healthy cells. Warburg observed that most cancer cells are programmed to increase glucose uptake in order to provide the necessary energy for their proliferative processes. Subsequently, much of the research into aberrant cancer metabolism has, thus far, been focused on targeting the glycolysis pathway, leaving the other metabolic pathways largely unexplored.¹ Recently, the contribution of fatty acid oxidation (FAO) to cancer cell function has gained more attention from the research community.^{2–4} It has been shown that there are specific cancer cell types; including prostate, ovarian and renal cell carcinoma, that are more reliant on fatty acids to satisfy their metabolic needs.⁵ It has also been demonstrated that following detachment of cancer cells from their extracellular matrix (i.e., the initial step of cancer metastasis), a metabolic switch towards increased fatty acid utilization for ATP generation can be seen even in the most glycolytic cancer cell types.² Finally, new research suggests that leukemia-initiating cancer (LIC) stem cells may be reliant on FAO for their maintenance and function, hinting at the

possibility of eradicating leukemia through the exhaustion of the chemo-resistant LIC pool via inhibition of FAO.⁶

Peroxisome proliferator-activated receptors (PPARs) are a group of DNA-binding transcription factors within which three distinct isoforms (i.e., α , β/δ and γ) have been identified.⁷ When activated by their respective endogenous ligands such as free fatty acids and eicosanoids, PPARs undergo a series of conformational changes to accommodate the recruitment of a suitable scaffolding co-activator protein and to facilitate their hetero-dimerization with 9-cis retinoic acid receptor (RXR). The resulting ternary complex then binds to the appropriate peroxisome proliferator response element (PPRE) on DNA and initiates the transcription and expression of its target genes. PPAR α is implicated primarily in the regulation of lipid metabolism and as such, its activation leads to an increase in uptake and catabolism of fatty acids. This phenotype is achieved via the up-regulation of genes involved in the binding and transport of fatty acids (i.e., CPT1/2, CACT, and others) for oxidative processing. Consequently, it follows that antagonism of the PPARa receptor, coupled with the knowledge that certain malignant cells rely on FAO, represents a novel paradigm to stop the proliferative and metastatic tendencies of these cancer cells.

Although numerous examples of potent and selective PPAR α agonists can be found in the literature; including those that have been approved clinically for the treatment of hypercholesterolemia, hypertriglyceridemia and other related metabolic disorders,⁸⁻¹² there are only a few isolated reports of confirmed

^{*} Corresponding author. Tel.: +1 858 354 9160. E-mail address: yaldabravo47@gmail.com (Y. Bravo).

Y. Bravo et al./Bioorg. Med. Chem. Lett. xxx (2014) xxx-xxx



Figure 1. Previously reported PPARa modulators.¹⁵

PPARα antagonists. These inhibitors (compounds **1** and **2**, Fig. 1) were reported to be discovered serendipitously, following independent SAR campaigns carried out on their respective PPARα agonist precursors (i.e., compounds **3** and **4**). The reported initial program goal was to identify a suitable replacement for their shared carboxylic acid warhead (highlighted in red). In both cases, replacement of the carboxylic acid group by either an inverse amide (i.e., GW6471, compound **1**) or an acyl sulfonamide (i.e., compound **2**) triggered an unanticipated agonist-to-antagonist switch that resulted from the lifting of the PPARα C-terminal AF-2 helix, which, in turn, favored the recruitment of a co-repressor peptide such as SMRT.

Both antagonists were found to dose-dependently inhibit the activation of PPARa-driven luciferase expression by GW7647 (a known PPAR α agonist) in a cell-based functional assay¹³ we employed to drive our SAR. Unfortunately, neither compound proved suitable for assessing whether the antagonism of PPAR α would be efficacious in our murine cancer models. Specifically, when mice were orally administered compound 1, no measurable drug levels could be detected in the mouse plasma regardless of when the blood was collected after dose. Although this lack of oral drug exposure could be mitigated via an alternative mode of administration (i.e., intra-peritoneal injection), compound 1's lack of solubility in conventional dosing vehicles necessitated that the compound be eventually formulated in neat DMSO. However, this vehicle choice complicated the interpretation of the resulting in vivo data. The DMSO vehicle arm was found to exhibit a significant anti-metastatic effect when compared to conventional vehicles such as 0.5% aqueous methocel or saline (data not shown).¹⁴ While compound 2 was found to possess physicochemical properties amenable for conventional formulation, we quickly discovered that this compound was very unstable in murine plasma and underwent an almost instantaneous enzyme-mediated hydrolytic cleavage to the acid; a potent PPARa agonist. As a result, we initiated separate SAR campaigns on both of these scaffolds with the aim of addressing their respective key liabilities that prohibited their profiling in vivo. This manuscript will focus on our efforts at improving the metabolic stability of compound **2** towards hydrolysis, while preserving its potency and selectivity against PPARα. Our effort around compound **1** will be disclosed separately in due course.

It was initially hypothesized that modification of the steric and/ or the electronic environment adjacent to the cleavage site in compound **2** might afford compounds that are more resistant towards hydrolysis. In this regard, we first evaluated the impact of the sulfonyl substituent on both stability in murine plasma (after a Table 1SAR of acyl sulfonamides



Compound	Х	R:	PPARa	% Parent compound remaining ^b			
			IC ₅₀ ª (nM)	30 min	60 min	24 h	
2	0	Ph	2.7 ± 0.9	6.7 ± 0.6	0.7 ± 0.1		
8	0	4-Me-Ph	2.8 ± 1.9	73.1 ± 1.4	63.7 ± 3.4		
9	0	4- ⁱ Pr-Ph	13 ± 0	89.5 ± 2.0	81.9 ± 3.3	2.7 ± 0.1	
10	0	4-CF ₃ -Ph	34 ± 18	102.7 ± 1.9	105.3 ± 3.1	53.7 ± 3.5	
11	0	3-Pyridyl	80 ± 18	82.4 ± 2.6	75.2 ± 1.1	0.2 ± 0.1	
12	0	Cyclohexyl	3 ± 0	60.5 ± 2.0	43.6 ± 2.0		
13	0	Cyclopentyl	4.6 ± 1.0	21.0 ± 0.8	7.0 ± 0.3		
14	0	Cyclopropyl	6.6 ± 3.0	1.7 ± 0.2	0.1 ± 0.04		
15	0	3-Furan	78 ± 0	75.4 ± 1.0	66.6 ± 2.0		
16	0	4-Pyran	42 ± 14	86.9 ± 2.9	86.5 ± 3.3		
17	С	Ph	60 ± 10	80.0 ± 2.8	62.3 ± 2.5		

^a Values are the mean of at least three experiments.

^b For experimental procedure see Ref. 20.

30 min and 60 min incubation period) and potency against PPAR α in our luciferase assay (Table 1). These compounds were readily accessed through EDC-mediated condensation of acid **5** and sulfonamide **7** in the presence of DMAP.¹⁶ Sulfonamides that were not commercially available were themselves synthesized from the corresponding sulfonyl chloride and liquid ammonia (Scheme 1).

The incorporation of an aliphatic group such as methyl (8) or isopropyl (9) at the para position of the terminal phenyl ring delivered PPAR α antagonists of comparable potency to compound **2**. Both modifications led to a significant improvement in plasma stability, with the larger isopropyl group being more resistant towards hydrolysis (82% of 9 remained after 60 min of incubation with murine plasma vs 64% of **8**). Unfortunately, the vast majority of acyl sulfonamide 9 (>97%) was still cleaved to PPARa agonist 5 after a 24 h incubation period. Although switching the isopropyl residue in **9** for its known electron-withdrawing isostere CF_3 (**10**) further improved the resulting compound's resistance towards hydrolytic cleavage at a cost of only a small drop in PPAR α potency, compound **10** could not survive the more rigorous 24 h incubation protocol unscathed. Replacement of the terminal benzene ring in compound 2 with a heteroaromatic plate such as 3-pyridyl (11) proved to be highly detrimental for PPARa antagonism. On the other hand, reducing it to the fully saturated cyclohexane (12) was tolerated in terms of potency against PPARa. These observations combine to reveal that the phenyl group in **2** occupies a large, hydrophobic pocket in the PPARa ligand binding domain but is



Scheme 1. Reagents and conditions: (a) liquid NH₃, DCM, -78 °C, 90–95%. (b) EDC, DMAP, DCM, 40–80%.

2

itself not participating in any stabilizing, π -stacking interactions. It is important to stress that while both of these acyl sulfonamides (11 and 12) were significantly more hydrolytically stable than 2, these alterations again failed to completely address compound 2's cleavage liability. Other aliphatic sulfonyl substituents were subsequently explored, including those cyclic (13 and 14) and heterocyclic (i.e., 15 and 16) in nature. As expected, an increase in the steric bulk around the site of cleavage was accompanied by an attendant improvement in plasma stability (compare cyclohexyl 12 > cyclopentyl 13 > cyclopropyl 14, and 4-pyran 16 > 3-furan 15). Analogues containing a heteroatom were consistently more resistant towards enzyme-mediated cleavage, but they were also found to be less effective at antagonizing PPAR α than those without (compare 13 vs 15, and 12 vs 16). Regardless, none of these alterations were able to deliver a potent, hydrolytically-stable PPAR α antagonist.

It has been previously observed that acyl sulfonamides bearing an oxygen atom β to the carbonyl group are much more susceptible

towards enzyme hydrolysis than those without¹⁷ and this property has been exploited for the design of pro-drugs that facilitate the delivery of non-orally bioavailable sulfonamide drugs. The synthesis of acyl sulfonamide 17 was carried out as detailed in Scheme 2. Briefly, palladium-catalyzed carbonylation of aryl bromide 18 was best carried out with triethylsilane as the reducing agent. The resulting benzaldehyde 19 then readily underwent Reformatsky reaction with the zincate generated in situ from commercially available ethyl α -bromoisobutyrate to deliver alcohol 20 in quantitative yield. Subsequent ionic de-oxygenation with triethylsilane and boron trifluoride etherate proceeded without incident and gave diester 21 in 67% yield. Selective saponification of the more sterically accessible methyl ester could be achieved with potassium trimethylsilanoate and the resulting carboxylic acid was then coupled with known hydrazine 22 in the presence of HATU and Hunig's base. When heated with an excess of CSA in ethyl acetate. hvdrazine carboxamide 23 underwent cvclization to afford. after hydrolysis with aqueous lithium hydroxide, triazalone 24. Finally,



Scheme 2. Reagents and conditions: (a) Pd(PPh_3)₂Cl₂, Na₂CO₃, Et₃SiH, CO, DMF, 80 °C, 53%. (b) Zn, I₂, BrC(Me)₂CO₂Et, sonication, rt, >99%. (c) BF₃·OEt₂, Et₃SiH, DCM, 0 °C \rightarrow rt, 67%. (d) KOTMS, THF, rt, 92%. (e) compound **22**, HATU, Hunig's base, DMF, rt, 64%. (f) CSA, EtOAc, 80 °C, 75%. (g) 2N aq LiOH, THF, MeOH, rt, 95%. (h) EDC, DMAP, DCM, 88%.



Scheme 3. Reagents and conditions: (a) CDI, THF, rt then NH₂NH₂·H₂O, rt, 95%. (b) EtN = C=O, DCM, rt, >99%. (c) KOH, MeOH, reflux, 85%. (d) 4-*tert*-butylbenzyl bromide, K₂CO₃, DMF, 45 °C, 93%. (e) Pd(dppf)Cl₂, KOAc, bis(pinacolato)diboron, dioxane, 85 °C, 82%. (f) Ar-Br, K₂CO₃, Pd(PPh₃)₄, DME, H₂O, 85 °C, 65–95%. (g) RSO₂Cl, pyridine, rt, 60–90%.

Please cite this article in press as: Bravo, Y.; et al. Bioorg. Med. Chem. Lett. (2014), http://dx.doi.org/10.1016/j.bmcl.2014.03.090

4

ARTICLE IN PRESS

Y. Bravo et al./Bioorg. Med. Chem. Lett. xxx (2014) xxx-xxx

using the conditions described earlier in Scheme 1, the requisite acyl sulfonamide **17** could be cleanly isolated in 88% yield. Unfortunately, although compound **17** was found to be much more stable towards hydrolysis than compound **2**, the removal of the oxygen linker did not completely protect **17** from enzyme-mediated hydrolysis.

Concurrent with our campaign to stabilize the acyl sulfonamide handle, we evaluated whether this labile motif in compound 2 could be replaced entirely without jeopardizing its PPARa antagonism. It was hypothesized that the fibrate core present in **2** could be effectively mimicked by a six-membered aryl ring.¹⁸ The requisite biaryl sulfonamide analogues were synthesized as shown in Scheme 3. Commercially available butanoic acid 25, after its initial conversion to the corresponding acyl imidazole with CDI, was coupled with hydrazine to give acylhydrazide 26. Subsequent treatment with ethyl isocvanate afforded hydrazine carboxamide **27**. which was then cyclized to triazalone **28** with KOH in refluxing methanol. This material was first alkylated with 4-tert-butylbenzyl bromide and then further transformed to pinacol boronate 29 under standard Suzuki-Miyaura borylation conditions. From this versatile intermediate, a variety of amino-aryl and amino-heteroaryl groups could then be appended via the Suzuki cross-coupling reaction. Finally, the desired biaryl sulfonamides (Table 2, compounds **31–39**) were obtained by treatment of aniline **30** with a selection of commercially available sulfonyl chlorides using pyridine as solvent.

Although replacement of the fibrate linker in compound **2** with a simple benzene spacer (**31**) led to a 200-fold drop in its ability to antagonize PPAR α , moving the sulfonamide group from the *meta* to

Table 2

SAR of biaryl triazalones



Compound	X ¹ / _{IJ} NHSO₂R	PPAR α IC ₅₀ ^a (nM)	$\%$ Parent compound remaining $^{\rm b}$ 24 h	Solubility in pH = 7.4 aqueous buffer (μM)
31	NHSO2Ph	386 ± 250	136 ± 8.4	0.037
32	NHSO ₂ Ph	21 ± 16	124.7 ± 4.6	0.018
33	NHSO ₂ Ph	77 ± 35	140.6 ± 2.4	1.5
34	NHSO ₂ Ph	122 ± 81	128.7 ± 2.3	0.014
35	NHSO ₂ Me	113 ± 63	104.8 ± 3.6	0.423
36	NHSO ₂ Bn	311 ± 66	133.9 ± 2.3	0.012
37	N SO OME	4800 ± 3492	119.2 ± 2.8	0.082
38	NHSO ₂ Ph	114 ± 57	119.4 ± 1.4	0.091
39	NHSO ₂ Ph	218 ± 153	131.2 ± 4.2	4.9

^a Values are the mean of at least three experiments.

^b For experimental procedure see Ref. 20.

^c For experimental procedure see Ref. 21.

Table 3

Pharmacokinetic profiles of compounds **33** and **37**

		33	37
CD1 mouse (I.P.)	Dose Vehicle Plasma AUC (h µg/mL) C _{max} (µM) C _{trough} (µM)	30 mpk Saline 3.4 1.2 1.0	30 mpk 98% saline/2% tween 23 7.2 0.29

the para position (32) recovered much of the initial loss in potency. However, compound **32** was found to be poorly soluble in a variety of aqueous formulations (0.018 μ M in pH = 7.4 aqueous buffer, Table 2) and could only be formulated in PEG400. As a consequence of its poor physiochemical properties, no bioavailability was achieved upon oral administration of compound 32 to rodents. In order to improve the solubility of this series, installation of heteroatoms were explored and 2-pyridyl was found to be optimal in terms of both PPARa potency and aqueous solubility (compare 32 vs 33 vs 34). Installation of the pyridine led to >80-fold increase in aqueous solubility. When increasing the nitrogen count at the terminal aromatic ring further to either a pyrimidine (38) or a pyrazine (39) potency was lost against PPARa. Sulfonamide 33 was also tested to see whether it recruits SMRT corepressor in a similar fashion to GW6471. Indeed, using a commercially available kit where the GST-tagged PPARα ligand binding domain is labeled with the Tb-anti-GST antibody and the SMRT co-repressor peptide is labeled with fluorescein,¹⁹ we witnessed a dose dependent quenching of the fluorescence signal with an EC₅₀ of 1.4μ M.

Y. Bravo et al./Bioorg. Med. Chem. Lett. xxx (2014) xxx-xxx

Table 4

Nuclear hormone receptor counter screen of compounds 33 and 37

	PPAR α antagonist IC ₅₀ ^a (μ M)	PPARα agonist EC ₅₀ ^{a,b} (μM)	PPAR IC ₅₀ ^{a,c} (μ M)	PPAR $\gamma IC_{50}^{a,c}$ (μM)	$ER\beta IC_{50}^{a,c} (\mu M)$	$\text{GR IC}_{50}{}^{\text{a,c}}\left(\mu M\right)$	$TR\beta IC_{50}^{a,c}(\mu M)$
33	0.077 ± 0.035	>100	6.0 ± 3.2	15 ± 19	15.2 ± 6.9	32.5 ± 19.3	>100
37	4.8 ± 3.5		>100	>100	34.8 ± 16.1	32 ± 19	>69

^a Values are the mean of at least three separate experiments.

^b For experimental procedure see Ref. 13.

^c For experimental procedure see Ref. 25.



Figure 2. Plasma FGF21 concentration after dosing of vehicle control versus **33** given at 30 mg/kg (IP) for 4 days * p < 0.05, t test.

With sulfonamide **33** as the reference point, we next examined the impact of the size of the sulfonyl substituent (R in Table 2) on PPAR α potency. Changing from phenyl to either methyl (**35**) or benzyl (**36**) proved to be detrimental. In fact, increasing the steric demand even further to sulfonamide **37** delivered an extremely weak PPAR α antagonist that would be suitable as a structurallysimilar, negative control to compound **33** in our proof-of-concept experiments. Therefore both compounds **33** and **37** were selected to be profiled further in vivo to assess whether sufficient drug concentrations could be achieved in mouse with conventional aqueous formulations (Table 3).

Gratifyingly, when the sodium salt of compound 33 was administered IP as a saline solution at a dose of 10 mpk in mouse, we observed good drug coverage across 24 h (i.e. C_{trough} = 1.0 μM) with a calculated plasma AUC of 3.4 h µg/mL. Furthermore with compound 33, the maximum plasma concentration was achieved 1 h post dose and found to be 1.2 µM. Similarly the sodium salt of compound 37, the proposed negative control to 33, also gave sufficient drug level in mouse using conventional formulation for the necessary in vivo proof-of-concept experiments. Our initial program goal was to design a molecule suitable for assessing whether PPAR α antagonism would be able to slow down or completely stop the proliferation of cancer cells both in vitro and in vivo. Therefore, it was prudent to assess our chosen antagonist and negative control's affinity for other nuclear hormone receptors; including estrogen receptor beta (ER β) and glucocorticoid receptor (GR), both of which have been implicated in the proliferation prostate²², breast²³ and renal²⁴ cancer cells (Table 4). As illustrated, compounds 33 and 37 exhibited minimal affinity for this small panel of nuclear hormone receptors.

Having identified a selective antagonist of the PPAR α nuclear hormone receptor with appreciable exposure after IP dosing we utilized a mouse model to assess compound **33**'s ability to inhibit PPAR α target genes in vivo. *Fgf21* (Fibroblast growth factor 21) is a known PPAR α target gene, induced by either fasting or administration of a fibrate agonist.²⁶ After 4 consecutive days dosing of compound **33** at 30 mg/kg (IP) to fasted mice,²⁷ we measured a significant decrease in plasma FGF21 in the treated animals versus control (Fig. 2) thus successfully demonstrating functional antagonism of the PPAR α receptor in vivo.²⁸

In summary, we have designed and synthesized the first series of potent and selective PPAR α antagonists that are suitable for in vivo proof-of-concept experiments. Optimizing the labile acyl sulfonamide linker in compound **2**, we arrived at biaryl sulfonamide **33**. Understanding of the binding tendencies of this novel series of compounds to PPAR α has facilitated the identification of compound **37** as a structurally similar, negative control. Both of these compounds will serve as invaluable tools to improve our understanding of the relationship between FAO and cancer proliferation and survival. Indeed, these data will be disclosed shortly.²⁹

Acknowledgments

We would like to thank Jill M. Baccei, Brian Stearns and Yen Truong for their contributions and insightful comments.

References and notes

- 1. Vander Heiden, M.; Cantley, C. L.; Thompson, C. B. Science 2009, 324, 1029.
- Schafer, Z. T.; Grassian, A. R.; Song, L.; Jiang, Z.; Gerhart-Hines, Z.; Irie, H. Y.; Gao, S.; Puigserver, P.; Brugge, J. S. Nature 2009, 461, 109.
- Samudio, I.; Harmancey, R.; Fiegl, M.; Kantarjian, H.; Konopleva, M.; Korchin, B.; Kaluarachchi, K.; Bornmann, W.; Duvvuri, S.; Taegtmeyer, H.; Andreeff, M. J. Clin. Invest. 2010, 120, 142.
- 4. Huqi, A. Heart Metab. 2011, 51, 27.
- Caro, P.; Kishan, A. U.; Norberg, E.; Stanley, I. A.; Chapuy, B.; Ficarro, S. B.; Plak, K.; Tondera, D.; Gounarldes, J.; Yin, H.; Zhou, F.; Green, M. R.; Chen, L.; Montl, S.; Marto, J. A.; Shipp, M. A.; Danial, N. N. *Cancer Cell* **2012**, *22*, 547.
- Ito, K.; Carracedo, A.; Weiss, D.; Arai, F.; Alau, U.; Avigan, D. E.; Schafer, Z. T.; Evans, R. M.; Suda, T.; Lee, C. H.; Pandolfi, P. P. *Nature Med.* 2012, *18*, 1350.
- 7. Berger, J.; Moller, D. E. Annu. Rev. Med. 2002, 53, 409. and references cited therein.
- Garcia-Casimiro, A.; Bigge, C. F.; Davis, J. A.; Padaline, T.; Pulaski, J.; Ohren, J. F.; McConnell, P.; Kane, C. D.; Royer, L. J.; Stevens, K. A.; Auerbach, B. J.; Collard, W. T.; McGregor, C.; Fakhoury, S. A.; Schaum, R. P.; Zhou, H. *Bioorg. Med. Chem.* 2008, 16, 4883.
- 9. Li, A. C.; Palinski, W. Annu. Rev. Pharmacol. Toxicol. 2006, 46, 1.
- Guerre-Millo, M.; Gervois, P.; Raspe, E.; Madsen, L.; Poulain, P.; Derudas, B.; Herbert, J.; Winegar, D. A.; Wilson, T. M.; Fruchart, J.; Berge, R. K.; Staels, B. J. Biol. Chem. 2000, 275, 16638.
- 11. Ye, J.; Doyle, P.; Iglesias, M.; Watson, D.; Cooney, G.; Kraegen, E. Diabetes 2001, 50, 411.
- Kobayashi, M.; Shigeta, Y.; Hirata, Y.; Omori, Y.; Sakamoto, N.; Nambu, S.; Baba, S. Diabetes Care 1988, 11, 495.
- 13. The screening of test compounds for activities against human PPAR α receptors were analyzed using a commercial kit, Human PPAR α Reporter Assay System (Indigo Biosciences, Cat. #IB00111). PPAR α Reporter cells were dispensed into 96-well assay plates and test compounds were then immediately added to the wells. For the antagonist assay, PPAR α agonist GW7647 (20 nM) was also added. Following an overnight incubation, the treatment media was discarded and Luciferase Detection Reagent (LDR) added. The luminescence intensity of light emission from the ensuing luciferase reaction is directly proportional to the relative level of PPAR α activation in the reporter cells. Luminescence was read using a Molecular Devices FlexStation 3.
- Lampugnani, M. G.; Pedenovi, M.; Niewiarowski, A.; Casali, B.; Donati, M. B.; Corbascio, G. C.; Marchisio, P. C. *Exp. Cell. Res.* 1987, 172, 385.
- (a) Xu, H. E.; Stanley, T. B.; Montana, V. G.; Lambert, M. H.; Shearer, B. G.; Cobb, J. E.; McKee, D. D.; Galardi, C. M.; Plunket, K. D.; Nolte, R. T.; Parks, D. J.; Moore, J. T.; Kliewer, S. A.; Willson, T. M.; Stimmel, J. B. *Nature* **2002**, *415*, 813; (b) Etgen, G. J.; Mantio, N. *Curr. Top. Med. Chem.* **2003**, *3*, 1649.
- 16. Xu, Y.; Mayhugh, D.; Saeed, A.; Wang, X.; Thompson, R. C.; Dominianni, S. J.; Kauffman, R. F.; Singh, J.; Bean, J. S.; Bensch, W. R.; Barr, R. J.; Osborne, J.;

Montrose-Rafizadeh, C.; Zink, R. W.; Yumibe, N. P.; Huang, N.; Luffer-Atlas, D.; Rungta, D.; Maise, D. E.; Mantlo, N. B. J. Med. Chem. **2003**, 46, 5121.

- 17. Larsen, J. D.; Bundgaard, H. Int. J. Pharm. 1987, 37, 87.
- 18. Analysis of the various known PPARα crystal structures in the RCSB protein data bank containing a fibrate-derived compound in their respective ligand binding domain (i.e. PDB: 1KKQ, ligand = GW6471, PDB: 2P54, ligand = GW735, and PDB: 4BCR, ligand = WY14643) revealed that the distance between the phenyl ring and the carbon in question ranged from 3.8 to 4.1 angstoms. Similarly, using the known biphenyl crystal structure (i.e., PDB: 2XRX), the distance between the phenyl ring on the left and the *meta* and *para* carbons of the phenyl ring on the right are 3.8 and 4.3 angstroms, respectively.



This would suggest that replacement of a fibrate moiety with an aromatic ring is feasible.

- 19. LanthaScreen TR-FRET Coactivator Assays, purchased from Invitrogen (PV4684), is used to identify antagonists of PPAR α following the manufacturer's instructions except that Fluorescein-SMRT ID2 corepressor peptide (Invitrogen, PV4423) was substituted for the Fluorescein-PGC1a peptide provided in the coactivator assay kit. Briefly, increasing concentrations of compound were added to LBD and corepressor peptide solutions with 10 nM of GW7647 (EC₈₀ of the PPAR α agonist, measured in this assay). After a 2 h incubation at room temperature, the 520/490 TR-FRET ratio was measured using a fluorescence microplate reader (Flex Station 3) excitation 340 nm, emission 495 nm, and emission 520 nm.
- 20. A 30 mM DMSO stock solution of test compound is diluted to $100 \,\mu$ M in acetonitrile. 5 μ L of this $100 \,\mu$ M solution is spiked into $495 \,\mu$ L of thawed mouse plasma (Bioreclamation Cat# MSEPLEDTA-F) yielding 1 μ M of test compound in plasma. Plasma vial is vortexed and a 50 μ L aliquot is immediately precipitated in 200 μ L acetonitrile containing 5 ng/mL buspirone. This is time zero (t = 0) test compound concentration. Plasma vial is incubated at 37 °C in water bath and additional 50 μ L plasma aliquots are

taken at 30 and 60 min time points and precipitated in the same fashion as time = 0. Precipitated samples are shaken for 5 min and centrifuged at 4000g for 10 min. 150 μ L of supernatant is collected and analyzed for test compound using LC-MS/MS. Percent test compound remaining is calculated by dividing concentration at specific time point by concentration at time equals zero and then multiplying by 100.

- 21. The test compounds and controls (10 mM in DMSO, 10 μ L/well) were added to a pH = 7.4 buffer (490 μ L/well) on a 96-well plate. The samples were then briefly vortexed for 2 min before it was shaken on an orbital shaker for 24 h at room temperature. 200 μ L of each sample solution was then transferred onto a MultiScreen filter plate equipped with a polycarbonate membrane and filtered using a Millipore vacuum manifold. The filtrate thus obtained was then analyzed and quantified by LCMSMS, using the calibration curve constructed from the corresponding test compound.
- 22. Signoretti, S.; Loda, M. Am. J. Pathol. 2001, 159, 13. and references cited therein.
- 23. Pan, D.; Kocherginksy, M.; Conzen, S. D. Cancer Res. 2011, 71, 6360.
- 24. Mink, K. J.; Jang, J. H.; Lee, J. T.; Choi, K. S.; Kwon, T. K. J. Mol. Med. 2012, 90, 309.
- 25. Screening of test compounds in Human PPARδ, PPARγ and the nuclear hormone receptors, estrogen receptor beta (ERβ), glucocorticoid receptor (GR) and thyroid hormone receptor beta (TRβ) were analyzed using reporter assay system kits from Indigo Biosciences according to the manufacturer's instructions.
- Badman, M. K.; Pissios, P.; Kennedy, A. R.; Koukos, G.; Flier, J. S.; Maratos-Flier, E. Cell Metab. 2007, 5, 426.
- 27. Mice (C57Bl/6) received an IP injection of **33** or vehicle control (saline) daily for four days. Starting on day 2 of antagonist treatment, mice were fasted for 48 h. 4 h after the final dose, blood was collected into EDTA tubes via terminal cardiac puncture. Plasma was isolated and analyzed for FGF21 using an ELISA (R&D Systems, Minneapolis, MN) and **33** concentration using LC/MS methods.
- Stebbins, K. J.; Broadhead, A. R.; Cabrera, G.; Correa, L. D.; Messmer, D.; Bundey, R.; Baccei, C.; Bravo, Y.; Chen, A.; Stock, N. S.; Prasit, P.; Lorrain, D. S. J. Pharm. Exp. Ther. 2014 (manuscript submitted).
- Messmer, D.; Lorrain, K.; Bravo, Y.; Stock, N.; Bundey, R.A.; Correa, L.; Poon, M.; Stebbins, K.; Cabrera, G.; Chen, A.; Jacintho, J.D.; Spaner, D.; Prasit, P.; Lorraine, D. Abstract of Papers, 54th Meeting of the American Society of Hematology, Atlanta, GA, 2011; Abstract 3879.