



Abstract

Vitamin A (retinol) and its metabolites play many physiological roles including cell differentiation, cell proliferation, energy homeostasis, circadian rhythm and immune response. These compounds are known to act through Retinoid Acid Receptors (RARs), retinoid-related Orphan Receptors (RORs) and Retinoid X Receptors (RXRs). These receptors are also important drug targets, although the specificity of many retinoid-like compounds for the retinoid receptors has not been carefully explored due to the lack of robust and selective assays. In the present work we examined 25 retinoid-like compounds for their ability to regulate RAR α , β , γ , RXR α , β , γ and ROR α , β , γ . Compounds were tested for agonistic as well as antagonistic activities towards these receptors. Additionally, cytotoxicity analyses were performed to confirm true antagonistic activity. Many of these compounds show a broad range of receptor activity, while others show moderate selectivity. Only a few of these test compounds show significant antagonistic activity to these retinoid receptors, illustrating the need to screen more broadly for drug candidates with improved specificity in inhibiting the human RXRs, RARs and RORs, as well as selectivity in targeting one receptor sub-type (α , β , or γ) over the others.

Introduction

Retinoic Acid Receptors (RARs), and Retinoid-related Orphan Receptors (RORs) belong to Thyroid hormone receptor-like subfamily, whereas the Retinoid X Receptors (RXRs) comprise their own subfamily. Each of the three receptor groups are comprised of three distinct variants (α , β and γ). Furthermore, each group shows a different mode of action: RARs form heterodimers with RXRs, RORs act as monomers, and RXRs partner with all of the other non-steroid hormone nuclear receptors to regulate gene transcription.

Dysfunctions of these receptors are associated with many disease conditions. For example, RARs are associated with cancer, neurodegenerative diseases, and schizophrenia; RORs are associated with metabolic disorders, inflammation, atherosclerosis and autoimmune diseases; and RXRs are associated with cancer. As such, all three groups of retinoid receptors present important targets for the development of new drugs that, ideally, would be *specific* to individual retinoid receptor groups and *selective* in targeting the activities of either the α , β or γ receptor variants.

The current study analyzes a library of 25 retinoid-like compounds for *specific* and *selective* bioactivities using commercially available cell-based assay systems for human RAR α , β and γ , RXR α , β and γ , and ROR γ (INDIGO Biosciences Inc; www.indigobiosciences.com). ROR α and β assays were developed for this study.

Materials and Methods

Test compounds- A portion of the test compounds analyzed in this study comprise the Retinoid Library offered by Enzo Life Sciences Inc.: Puerarin, ec23, ATRA, 9CRA, 13CRA, 4 OH-phenylretinamide, AM 580, TTNPB, Methoprene acid, Methoprene, Acitretin, 4-OH retinoic acid, RO 41-5253, adapalene, and AC-55469. DHA, EPA, and Ursolic Acid (Sigma-Aldrich Corp.), TO901317, BMS961, ER50891, HX531, HX630, and MM11253 (Tocris Biosciences), and SR1001 (Cayman Chemicals) were also tested.

Retinoid Receptor Assays- Human RAR α , β and γ , RXR α , β and γ , and ROR γ Reporter Assay Systems are commercially available in kit formats from INDIGO Biosciences Inc. INDIGO's Retinoid Receptor Assays utilize division competent proprietary cells engineered to provide constitutive, high-level expression of the respective retinoid receptors, and include the luciferase reporter gene functionally linked to a responsive promoter. Luciferase gene expression occurs after ligand-bound receptor undergoes nuclear translocation, DNA binding, recruitment and assembly of the co-activators and accessory factors required to form a functional transcription complex, culminating in expression of the target gene. The Human ROR α and β assays were developed for this study. The scheme for performing agonist-mode retinoid receptor assays is depicted in **Figure 1**. Antagonist-mode retinoid receptor assays were performed in combination with INDIGO's Live Cell Multiplex (LCM) Assay, as depicted in **Figure 2**.

Figure 1. Agonist-mode Nuclear Receptor assay scheme.

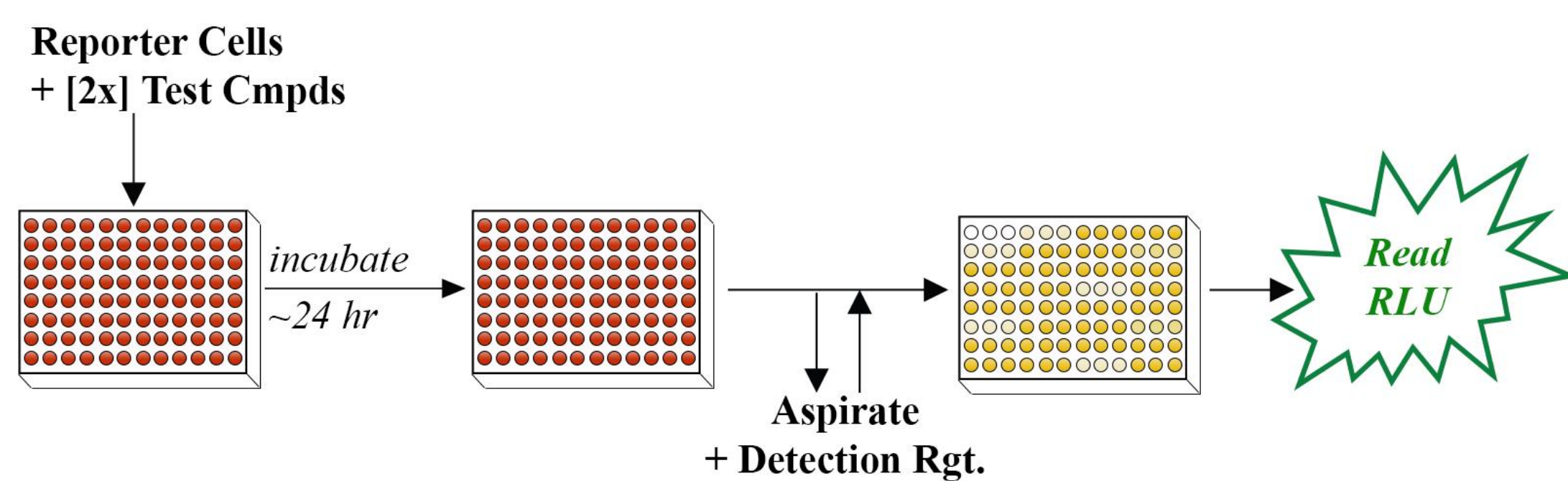
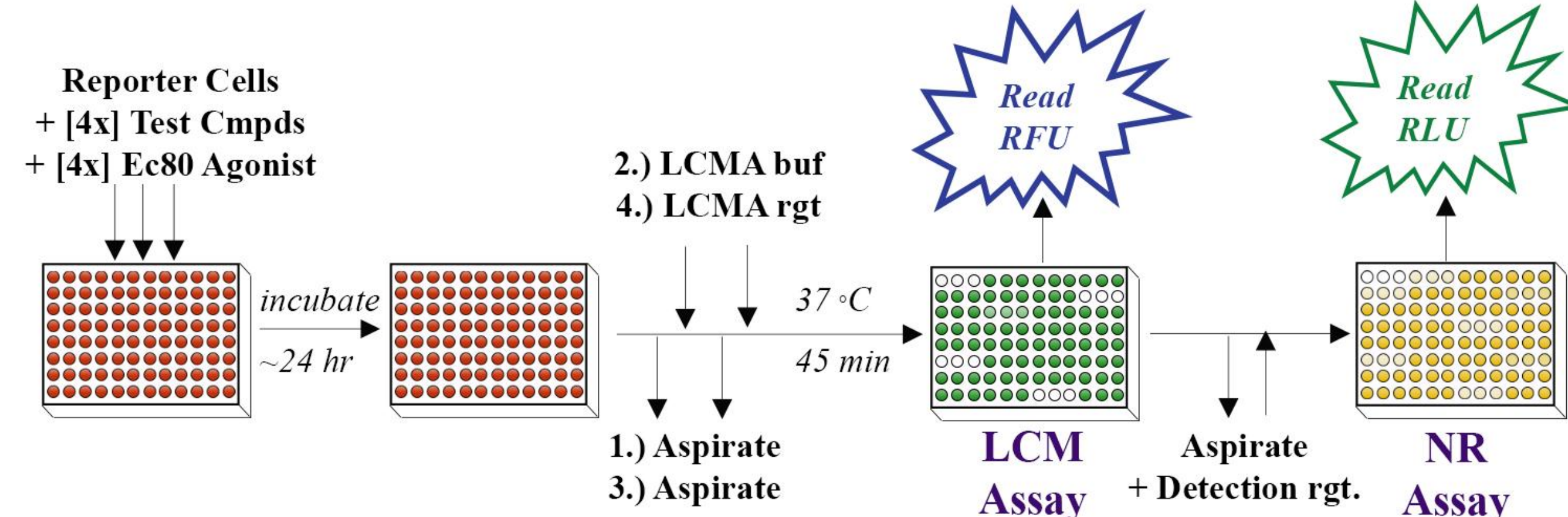


Figure 2. Antagonist Nuclear Receptor assay and integrated Live Cell Multiplex (LCM) assay.



Primary Screen

A panel of 25 'retinoid like' compounds were screened at 3 concentrations against each of the RARs, RORs and RXRs to assess their bioactivity and selectivity. Test compounds were screened at: + 5 μ M, + 1.67 μ M, and + 0.55 μ M (except HX531 + 2.5 μ M, + 0.83 μ M, and + 0.27 μ M); n=3

Figure 3. Human RAR alpha Assays.

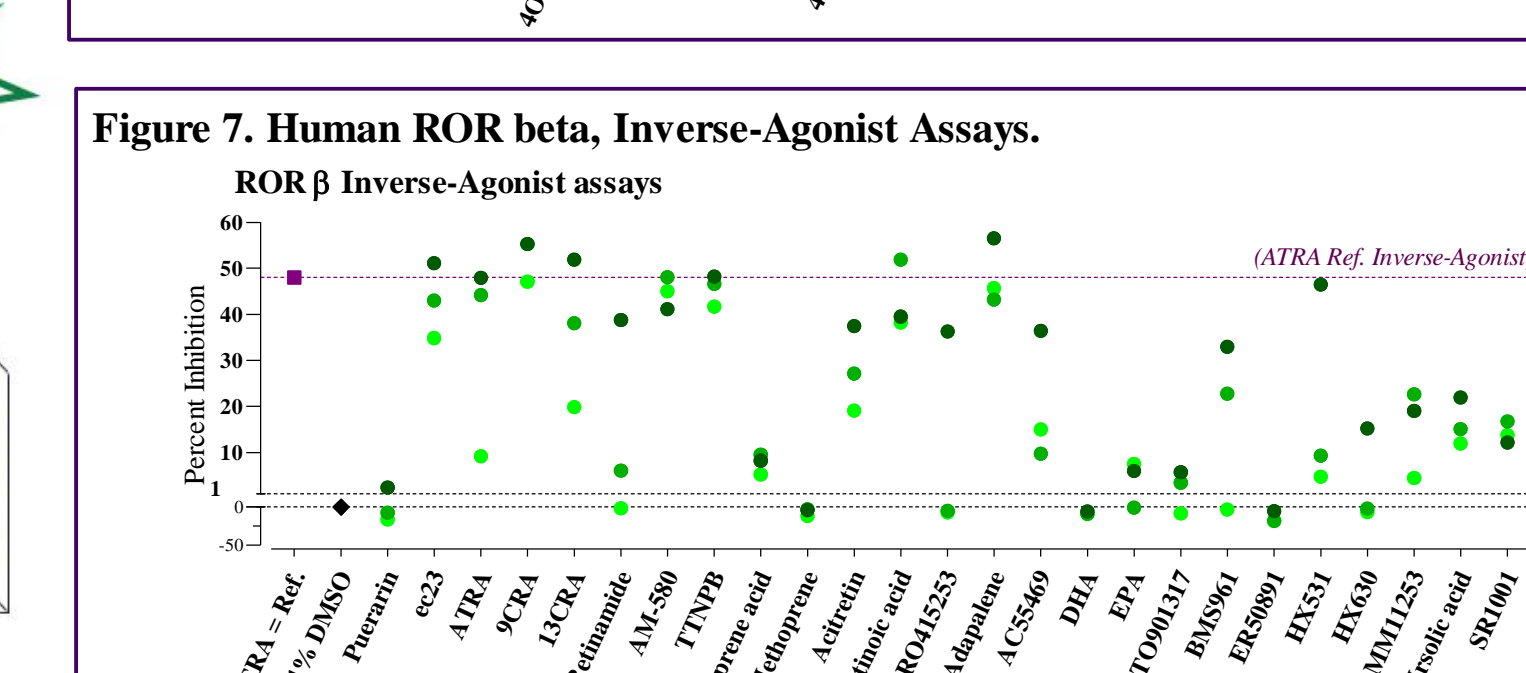
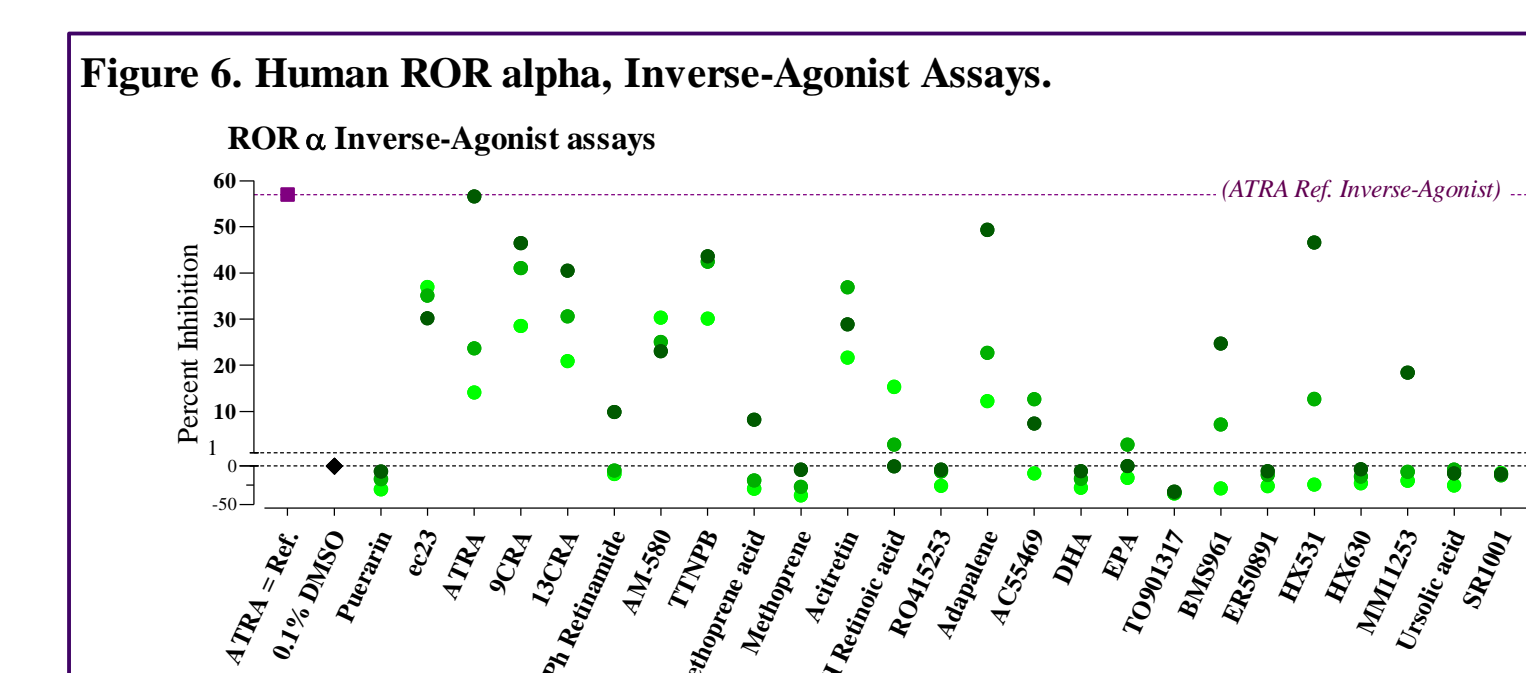
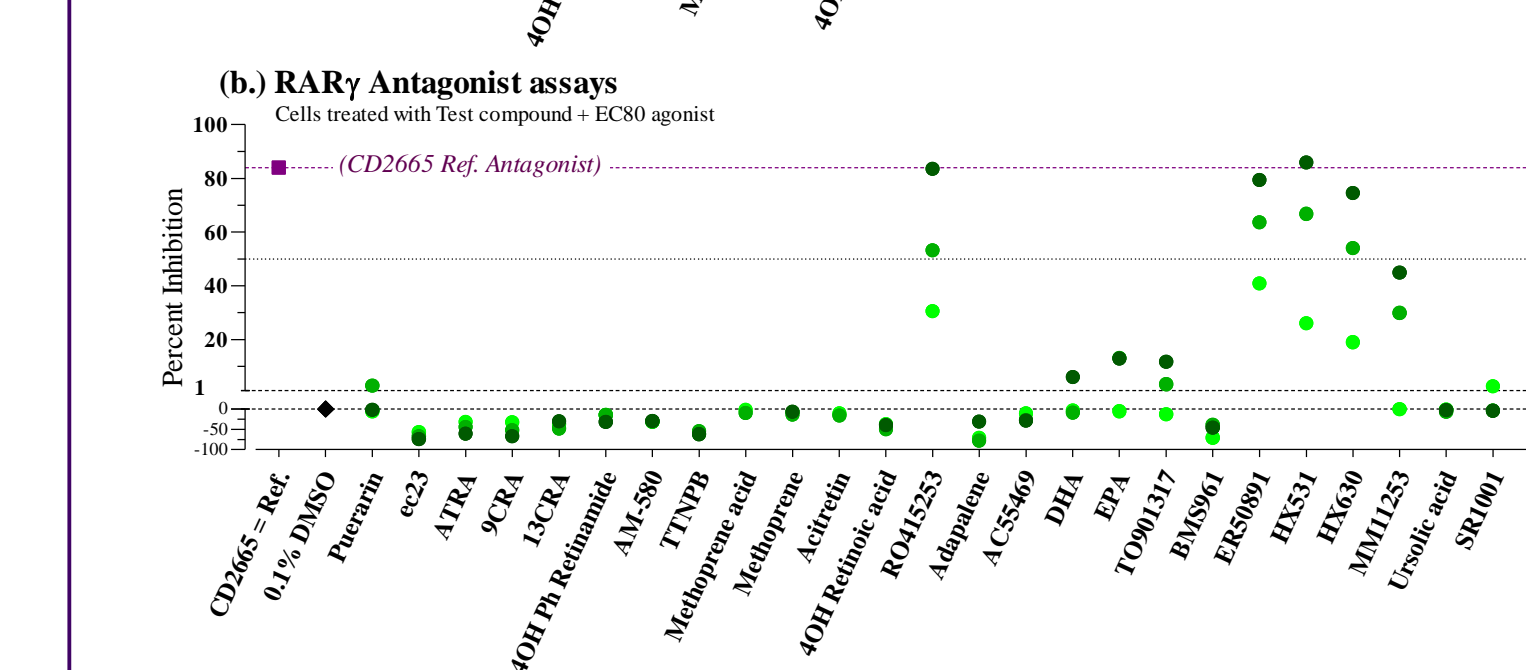
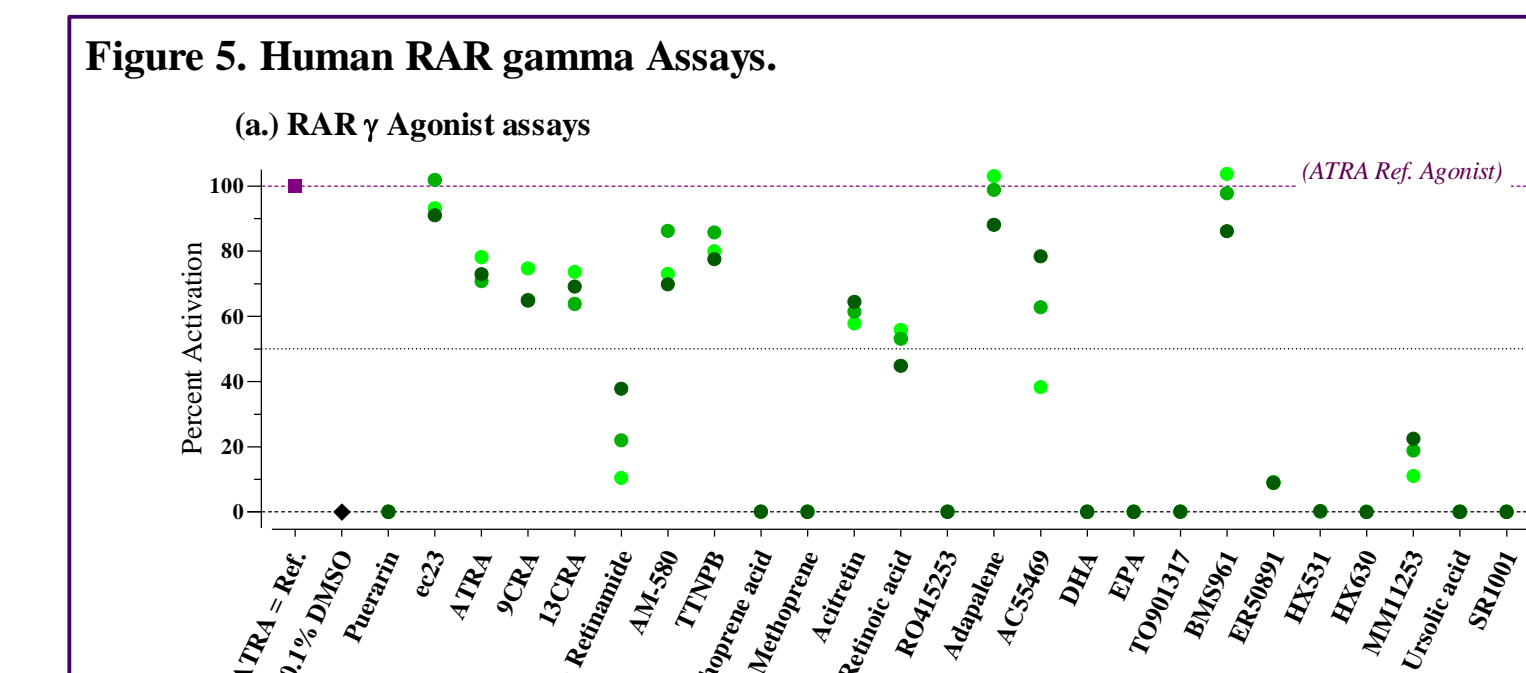
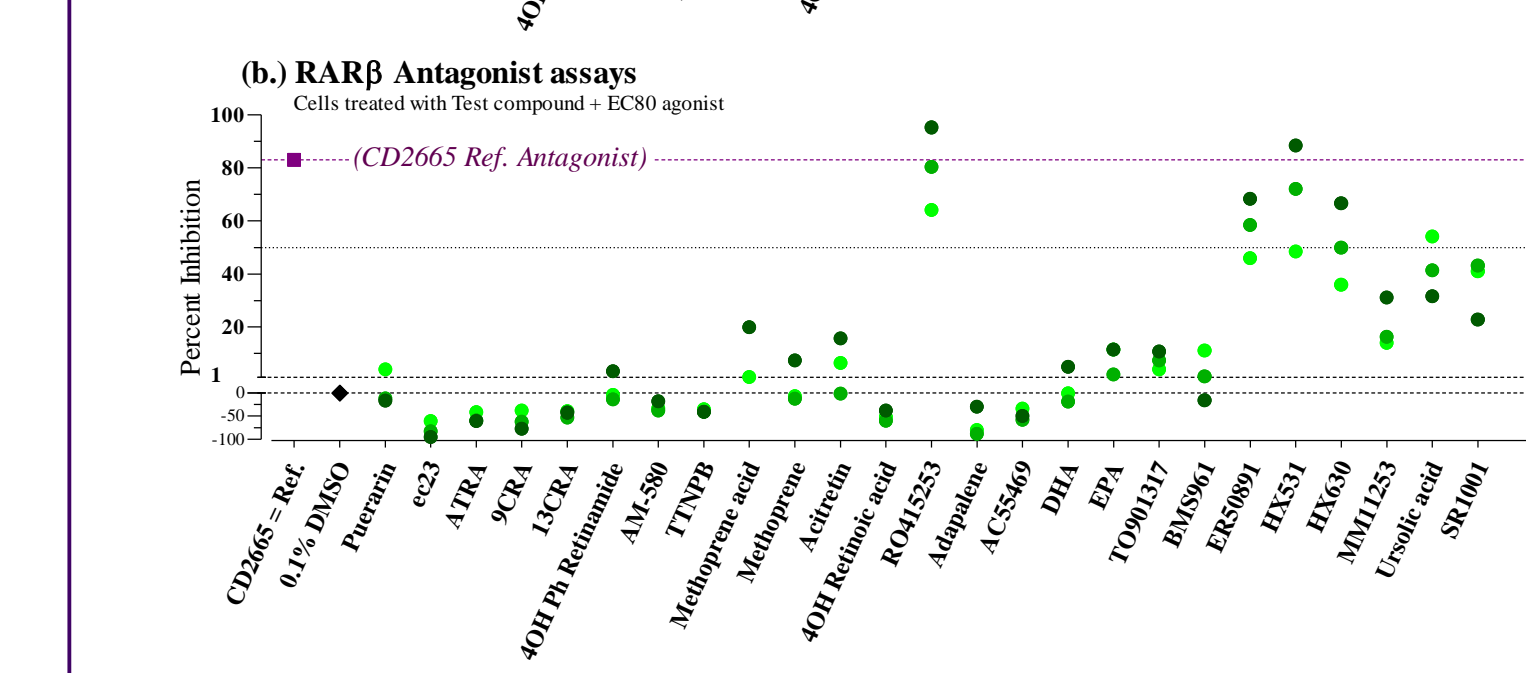
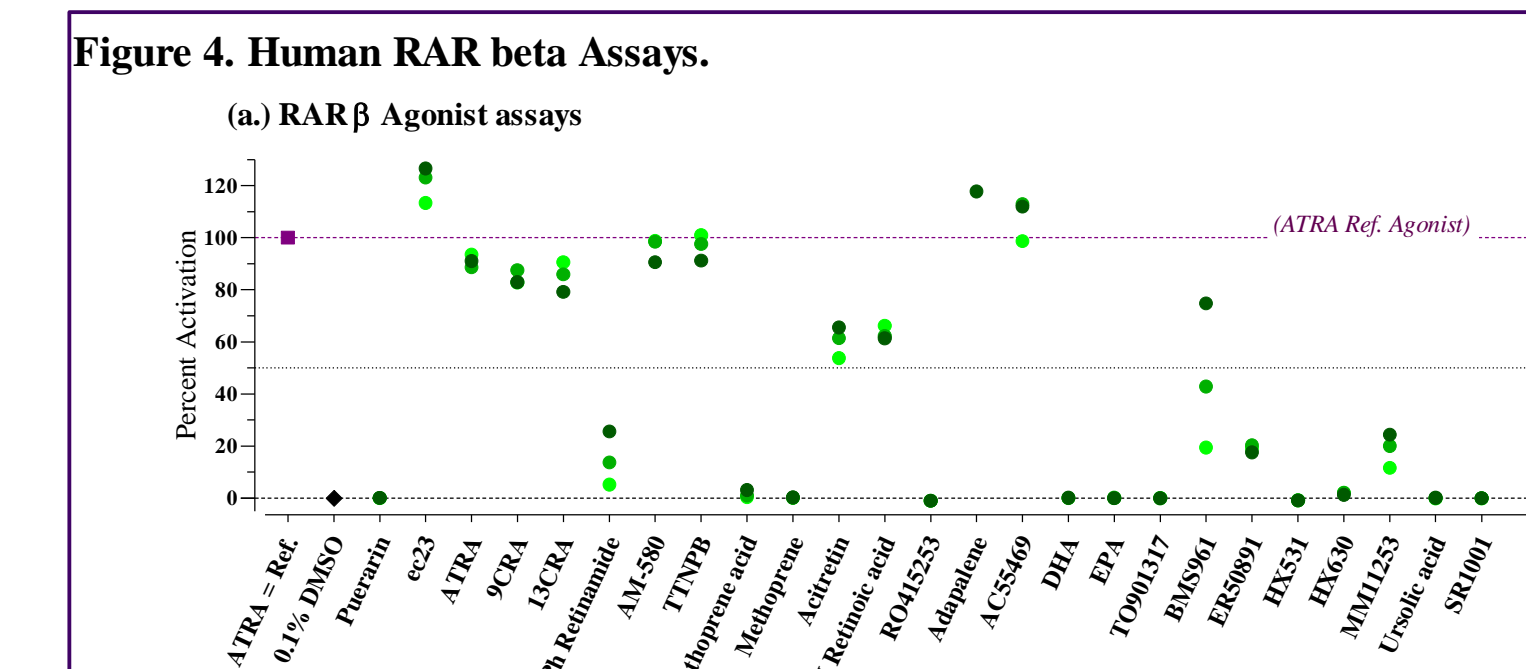
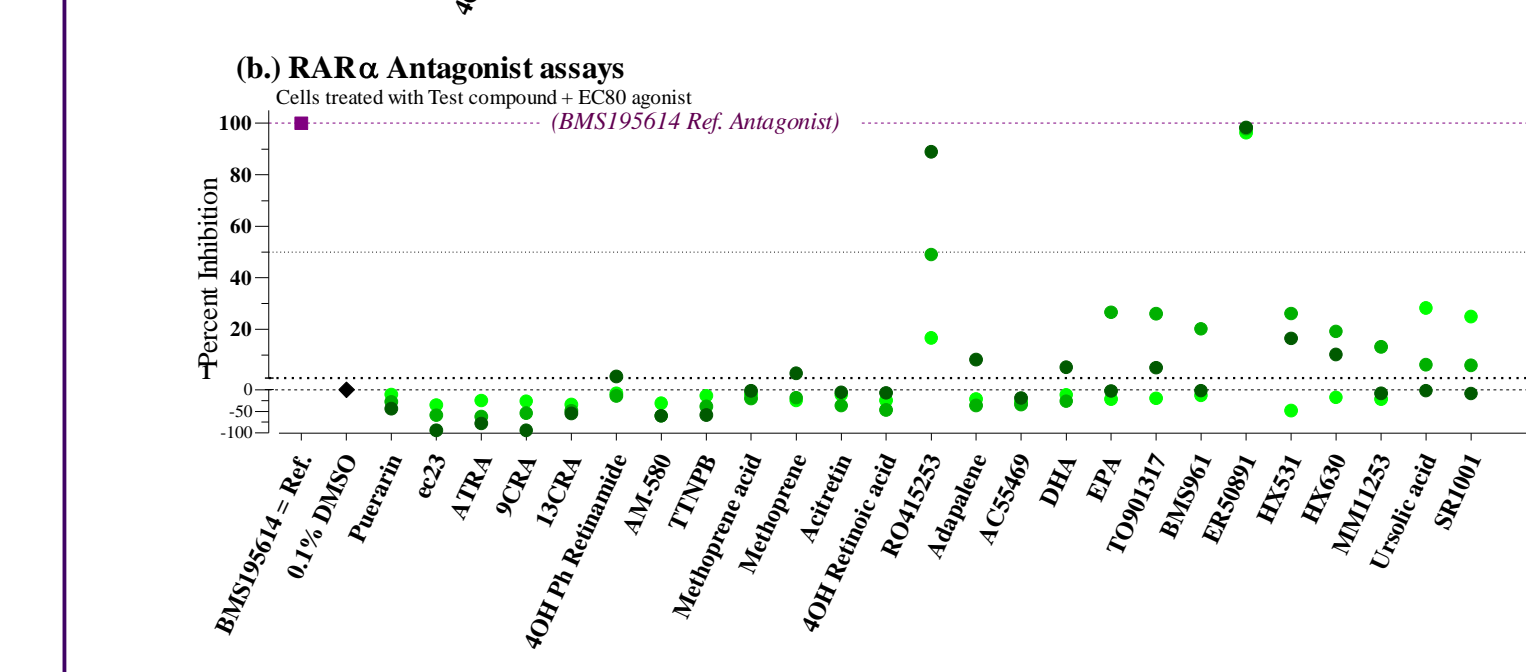
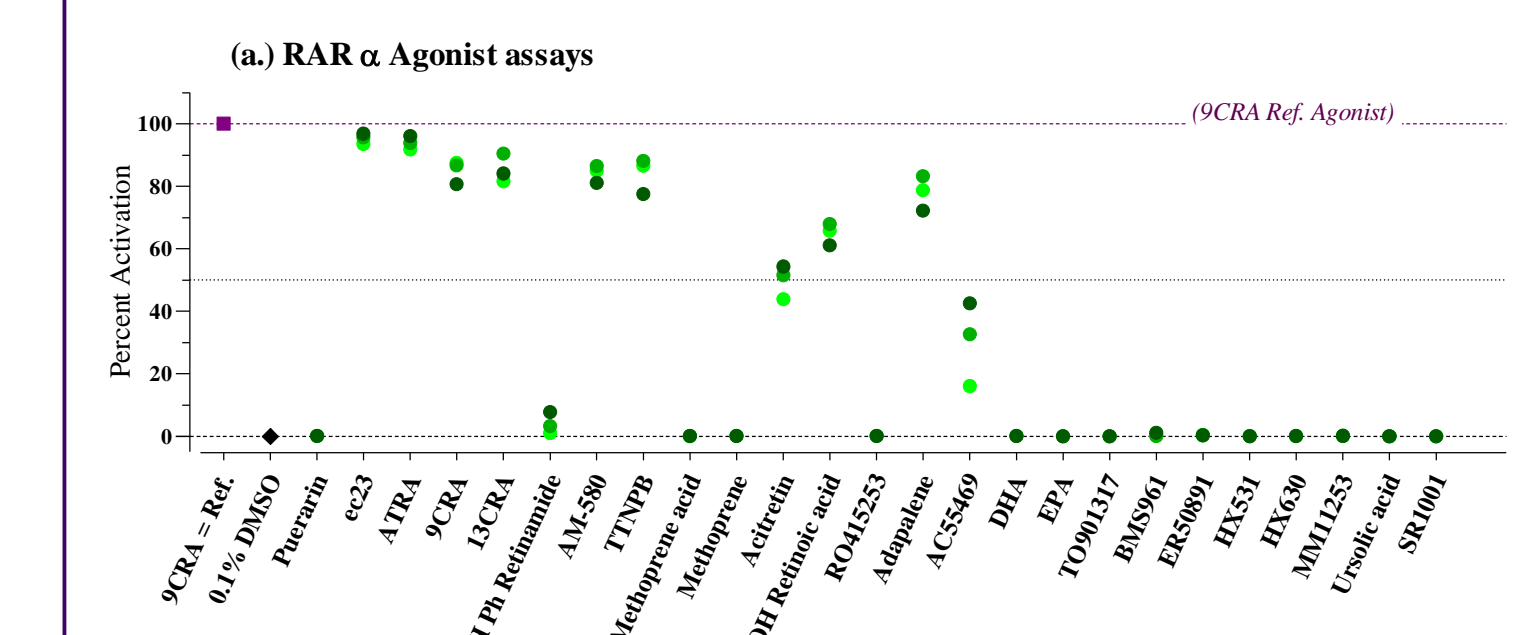
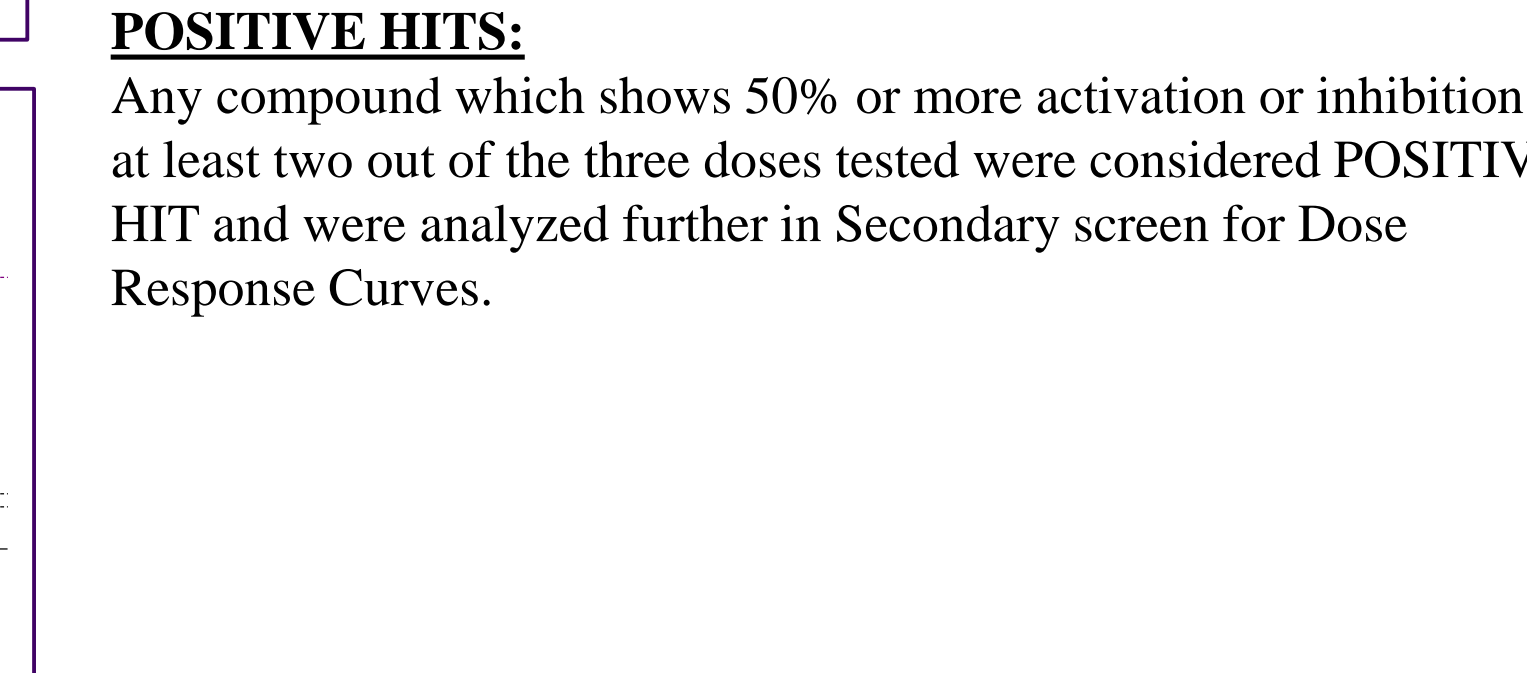
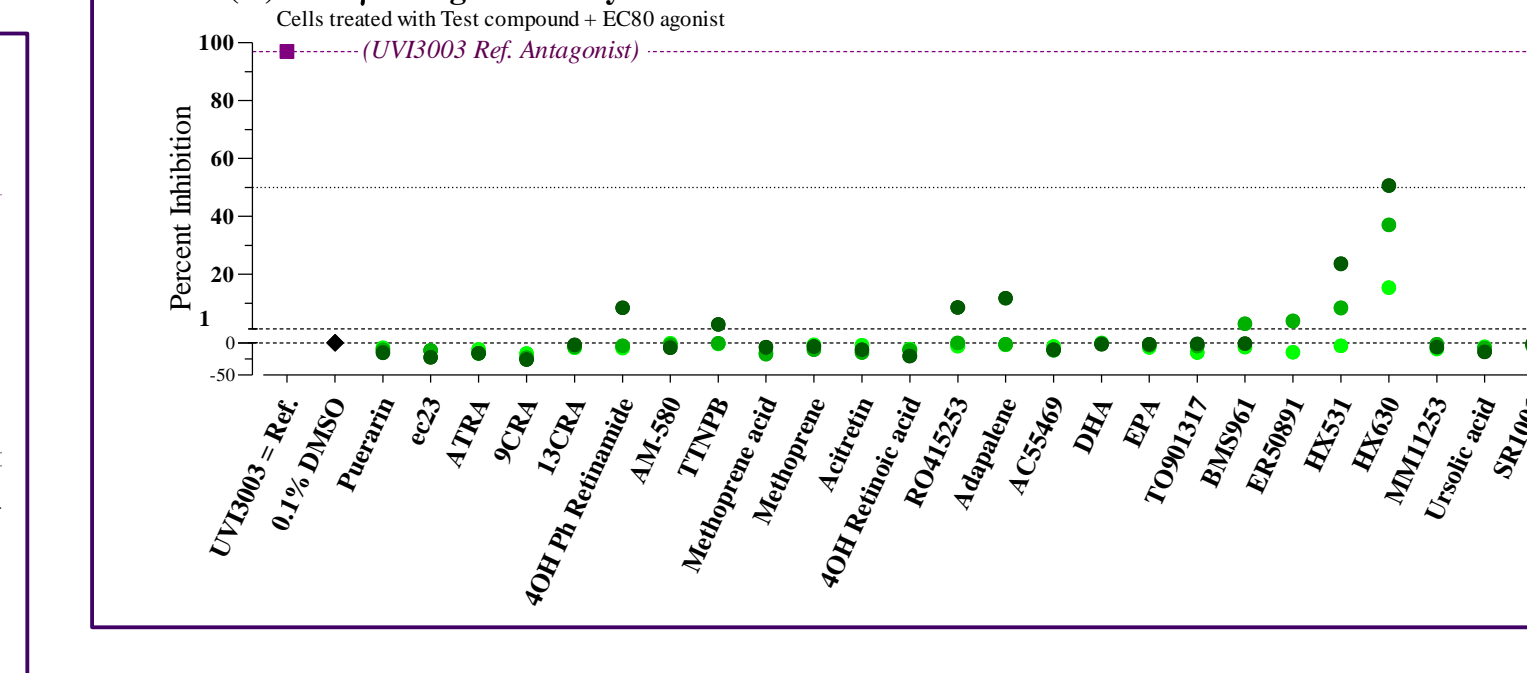
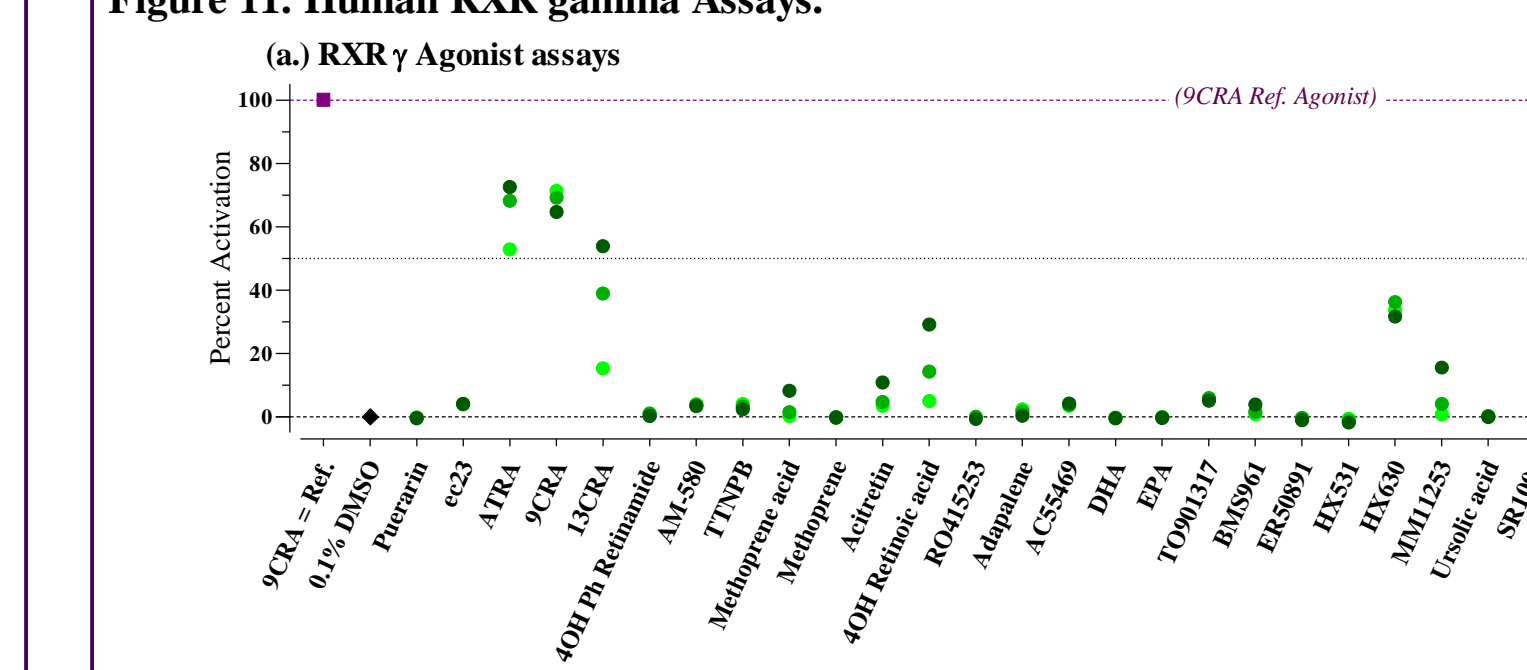
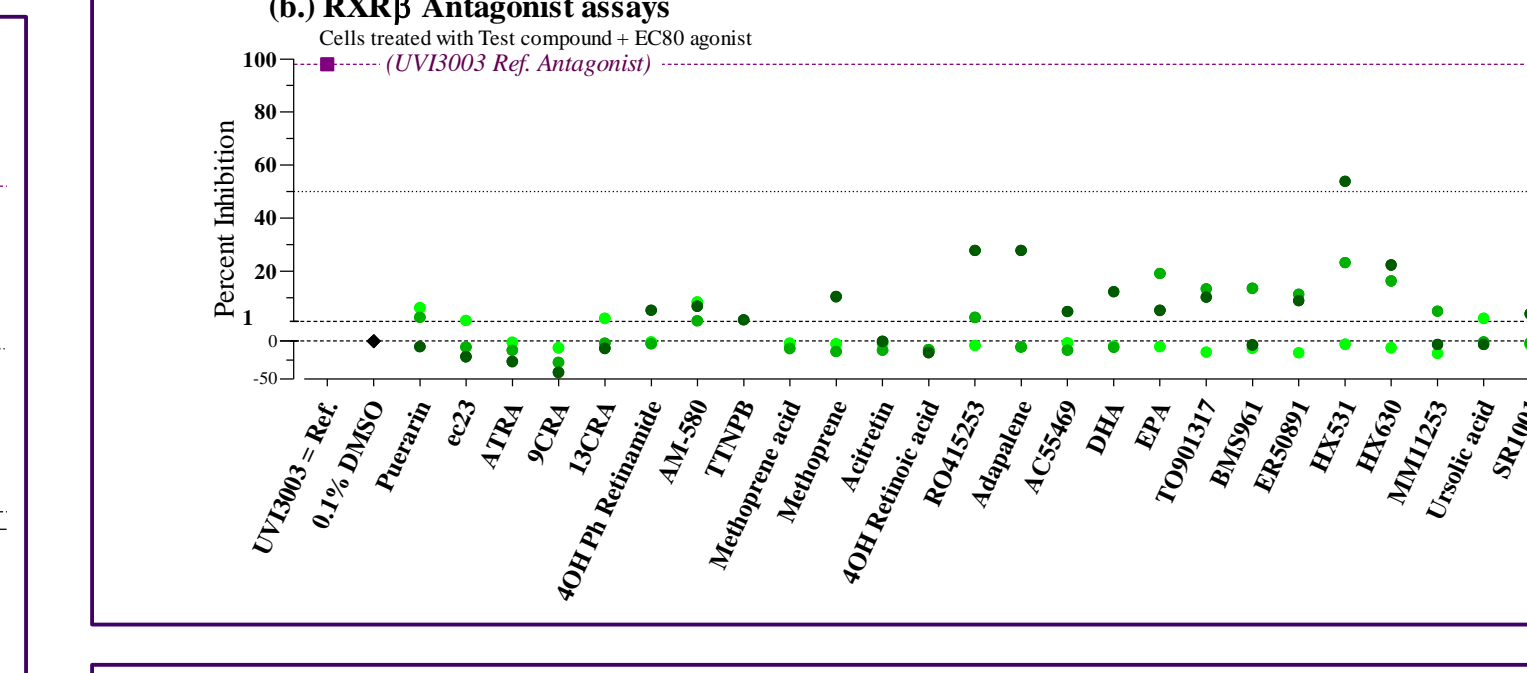
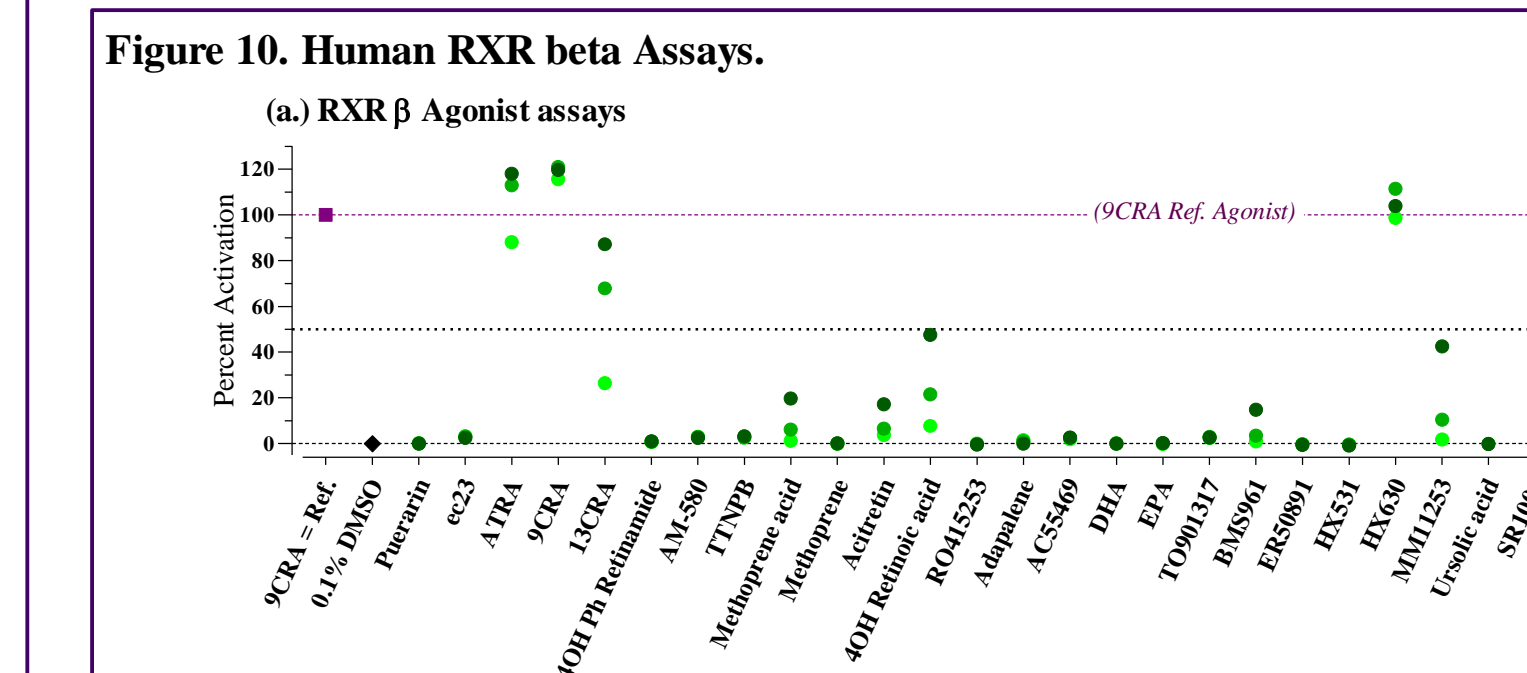
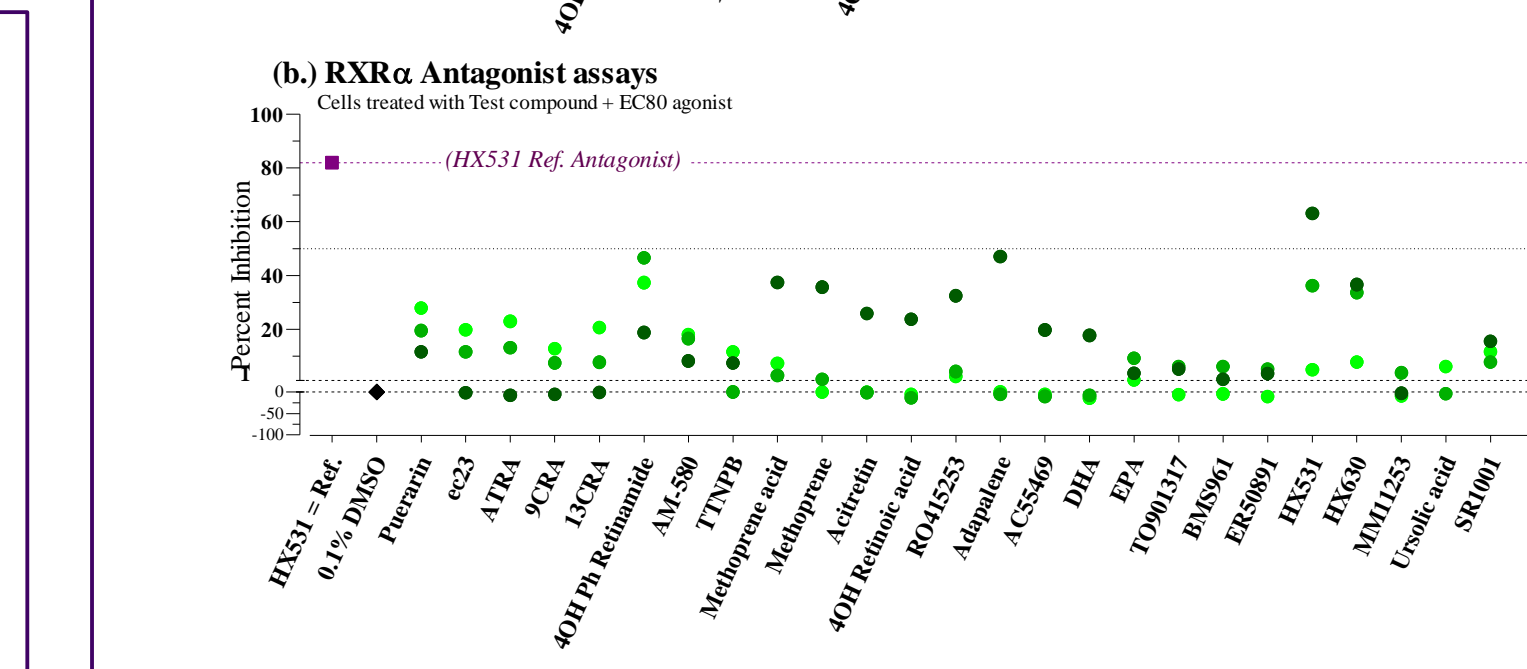
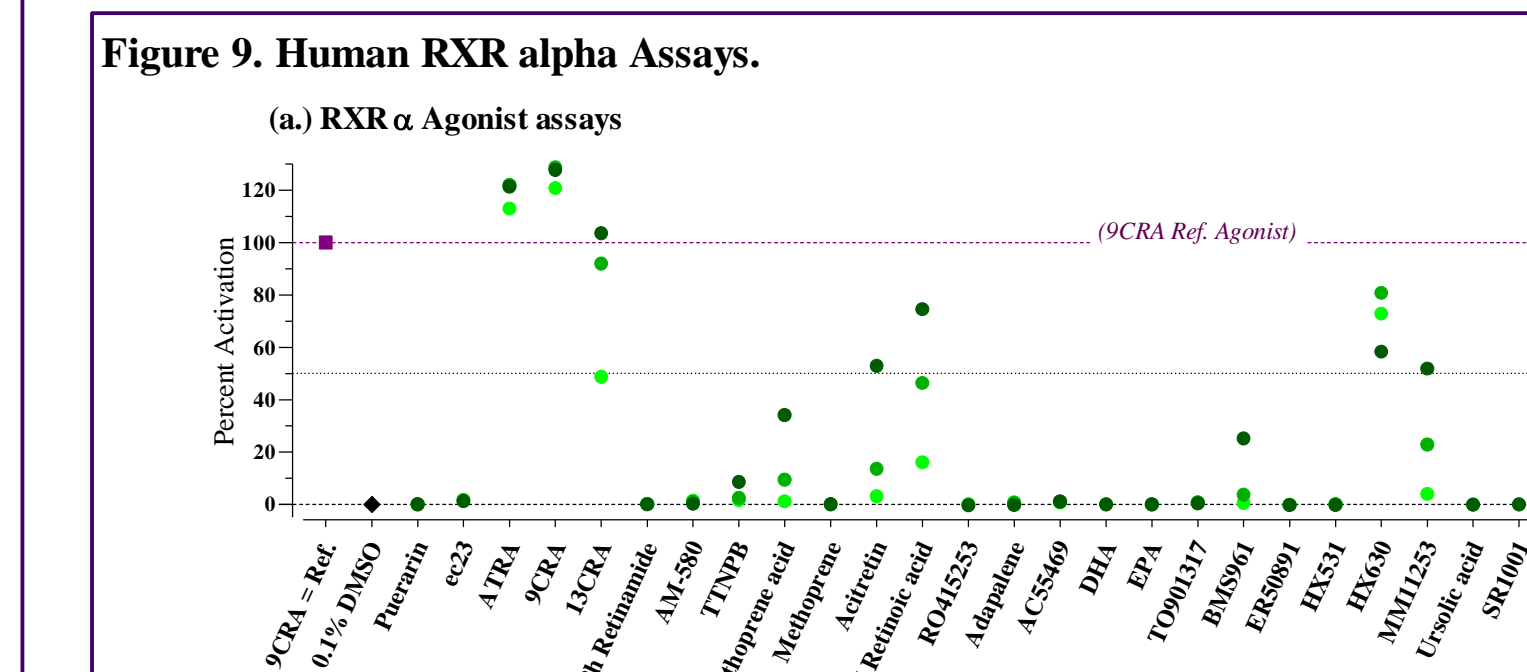
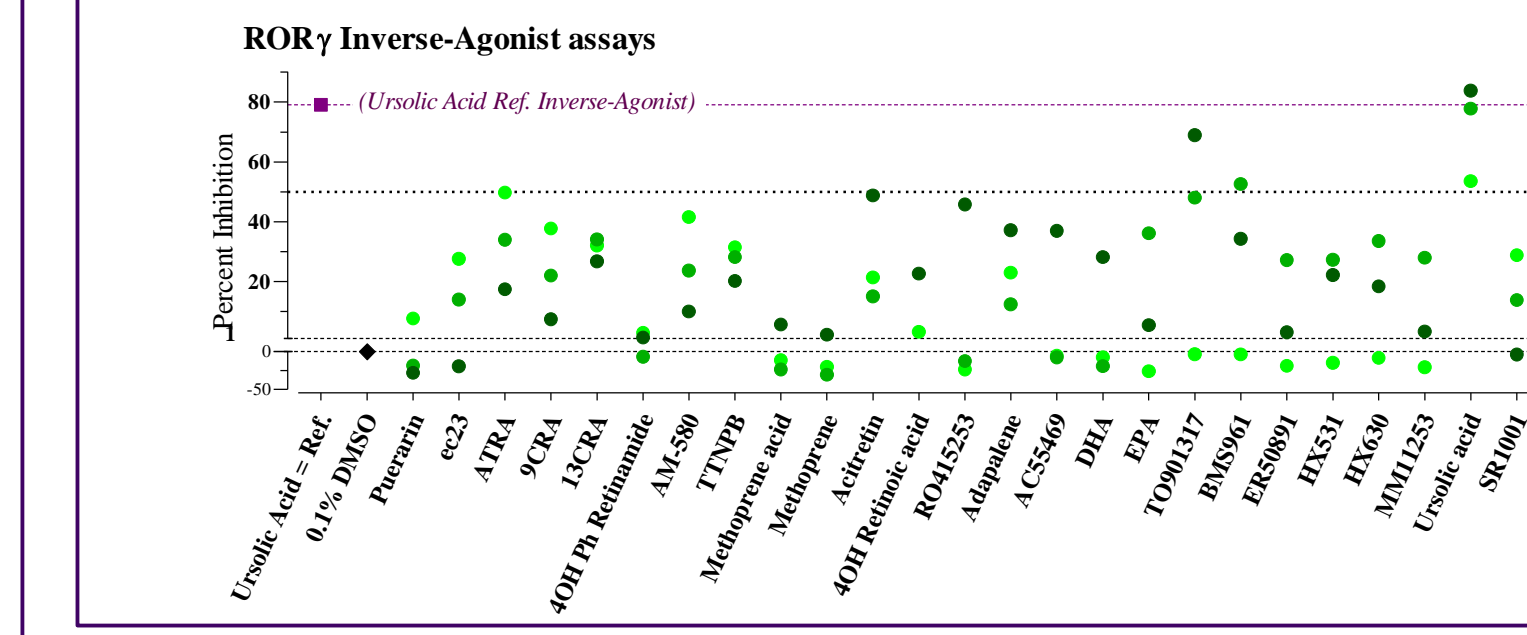


Figure 8. Human ROR gamma, Inverse-Agonist Assays.



Positive compounds in AGONIST mode

Compound	[nM]	Percent (%) ACTIVATION					
		RAR α	RAR β	RAR γ	RXR α	RXR β	RXR γ
ec23	555	94	113	93	17	33	39
	1,667	96	123	102	17	30	39
	5,000	97	127	91	12	26	40
ATRA	555	92	94	78	113	88	53
	1,667	94	89	71	122	113	68
	5,000	96	91	73	121	118	72
9 cis-Retinoic Acid	555	88	83	75	121	116	71
	1,667	87	88	65	129	121	69
	5,000	81	83	65	128	120	65
13 cis-Retinoic Acid	555	82	91	74	49	26	15
	1,667	90	86	64	92	68	39
	5,000	84	79	69	104	87	54
AM-580	555	85	99	73	13	29	40
	1,667	87	98	86	10	29	37
	5,000	81	91	70	0	25	33
TTNPB	1,667	88	98	86	24	27	32
	5,000	77	91	78	8.6	3.3	2.4

Compound	[nM]	Percent (%) ACTIVATION					
		RAR α	RAR β	RAR γ	RXR α	RXR β	RXR γ
Acitretin	555	44	54	58	3.1	3.7	3.5
	1,667	52	61	61	14	6.6	4.7
	5,000	54	66	64	53	17.1	11
4-OH Retinoic Acid	555	68	66	66	56	7.7	5.0
	1,667	68	62	53	46	22	14
	5,000	61	61	45	75	47	29
Adapalene	555	79	131	103	0	1.6	2.3
	1,667	83	130	99	0	1.0	1.6
	5,000	72	118	88	0	0	0
AC-55469	555	16	19	104	0	1.0	0
	1,667	33	113	63	1.0	2.4	3.9
	5,000	43	112	78	1.1	2.7	4.2
BMS961	555	0	49	104	0	1.0	0
	1,667	0	13	98	3.6	3.5	1.6
	5,000	1.1	75	86	25	15	3.9
HX630	555	0	2.1	0	73	99	34
	1,667	0	2.1	0	81	111	36
	5,000	0	1.3	0	58	104	32

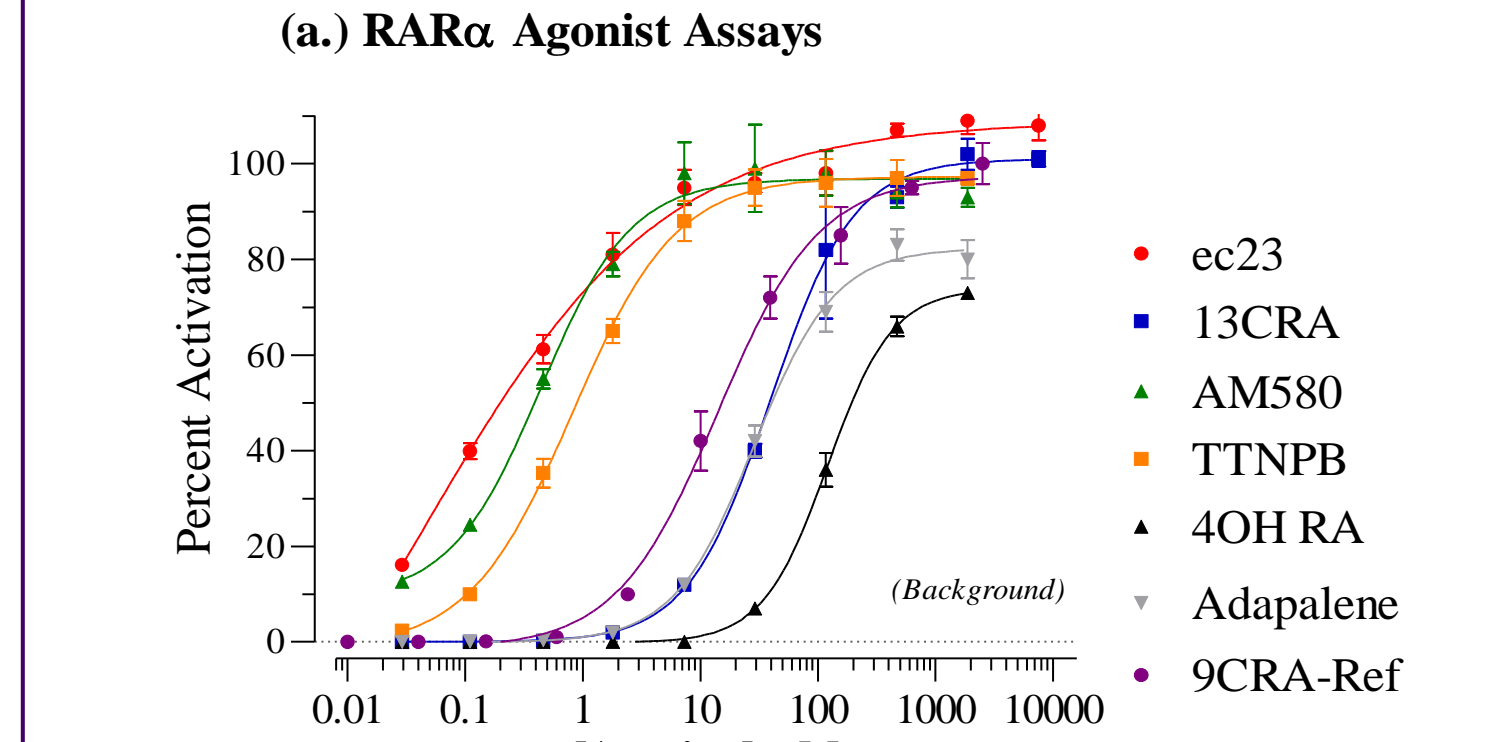
Positive compounds in ANTAGONIST mode

Compound	[nM]	Percent (%) INHIBITION								
		RAR α	RAR β	RAR γ	ROR α	ROR β	ROR γ	RXR α	RXR β	RXR γ
trans-Retinoic Acid	555	0	0	0	14	9	50	23	0	0
	1,667	0	0	0	24	44	34	13	0	0
	5,000	0	0	0	27	48	18	0	0	0
RO41-5253	555	17	64	31	0	0	0	2.4	0	0
	1,667	49	80	53	0	0	0	4.3	2.5	0
	5,000	89	95	84	0	0	0	33	28	8.5
ER50891	555	96	96	41	0	0	0	0	0	0
	1,667	98	98	64	0	0	0	27	5.2	11
	5,000	98	68	79	0	0	0	3.2	3.6	8.8
HX531	555	0	49	26	0	0	4.7	0	5.0	0
	1,667	26	72	67	13	9.3	27	36	23	8.3
	5,000	16	88	86	47	46	22	63	54	24
HX630	555	0	36	19	0	0	0	7.9	0	15
	1,667	19	50	54	0	0	34	34	16	37
	5,000	10	67	75	0	15	18	37	22	51
Ursolic Acid	555	28	54	0	0	12	54	6.1	2.2	0
	1,667	6.2	41	0	0	15	78	0	0	0
	5,000	0	32	0	0	22	84	0.7	0	0

Secondary Screen

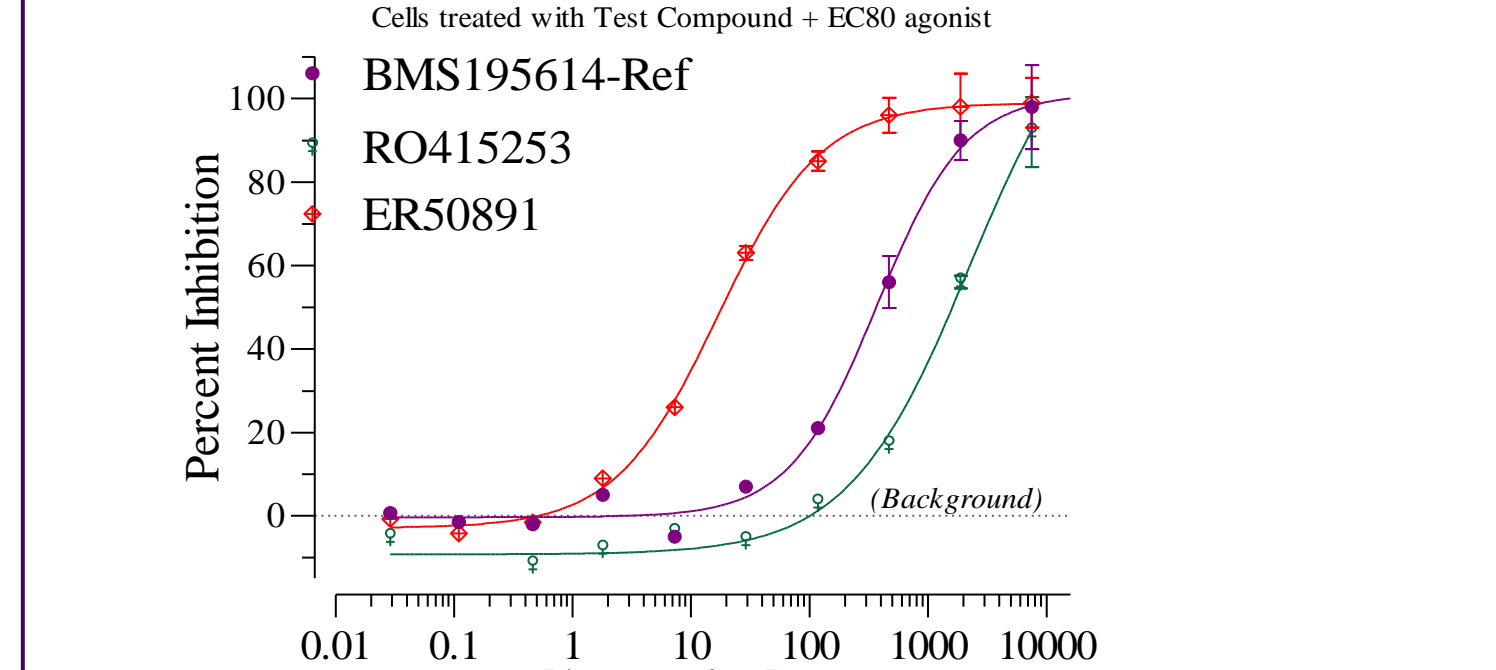
Test compounds showing positive hits in 1^o screen were tested at concentrations starting at 10 μ M (serial dilutions of 1:3). The reference compounds in each assay are denoted by suffix "Ref". Reference compounds were tested at same concentrations as the test compounds. Results from the Live Cell Multiplex Assay demonstrated that these test compounds did not show cytotoxicity at the highest dose tested in antagonist assays, indicating that the positive hits in antagonist mode are not 'False Positives' (results not shown).

Figure 12. Human RAR alpha dose-response assays.



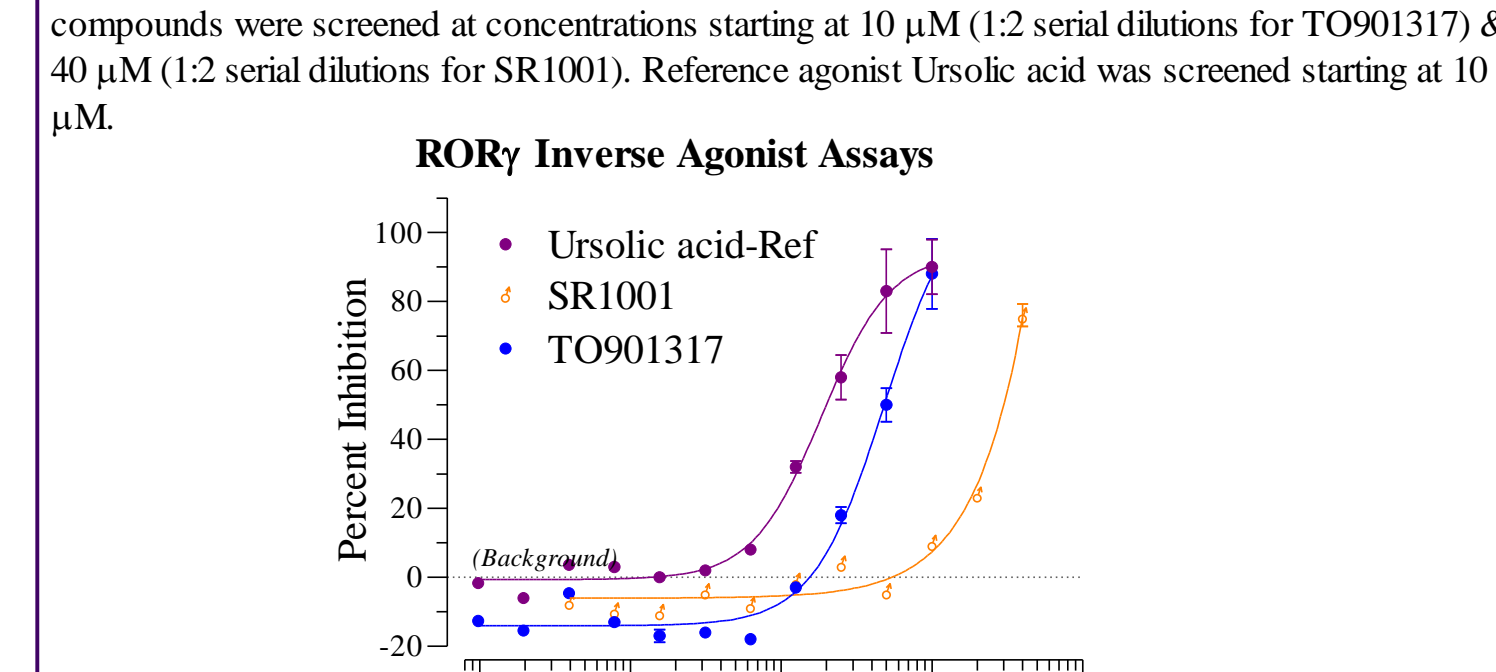
	ec23	13CRA	AM580	TTNPB	4OH RA	Adapalene	9CRA-Ref
HillSlope	0.4218	1.235	1.132	0.9860	1.377	1.268	1.008
EC50	0.04136	39.34	0.4384	0.8181	120.6	28.66	14.08
Span	172.7	101.0	87.70	98.62	74.07	82.44	98.67

Figure 13. Human RAR beta dose-response assays.



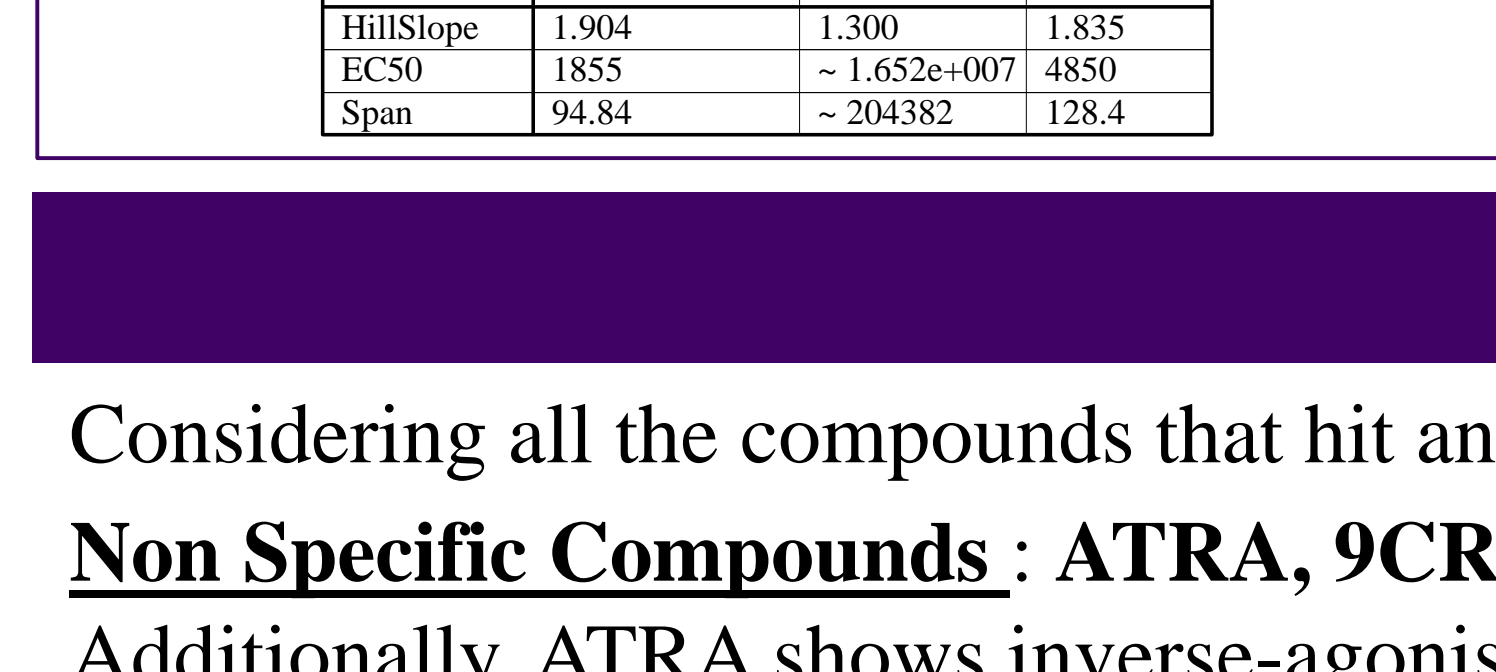
	BMS195614-Ref	RO415253	ER50891
HillSlope	1.176	0.6787	0.9728
EC50	370.8	2243	16.89
Span	101.6	138.2	102.0

Figure 14. Human RAR gamma dose-response assays.



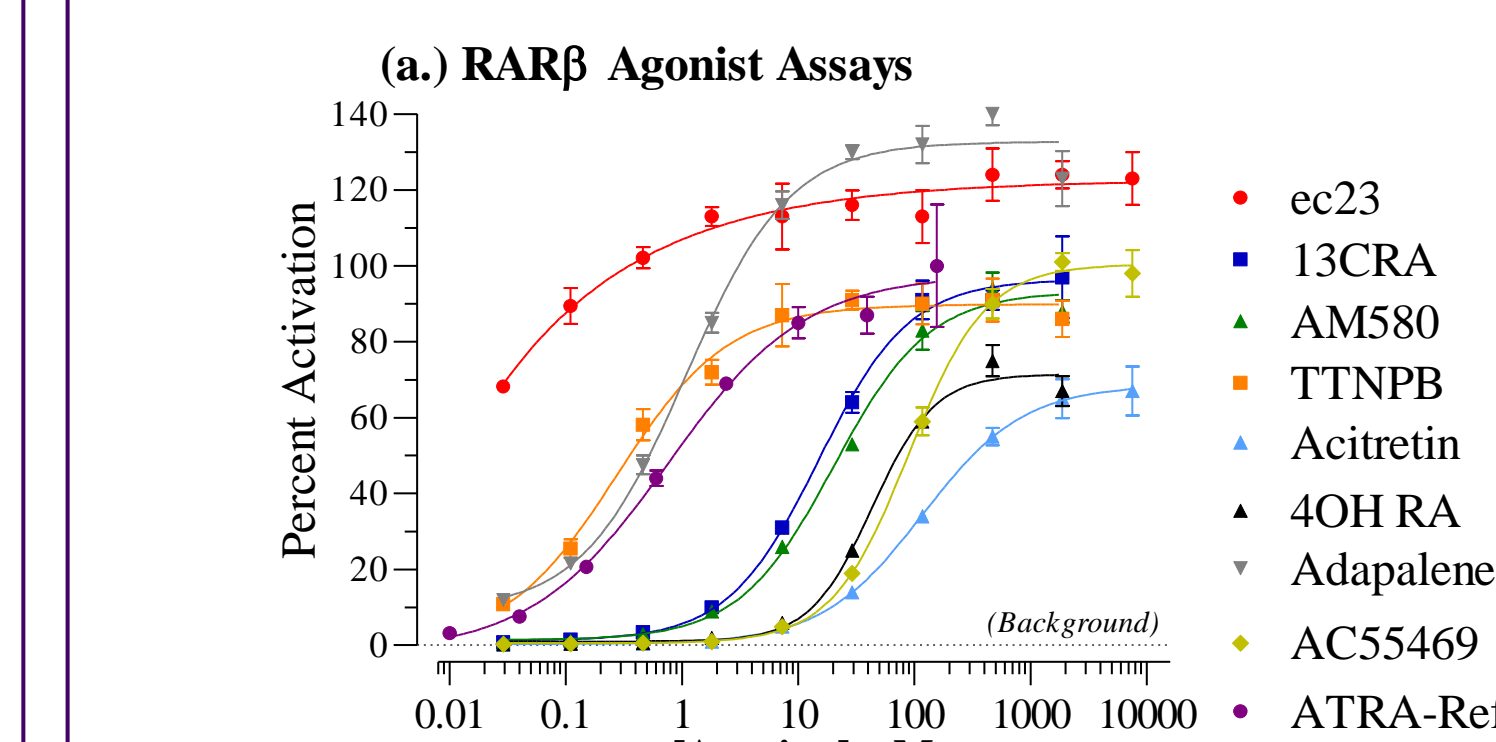
	C2665-Ref	RO415253	ER50891	HX531	HX630
HillSlope	0.7568	0.6787	0.9728	0.5377	0.5083
EC50	917.5	1222	1201	14102	~4.035e+010
Span	93.62	135.2	96.82	304.9	~197225

Figure 15. Human ROR gamma inverse-agonist dose-response assays.



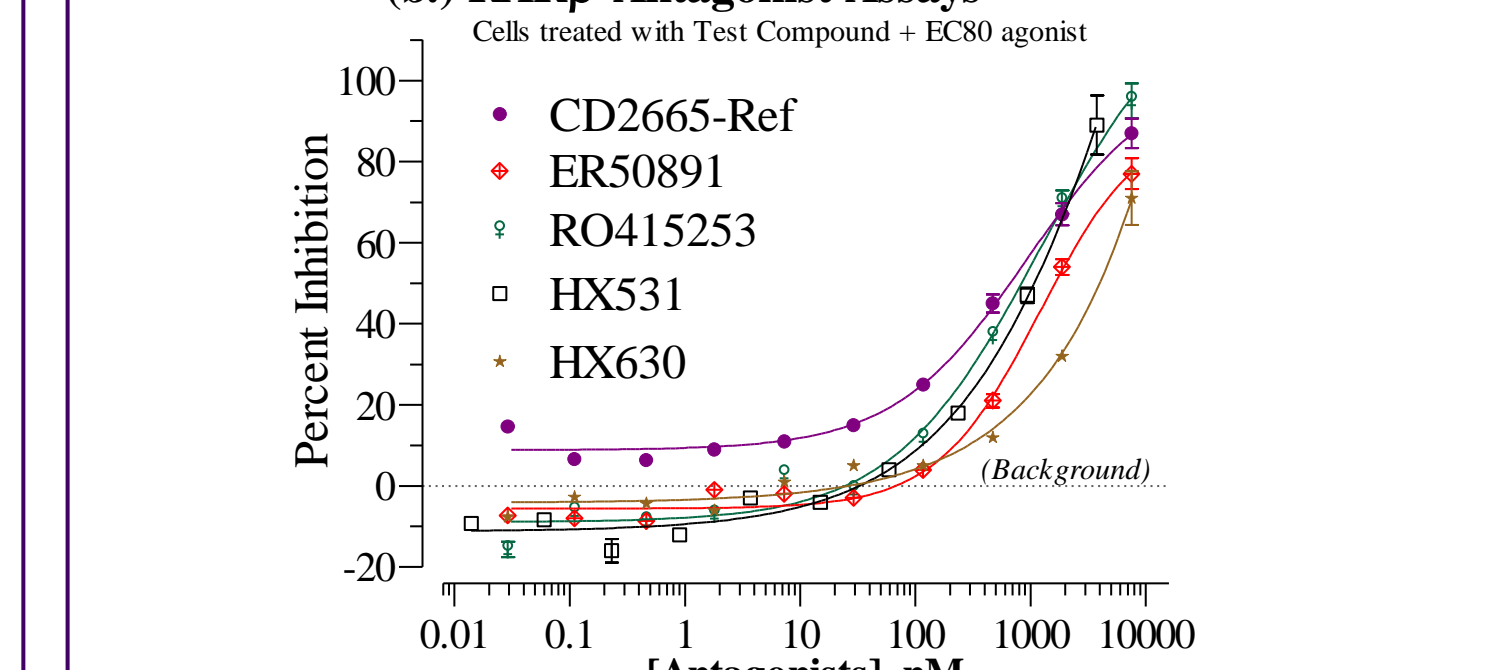
	Ursolic acid-Ref	SR1001	TO901317
HillSlope	1.904	1.300	1.835
EC50	1855	~1.652e+007	4850
Span	94.84	~204382	128.4

Figure 16. Human RXR alpha agonist dose-response assays.



	13CRA	9CRA-Ref	HX630
HillSlope	1.561	1.301	1.335
EC50	570.8	19.18	52.52
Span	74.95	98.66	63.22

Figure 17. Human RXR beta agonist dose-response assays.



	13CRA	9CRA-Ref	HX630
HillSlope	1.947	1.616	1.494
EC50	1142	41.77	130.3
Span	64.12	97.18	82.07

Conclusions

Considering all the compounds that hit any receptor at percent activation or inhibition of more than 50%, they can be classified as:

Non Specific Compounds: ATRA, 9CRA, and 13CRA are not specific agonists as they show activity against all 6 receptors. Additionally, ATRA shows inverse-agonistic activity against ROR α and ROR γ .

Moderately Specific Compounds: 4-hydroxy retinoic acid show agonistic activity against all three RARs and RXR α . HX630 shows agonistic activity against RXR β and RXR γ and antagonistic activity against RAR α , RAR β and RXR γ . HX531 shows antagonistic activity against RAR α , RAR β and RXR α , RXR β .

Specific Compounds: ec23, AM580, TTNPB, Acitretin, and Adapalene show agonistic activity against RARs, whereas AC55469 and BMS961 show agonistic activity against RAR β and RAR γ with little or no activity against RAR α , respectively. RO415253 and ER50891 show antagonistic activity only against the RARs. Ursolic acid shows inverse-agonistic (or antagonistic activity) against ROR γ and RAR β .

POSITIVE HITS:

Any compound which shows 50% or more activation or inhibition in at least two out of the three doses tested were considered POSITIVE HIT and were analyzed further in Secondary screen for Dose Response Curves.