JBC Papers in Press. Published on October 16, 2014 as Manuscript M114.610501

The latest version is at http://www.jbc.org/cgi/doi/10.1074/jbc.M114.610501

 β -apo-13-Carotenone as an RXR antagonist

 $\label{eq:based} \begin{array}{l} \beta \text{-apo-13-Carotenone Regulates Retinoid X Receptor Transcriptional Activity Through} \\ \text{Tetramerization of the Receptor*} \end{array}$

Jian Sun¹, Sureshbabu Narayanasamy^{1, 2}, Robert W. Curley Jr.², Earl H. Harrison¹

From the ¹Department of Human Sciences and ²College of Pharmacy, The Ohio State University, Columbus, OH 43210 USA

*Running title: β -apo-13-Carotenone as an RXR antagonist

To whom correspondence should be addressed: Earl H. Harrison, 314 Campbell Hall, 1787 Neil Ave., Columbus, OH, 43210 USA, Tel.: (614) 292-8189; Fax: (614) 292-4339; Email:harrison.304@osu.edu.

KEYWORDS: retinoid receptors; tetramerization; apocarotenoid

β-apo-13-Carotenone, **Background:** а naturally-occurring apocarotenoid, functions as an antagonist of the retinoid X receptor (RXR). β-apo-13-Carotenone **Results:** inhibits transactivation of RXRα but does not interfere with coactivator binding to the receptor like the antagonist known UVI3003. **Conclusion:** β-apo-13-Carotenone induces the formation of a transcriptionally silent RXR tetramer.

Significance: β -apo-13-Carotenone is a naturally-occurring rexinoid with a novel mechanism of antagonism.

ABSTRACT

Retinoid X receptor (RXRa) is activated by 9-cis-retinoic acid (9cRA) and regulates transcription as a homodimer or as a heterodimer with other nuclear receptors. We have previously demonstrated that β -apo-13carotenone, an eccentric cleavage product of βcarotene, antagonizes the activation of RXRa by 9cRA in mammalian cells overexpressing this receptor. However. the molecular mechanism of β -apo-13-carotenone's modulation on RXRa transcriptional activity is not understood and is the subject of this report. We performed transactivation assays using full length RXRa and reporter gene constructs (RXRE-Luc) transfected into COS-7 cells and luciferase activity was examined. β-apo-13-Carotenone was compared with the RXRa antagonist UVI 3003. The results showed that

both β-apo-13-carotenone and UVI 3003 shifted the dose-dependent RXRα activation by 9cRA. In contrast, results of assays using a Gal4-DBD:RXRa-LBD hvbrid receptor reporter cell assay that detects 9cRA-induced coactivator binding to the ligand binding demonstrated that domain UVI3003 significantly inhibited 9cRA-induced coactivator binding to RXRaLBD, but β-apo-13-carotenone did not. However, both β-apo-13-carotenone and UVI 3003 inhibited 9-cRA induction of caspase 9 gene expression in the mammary carcinoma cell line MCF7. In order to resolve this apparent contradiction we investigated the effect of β -apo-13-carotenone on the oligomeric state of purified recombinant RXR α LBD. β -apo-13-carotenone induces tetramerization of the RXRaLBD whereas UVI3003 had no effect on the oligomeric state. These observations suggest that β -apo-13carotenone regulates RXRa transcriptional activity through inducing the formation of the "transcriptionally silent" RXRα tetramer.

Retinoid X receptors (RXR α , RXR β and RXR γ) are members of nuclear receptor (NR) family and play a central role in NR-regulated signaling pathways. RXRs are involved in biological processes, including cell growth and differentiation, metabolism, morphogenesis, and embryogenic development (1 – 7) The active form of RXR is a dimer or heterodimer (8, 9). Besides the RXR homodimer, RXR also forms heterodimers with other NR family members, including retinoic acid receptors (RARs), the vitamin D receptor (VDR), peroxisome proliferator-activated receptors (PPARs), the farnesoid X receptor (FXR) and the liver X receptors (LXRs) (10, 11). RXR naturally forms into tetramers which are transcriptionally inactive (12).

RXRs are primarily made up of two modular domains: a central DNA binding domain (DBD) and a carboxy-terminal ligand binding domain (LBD). In addition to its role in binding of ligands, the LBD contains dimerization motifs and an activation function 2 (AF-2) domain (13, 14). Ligand free RXR represses transcription of target genes through interaction with corepressor proteins. Ligand binding induces a conformational change of the AF-2 helix that releases corepressor protein and allows recruiting of coactivator complexes. compounds synthesized Numerous as antagonists, such as UVI 3003, target the AF-2 helix (13).

Ligand free RXR tends to associate into homotetramers both in solution and when bound to DNA. However, RXR tetramers rapidly dissociate into active dimers upon binding of an agonist such as 9-cis retinoic acid (9cRA). RXR heterodimers bind in regulatory regions of their target genes by associating with response elements (REs). RXR homodimers bind to a retinoid DNA response element (RXRE). Activation of DNA-bound dimers by ligands promotes the recruiting of transcriptional coactivators to the promoters of target genes and enhances transcription rate. In vitro studies have indicated that full-length RXR self-associates into tetramers and the LBD alone is sufficient to mediate tetramer formation with 3-5 nM affinity between the dimers (12, 15). Studies have substantiated the existence in vivo of an RXR tetramer (12, 15). It also has been shown that the RXR tetramer is transcriptionally silent based on the correlation between the transcriptional activity of RXR mutants and their ability to form tetramers (16).

The vitamin A metabolite 9cRA is a ligand of RXR (17). Binding of 9cRA as an agonist induces the dissociation of the tetramer into dimer, which is the first step for RXR activation (18.19). Carotenoids are polyisoprenoids that are biosynthesized in plants, fungi and bacteria. Approximately 50 -60 carotenoids that contain at least one unsubstituted *B*-ionone ring and the correct number and position of methyl groups in the polyene chain exhibit provitamin A activity (20, 21). Dietary provitamin A, β -carotene, can be metabolized in mammals through two pathways (22). β-Carotene oxygenase 1 (BCO1) catalyzes the cleavage of the 15,15' double bond resulting in two retinaldehyde molecules, and the eccentric cleavage takes place at double bonds other than the central 15.15' double bond to produce β -apocarotenoids with different chain lengths. β-Apocarotenoids have been detected in foods (23) and the blood of both humans (24) and animals (25). Recently, β -apo-8'-carotenal was detected in plasma after ingestion of β -carotene by a healthy human subject (24).Our previous studies demonstrated β-apo-13-carotenone that functioned as an antagonist in transactivation assay using full length RXR α (26) and the retinoic acid receptors (RAR) α , β , and γ (27). We have reported that β -apo-13-carotenone competes for 9cRA binding to RXRa with an affinity (7-8 nM) identical to 9cRA itself (27). However, the molecular mechanism of β -apo-13-carotenone's modulation of transcriptional activity is not understood yet. This study focused on the mechanism by which β -apo-13antagonizes 9cRA carotenone induced activation of RXRa. Our results show that Bapo-13-carotenone induces formation of the RXRα transcriptionally-silent tetramer but does not inhibit coactivator recruitment to the isolated LBD.

EXPERIMENTAL PROCEDURES

Materials

COS-7 African green monkey kidney cells and MCF7 mammary carcinoma cells from ATCC (Rockville, MD) were cultured in

 β -apo-13-Carotenone as an RXR antagonist

DMEM supplemented with 10% FBS. Cells were maintained at 37° C with 10% CO₂. 9-cis Retinoic acid and UVI3003 were purchased from Santa Cruz. β -apo-13-carotenone was synthesized as described previously (27). All other chemicals were from Sigma-Aldrich.

Nuclear Receptor Reporter Cell Assay with Full Length hRXRa

COS-7 cells were cultured in 96-well plates overnight. cDNA of full length human RXRa in pSV sport vector (Addgene) was cotransfected with Renilla (pRL-tk) and Firefly luciferase (RXRE-Luc) reporter constructs into COS-7 cells in serum free DMEM with XtremeGENE 9 DNA (Roche). Twenty four hours after transfection, COS-7 cells in DMEM with 10% charcoal stripped FBS were then treated with 9cRA in presence or absence of β-apo-13-carotenone or UVI 3003 for an additional 24 hours. Cell lysates were used in the dual luciferase assay (Promega) to determine the activation of hRXRa by 9cRA and the inhibition by β -apo-13-carotenone and UVI 3003. For each experiment, the firefly luciferase (experimental reporter) activity was normalized to Renilla luciferase (control reporter).

Human RXRα-LBD Reporter Cell Assay

Reporter cells expressing human RXRaLBD fused to the GAL-4 DBD (Indigo Biosciences, State College PA) were treated according to the manufacturer's protocol. Reporter cells were incubated with 0, 0.32, 1.6, 8, 40, 200, 1000, and 5000 nM 9cRA for 24 hours at 37 ° C in the presence or absence of fixed concentrations of either β -apo-13carotenone or UVI 3003. Luminescence was detected with Glomax96 luminometer (Promega).

Quantitative Real-Time PCR

Breast cancer cells MCF7 were cultured in 6-well plates and starved for 24 hours in serum free DMEM. MCF7 cells were then treated with ligands in serum free medium for 4 hours. Total RNA was isolated using NucleoSpin RNA II (Macherey-Nagel). Two micrograms of RNA was reversed-transcribed into cDNA using the high-capacity cDNA reverse transcription kit (Applied Biosystems). Quantitative real-time PCR analysis was performed in quadruplicates with TaqMan Chemistry and probed for caspase 9 (Hs00154260-m1) (Applied Biosystems). Human GAPDH (4333764F) was used as housekeeping gene. The comparative C_t method ($\Delta\Delta C_t$) was used to analyze results.

mRXRaLBD Expression in E. Coli and Protein Purification

mouse RXRa LBD N-His-tagged (pET15b) was transformed into BL21(DE3). The E coli culture was grown at 37°C to OD600 of 0.6. After induction with 0.5 mM IPTG, cells was incubated for another $2 \sim 4$ hours at 25 ° C. Cells were harvested and lysed in lysis buffer (20 mM Tris, 500 mM NaCl, 5 mM imidazole and 3 mM DTT, pH8.0). The supernatant was loaded onto a HisPur Ni-NTA affinity column followed by extensive washing with 20 mM imidazole in lysis buffer. HismRXRaLBD was eluted with 500 mM imidazole in lysis buffer. The concentrated protein peak fraction was then applied to Hiload Superdex 200 gel filtration column for isolation of mRXRaLBD dimer and tetramer. The gel filtration column was calibrated with protein standards of 13.7, 25, 43, 67, 158, 232, and 440 kDa and blue dextran 2000 in order to molecular weights of the confirm the mRXRaLBD dimer and tetramer. Protein concentration was determined with the Bradford reagent (Bio-Rad) with bovine serum albumin as a standard. The purity of protein was assessed by SDS-PAGE and Coomassie Blue staining.

Gel filtration Chromatography for Detection of mRXRaLBD Dimer and Tetramer

Purified mRXR α LBD dimer (50 μ M in monomer concentration) was incubated with β apo-13-carotenone or UVI 3003 in various concentrations on ice for 3 hours or overnight. Tween 40 was added in order to increase ligand solubility in the aqueous buffer. In other experiments, mRXR α LBD tetramer (50 μ M in monomer concentration) was firstly saturated with 100 μ M β -apo-13-carotenone then was incubated with 9cRA. After ligand treatment mRXR α LBD was subjected to gel filtration chromatography on a Superdex 200 HR column controlled by an AKTA FPLC system (GE). The running buffer contained 20mM Tris, 150mM NaCl, pH7.5 at 4°C. Protein chromatograms were monitored at 280nm. As above, this gel filtration column was also calibrated with proteins of known molecular weight to confirm the retention volumes of the mRXR α LBD dimers and tetramers.

RESULTS

 β -apo-13-Carotenone and UVI 3003 Antagonize 9cRA Induced Transactivation of Full Length RXR α

To investigate the effect of β -apo-13carotenone and UVI 3003 on the functional RXRα. **RXRE-luciferase** role of an receptor/reporter transactivation assay was performed. Full length hRXRa was transiently co-transfected in COS-7 cells with two reporter plasmids. а firefly luciferase reporter containing RXRE from CRBP-II and Renilla luciferase as an internal control. In transfected cells 9cRA induced luciferase activity in a dose-dependent manner over a concentration range of 5 x 10^{-5} M (50 μ M) – 3.2 x 10^{-10} M (0.32 nM), as shown in Figure 1. To determine the antagonist function of β -apo-13-carotenone and UVI 3003, cells were treated with 9cRA in a presence of β -apo-13-carotenone or UVI 3003 at a constant concentration of 200 nM. We observed a shift in the 9cRA dose-response curve induced by both β -apo-13-carotenone and the known antagonist UVI 3003. β-apo-13-Carotenone alone did not induce the activation of RXR α (data not shown). This suggests that β-apo-13-carotenone antagonizes 9cRA activation of full length hRXRa with a similar efficiency as the known antagonist UVI 3003.

β-apo-13-Carotenone and UVI 3003 Inhibit 9cRA Induction of an Endogenous RXRα Responsive Gene

It was previously reported that caspase 9 and Btg2 are direct target genes for RAR-RXR heterodimers (28 - 30). Thus, we asked whether β-apo-13-carotenone and UVI 3003 would inhibit the 9cRA-induced transcription of the endogenous gene caspase 9. For these experiments MCF-7 cells were serum-starved for 24 hours and followed by incubation with 9cRA, β-apo-13-carotenone or UVI 3003 at concentrations of 200 nM for 4 hours. As shown in Figure 2, 9cRA upregulated the expression of mRNA for caspase 9 3-4 fold. Both β-apo-13-carotenone or UVI 3003 inhibited 9cRA-induced gene expression. These data support the results shown above that β -apo-13-carotenone and UVI 3003 antagonize 9cRA induced transactivation of RXRα in transactivation assays.

UVI 3003 Inhibits 9cRA-induced Coactivator Binding to the RXR α Ligand Binding Domain but β -apo-13-Carotenone Does Not

To further characterize the mechanisms of β -apo-13-carotenone and UVI 3003 as antagonists for RXR α , we used cells that stably express a fusion protein containing the Gal4-DBD linked to the ligand binding domain (LBD) of RXRa. The luciferase reporter gene utilized in these assays contains the Gal4 upstream activation sequence (UAS) linked to the luciferase reporter gene. 9cRA activated transcription. however β-apo-13the carotenone alone did not activate the RXRa reporter assay (Figure 3A). Although we tested co-treatment of β -apo-13-carotenone at several different concentrations (5, 10, 100, 200, 500, 1000 nM) with 9cRA, no marked shift of the 9cRA dose-response curve was observed (Figure 3B). In contrast, 200 nM or 500 nM UVI 3003 prominently shifted the 9cRA doseresponse curve, as shown in Figure 3 C. In this assay the activation of RXRa does not require the formation of RXRaLBD dimer. 9cRA binding to the ligand binding domain of RXRa provokes a conformational change of the AF-2 motif that produces a suitable binding surface for recruitment of coactivators. Previous structural studies have shown that the binding of antagonist UVI 3003 to LBD of RXRa disturbs the conformation of helix 12 (H12)

and leads to inhibition of coactivator recruitment (31). Strikingly, in the experiments reported here using the hybrid receptor, β -apo-13-carotenone had no effect on coactivator binding to the RXR α LBD.

β-apo-13-Carotenone Regulates RXRα Through Tetramerization of the Receptor

To further elucidate the mechanism regulation of underlying the RXRα transcriptional activity by β -apo-13-carotenone, we investigated the dimer-tetramer equilibrium of RXRaLBD after exposure to ligand. Mouse RXRaLBD was expressed in E. coli and purified to homogeneity as shown in Figure 4. Gel filtration chromatography on calibrated columns of Superdex 200 was used to isolate the mRXRaLBD dimer and tetramer used in the following experiments. Recombinant mouse RXRaLBD dimer 50 µM (calculated as monomer concentration) was incubated with increasing concentrations (100, 250, 500 µM) of β -apo-13-carotenone on ice for 3.5 hours or overnight. Gel filtration chromatography demonstrated *B*-apo-13-carotenone induced formation of RXRaLBD tetramer (Figure 5). Treatment with 500 µM of β-apo-13carotenone for 3.5 hours, in the molar ratio to monomer receptor of 5:1, led to 33% tetramer formation of the RXRaLBD, whereas if treatment was extended overnight 50% tetramer RXRaLBD formed. In contrast, the antagonist UVI 3003 did not induce tetramer formation at any of the tested concentrations (Figure 6) even if incubations were extended overnight (data not shown). Finally, β-apo-13carotenone saturated RXRaLBD tetramer, in $(\beta$ -apo-13-carotenone: molar ratio 2:1monomer RXRaLBD), was incubated with the agonist 9cRA. RXRaLBD tetramer dissociated to dimer with the addition of agonist 9cRA. At 50 µM, in an equal molar concentration to monomer RXRaLBD, 9cRA induced approximately 55% of dimer, whereas higher concentrations of 9cRA almost completely converted tetramer to dimer (Figure 7). The gel filtration chromatography results showed that β-apo-13-carotenone induced the tetramerization of RXRaLBD, which was reversed with addition of 9cRA. In contrast,

the antagonist UVI 3003 did not influence the tetramer-dimer equilibrium of $RXR\alpha LBD$.

DISCUSSION

In this study we characterize the activity of β -apo-13-carotenone as an antagonist to RXRa and reveal the mechanism of RXRa antagonism by β-apo-13-carotenone-induced tetramerization of the receptor. The comparison of experimental data of B-apo-13carotenone and UVI 3003 indicated that these RXRα antagonists use two distinct β-apo-13-Carotenone, mechanisms. а naturally-occurring β -apocarotenoid which can be obtained from either the diet directly or β-carotene eccentric cleavage from of functioned as an antagonist of RXRa. UVI 3003 is a selective antagonist of RXRa whose inhibitory effect results from an interference of its long side chains with L451 of H12 (31). Both β-apo-13-carotenone and UVI 3003 inhibited 9cRA induced transactivation of full length RXR α in a dual luciferase assay, whereas higher concentrations of 9cRA overcame the inhibition by the two antagonists. Both antagonists inhibited 9cRA induction of the expression of the caspase 9 gene in MCF7 cells.

The tetramerization of RXRa induced by β -apo-13-carotenone is supported by reporter cell based assays. The 'Gal4-DBD:RXRa-LBD' receptor expressed in the reporter cells will bind ligand, translocate to the nucleus, bind to the Gal4 UAS sequence on the reporter gene, recruit co-activator proteins, and lead to the transcription of luciferase. It is important to note that the whole process of transcription in these reporter cells does not require dimerization of RXRa. The conformational change of RXRaLBD due to ligand binding is sufficient to activate coactivator recruitment and subsequent luciferase transcription. β -apo-13-carotenone is inactive in this assay. In contrast, full length RXRα expressed in Cos-7 cells undergoes nuclear translocation, dimer formation upon agonist binding, binding to the RXRE. coactivator recruitment and luciferase

transcription. The distinctive difference in mechanism between the reporter cell assay with Gal4-DBD:RXR α -LBD and the transactivation assay with full length RXR α is that dimer formation is obligatory for the latter; and β -apo-carotenone is only effective as an antagonist in this assay. These observations suggest that β -apo-13-carotenone inhibits 9cRA induced RXR α transcription through the formation of the RXR α tetramer.

In an X-ray crystal study, atRA has been shown to bind to the transcriptionally silent tetrameric RXR α in a unique conformation (32). Previously, we showed using molecular modeling that when this bound atRA is computationally removed from the tetrameric RXR protein and redocked it assumes the identical position as in the crystal structure (26). In addition, when β -apo-13-carotenone is built in a similar conformation to this atRA and docked into this RXRa tetramer it occupies the same position and has the same conformation as the bound atRA. Alternatively, when we built β-apo-13-carotenone in a conformation similar to RXRa bound 9cRA (33), and attempted to dock this molecule into the dimeric RXRa it assumes a very different position than the agonist ligand. Thus we suggested that β -apo-13-carotenone should be capable of acting as an antagonist of RXR α by stabilizing the transcriptionally silent tetramer, but we had no direct biochemical evidence for that suggestion at that time.

The crystal structure study of the ligandbinding domain of the RXRa suggested that a cavity corresponded to the RXR ligandbinding site and that 9cRA binding triggered a conformational modification of H11 which lead to ligand-dependent transactivation by AF-2 (33). It has also been reported that the tetramerization domain is located in helix 11 at the RXRaLBD and tetramerization does not interfere with the function of helix12 (34). Thus β -apo-13-carotenone could cause tetramerization of RXRa by interacting with helix 11, and not affecting helix 12 (or coactivator binding). In contrast, inhibition of 9cRA induced RXRa transcription by UVI 3003 is due to the blockage of helix 12 as pointed out above (31).

RXR α tetramer formation induced by β apo-13-carotenone was confirmed biochemically in the present studies by the observations of gel filtration chromatography with purified recombinant mouse RXRa LBD. Comparison between the β -apo-13-carotenone and UVI 3003 treated dimeric RXRa LBD indicates that β -apo-13-carotenone regulates RXRa transcription through tetramerization, whereas inhibition by the antagonist UVI 3003 is due to interference with helix 12. The complete dissociation of the tetramer RXRa LBD saturated with β -apo-13-carotenone to dimer by 9cRA shows that tetramerization is reversible when the agonist is in sufficient concentration. Thus, the equilibrium of RXRa dimer and tetramer could be controlled by the availability of ligands. A model of these various effects of ligands on RXRa is shown in Fig. 8.

In summary, the present study revealed the mechanism of ligand-dependent regulation of RXRa transcriptional activity by the antagonist, β -apo-13-carotenone. The findings imply that tetramerization of RXR and factors that modulate the oligomer state may contribute to regulation of cellular signaling. βapo-13-Carotenone induced tetramerization could conserve RXRa as an inactive nuclear receptor pool which can rapidly supply dimeric or monomeric RXRa upon 9cRA generation. This may also suggest a ligand-dependent modulation controlling the availability of RXRa for the heterodimerization with other nuclear receptor partners engaged in multiple signaling pathways.

REFERENCES

1. Sakashita, A., Kizaki, M., Pakkala, S., Schiller, G., Tsuruoka, N., Tomosaki, R., Cameron, J.F., Dawson, M.I., Koeffler, H.P. (1993) 9-cis-retinoic acid: effects on normal and leukemic hematopoiesis in vitro. *Blood* **15**, 1009-1016

2. Robertson, K.A., Emami, B., Mueller, L., Collins, S.J. (1992) Multiple members of the retinoic acid receptor family are capable of mediating the granulocytic differentiation of HL-60 cells. *Mol. Cell Biol.* **12**, 3743-3749

3. Sanz, M.J., Albertos, F., Otero, E., Juez, M., Morcillo, E.J., Piqueras, L. (2012) Retinoid X receptor agonists impair arterial mononuclear cell recruitment through peroxisome proliferator-activated receptor- γ activation. *J. Immunol.* **189**, 411-424

4. Qian, L., Zolfaghari, R., Ross, A.C. (2010) Liver-specific cytochrome P450 CYP2C22 is a direct target of retinoic acid and a retinoic acid-metabolizing enzyme in rat liver. *J. Lipid Res.* **51**, 1787-1792

5. Thaller, C., Hofmann, C., Eichele, G. (1993) 9-cis-retinoic acid, a potent inducer of digit pattern duplications in the chick wing bud. *Development* **118**, 957-965

6. Pijnappel, W.W., Hendriks, H.F., Folkers, G.E., van den Brink, C.E., Dekker, E.J., Edelenbosch, C., van der Saag, P.T., Durston, A.J. (1993) The retinoid ligand 4-oxo-retinoic acid is a highly active modulator of positional specification. *Nature* **366**, 340-344

7. Carter, C.J., Farrar, N., Carlone, R.L., Spencer, G.E. (2010) Developmental expression of a molluscan RXR and evidence for its novel, nongenomic role in growth cone guidance. *Dev. Biol.* **343**, 124-137

8. Zhang, X.K., Lehmann, J., Hoffmann, B., Dawson, M.I., Cameron, J., Graupner, G., Hermann, T., Tran, P., Pfahl, M. (1992) Homodimer formation of retinoid X receptor induced by 9-cis retinoic acid. *Nature* **358**, 587-591

9. Zhang, X.K., Hoffmann, B., Tran, P.B., Graupner, G., Pfahl, M. (1992) Retinoid X receptor is an auxiliary protein for thyroid hormone and retinoic acid receptors. *Nature* **355**, 441-446

10. Lefebvre, P., Benomar, Y., Staels, B. (2010) Retinoid X receptors: common hetero-dimerization partners with distinct functions. *Trends Endocrinol. Metab.* **21**, 676-683

11. Evans, R.M., Mangelsdorf, D.J. (2014) Nuclear Receptors, **RXR**, and the Big Bang. *Cell* **157**, 255-266

12. Kersten, S., Kelleher, D., Chambon, P., Gronemeyer, H., and Noy, N. (1995) Retinoid X receptor alpha forms tetramers in solution. *Proc. Natl. Acad. Sci.* **92**, 8645-8649

13. Dawson, M. I., Xia, Z. (2012) The retinoid X receptors and their ligands. *Biochim. Biophy. Acta.* **1821**, 21-56

14. de Lera, A. R., Bourguet, W., Altucci, L., Gronemeyer, H. (2007) Design of selective nuclear receptor modulators : RAR and RXR as a case study. *Nature Rev.* **6**, 811-820

15. Kersten, S., Pan, L., Chambon, P., Gronemeyer, H., and Noy, N. (1995) Role of ligand in retinoid signaling. 9-cis-retinoic acid modulates the oligomeric state of the retinoid X receptor. *Biochem.* **34**, 13717-13721

16. Kersten, S., Dong, D., Lee, W., Reczek, P.R., and Noy, N. (1998) Auto-silencing by the retinoid X receptor. *J. Mol. Biol.* **284**, 21-32

17. Mangelsdorf, D.J., Borgmeyer, U., Heyman, R.A., Zhou, J.Y., Ong, E.S., Oro, A.E., Kakizuka, A., Evans, R.M. (1992) Characterization of three RXR genes that mediate the action of 9-cis retinoic acid. *Genes Dev.* **6**, 329-344

18. Heyman, R. A., Mangelsdorf, D. J., Dyck, J.A., Stein, R. B., Eichele, G., Evans, R. M., Thaller, C. (1992) 9-cis retinoic acid is a high affinity ligand for the retinoid X receptor. *Cell* **68**, 397-406

19. Levin, A. A, Sturzenbecker, L. J., Kazmer, S., Bosakowski, T., Huselton, C., Allenby, G., Speck, J., Kratzeisen, C., Rosenberger, M., Lovey, A., and Grippo, J. F. (1992) 9-cis retinoic acid stereoisomer binds and activates the nuclear receptor RXR alpha. *Nature* **355**, 359-361

20. Olson, J. A., and Krinsky, N. I. (1995) Introduction: the colorful, fascinating world of the carotenoids: important physiologic modulators. *FASEB J.* **9**, 1547-1550

21. Parker, R. S. (1996) Absorption, metabolism, and transport of carotenoids. FASEB J. 10, 542-551

22. Harrison, E. H. (2005) Mechanisms of digestion and absorption of dietary vitamin A. Ann. Rev. Nutr. 25, 87-103

23. Fleshman, M. K., Lester, G. E., Riedl, K. M., Kopec, R. E., Narayanasamy, S., Curley, R. W. Jr, Schwartz, S. J., Harrison, E. H. (2011) Carotene and novel apocarotenoid concentrations in orange-fleshed Cucumis melo melons: determinations of β -carotene bioaccessibility and bioavailability. *J. Agric. Food Chem.* **59**, 4448-4454

24. Ho, C. C., de Moura, F. F., Kim, S. H., Clifford, A. J. (2007) Excentral cleavage of beta-carotene in vivo in a healthy man. *Am. J. Clin. Nutr.* **85**, 770-777

25. Shmarakov, I, Fleshman, M. K., D'Ambrosio, D. N., Piantedosi, R., Riedl, K. M., Schwartz, S. J., Curley, R. W. Jr, von Lintig, J., Rubin, L. P., Harrison, E. H., Blaner, W. S. (2012) Hepatic stellate cells are an important cellular site for β -carotene conversion to retinoid. *Arch. Biochem. Biophys.* **504**, 3-10

26. Eroglu, A., Hruszkewycz, D. P., Curley, R. W. Jr, Harrison, E. H. (2010) The eccentric cleavage product of β -carotene, β -apo-13-carotenone, functions as an antagonist of RXR α . Arch. Biochem. Biophys. **504**, 11-16

27. Eroglu, A, Hruszkewycz, D. P, dela Sena C., Narayanasamy, S., Riedl, K. M., Kopec, R. E., Schwartz, S. J., Curley, R. W. Jr, Harrison, E. H. (2012) Naturally occurring eccentric cleavage products of provitamin A β -carotene function as antagonists of retinoic acid receptors. *J. Biol. Chem.* **287**, 15886-15895

28. Donato, L. J., Noy, N. (2005) Suppression of mammary carcinoma growth by retinoic acid: proapoptotic genes are targets for retinoic acid receptor and cellular retinoic acid-binding protein II signaling. *Cancer Res.* **65**, 8193-8199

29. Donato, L. J., Suh, J. H., Noy, N. (2007) Suppression of mammary carcinoma cell growth by retinoic acid: the cell cycle control gene Btg2 is a direct target for retinoic acid receptor signaling. *Cancer Res.* **67**, 609-615

30. Yasmin, R., Kannan-Thulasiraman, P., Kagechika, H., Dawson, M. I., Noy, N. (2010) Inhibition of mammary carcinoma cell growth by RXR is mediated by the receptor's oligomeric switch. *J. Mol. Biol.* **397**, 1121-1131

31. Nahoum, V., Pérez, E., Germain, P., Rodríguez-Barrios, F., Manzo, F., Kammerer, S., Lemaire, G., Hirsch, O., Royer, C.A., Gronemeyer, H., de Lera, A. R., Bourguet, W. (2007) Modulators of the structural dynamics of the retinoid X receptor to reveal receptor function. *Proc. Natl. Acad. Sci.* **104**, 17323-17328

32. Gampe, R. T., Montana, V. G., Lambert, M. H., Wisely, G. B., Milburn, M. V., and Xu, H. E. (2000) Structural basis for autorepression of retinoid X receptor by tetramer formation and the AF-2 helix. *Genes Dev.* **14**, 2229-2241

33. Bourguet, W., Ruff, M., Chambon, P., Gronemeyer, H., Moras, D. (1995) Crystal structure of the ligand-binding domain of the human nuclear receptor RXR-alpha. *Nature* **375**, 377-382

34. Kersten, S., Reczek, P. R., Noy, N. (1997) The tetramerization region of the retinoid X receptor is important for transcriptional activation by the receptor. *J. Biol. Chem.* **272**, 29759-29768

Acknowledgments-The authors would like to thank Dr. Noa Noy of Case Western Reserve University for the mRXRαLBD-pET-28a plasmid.

FOOTNOTES

This work was supported by NIH Grant R01-HL049879 (National Heart, Lung and Blood Institute and the NIH Office of Dietary Supplements).

The abbreviations used are: RXR-retinoid x receptor, 9cRA-9-cis-retinoic acid, LBD-ligand binding domain, DBD-DNA binding domain, RXRE –retinoid x response element, MCF-7-Michigan Cancer Foundation-7 (a human mammary cancer cell line),

Figure Legends

Figure 1. The antagonist effects of β -apo-13-carotenone and UVI 3003 on transactivation of full length RXR α . COS-7 cells were co-transfected with full length hRXR α , pRL-tk-luc(Renilla) and RXRE-luc(firefly). Twenty-four hours after transfection, cells were treated with 9cRA at concentrations of 0, 0.064, 0.32, 1.6, 8, 40, 200, 1000, 5000,25000, 50000 nM in presence or absence of 200 nM β -apo-13-carotenone or UVI 3003 for another 24 hours in 37 °C incubator. 9cRA alone (Blue), 9cRA plus β -apo-13-carotenone at 200 nM (Red), 9cRA plus UVI 3003 a 200 nM (Green). Luciferase activity is shown on the y-axis.

Figure 2. β -apo-13-carotenone and UVI 3003 inhibit 9cRA-induced expression of mRNA for caspase 9. MCF7 cells were cultured in 6-well plate and serum-starved for 24 hours before ligand treatment. Cells were incubated with 200 nM ligands (9cRA, β -apo-13-carotenone or UVI 3003) in serum-free medium for 4 hour. The comparative Ct method ($\Delta\Delta$ Ct) was used to analyze results. Data are presented as means ± SE, n=5. Antagonist treatment (c) significantly inhibits 9cRA induced Casp9 gene expression (b) at P < 0.006).

Figure 3. β -apo-13-carotenone does not antagonize 9cRA induced transactivation in cells expressing Gal4-DBD:RXR α -LBD and a UAS-driven luciferase reporter. Cells were incubated with 9cRA or β -apo-13-carotenone at concentrations of 0, 0.32, 1.6, 8, 40, 200, 1000, and 5000 nM for 24 hours at 37 ⁰ C (A); cells were incubated with these same concentrations of 9cRA in presence or absence of fixed concentrations of β -apo-13-carotenone (B) or UVI 3003(C).

Figure 4. Purification of recombinant mRXR α LBD. Histidine tagged mRXR α LBD (theoretical molecular mass for monomer = 28,821 Da) was expressed in *E. coli* strain BL21(DE3) and purified firstly with a Ni-NTA affinity column. The elution fraction 1 of the affinity column was then applied onto a gel filtration column of Hiload Superdex200 16x60 column to separate the mRXR α LBD tetramer and dimer. A, SDS-PAGE of the fractions from Ni-NTA affinity column. Lane 9 was the mRXR α LBD elution fraction collected for second step purification with gel filtration. B, Separation of mRXR α LBD tetramer and dimer with gel filtration column. mRXR α LBD tetramer was collected between the retention volume of 61.4 ml to 68.4 ml. mRXR α LBD dimer was collected between retention volume 72.4 ml to 84.4 ml. The gel filtration column was calibrated with standard proteins of known molecular weight as described in the text.

Figure 5. Gel filtration analysis for tetramerization of RXR α LBD induced by β -apo-13-carotenone. Purified mRXR α LBD dimer (50 μ M in monomer concentration) was incubated with β -apo-13-carotenone in concentrations of 0, 100, 250, 500 μ M on ice for 3 hours or overnight. Tween 40 of 0.75% (w/v) was added in order to increase ligand solubility in the aqueous buffer. Superdex 200 HR column and elution buffer were well equilibrated at 4 °C with elution buffer consisted of 20 mM Tris, 150 mM NaCl, and 3 mM DTT, pH 7.4. Protein chromatograms were monitored at 280nm.

Figure 6. UVI 3003 does not induce tetramerization of RXR α LBD. Purified mRXR α LBD dimer (50 μ M in monomer concentration) was incubated with UVI 3003 in concentrations of 0, 250, 500 μ M on ice for 3 hours. Superdex 200 HR column and elution buffer were well equilibrated at 4 °C with elution buffer consisted of 20 mM Tris, 150 mM NaCl, and 3 mM DTT, pH 7.4. Protein chromatograms were monitored at 280nm. Extending the incubation time with 500 μ M UVI 3003 overnight did not lead to tetramer formation.

Figure 7. 9cRA dissociates the tetramer of RXR α LBD induced by β -apo-13-carotenone. mRXR α LBD tetramer (50 μ M in monomer concentration) was firstly saturated with 100 μ M β -apo-13-carotenone then was incubated with 9cRA in concentrations of 0, 50, 100, 250 μ M on ice for 3 hours. Protein chromatograms were monitored at 280nm.

Figure 8. Model depicting the effects of ligands on the oligomeric state and transcriptional activity of $RXR\alpha$. See text for details.













С

Figure 4A



1	2	3	4	5	6	7	8	9	10
Maker	Whole	Pellet	Super-	Flow	Wash	Wash	Wash	Fraction	Fraction
	lysate		natant	through	1	2	3	1	2

Figure 4B



Figure 5





Figure 6





Figure 8





Metabolism:

β-apo-13-Carotenone Regulates Retinoid X Receptor Transcriptional Activity Through Tetramerization of the Receptor



Jian Sun, Sureshbabu Narayanasamy, Robert W. Curley, Jr. and Earl H. Harrison J. Biol. Chem. published online October 16, 2014

Access the most updated version of this article at doi: 10.1074/jbc.M114.610501

Find articles, minireviews, Reflections and Classics on similar topics on the JBC Affinity Sites.

Alerts:

- · When this article is cited
- When a correction for this article is posted

Click here to choose from all of JBC's e-mail alerts

This article cites 0 references, 0 of which can be accessed free at http://www.jbc.org/content/early/2014/10/16/jbc.M114.610501.full.html#ref-list-1