

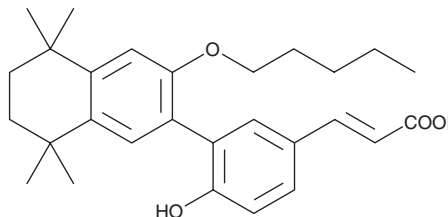
PRODUCT INFORMATION



UVI3003

Item No. 16026

CAS Registry No.: 847239-17-2
Formal Name: 3-[4-hydroxy-3-[5,6,7,8-tetrahydro-5,5,8,8-tetramethyl-3-(pentyloxy)-2-naphthalenyl]phenyl]-2-propenoic acid
MF: C₂₈H₃₆O₄
FW: 436.6
Purity: ≥90%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

UVI3003 is supplied as a crystalline solid. A stock solution may be made by dissolving the UVI3003 in the solvent of choice, which should be purged with an inert gas. UVI3003 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of UVI3003 in these solvents is approximately 2, 1, and 0.5 mg/ml, respectively.

Description

UVI3003 is a full antagonist of RXR that demonstrates potent, nanomolar binding affinity.¹ At 1 μM, UVI3003 does not affect the corepressor interaction capacity of the RARα subunit in the RAR-RXR heterodimer configuration.¹ This compound can be used to isolate the contribution of RXR to the function of RXR heterodimers.^{1,2}

References

1. Nahoum, V., Pérez, E., Germain, P., *et al.* Modulators of the structural dynamics of the retinoid X receptor to reveal receptor function. *Proc. Natl. Acad. Sci. USA* **104(44)**, 17323-17328 (2007).
2. le Maire, A., Grimaldi, M., Roecklin, D., *et al.* Activation of RXR-PPAR heterodimers by organotin environmental endocrine disruptors. *EMBO reports* **10(4)**, 367-373 (2009).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/10/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM