

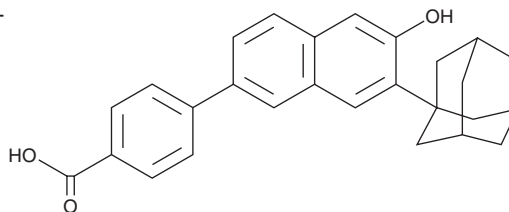
PRODUCT INFORMATION



CD1530

Item No. 19062

CAS Registry No.: 107430-66-0
Formal Name: 4-(6-hydroxy-7-tricyclo[3.3.1.1^{3,7}]dec-1-yl-2-naphthalenyl)-benzoic acid
MF: C₂₇H₂₆O₃
FW: 398.5
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 277, 325 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CD1530 is supplied as a crystalline solid. A stock solution may be made by dissolving the CD1530 in the solvent of choice, which should be purged with an inert gas. CD1530 is soluble at 1 molar equivalent NaOH to 50 mM. We do not recommend storing the aqueous solution for more than one day.

Description

CD1530 is a selective retinoic acid receptor (RAR γ) agonist (K_ds = 150, 1,500, and 2,750 nM for RAR γ , RAR β , and RAR α receptors, respectively).¹ It demonstrates transcriptional activity at RAR γ with an AC₅₀ value of 1.8 nM.¹ It has been used in combination with bexarotene (Item No. 11571), a retinoid X receptor agonist, to inhibit oral carcinogenesis induced by the carcinogen 4-nitroquinoline 1-oxide in a mouse model of human oral-cavity and esophageal squamous-cell carcinoma.²

References

1. Bernard, B.A., Bernardon, J.-M., Delescluse, C., *et al.* Identification of synthetic retinoids with selectivity for human nuclear retinoic acid receptor γ . *Biochem. Biophys. Res. Commun.* **186(2)**, 977-983 (1992).
2. Tang, X.-H., Osei-Sarfo, K., Urvalek, A.M., *et al.* Combination of bexarotene and the retinoid CD1530 reduces murine oral-cavity carcinogenesis induced by the carcinogen 4-nitroquinoline 1-oxide. *Proc. Natl. Acad. Sci. USA* **111(24)**, 8907-8912 (2014).

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, 11/11/2021

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