

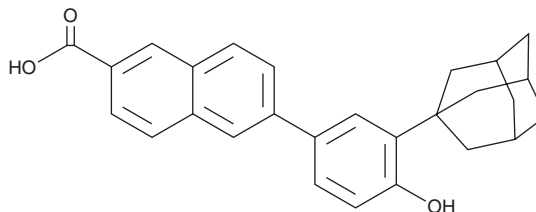
PRODUCT INFORMATION



CD437

Item No. 71610

CAS Registry No.: 125316-60-1
Formal Name: 6-(4-hydroxy-3-tricyclo[3.3.1.1^{3,7}]-naphthalenecarboxylic acid
Synonym: AHPN
MF: C₂₇H₂₆O₃
FW: 398.5
Purity: ≥95%
UV/Vis.: λ_{max}: 234, 274, 325 nm
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

CD437 is supplied as a crystalline solid. A stock solution may be made by dissolving the CD437 in the solvent of choice, which should be purged with an inert gas. CD437 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CD437 in these solvents is approximately 1, 16, and 10 mg/ml.

CD437 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CD437 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CD437 has a solubility of approximately 0.05 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CD437 is the prototypical adamantyl arotinoid of the retinoid-related molecule family that acts as a selective agonist of retinoic acid receptor (RAR)γ (K_ds = 6.5 μM, 2.5 μM, and 77 nM for RARα, β, and γ, respectively).^{1,2} CD437 is cytotoxic to the acute myeloid leukemia cell line NB4 (EC₅₀ = 0.17 μM), inducing DNA damage by producing DNA double strand breaks.³ CD437 also induces cell cycle arrest and apoptosis in various other cancer cells including melanoma, breast, non-small lung, and prostate cancer cells through RAR-dependent and -independent signaling pathways.^{4,5}

References

1. Pérez-Rodríguez, S., Ortiz, M.A., Pereira, R., *et al.* Highly twisted adamantyl arotinoids: Synthesis, antiproliferative effects and RXR transactivation profiles. *Eur. J. Med. Chem.* **44(6)**, 2434-2446 (2009).
2. 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).
3. Valli, C., Paroni, G., Di Francesco, A.M., *et al.* Atypical retinoids ST1926 and CD437 are S-phase-specific agents causing DNA double-strand breaks: Significance for the cytotoxic and antiproliferative activity. *Mol. Cancer Ther.* **7(9)**, 2941-2954 (2008).
4. Fontana, J.A. and Rishi, A.K. Classical and novel retinoids: Their targets in cancer therapy. *Leukemia* **16(4)**, 463-472 (2002).
5. Jin, F., Liu, X., Zhou, Z., *et al.* Activation of nuclear factor-κB contributes to induction of death receptors and apoptosis by the synthetic retinoid CD437 in DU145 human prostate cancer cells. *Cancer Res.* **65(14)**, 6354-6363 (2005).

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, 11/21/2022

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