

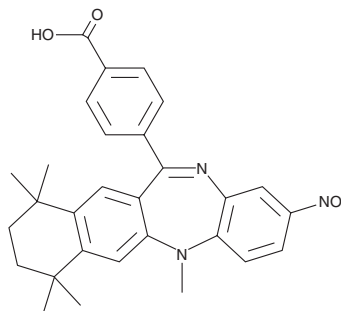
PRODUCT INFORMATION



HX 531

Item No. 20762

CAS Registry No.: 188844-34-0
Formal Name: 4-(7,8,9,10-tetrahydro-5,7,7,10,10-pentamethyl-2-nitro-5H-benzo[b]naphtho[2,3-e][1,4]diazepin-12-yl)-benzoic acid
MF: C₂₉H₂₉N₃O₄
FW: 483.6
Purity: ≥95%
UV/Vis.: λ_{max}: 264 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

HX 531 is supplied as a crystalline solid. A stock solution may be made by dissolving the HX 531 in the solvent of choice, which should be purged with an inert gas. HX 531 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HX 531 in these solvents is approximately 0.12, 10, and 15 mg/ml, respectively.

HX 531 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HX 531 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HX 531 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HX 531 is an antagonist of the retinoid X receptor (RXR) with an IC₅₀ value of 18 nM that inhibits 9-*cis* retinoic acid (Item No. 14587)-induced transactivation of RXR.¹ It has been shown to reduce triglyceride content in white adipose tissue, skeletal muscle, and the liver of mice on a high fat diet.²

References

1. Konta, T., Xu, Q., Furusu, A., *et al.* Selective roles of retinoic acid receptor and retinoid X receptor in the suppression of apoptosis by all-*trans*-retinoic acid. *J. Biol. Chem.* **276**(16), 12697-12701 (2001).
2. Yamauchi, T., Waki, H., Kamon, J., *et al.* Inhibition of RXR and PPARγ ameliorates diet-induced obesity and type 2 diabetes *J. Clin. Invest.* **108**(7), 1001-1013 (2001).

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, 11/29/2022

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