

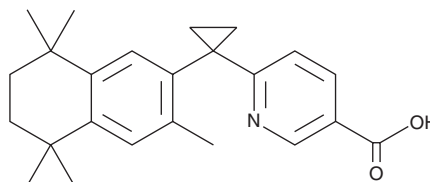
PRODUCT INFORMATION



LG 100268

Item No. 21606

CAS Registry No.: 153559-76-3
Formal Name: 6-[1-(5,6,7,8-tetrahydro-3,5,5,8,8-pentamethyl-2-naphthalenyl)cyclopropyl]-3-pyridinecarboxylic acid
Synonyms: AGN 192620, CD3127
MF: C₂₄H₂₉NO₂
FW: 363.5
Purity: ≥98%
UV/Vis.: λ_{max}: 271, 318 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

LG 100268 is supplied as a crystalline solid. A stock solution may be made by dissolving the LG 100268 in the solvent of choice. LG 100268 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of LG 100268 in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

LG 100268 is an agonist of retinoid X receptors (RXRs; $K_d = 3 \text{ nM}$).¹ It is selective for RXRs over retinoic acid receptors (RARs; $K_d = >10,000 \text{ nM}$).² *In vitro*, LG 100268 (1 μM) induces apoptosis in HL-60 human leukemia cells when used in combination with TTNPB (Item No. 16144).¹ LG 100268 (100, 300, and 1,000 nM) dose-dependently inhibits COX-2 expression and increases ATP-binding cassette transporter (ABCA1) levels in macrophage-like RAW 264.7 cells.³ *In vivo*, it inhibits vinyl carbamate-induced lung tumor growth in mice when administered at a dose of 40 mg/kg.

References

- Boehm, M.F., Zhang, L., Zhi, L., *et al.* Design and synthesis of potent retinoid X receptor selective ligands that induce apoptosis in leukemia cells. *J. Med. Chem.* **38**(16), 3146-3155 (1995).
- Bissonnette, R.P., Brunner, T., Lazarchik, S.B., *et al.* 9-*cis* retinoic acid inhibition of activation-induced apoptosis is mediated via regulation of fas ligand and requires retinoic acid receptor and retinoid X receptor activation. *Mol. Cell. Biol.* **15**(10), 5576-5585 (1995).
- Cao, M., Royce, D.B., Risingsong, R., *et al.* The rexinoids LG100268 and LG101506 inhibit inflammation and suppress lung carcinogenesis in A/J mice. *Cancer Prev. Res. (Phila)*. **9**(1), 105-114 (2016).

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, 10/24/2022

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