

PRODUCT INFORMATION

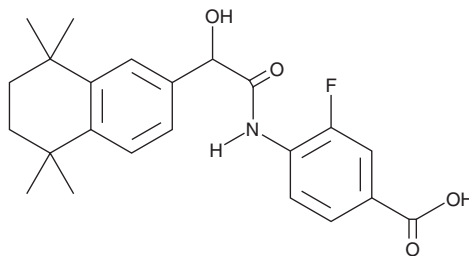


BMS 189961

Item No. 29380

CAS Registry No.: 185629-22-5
Formal Name: 3-fluoro-4-[[2-hydroxy-2-(5,6,7,8-tetrahydro-5,5,8,8-tetramethyl-2-naphthalenyl)acetyl]amino]benzoic acid

Synonym: BMS 961
MF: C₂₃H₂₆FNO₄
FW: 399.5
Purity: ≥98%
UV/Vis.: λ_{max}: 268 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

BMS 189961 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 189961 in the solvent of choice, which should be purged with an inert gas. BMS 189961 is soluble in the organic solvent DMSO.

Description

BMS 189961 is an agonist of retinoic acid receptor γ (RAR γ ; EC₅₀ = 30 nM in a transactivation assay).¹ It is selective for RAR γ over RAR β (EC₅₀ = 1,000 nM). *In vitro*, BMS 189961 decreases long bone growth and digit formation of isolated mouse embryonic limb buds when used at concentrations of 0.01 and 0.1 μ M.² Topical administration of BMS 189661 (40 nmol) increases the skin expression of retinoid metabolism-related genes, including *Rdh10*, *Crabp2*, and *Krt4*, as well as epidermal barrier genes, including *Spink5*, *Klk5*, and *Klk7*, in mice.³

References

1. Klaholz, B.P., Mitschler, A., Belema, M., *et al.* Enantiomer discrimination illustrated by high-resolution crystal structures of the human nuclear receptor hRAR γ . *Proc. Natl. Acad. Sci. USA* **97(12)**, 6322-6327 (2000).
2. Galdones, E. and Hales, B.F. Retinoic acid receptor gamma-induced misregulation of chondrogenesis in the murine limb bud *in vitro*. *Toxicol. Sci.* **106(1)**, 223-232 (2008).
3. Gericke, J., Ittensohn, J., Mihály, J., *et al.* Regulation of retinoid-mediated signaling involved in skin homeostasis by RAR and RXR agonists/antagonists in mouse skin. *PLoS One* **8(4)**, e62643 (2013).

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, 10/04/2022

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