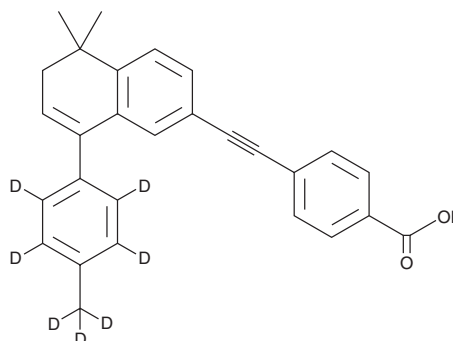


PRODUCT INFORMATION



AGN 193109-d₇
Item No. 23976

CAS Registry No.: 1216429-25-2
Formal Name: 4-[2-[5,6-dihydro-5,5-dimethyl-8-(4-methylphenyl-d₇)-2-naphthalenyl]ethynyl]-benzoic acid
MF: C₂₈H₁₇D₇O₂
FW: 399.5
Chemical Purity: ≥98% (AGN 193109)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
Supplied as: A 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

AGN 193109-d₇ is intended for use as an internal standard for the quantification of AGN 193109 (Item No. 23975) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

AGN 193109-d₇ is supplied as a solid. A stock solution may be made by dissolving the AGN 193109-d₇ in the solvent of choice, which should be purged with an inert gas. AGN 193109-d₇ is slightly soluble in DMSO and methanol.

Description

AGN 193109 is an antagonist of retinoic acid receptors (RARs).¹ AGN 193109 decreases expression of cytokeratin K5-8, 13, 14, 16, 17, and 19 genes, markers of retinoid action in ECE16-1 cells, when co-administered with TTNPB (Item No. 16144) but not when used alone. *In vivo*, AGN 193109 induces cleft palate or frontonasal dysplasia and eye malformations in fetuses of pregnant mice following a single oral dose of 1 mg/kg.²

References

1. Agarwal, C., Chandraratna, A.S., Johnson, A.L., *et al.* AGN193109 is a highly effective antagonist of retinoid action in human ectocervical epithelial cells. *J. Biol. Chem.* **271(21)**, 12209-12212 (1996).
2. Kochhar, D.M., Jiang, H., Penner, J.D., *et al.* The use of a retinoid receptor antagonist in a new model to study vitamin A-dependent developmental events. *Int. J. Dev. Biol.* **42(4)**, 601-608 (1998).

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, 11/16/2022

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