

PRODUCT INFORMATION



13-*cis*-Retinoic Acid-d₅

Item No. 21855

Formal Name: 9-(6,6-dimethyl-2-(methyl-d₃)cyclohex-1-en-1-yl-3,3-d₂)-3,7-dimethylnona-2Z,4E,6E,8E-tetraenoic acid

Synonyms: Accutane-d₅, Isotretinoin-d₅, 13-*cis*-Vitamin A Acid-d₅

MF: C₂₀H₂₃D₅O₂

FW: 305.5

Chemical Purity: 90% (13-*cis*-Retinoic Acid)

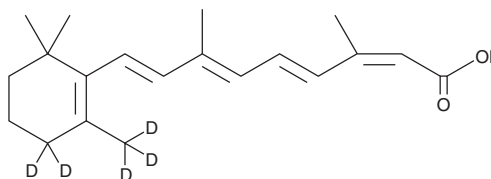
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀

Supplied as: A solid

Storage: -80°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

13-*cis*-Retinoic acid-d₅ is intended for use as an internal standard for the quantification of 13-*cis*-retinoic acid 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).

13-*cis*-Retinoic acid-d₅ is supplied as a solid. A stock solution may be made by dissolving the 13-*cis*-retinoic acid-d₅ in the solvent of choice, which should be purged with an inert gas. 13-*cis*-retinoic acid-d₅ is soluble in methanol.

Description

13-*cis*-Retinoic acid is a naturally occurring retinoid that inhibits retinol oxidation and dehydrogenation of 17β-hydroxy steroids.¹

Reference

1. Blaner, W.S. Cellular metabolism and actions of 13-*cis*-retinoic acid. *J. Am. Acad. Dermatol.* **45(5)**, S129-S135 (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, 11/10/2022

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