

PRODUCT INFORMATION

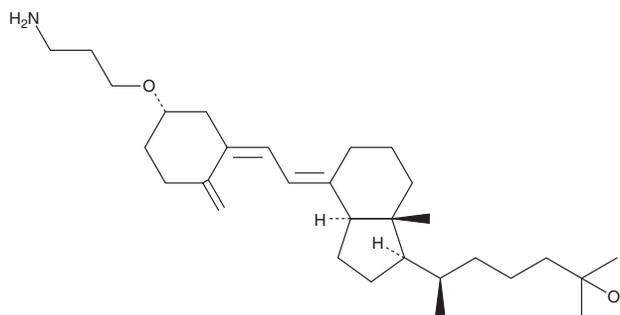


25-hydroxy Vitamin D₃ 3,3'-aminopropyl ether

Item No. 23122

CAS Registry No.: 163018-26-6
Formal Name: (εR,1R,3aS,4E,7aR)-4-[(2Z)-2-[(5S)-5-(3-aminopropoxy)-2-methylenecyclohexylidene]ethylidene]octahydro-α,α,ε,7a-tetramethyl-1H-indene-1-pentanol

MF: C₃₀H₅₁NO₂
FW: 457.7
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 265 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

25-hydroxy Vitamin D₃ 3,3'-aminopropyl ether is supplied as a crystalline solid. A stock solution may be made by dissolving the 25-hydroxy vitamin D₃ 3,3'-aminopropyl ether in the solvent of choice, which should be purged with an inert gas. 25-hydroxy Vitamin D₃ 3,3'-aminopropyl ether is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 25-hydroxy vitamin D₃ 3,3'-aminopropyl ether in these solvents is approximately 20, 10, and 25 mg/ml, respectively.

25-hydroxy Vitamin D₃ 3,3'-aminopropyl ether is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 25-hydroxy vitamin D₃ 3,3'-aminopropyl ether should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 25-hydroxy Vitamin D₃ 3,3'-aminopropyl ether has a solubility of approximately 0.3 mg/ml in a 1:12 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

25-hydroxy Vitamin D₃ 3,3'-aminopropyl ether is an analog of vitamin D₃ (Item No. 11792).¹ It contains a highly reactive amine group that can be coupled to a sepharose matrix for affinity purification of vitamin D binding protein. 25-hydroxy Vitamin D₃ 3,3'-aminopropyl ether also inhibits 25-hydroxy vitamin D₃-1α-hydroxylase.

Reference

1. Roy, A. and Ray, R. Aminopropylation of vitamin D hormone (1 α,25-dihydroxyvitamin D₃), its biological precursors, and other steroidal alcohols: An anchoring moiety for affinity studies of sterols. *Steroids* **60**(8), 530-533 (1995).

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, 11/15/2022

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