

# PRODUCT INFORMATION



## 3-*epi*-25-hydroxy Vitamin D<sub>3</sub>

Item No. 34200

**CAS Registry No.:** 73809-05-9  
**Formal Name:** (3aS)-octahydro-4E-[(2Z)-2-[(5R)-5-hydroxy-2-methylenecyclohexylidene]ethylidene]-α,α,εR,7aR-tetramethyl-1H-indene-1R-pentanol

**Synonyms:** 3-*epi*-25(OH)D<sub>3</sub>,  
3-*epi*-25-hydroxy Cholecalciferol

**MF:** C<sub>27</sub>H<sub>44</sub>O<sub>2</sub>

**FW:** 400.6

**Purity:** ≥95%

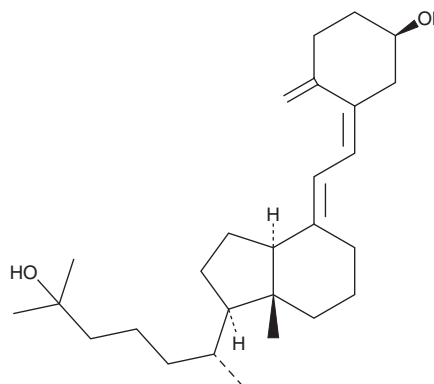
**Supplied as:** A solid

**Storage:** -80°C

**Stability:** ≥2 years

**Special Conditions:** Light sensitive

**Item Origin:** Synthetic



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

### Laboratory Procedures

3-*epi*-25-hydroxy Vitamin D<sub>3</sub> is supplied as a solid. A stock solution may be made by dissolving the 3-*epi*-25-hydroxy vitamin D<sub>3</sub> in the solvent of choice, which should be purged with an inert gas. 3-*epi*-25-hydroxy Vitamin D<sub>3</sub> is slightly soluble in chloroform and ethyl acetate.

### Description

3-*epi*-25-hydroxy Vitamin D<sub>3</sub> is the C-3 epimer of 25-hydroxy vitamin D<sub>3</sub> (Item No. 9000683).<sup>1</sup> Dietary administration of 3-*epi*-25-hydroxy vitamin D<sub>3</sub> (0.5 and 1 IU/g) decreases levels of serum parathyroid hormone (PTH) in male, but not female, weanling rats.

### Reference

1. Djekic-Ivankovic, M., Lavery, P., Agellon, S., *et al.* The C-3α epimer of 25-hydroxycholecalciferol from endogenous and exogenous sources supports normal growth and bone mineral density in weanling rats. *J. Nutr.* **147**(2), 141-151 (2017).

#### 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, 06/29/2021

#### 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