

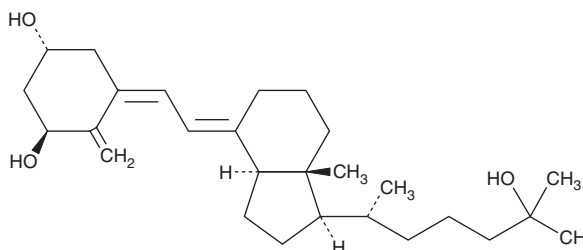
PRODUCT INFORMATION



Calcitriol

Item No. 71820

CAS Registry No.: 32222-06-3
Formal Name: 9,10-secocholesta-5Z,7E,10(19)-triene-1 α ,3 β ,25-triol
Synonym: 1 α ,25-dihydroxy Vitamin D₃
MF: C₂₇H₄₄O₃
FW: 416.6
Purity: \geq 97%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



Special Conditions: Product is light sensitive and air sensitive. Handle under inert conditions.

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

Laboratory Procedures

Calcitriol is supplied as a crystalline solid. A stock solution may be made by dissolving the calcitriol in the solvent of choice, which should be purged with an inert gas. Calcitriol is soluble in organic solvents such as ethanol, methanol, and DMSO. The solubility of calcitriol in ethanol is approximately 1 mg/ml and approximately 50 mg/ml in methanol and DMSO.

Description

Calcitriol is synthesized from 7-dehydro cholesterol in humans *via* a non-enzymatic photochemical reaction with 290-310 nm UV light in the skin.¹ Hydroxylation of the resulting cholecalciferol in the liver produces 25-hydroxy vitamin D₃, the principal circulating form of vitamin D. A second, tightly regulated hydroxylation in the kidney produces calcitriol. Plasma calcitriol levels range from 10-70 pg/ml and are influenced by numerous dietary and hormonal factors.² The main physiologic effects of calcitriol are to increase the absorption of calcium at the level of the intestinal epithelium, and to increase the mineralization of bone *via* the direct stimulation of osteoblasts.³

References

1. Kumar, R. Metabolism of 1,25-dihydroxyvitamin D₃. *Physiol. Rev.* **64**(2), 478-505 (1984).
2. Bikle, D.D., Gee, E., Halloran, B., *et al.* Free 1,25-dihydroxyvitamin D levels in serum from normal subjects, pregnant subjects, and subjects with liver disease. *J. Clin. Invest.* **74**(6), 1966-1971 (1984).
3. Portale, A.A., Halloran, B.P., and Morris, R.C., Jr. Physiologic regulation of the serum concentration of 1,25-dihydroxyvitamin D by phosphorus in normal men. *J. Clin. Invest.* **83**(5), 1494-1499 (1989).

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, 10/30/2020

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