

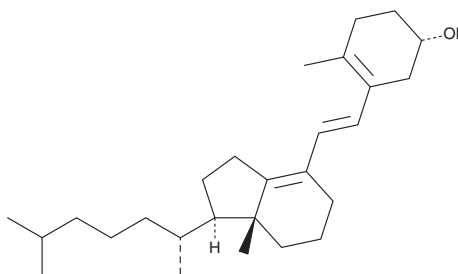
PRODUCT INFORMATION



Isotachysterol 3

Item No. 35751

CAS Registry No.: 22350-43-2
Formal Name: (3 β ,6E)-9,10-secocholesta-5(10),6,8(14)-trien-3-ol
MF: C₂₇H₄₄O
FW: 384.6
Purity: \geq 98%
Supplied as: A solution in chloroform
Storage: -20°C
Stability: \geq 2 years
Item Origin: Synthetic



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

Laboratory Procedures

Isotachysterol 3 is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. Isotachysterol 3 is slightly soluble in acetone, methanol, and ethyl acetate.

Description

Isotachysterol 3 is an impurity in the preparation of vitamin D₃ (Item No. 17890).¹ It stimulates intestinal calcium transport and bone calcium mobilization in control, as well as in anephric, rats fed a low-calcium and vitamin D-deficient diet when administered at a dose of 5 μ g/animal.²

References

1. Tartivita, K.A., Sciarello, J.P., and Rudy, B.C. High-performance liquid chromatographic analysis of vitamins I: Quantitation of cholecalciferol or ergocalciferol in presence of photochemical isomers of the provitamin and application to cholecalciferol resins. *J. Pharm. Sci.* **65(7)**, 1024-1027 (1976).
2. Holick, M.F., DeLuca, H.F., Kasten, P.M., *et al.* Isotachysterol₃ and 25-hydroxyisotachysterol₃: Analogs of 1,25-dihydroxvitamin D₃. *Science* **180(4089)**, 964-966 (1973).

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, 12/13/2022

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