

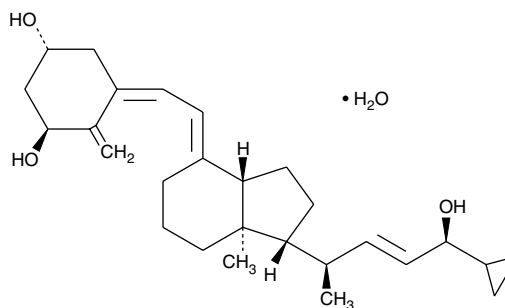
# Product Information



## Calcipotriol (hydrate)

Item No. 10009599

**CAS Registry No.:** 112828-00-9  
**Formal Name:** 24-cyclopropyl-(1 $\alpha$ ,3 $\beta$ ,5Z,7E,22E,24S)-9,10-secochola-5,7,10(19),22-tetraene-1,3,24-triol, monohydrate  
**Synonym:** Calcipotriene  
**MF:** C<sub>27</sub>H<sub>40</sub>O<sub>3</sub> • H<sub>2</sub>O  
**FW:** 412.6  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that calcipotriol (hydrate) be stored as supplied at -20°C. It will be stable for at least two years.

Calcipotriol (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the calcipotriol (hydrate) in an organic solvent purged with an inert gas. Calcipotriol (hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of calcipotriol (hydrate) in these solvents is approximately 50 mg/ml.

Calcipotriol (hydrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Calcipotriol (hydrate) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Calcipotriol (hydrate) has a solubility of approximately 0.15 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Calcipotriol (hydrate) is a low-calcemic vitamin D receptor (VDR) agonist. Calcipotriol (hydrate) is about 200 times less potent in its effects on calcium metabolism than vitamin D (1,25[OH]D<sub>3</sub>).<sup>1</sup> Binding of calcipotriol (hydrate) to the VDR increases AP-1, a transcription factor important for keratinocyte differentiation, and reduces expression of JunB protein, a transcriptional activator in the inflammatory response.<sup>2</sup> Calcipotriol (hydrate) also induces expression of thymic stromal lymphopoietin, which triggers T-cell differentiation in keratinocytes.<sup>3</sup>

### References

1. Kragballe, K. Vitamin D analogues in the treatment of psoriasis. *J. Cell. Biochem.* **49**, 46-52 (1992).
2. Johansen, C., Kragballe, K., Rasmussen, M., *et al.* Activator protein 1 DNA binding activity is decreased in lesional psoriatic skin compared with nonlesional psoriatic skin. *Br. J. Dermatol.* **151**, 600-607 (2004).
3. Li, M., Hener, P., Zhang, Z., *et al.* Topical vitamin D<sub>3</sub> and low-calcemic analogs induce thymic stromal lymphopoietin in mouse keratinocytes and trigger an atopic dermatitis. *Proc. Natl. Acad. Sci. USA* **103(31)**, 11736-11741 (2006).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10009599](http://www.caymanchem.com/catalog/10009599)

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 02/09/2012