

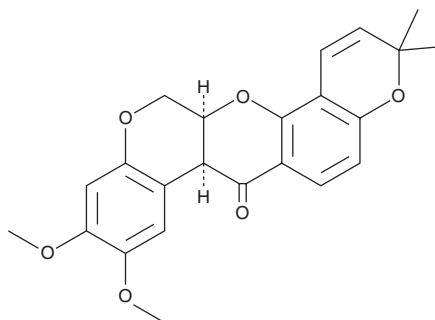
Product Information



(-)-Deguelin

Item No. 10010706

CAS Registry No.: 522-17-8
Formal Name: 13,13aS-dihydro-9,10-dimethoxy-3,3-dimethyl-3H-[1]benzopyrano[3,4-b]pyrano[2,3-h][1]benzopyran-7(7aS)-one
Synonym: (-)-*cis*-Deguelin
MF: C₂₃H₂₂O₆
FW: 394.06
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

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

(-)-Deguelin is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-deguelin in an organic solvent purged with an inert gas. (-)-Deguelin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (-)-deguelin in these solvents is approximately 25 mg/ml.

(-)-Deguelin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (-)-deguelin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (-)-Deguelin has a solubility of approximately 0.33 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Rotenoids, deriving from the flavonoid family of compounds, act as chemopreventive agents that inhibit NADH:ubiquinone oxidoreductase activity and suppress phorbol ester-induced ornithine decarboxylase (ODC) activity.¹ The rotenoid compound deguelin, originally isolated from the bark of *M. sericea* (*Leguminosae*) to be used as an insecticide and fish poison, has chemopreventive and chemosensitizing effects in models of skin, mammary, colon, and lung carcinogenesis.²⁻⁴ Deguelin inhibits cell growth (IC₅₀ = <10⁻⁸ M), blocks PI3K/Akt activation, suppresses COX-2 expression, and induces apoptosis in premalignant and squamous human bronchial epithelial (HBE) cells without affecting normal HBE cells.^{2,5} At 6 mg/kg, deguelin induces Parkinson's disease-like symptoms in rats after six days of subcutaneous infusion and therefore may also serve as a useful model for Parkinson's disease research.⁶

References

1. Fang, N. and Casida, J.E. *Proc. Natl. Acad. Sci. USA* **95**, 3380-3384 (1998).
2. Chun, K.-H., Kosmeder, J.W.II., Sun, S., *et al. Journal of the National Cancer Institute* **95**(4), 291-302 (2003).
3. Lee, H.-Y., Oh, S.-H., Woo, J.K., *et al. Journal of the National Cancer Institute* **97**(22), 1695-1699 (2005).
4. Peng, X.-H., Karna, P., O'Regan, R.M., *et al. Mol. Pharmacol.* **71**, 101-111 (2007).
5. Lee, H.-Y., Suh, Y.-A., Kosmeder, J.W., *et al. Clinical Cancer Research* **10**, 1074-1079 (2004).
6. Caboni, P., Sherer, T.B., Zhang, N., *et al. Chem. Res Toxicol.* **17**, 1540-1548 (2004).

Related Product

Epoxomicin - Item No. 10007806

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

Web

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, 02/18/2011