

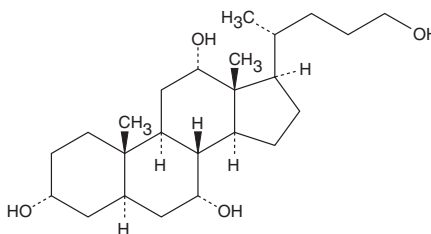
Product Information



Petromyzonol

Catalog No. 98250

CAS Registry No.: 28979-29-5
Formal Name: 3 α ,7 α ,12 α ,24-tetrahydroxy-5 α -cholane
MF: C₂₄H₄₂O₄
FW: 394.6
Purity: \geq 98%
Stability: \geq 1 year at room temperature
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that petromyzonol be stored as supplied at room temperature. It should be stable for at least one year.

Petromyzonol is supplied as a crystalline solid. A stock solution may be made by dissolving the petromyzonol in the solvent of choice. Petromyzonol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of petromyzonol in these solvents is approximately 1.8 mg/ml.

Petromyzonol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, petromyzonol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Petromyzonol has a solubility of approximately 0.05 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Petromyzonol is a tetrahydroxy stearyl produced in the bile of sea lamprey larvae from the bile acid precursor allocholic acid. While the adult sea lamprey is relatively insensitive to petromyzonol itself, the C-24 sulfate ester (petromyzonol sulfate) is a pheromone and a spawning chemoattractant which can be detected at very low concentrations by lamprey olfactory chemoreceptors. Petromyzonol, petromyzonol sulfate, and allocholic acid are all found in water samples from fresh water streams bearing larval lamprey populations.¹

References

1. Li, W., Sorensen, P.W., and Gallaher, D.D. The olfactory system of migratory adult sea lamprey (*Petromyzon marinus*) is specifically and acutely sensitive to unique bile acids released by conspecific larvae. *J. Gen. Physiol.* **105**, 569-587 (1995).

Related Products

3-keto Petromyzonol - Cat. No. 10007055

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, 11/17/2009