

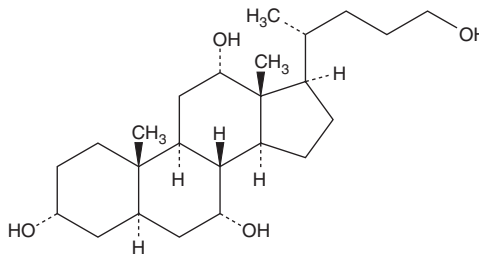
PRODUCT INFORMATION



Petromyzonol

Item 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%
Supplied as: A crystalline solid
Storage: Room temperature
Stability: \geq 4 years



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

Laboratory Procedures

Petromyzonol is supplied as a crystalline solid. A stock solution may be made by dissolving the petromyzonol in the solvent of choice, which should be purged with an inert gas. Petromyzonol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of petromyzonol in ethanol and DMF is approximately 2 mg/ml and approximately 1 mg/ml in DMSO.

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 50 μ g/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.

Description

Petromyzonol is a tetrahydroxy stearyl which serves as the primary spawning pheromone in adult sea lamprey. It is 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

Reference

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**(5), 569-587 (1995).

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, 01/11/2022

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