

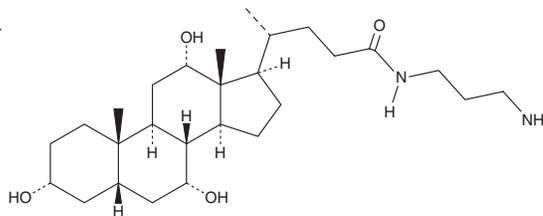
PRODUCT INFORMATION



Cholyl-1,3-diaminopropane

Item No. 40898

CAS Registry No.: 136533-45-4
Formal Name: (5 β)-N-(3-aminopropyl)-3 α ,7 α ,12 α -trihydroxy-cholan-24-amide
MF: C₂₇H₄₈N₂O₄
FW: 464.7
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Cholyl-1,3-diaminopropane is supplied as a solid. A stock solution may be made by dissolving the cholyl-1,3-diaminopropane in the solvent of choice, which should be purged with an inert gas. Cholyl-1,3-diaminopropane is soluble (\geq 10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of cholyl-1,3-diaminopropane can be prepared by directly dissolving the solid in aqueous buffers. Cholyl-1,3-diaminopropane is soluble (\geq 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Cholyl-1,3-diaminopropane is a polyamine bile amidate.¹ It has been used as a standard for the detection of endogenous bile acid amidates in mammalian intestinal and fecal samples.

Reference

1. Mohanty, I., Mannocho-Russo, H., Schweer, J.V., *et al.* The underappreciated diversity of bile acid modifications. *Cell* **187**(7), 1801-1818 (2024).

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, 06/07/2024

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