

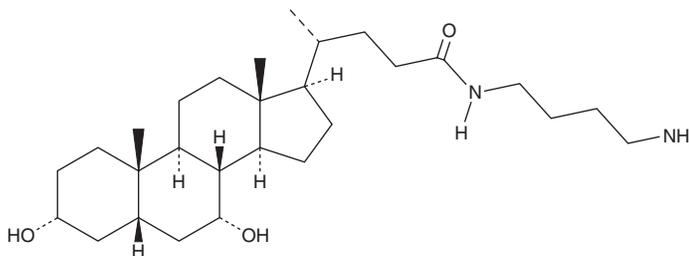
PRODUCT INFORMATION



Chenodeoxycholyl-putrescine

Item No. 40896

CAS Registry No.: 142970-48-7
Formal Name: (5β)-N-(4-aminobutyl)-3α,7α-dihydroxy-cholan-24-amide
Synonym: GCDCA-putrescine
MF: C₂₈H₅₀N₂O₃
FW: 462.7
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Chenodeoxycholyl-putrescine is supplied as a crystalline solid. A stock solution may be made by dissolving the chenodeoxycholyl-putrescine in the solvent of choice, which should be purged with an inert gas. Chenodeoxycholyl-putrescine is sparingly soluble (1-10 mg/ml) in organic solvents ethanol and 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 chenodeoxycholyl-putrescine can be prepared by directly dissolving the crystalline solid in aqueous buffers. Chenodeoxycholyl-putrescine is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Chenodeoxycholyl-putrescine is a polyamine bile amidate.¹ It has been found in lion, but not mouse, fecal samples. Chenodeoxycholyl-putrescine has been used as a standard in the detection of chenodeoxycholyl-putrescine 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, 06/11/2024

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