

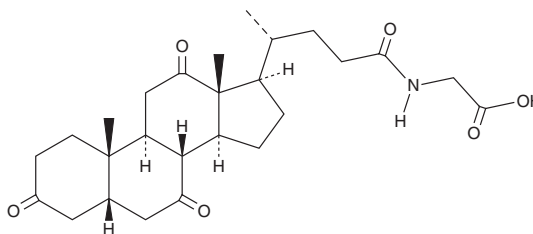
PRODUCT INFORMATION



Glycodehydrocholic Acid

Item No. 29892

CAS Registry No.: 3415-45-0
Formal Name: N-[(5 β)-3,7,12,24-tetraoxocholan-24-yl]-glycine
Synonym: Glycine Dehydrocholate
MF: C₂₆H₃₇NO₆
FW: 459.6
Purity: \geq 95%
Supplied as: A crystalline 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

Glycodehydrocholic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the glycodehydrocholic acid in the solvent of choice, which should be purged with an inert gas. Glycodehydrocholic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of glycodehydrocholic acid in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

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 glycodehydrocholic acid can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of glycodehydrocholic acid in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Glycodehydrocholic acid is a glycine-conjugated form of the synthetic bile acid dehydrocholic acid (Item No. 29176).¹

Reference

1. Beher, W.T., Stradnieks, S., Beher, G.R., *et al.* The hydrolysis of bile acid conjugates. *Steroids* **32**(3), 355-363 (1978).

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, 11/23/2022

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