

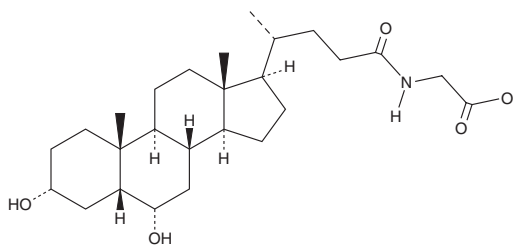
PRODUCT INFORMATION



Glycohyodeoxycholic Acid

Item No. 22643

CAS Registry No.: 13042-33-6
Formal Name: N-[(3 α ,5 β ,6 α)-3,6-dihydroxy-24-oxocholan-24-yl]-glycine
Synonyms: GHDCa, Glycine hyodeoxycholate
MF: C₂₆H₄₃NO₅
FW: 449.6
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

Glycohyodeoxycholic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the glycohyodeoxycholic acid in the solvent of choice. Glycohyodeoxycholic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of glycohyodeoxycholic acid in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

Glycohyodeoxycholic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, glycohyodeoxycholic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Glycohyodeoxycholic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Glycohyodeoxycholic acid is a major metabolite of the secondary bile acid hyodeoxycholic acid (Item No. 20294) in humans.¹ Glycohyodeoxycholic acid supplementation in prairie dogs fed a lithogenic diet decreases the frequency of cholesterol crystals in the gallbladder and inhibits the activity of cholesterol 7 α -hydroxylase.²

References

1. Sacquet, E., Parquet, M., Riottot, M., *et al.* Intestinal absorption, excretion, and biotransformation of hyodeoxycholic acid in man. *J. Lipid Res.* **24(5)**, 604-613 (1983).
2. McSherry, C.K., Mosbach, E.H., Cohen, B.I., *et al.* Hyodeoxycholic acid: A new approach to gallstone prevention. *Am. J. Surg.* **149(1)**, 126-132 (1985).

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, 11/11/205

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