

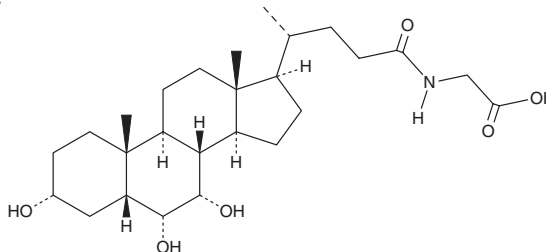
PRODUCT INFORMATION



Glycohyocholic Acid

Item No. 22670

CAS Registry No.: 32747-08-3
Formal Name: N-[(3 α ,5 β ,6 α ,7 α)-3,6,7-trihydroxy-24-oxocholan-24-yl]-glycine
Synonyms: GHCA, Glycine Hyocholate, Glycohyocholate
MF: C₂₆H₄₃NO₆
FW: 465.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 202 nm
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

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

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

Description

GHCA is a glycine-conjugated form of the primary bile acid hyocholic acid (Item No. 20293).¹ GHCA is upregulated 6.6-fold in the serum of patients with cirrhosis induced by hepatitis C virus (HCV) compared to healthy controls.² Plasma levels of GHCA increase in patients who are no longer diabetic following gastric bypass surgery.³

References

1. Guo, L., Milburn, M.V., Ryals, J.A., *et al.* Plasma metabolomic profiles enhance precision medicine for volunteers of normal health. *Proc. Nat. Acad. Sci. USA* **112(35)**, E4901-E4910 (2015).
2. Fitian, A.I., Nelson, D.R., Liu, C., *et al.* Integrated metabolomic profiling of hepatocellular carcinoma in hepatitis C cirrhosis through GC/MS and UPLC/MS-MS. *Liver Int.* **34(9)**, 1428-1444 (2014).
3. Yu, H., Ni, Y., Bao, Y., *et al.* Chenodeoxycholic acid as a potential prognostic marker for roux-en-Y gastric bypass in Chinese obese patients. *J. Clin. Endocrinol. Metab.* **100(11)**, 4222-4230 (2015).

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