

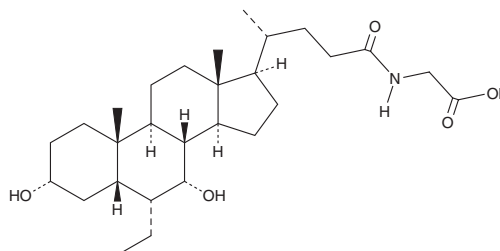
PRODUCT INFORMATION



Glyco-Obeticholic Acid

Item No. 28242

CAS Registry No.: 863239-60-5
Formal Name: N-[(3 α ,5 β ,6 α ,7 α)-6-ethyl-3,7-dihydroxy-24-oxocholan-24-yl]-glycine
Synonym: Obeticholic Acid Glycine Conjugate
MF: C₂₈H₄₇NO₅
FW: 477.7
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Glyco-obeticholic acid is supplied as a solid. A stock solution may be made by dissolving the glyco-obeticholic acid in the solvent of choice, which should be purged with an inert gas. Glyco-obeticholic acid is slightly soluble in DMSO and methanol.

Description

Glyco-obeticholic acid is an active metabolite of obeticholic acid (6-ECDCA; Item No. 11031), which is a farnesoid X receptor (FXR) agonist and semisynthetic derivative of chenodeoxycholic acid (Item No. 10011286).¹ Glyco-obeticholic acid is formed from obeticholic acid by glycine conjugation in the liver but can be reconverted back to obeticholic acid by microorganism-mediated deconjugation in the ileum and colon. It has been used as a precursor in the synthesis of bile acid analogs as agonists of the farnesoid X receptor (FXR) and TGR5.²

References

1. Markham, A. and Keam, S.J. Obeticholic Acid: First global approval. *Drugs* **76**(12), 1221-1226 (2016).
2. Or, Y.S., Wang, G., Shen, R., *et al.* Bile acid analogs and FXR/TGR5 agonists and methods of use thereof. *Enanta Pharmaceuticals, Inc.* **WO 2016/073767 A1** (2016).

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, 12/07/2022

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