

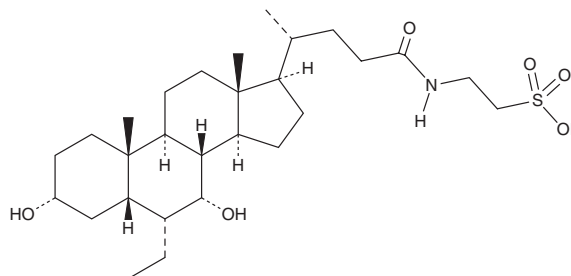
PRODUCT INFORMATION



Tauro-Obeticholic Acid

Item No. 28243

CAS Registry No.: 863239-61-6
Formal Name: 2-[[[(3 α ,5 β ,6 α ,7 α)-6-ethyl-3,7-dihydroxy-24-oxocholan-24-yl]amino]-ethanesulfonic acid
Synonyms: Obeticholic Acid Taurine Conjugate, T-ECDCA, Tauro-OCA, Tauro-6-Ethyl-CDCA, Tauro 6-Ethylchenodeoxycholic Acid
MF: C₂₈H₄₉NO₆S
FW: 527.8
Purity: $\geq 95\%$
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

Tauro-obeticholic acid is supplied as a solid. A stock solution may be made by dissolving the tauro-obeticholic acid in the solvent of choice, which should be purged with an inert gas. Tauro-obeticholic acid is soluble in organic solvents such as methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Tauro-Obeticholic acid is an active metabolite of obeticholic acid, which is a farnesoid X receptor (FXR) agonist and semisynthetic derivative of chenodeoxycholic acid (Item No. 10011286).¹ Tauro-obeticholic acid is formed from obeticholic acid by taurine conjugation in the liver but can be reconverted back to obeticholic acid by intestinal microflora.

Reference

1. Markham, A. and Keam, S.J. Obeticholic acid: First global approval. *Drugs* **76(12)**, 1221-1226 (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, 07/13/2023

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