

PRODUCT INFORMATION



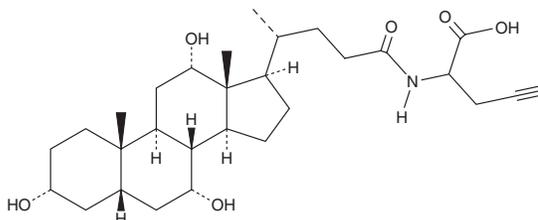
Clickable Cholic Acid

Item No. 40735

Formal Name: 2-((R)-4-((3R,5S,7R,8R,9S,10S,12S,13R,14S,17R)-3,7,12-trihydroxy-10,13-dimethylhexadecahydro-1H-cyclopenta[a]phenanthren-17-yl)pentanamido)pent-4-ynoic acid

Synonyms: Cholalic Acid, Cholalin, ClickTag™ Cholic Acid

MF: C₂₉H₄₅NO₆
FW: 503.7
Purity: ≥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

Clickable cholic acid is supplied as a solid. A stock solution may be made by dissolving the clickable cholic acid in the solvent of choice, which should be purged with an inert gas. Clickable cholic acid is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Clickable cholic acid is a clickable form of the primary bile acid cholic acid (Item No. 20250). It has been used in the synthesis of bile acid-nucleoside conjugates with anticancer and antimycobacterial activities.¹

Reference

1. Agarwal, D.S., Siva Krishna, V., Sriram, D., *et al.* Clickable conjugates of bile acids and nucleosides: Synthesis, characterization, *in vitro* anticancer and antituberculosis studies. *Steroids* **139**, 35-44 (2018).

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, 06/11/2024

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