

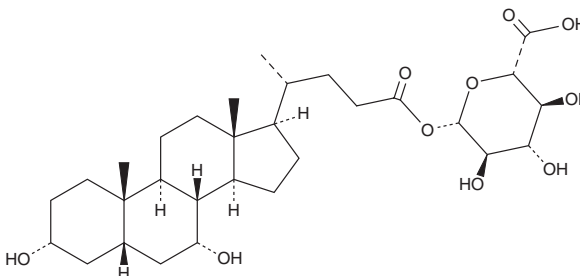
PRODUCT INFORMATION



Chenodeoxycholic Acid 24-Acyl-β-D-Glucuronide

Item No. 33227

CAS Registry No.: 208038-27-1
Formal Name: 1-[(3α,5β,7α)-3,7-dihydroxycholan-24-oate]β-D-glucopyranuronic acid
Synonyms: CDCA-24G, CDCA-24-glucuronide
MF: C₃₀H₄₈O₁₀
FW: 568.7
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

Chenodeoxycholic acid 24-acyl-β-D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the chenodeoxycholic acid 24-acyl-β-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Chenodeoxycholic acid 24-acyl-β-D-glucuronide is slightly soluble in DMSO.

Description

Chenodeoxycholic acid 24-acyl-β-D-glucuronide (CDCA-24G) is a metabolite of CDCA (Item No. 10011286).^{1,2} It is formed from CDCA by the UDP-glucuronosyltransferase (UGT) isoform UGT1A3.¹

References

1. Caron, P., Trottier, J., Verreault, M., *et al.* Enzymatic production of bile acid glucuronides used as analytical standards for liquid chromatography-mass spectrometry analyses. *Mol. Pharm.* **3(3)**, 293-302 (2006).
2. Takehara, I., Terashima, H., Nakayama, T., *et al.* Investigation of glycochenodeoxycholate sulfate and chenodeoxycholate glucuronide as surrogate endogenous probes for drug interaction studies of OATP1B1 and OATP1B3 in healthy Japanese volunteers. *Pharm. Res.* **34(8)**, 1601-1614 (2017).

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, 12/06/2022

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