

PRODUCT INFORMATION



Chenodeoxycholic Acid 3-Glucuronide

Item No. 36289

CAS Registry No.: 58814-71-4

Formal Name: β -D-glucopyranosiduronic acid,
(3 α ,5 β ,7 α)-23-carboxy-7-hydroxy-
24-norcholan-3-yl

Synonyms: CDCA-3G, CDCA-3-Glucuronide

MF: C₃₀H₄₈O₁₀

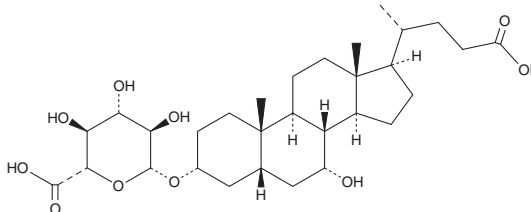
FW: 568.7

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

Chenodeoxycholic acid 3-glucuronide (CDCA-3G) is supplied as a solid. A stock solution may be made by dissolving the CDCA-3G in the solvent of choice, which should be purged with an inert gas. CDCA-3G is slightly soluble in methanol and DMSO.

Description

CDCA-3G is a metabolite of CDCA (Item No. 10011286).^{1,2} It is formed from CDCA in human liver microsomes.

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, 12/19/2022

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