

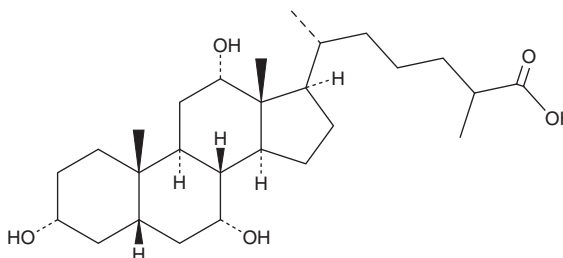
PRODUCT INFORMATION



Trihydroxycholestanoic Acid

Item No. 26098

CAS Registry No.: 547-98-8
Formal Name: (5 β)-3 α ,7 α ,12 α -trihydroxy-cholestan-26-oic acid
Synonym: Trihydroxycoprostanic Acid
MF: C₂₇H₄₆O₅
FW: 450.7
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

Trihydroxycholestanoic acid is supplied as a solid. A stock solution may be made by dissolving the trihydroxycholestanoic acid in the solvent of choice, which should be purged with an inert gas. Trihydroxycholestanoic acid is slightly soluble in chloroform (warmed), ethanol, and DMSO.

Description

Trihydroxycholestanoic acid is an intermediate in the biosynthesis of cholic acid (Item No. 20250).¹ Elevated plasma levels of trihydroxycholestanoic acid have been found in patients with Zellweger syndrome, a neurological disorder characterized by mutations in *PEX* genes which result in defects in peroxisome formation.^{2,3}

References

1. Keane, M.H., Overmars, H., Wikander, T.M., *et al.* Bile acid treatment alters hepatic disease and bile acid transport in peroxisome-deficient *PEX2* Zellweger mice. *Hepatology* **45**(4), 982-997 (2007).
2. Ferdinandusse, S., Overmars, H., Denis, S., *et al.* Plasma analysis of di- and trihydroxycholestanoic acid diastereoisomers in peroxisomal α -methylacyl-CoA racemase deficiency. *J. Lipid Res.* **42**(1), 137-141 (2001).
3. Klouwer, F.C.C., Berendse, K., Ferdinandusse, S., *et al.* Zellweger spectrum disorders: Clinical overview and management approach. *Orphanet J. Rare Dis.* **10**, 151 (2015).

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, 12/01/2022

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