

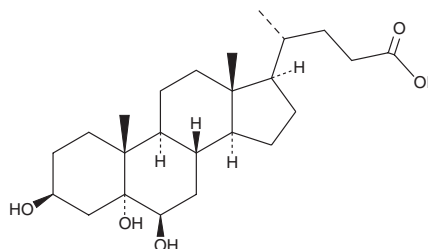
PRODUCT INFORMATION



3 β ,5 α ,6 β -Trihydroxycholan-24-oic Acid

Item No. 19636

CAS Registry No.: 16030-92-5
Formal Name: 3 β ,5 α ,6 β -trihydroxy-cholan-24-oic acid
MF: C₂₄H₄₀O₅
FW: 408.6
Purity: \geq 90%
Supplied as: A solid
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

3 β ,5 α ,6 β -Trihydroxycholan-24-oic acid is supplied as a solid. A stock solution may be made by dissolving the 3 β ,5 α ,6 β -trihydroxycholan-24-oic acid in the solvent of choice, which should be purged with an inert gas. 3 β ,5 α ,6 β -Trihydroxycholan-24-oic acid is soluble in methanol and DMSO.

Description

3 β ,5 α ,6 β -Trihydroxycholan-24-oic acid is a metabolite of 5 α ,6 β -dihydroxycholestanol (Item No. 25538).^{1,2} It is increased in dried blood spots from patients with Niemann-Pick disease type C, a neurodegenerative cholesterol-sphingolipid lysosomal storage disorder.²

References

1. Kikuchi, S., Imai, Y., Suzuoki, Z., *et al.* Biologic studies of cholestane-3 β ,5 α ,6 β -triol and its derivatives. III. The metabolic fate and metabolites of cholestane-3 β ,5 α ,6 β -triol in animals. *J. Pharmacol. Exp. Ther.* **159**(2), 399-408 (1968).
2. Jiang, X., Sidhu, R.S., Mydock-McGrane, L., *et al.* Development of a bile acid-based newborn screen for Niemann-Pick disease type C. *Sci. Transl. Med.* **8**(337), 337ra63 (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, 07/07/2021

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