

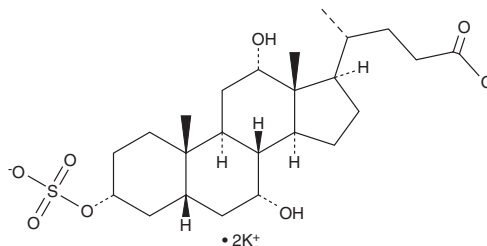
PRODUCT INFORMATION



Cholic Acid 3-sulfate (potassium salt)

Item No. 38019

Formal Name: (3 α ,5 β ,7 α ,12 α)-7,12-dihydroxy-3-(sulfooxy)-
cholan-24-oic acid, dipotassium salt
Synonyms: 3-Sulfocholate, 3-Sulfocholic Acid
MF: C₂₄H₃₈O₈S • 2K
FW: 564.8
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



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

Laboratory Procedures

Cholic acid 3-sulfate (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the cholic acid 3-sulfate (potassium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Cholic acid 3-sulfate is a metabolite of the primary bile acid cholic acid (Item No. 20250).^{1,2} Plasma levels of cholic acid 3-sulfate are increased in patients with intrahepatic cholestasis and mutations in the gene encoding the bile salt export pump (BSEP), *ABCB11*.³

References

1. Alnouti, Y. Bile acid sulfation: A pathway of bile acid elimination and detoxification. *Toxicol. Sci.* **108**(2), 225-246 (2009).
2. Goto, J., Kato, H., Saruta, Y., *et al.* Studies on steroids. CLXX. Separation and determination of bile acid 3-sulfates in human bile by high-performance liquid chromatography. *J. Chromatogr.* **226**(1), 13-24 (1981).
3. Liu, T., Wang, R.-X., Han, J., *et al.* Comprehensive bile acid profiling in hereditary intrahepatic cholestasis: Genetic and clinical correlations. *Liver Int.* **38**(9), 1676-1685 (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, 11/06/2024

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