

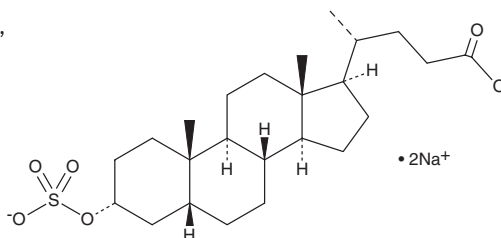
PRODUCT INFORMATION



Lithocholic Acid 3-sulfate (sodium salt)

Item No. 20676

CAS Registry No.: 64936-81-8
Formal Name: (3 α ,5 β)-3-(sulfooxy)-cholan-24-oic acid, disodium salt
Synonyms: LCA 3-sulfate, Lithocholate 3-sulfate
MF: C₂₄H₃₈O₆S • 2Na
FW: 500.6
Purity: \geq 98%
Supplied as: A crystalline 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

Lithocholic acid 3-sulfate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the lithocholic acid 3-sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Lithocholic acid 3-sulfate (sodium salt) is soluble in the organic solvent methanol. It is also soluble in water. The solubility of lithocholic acid 3-sulfate (sodium salt) in methanol and water is approximately 1 and 5 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Lithocholic acid 3-sulfate is a metabolite of the secondary bile acid lithocholic acid (Item No. 20253).¹ It is water soluble and the primary form of lithocholic acid in the bile duct. Lithocholic acid 3-sulfate forms a complex with calcium *in vitro*.² Biliary secretion of lithocholic acid 3-sulfate is increased in a rat model of diabetes induced by streptozotocin (Item No. 13104).³

References

1. Borriello, S.P. and Owen, R.W. The metabolism of lithocholic acid and lithocholic acid-3- α -sulfate by human fecal bacteria. *Lipids* **17(7)**, 477-482 (1982).
2. Oelberg, D.G., Dubinsky, W.P., Adcock, E.W., *et al.* Calcium binding by lithocholic acid derivatives. *Am. J. Physiol.* **247(1 Pt 1)**, G112-G115 (1984).
3. Kirkpatrick, R.B. and Kraft, B.G. Effect of streptozotocin-induced diabetes on bile acid sulfation in male rat liver. *Am. J. Physiol.* **247(3 Pt 1)**, G226-G230 (1984).

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, 04/28/2022

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