

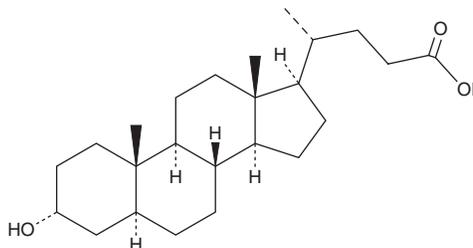
PRODUCT INFORMATION



Allolithocholic Acid

Item No. 38158

| | |
|--------------------------|--|
| CAS Registry No.: | 2276-94-0 |
| Formal Name: | (3 α ,5 α)-3-hydroxy-cholan-24-oic acid |
| Synonyms: | 3 α -hydroxy-5 α -Cholanoic Acid, <i>allo</i> -LCA |
| MF: | C ₂₄ H ₄₀ O ₃ |
| FW: | 376.6 |
| Purity: | ≥98% |
| 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

Allolithocholic acid is supplied as a solid. A stock solution may be made by dissolving the allolithocholic acid in the solvent of choice, which should be purged with an inert gas. Allolithocholic acid is slightly soluble in ethanol, methanol, and DMSO.

Description

Allolithocholic acid is a secondary allomonohydroxy bile acid.¹ It activates large-conductance calcium-activated potassium channels (BK/K_{CaS}; EC₅₀ = 44.21 μ M in *Xenopus* oocytes expressing the rat channel). Fecal levels of allolithocholic acid are increased in patients with colorectal cancer.² Allolithocholic acid has also been found in the urine of infants with biliary atresia.³

References

1. Bukiya, A.N., McMillan, J., Parrill, A.L., *et al.* Structural determinants of monohydroxylated bile acids to activate β_1 subunit-containing BK channels. *J. Lipid Res.* **49(11)**, 2441-2451 (2008).
2. Tadano, T., Kanoh, M., Kondoh, H., *et al.* Kinetic analysis of bile acids in the feces of colorectal cancer patients by gas chromatography-mass spectrometry (GC-MS). *Rinsho Byori.* **55(5)**, 417-427 (2007).
3. Makino, I., Sjövall, J., Norman, A., *et al.* Excretion of 3 β -hydroxy-5-cholenoic and 3 α -hydroxy-5 α -cholanoic acids in urine of infants with biliary atresia. *FEBS Lett.* **15(2)**, 161-164 (1971).

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, 08/22/2024

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