

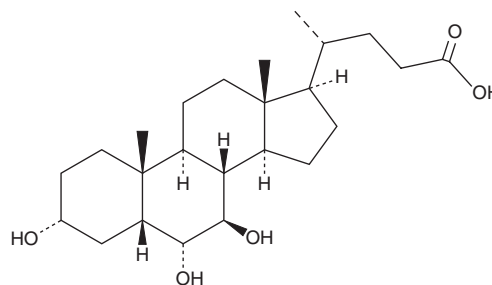
PRODUCT INFORMATION



ω -Muricholic Acid

Item No. 20292

CAS Registry No.: 6830-03-1
Formal Name: (3 α ,5 β ,6 α ,7 β)-3,6,7-trihydroxy-
cholan-24-oic acid
Synonym: ω -MCA
MF: C₂₄H₄₀O₅
FW: 408.6
Purity: \geq 95%
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

ω -Muricholic acid (ω -MCA) is supplied as a crystalline solid. A stock solution may be made by dissolving the ω -MCA in the solvent of choice. ω -MCA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ω -MCA in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

ω -MCA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ω -MCA should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ω -MCA has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ω -MCA is a murine-specific secondary bile acid.^{1,2} It is formed *via* enzymatic conversion of β -MCA (Item No. 20287) by a variety of intestinal microorganisms, including *Clostridium*, in rat and mouse gut.

References

1. Eyssen, H., De Pauw, G., Stragier, J., *et al.* Cooperative formation of ω -muricholic acid by intestinal microorganisms. *Appl. Environ. Microbiol.* **45(1)**, 141-147 (1983).
2. Sacquet, E.C., Raibaud, P.M., Mejean, C., *et al.* Bacterial formation of ω -muricholic acid in rats. *Appl. Environ. Microbiol.* **37(6)**, 1127-1131 (1979).

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, 09/25/2018

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