

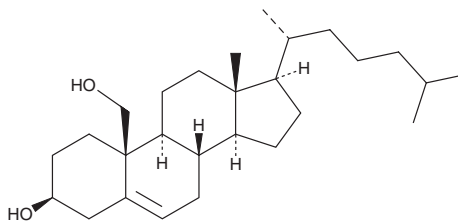
PRODUCT INFORMATION



19-hydroxy Cholesterol

Item No. 15209

CAS Registry No.: 561-63-7
Formal Name: cholest-5-ene-3 β ,19-diol
MF: C₂₇H₄₆O₂
FW: 402.7
Purity: $\geq 98\%$
Supplied as: A crystalline 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

19-hydroxy Cholesterol is supplied as a crystalline solid. A stock solution may be made by dissolving the 19-hydroxy cholesterol in the solvent of choice. 19-hydroxy Cholesterol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 19-hydroxy cholesterol in these solvents is approximately 20, 0.1, and 2 mg/ml, respectively.

19-hydroxy Cholesterol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 19-hydroxy cholesterol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 19-hydroxy Cholesterol has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

19-hydroxy Cholesterol is formed during metabolic oxidation of cholesterol. It has been used as an internal standard for the quantitative determination of sterols by mass spectroscopic analysis.¹

References

1. Yamaga, N., Ogura, Y., Islam, A., *et al.* Utility of 19-hydroxycholesterol as an internal standard compound for the quantitative determination of sterols using capillary gas chromatograph. *Yonago Acta Med.* **45**, 27-33 (2002).

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, 10/31/2022

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