

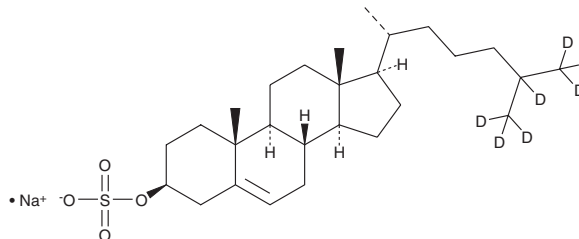
PRODUCT INFORMATION



Cholesterol Sulfate-d₇ (sodium salt)

Item No. 45740

CAS Registry No.: 2687961-01-7
Formal Name: (3β)-3-(hydrogen sulfate), cholest-5-en-25,26,26,26,27,27,27-d₇-3-ol, monosodium salt
Synonyms: SCS-d₇, Sodium Cholesteryl Sulfate-d₇
MF: C₂₇H₃₈D₇O₄S • Na
FW: 495.7
Chemical Purity: ≥98% (cholesterol sulfate)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
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

Cholesterol sulfate-d₇ (sodium salt) is intended for use as an internal standard for the quantification of cholesterol sulfate (Item No. 15106) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Cholesterol sulfate-d₇ (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the cholesterol sulfate-d₇ (sodium salt) in the solvent of choice, which should be purged with an inert gas. Cholesterol sulfate-d₇ (sodium salt) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of cholesterol sulfate-d₇ (sodium salt) in these solvents is approximately 10 mg/ml and 20 mg/ml, respectively.

Description

Cholesterol sulfate is an endogenous component of cell membranes, where it serves as a substrate for steroid synthesis, influences lipid metabolism, stabilizes cell membranes, regulates the activity of functional proteins, and plays a role in keratinocyte differentiation and development of the epidermal barrier.¹

Reference

1. Strott, C.A. and Higashi, Y. Cholesterol sulfate in human physiology: What's it all about? *J. Lipid Res.* **44**(7), 1268-1278 (2003).

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, 06/02/2026

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