

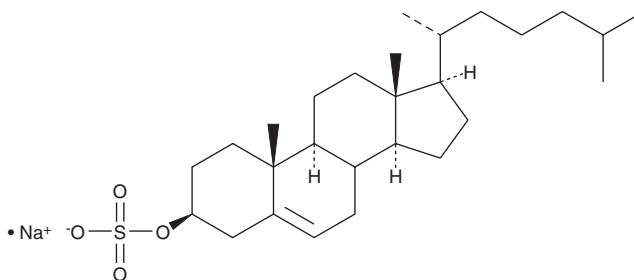
PRODUCT INFORMATION



Cholesterol Sulfate (sodium salt)

Item No. 15106

CAS Registry No.: 2864-50-8
Formal Name: (3 β)-3-(hydrogen sulfate)-cholest-5-en-3-ol, monosodium salt
Synonyms: SCS, Sodium Cholesteryl Sulfate
MF: C₂₇H₄₅O₄S • Na
FW: 488.7
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Cholesterol sulfate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the cholesterol sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Cholesterol sulfate (sodium salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of cholesterol sulfate (sodium salt) in these solvents is approximately 10 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 This synthetic formulation can be used to form nonconventional liposomes in order to study cellular membrane properties.²

References

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).
2. Méléard, P., Gerbeaud, C., Pott, T., *et al.* Bending elasticities of model membranes: Influences of temperature and sterol content. *Biophys. J.* **72**(6), 2616-2629 (1997).

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, 12/12/2022

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