

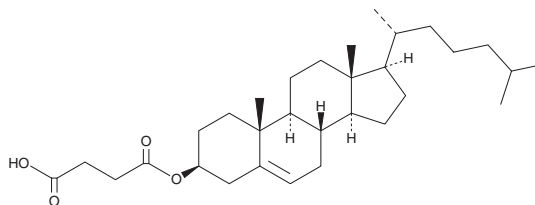
PRODUCT INFORMATION



Cholesteryl Hemisuccinate

Item No. 25698

CAS Registry No.: 1510-21-0
Formal Name: (3 β)-cholest-5-en-3-ol 3-(hydrogen butanedioate)
Synonym: Cholesteryl hydrogen succinate
MF: C₃₁H₅₀O₄
FW: 486.7
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

Cholesteryl hemisuccinate is supplied as a crystalline solid. A stock solution may be made by dissolving the cholesteryl hemisuccinate in the solvent of choice, which should be purged with an inert gas. Cholesteryl hemisuccinate is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

Cholesteryl hemisuccinate is a cholesterol ester with anticancer activity.¹ It inhibits the growth of murine C1498 myeloid and L1210 lymphocytic leukemia cells when used at concentrations of 50 and 150 μ M, respectively.¹ Cholesteryl hemisuccinate acts as an ionizable anionic detergent and is commonly used to stabilize unilamellar vesicles and liposomes.² It has also been used as an emulsifying agent in various vesicular drug delivery systems for anticancer drugs, antibiotics, and oligonucleotides and to solubilize various proteins including chemokine receptor 1 as well as erythrocyte ghosts.³⁻⁵

References

1. Fariss, M.W., Fortuna, M.B., Everett, C.K., *et al.* The selective antiproliferative effects of α -tocopheryl hemisuccinate and cholesteryl hemisuccinate on murine leukemia cells result from the action of the intact compounds. *Cancer Res.* **54(13)**, 3346-3351 (1994).
2. Ding, W.X., Qi, X.R., Li, P., *et al.* Cholesteryl hemisuccinate as a membrane stabilizer in dipalmitoylphosphatidylcholine liposomes containing saikosaponin-d. *Int. J. Pharm.* **300(1-2)**, 38-47 (2005).
3. Simoes, S., Moreira, J.N., Fonseca, C., *et al.* On the formulation of pH-sensitive liposomes with long circulation times. *Adv. Drug Deliv. Rev.* **56(7)**, 947-965 (2004).
4. Allen, S.J., Ribeiro, S., Horuk, R., *et al.* Expression, purification and *in vitro* functional reconstitution of the chemokine receptor CCR1. *Protein Expr. Purif.* **66(1)**, 73-81 (2009).
5. Grodecka, M., Bertrand, O., Karolak, E., *et al.* One-step immunopurification and lectinochemical characterization of the Duffy atypical chemokine receptor from human erythrocytes. *Glycoconj. J.* **29(2-3)**, 93-105 (2012).

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, 12/13/2022

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