

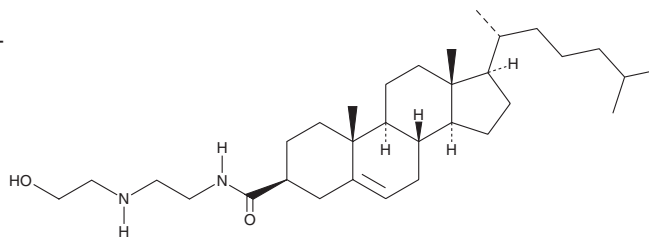
PRODUCT INFORMATION



OH-Chol

Item No. 26586

CAS Registry No.: 191173-82-7
Formal Name: N-[2-[(2-hydroxyethyl)amino]ethyl]-(3 β)-cholest-5-ene-3-carboxamide
MF: C₃₂H₅₆N₂O₂
FW: 500.8
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

OH-Chol is supplied as a crystalline solid. A stock solution may be made by dissolving the OH-chol in the solvent of choice, which should be purged with an inert gas. OH-Chol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of OH-chol in these solvents is approximately 10 mg/ml.

OH-Chol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, OH-chol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. OH-Chol has a solubility of approximately 0.14 mg/ml in a 1:6 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

OH-Chol is a cationic cholesterol derivative.¹ OH-Chol, as a component of lipoplexes with DOPE (Item No. 15091), has been used for siRNA delivery and gene silencing in MCF-7 cells, as well as in mice *via* intravenous injection, resulting in lipoplex accumulation in the liver. It has also been used in cationic nanoparticles in combination with Tween 80 to transfect pDNA and siRNA into PC3 mouse xenografts *via* intratumoral injection and with Tween 80 and folate-PEG₂₀₀₀-DSPE in a KB mouse xenograft model for intratumoral gene delivery.²

References

1. Hattori, Y., Nakamura, M., Takeuchi, N., *et al.* Effect of cationic lipid in cationic liposomes on siRNA delivery into the lung by intravenous injection of cationic lipoplex. *J. Drug. Target* **27(2)**, 217-227 (2019).
2. Hattori, Y. Development of non-viral vector for cancer gene therapy. *Yakugaku Zasshi* **130(7)**, 917-923 (2010).

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, 10/14/2019

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