PRODUCT INFORMATION



OH-C-Chol

Item No. 26587

CAS Registry No.: 496801-51-5

Formal Name: (3β)-cholest-5-en-3-ol 3-[N-[2-[(2-

hydroxyethyl)aminolethyl]carbamatel

MF: $C_{32}H_{56}N_2O_3$ FW: 516.8 **Purity:** ≥95%

Supplied as: A crystalline solid

Storage: -20°C Stability: ≥2 years

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

Laboratory Procedures

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

OH-C-Chol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, OH-C-chol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. OH-C-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-C-Chol is a cationic cholesterol derivative. OH-C-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.

Reference

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).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

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, 02/18/2019

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