

PRODUCT INFORMATION



MHAPC-Chol

Item No. 26585

CAS Registry No.: 1027801-74-6
Formal Name: (3 β)-cholest-5-en-3-ol, 3-[N-

[3-[(2-hydroxyethyl)methyl
amino]propyl]carbamate]

MF: C₃₄H₆₀N₂O₃

FW: 544.9

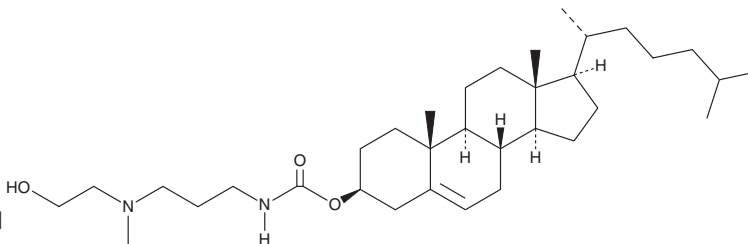
Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 219 nm

Supplied as: A 5 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥ 2 years



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

Laboratory Procedures

MHAPC-Chol is supplied as a 5 mg/ml solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. MHAPC-Chol is soluble in dimethyl formamide at a concentration of 10 mg/ml.

MHAPC-Chol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of MHAPC-chol should be diluted with the aqueous buffer of choice. MHAPC-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

MHAPC-Chol is a cationic cholesterol.¹ MHAPC-Chol, as part of a lipoplex with DOPE (Item No. 15091), has been used for siRNA delivery and gene silencing in MCF-7 cells in a luciferase assay without affecting cell viability. It has also been used to deliver siRNA into mice *via* intravenous injection, resulting in MHAPC-chol 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.

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, 11/24/2025

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