

PRODUCT INFORMATION

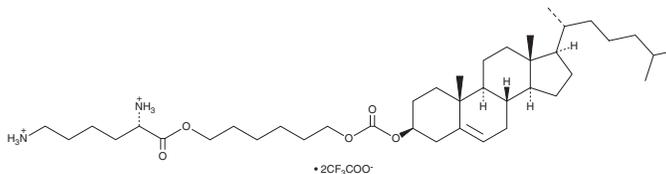


Cho-es-Lys

Item No. 37337

CAS Registry No.: 1356250-80-0
Formal Name: L-lysine, 6-[[[(3 β)-cholest-5-en-3-yloxy]carbonyl]oxy]hexyl ester, 2,2,2-trifluoroacetate

Synonym: CEL
MF: C₄₀H₇₂N₂O₅ • 2CF₃COO
FW: 887.1
Purity: ≥95%
Supplied as: A 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

Cho-es-Lys is supplied as a solid. A stock solution may be made by dissolving the cho-es-Lys in the solvent of choice, which should be purged with an inert gas. Cho-es-Lys is soluble in the organic solvent ethanol at a concentration of approximately 2 mg/ml.

Description

Cho-es-Lys is a cationic cholesterol derivative composed of cholesterol (Item No. 9003100) conjugated to an L-lysine head group containing two primary amines (pK_a s = 10.48 and 9.46) via a hexanedioyl linker.^{1,2} It has been used for transfection of Cy3-pDNA in COS-7 cells. Cho-es-Lys nanocomplexes containing luciferase siRNA reduce luciferase levels *in vitro*.²

References

- Sheng, R., Luo, T., Li, H., *et al.* Cholesterol-based cationic lipids for gene delivery: Contribution of molecular structure factors to physico-chemical and biological properties. *Colloids Surf. B Biointerfaces* **116**, 32-40 (2014).
- Zhu, Z., Zhang, L., Sheng, R., *et al.* Microfluidic-based cationic cholesterol lipid siRNA delivery nanosystem: Highly efficient *in vitro* gene silencing and the intracellular behavior. *Int. J. Mol. Sci.* **23(7)**, 3999 (2022).

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, 05/02/2023

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