

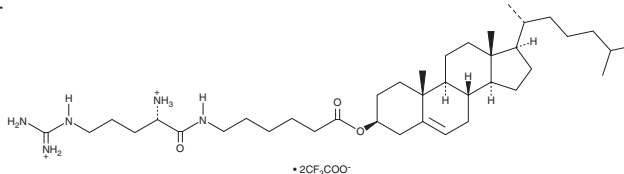
PRODUCT INFORMATION



Cho-Arg (trifluoroacetate salt)

Item No. 25943

CAS Registry No.: 1609010-56-1
Formal Name: cholest-5-en-3-ol (3 β)-3-[6-[[[(2S)-2-amino-5-[(aminoiminomethyl)amino]-1-oxopentyl]amino]hexanoate], 2,2,2-trifluoroacetate (1:2)
MF: C₃₉H₇₁N₅O₃ • 2CF₃COO
FW: 884.1
Purity: ≥95%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

Cho-Arg (trifluoroacetate salt) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of Cho-Arg (trifluoroacetate salt) in these solvents is approximately 50 mg/ml.

Cho-Arg (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of Cho-Arg (trifluoroacetate salt) should be diluted with the aqueous buffer of choice. Cho-Arg (trifluoroacetate salt) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method.

Description

Cho-Arg is a steroid-based cationic lipid that contains a cholesterol skeleton coupled to an L-arginine head group.¹ It forms a complex with plasmid DNA and decreases plasmid DNA migration in an agarose-gel retardant assay at charge ratios greater than or equal to 4. Cho-Arg facilitates transfection of plasmid DNA into H1299 and HeLa cells in the presence or absence of fetal bovine serum, an effect that is reversed by the lipid raft-mediated endocytosis inhibitor methyl- β -cyclodextrin (Item No. 21633) and the caveolae-mediated endocytosis inhibitor genistein (Item No. 10005167). It is cytotoxic to H1299 cells (IC₅₀ = 88.5 μ g/ml).

Reference

1. Sheng, R., Wang, Z., Luo, T., *et al.* Skeleton-controlled pDNA delivery of renewable steroid-based cationic lipids, the endocytosis pathway analysis and intracellular localization. *Int. J. Mol. Sci.* **19**(2), e369 (2018).

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, 03/25/2026

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