

PRODUCT INFORMATION



DOSEH (chloride)

Item No. 41601

Formal Name: (2-hydroxyethyl)dioctadecylsulfonium, monochloride

MF: C₃₈H₇₉OS • Cl

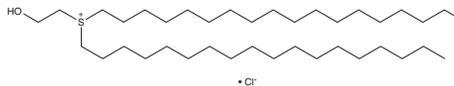
FW: 619.6

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

DOSEH (chloride) is supplied as a solid. A stock solution may be made by dissolving the DOSEH (chloride) in the solvent of choice, which should be purged with an inert gas. DOSEH (chloride) is slightly soluble (0.1-1 mg/ml) in acetonitrile and chloroform.

Description

DOSEH is a cationic sulfonium lipid that has been used in the generation of sulfonium lipid nanoparticles (sLNPs) for the delivery of mRNA *in vivo*.¹ LNPs containing DOSEH and encapsulating an mRNA reporter accumulate primarily in the lungs after intravenous administration in mice. DOSEH-containing LNPs encapsulating Cre mRNA induce the expression of tdTomato in the lungs of Ai14 mice engineered to express tdTomato upon Cre-mediated recombination.

Reference

1. Popoola, D.O., Cao, Z., Men, Y., *et al.* Lung-specific mRNA delivery enabled by sulfonium lipid nanoparticles. *Nano Lett.* **24(26)**, 8080-8088 (2024).

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, 10/24/2024

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