

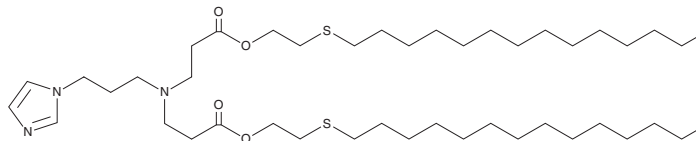
PRODUCT INFORMATION



93-O17S

Item No. 34367

CAS Registry No.: 2227008-67-3
Formal Name: bis(2-(tetradecylthio)ethyl)
3,3'-((3-(1H-imidazol-1-yl)
propyl)azanediyl)dipropionate
MF: C₄₄H₈₃N₃O₄S₂
FW: 782.3
Purity: ≥95%
Supplied as: A 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

93-O17S 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. 93-O17S is soluble in chloroform at a concentration of approximately 100 mg/ml.

Description

93-O17S is a chalcogen-containing ionizable cationic lipidoid.¹ It has been used in the synthesis of lipid nanoparticles (LNPs) for the delivery of Cre recombinase and ribonucleoproteins for genome editing in mice. LNPs containing 93-O17S have been used for the intratumoral delivery of cGAMP to enhance cross-presentation of tumor antigens and stimulation of interferon genes (STING) activation in a B16/F10 murine melanoma model.²

References

1. Li, Y., Yang, T., Yu, Y., *et al.* Combinatorial library of chalcogen-containing lipidoids for intracellular delivery of genome-editing proteins. *Biomaterials* **178**, 652-662 (2018).
2. Chen, J., Qiu, M., Ye, Z., *et al.* In situ cancer vaccination using lipidoid nanoparticles. *Sci. Adv.* **7(19)**, eabf1244 (2021).

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, 11/10/2022

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