

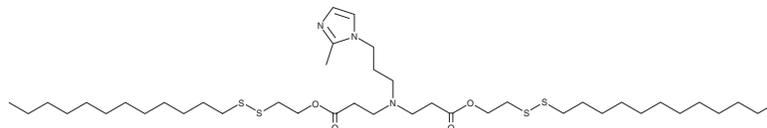
PRODUCT INFORMATION



9322-O16B

Item No. 40208

CAS Registry No.: 2909427-23-0
Formal Name: bis(2-(dodecylsulfaneyl)ethyl)
3,3'-((3-(2-methyl-1H-imidazol-1-yl)propyl)
azanediyl)dipropionate
MF: C₄₁H₇₇N₃O₄S₄
FW: 804.3
Purity: ≥95%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

9322-O16B 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. 9322-O16B is soluble in chloroform.

Description

9322-O16B is an ionizable cationic lipid.¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vitro*. LNPs containing 9322-O16B and encapsulating mRNA encoding chimeric antigen receptors (CARs) have been used in the generation of anti-CD19 CAR macrophages with anticancer activity.

Reference

1. Ye, Z., Chen, J., Zhao, X., *et al.* In vitro engineering chimeric antigen receptor macrophages and T cells by lipid nanoparticle-mediated mRNA delivery. *ACS Biomater. Sci. Eng.* **8(2)**, 722-733 (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, 11/18/2025

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