

PRODUCT INFORMATION



Lipid 8

Item No. 38746

CAS Registry No.: 2226547-25-5
Formal Name: 4-(dimethylamino)-butanoic acid,
2-[di-(9Z,12Z)-9,12-octadecadien-
1-ylamino]ethyl ester

MF: C₄₄H₈₂N₂O₂

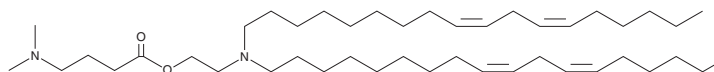
FW: 671.1

Purity: ≥95%

Supplied as: A 25 mg/ml solution in methyl acetate

Storage: -20°C

Stability: ≥2 years



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

Description

Lipid 8 is an ionizable cationic lipid that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).¹ LNPs containing lipid 8 and encapsulating polo-like kinase 1 (Plk1) siRNA induce apoptosis in U266 multiple myeloma cells.¹

Reference

1. Ramishetti, S., Hazan-Halevy, I., Chatterjee, S., *et al.* A combinatorial library of lipid nanoparticles for RNA delivery to leukocytes. *Adv. Mater.* **32(12)**, e1906128 (2020).

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, 11/11/2025

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