

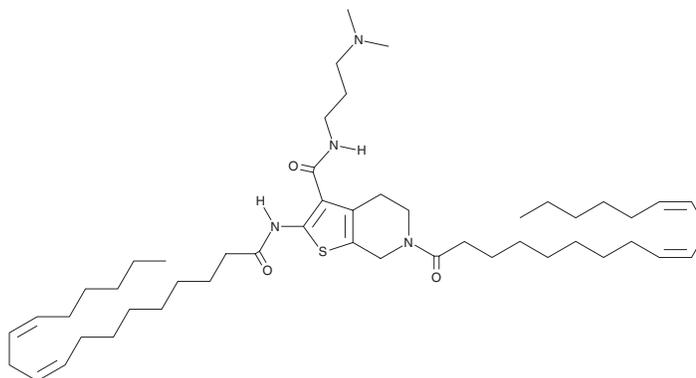
PRODUCT INFORMATION



Lipid 29d

Item No. 40913

CAS Registry No.: 2921586-67-4
Formal Name: N-[3-(dimethylamino)propyl]-4,5,6,7-tetrahydro-6-[(9Z,12Z)-1-oxo-9,12-octadecadien-1-yl]-2-[[[(9Z,12Z)-1-oxo-9,12-octadecadien-1-yl]amino]-thieno[2,3-c]pyridine-3-carboxamide
MF: C₄₉H₈₂N₄O₃S
FW: 807.3
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.

Description

Lipid 29d is an unsaturated ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ Intravenous administration of LNPs containing lipid 29d and encapsulating mRNA encoding a luciferase reporter induce bioluminescence in the lungs and spleen of mice.

Reference

1. Eygeris, Y., Gupta, M., Kim, J., *et al.* Thiophene-based lipids for mRNA delivery to pulmonary and retinal tissues. *Proc. Natl. Acad. Sci. USA* **121(11)**, e2307813120 (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, 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