

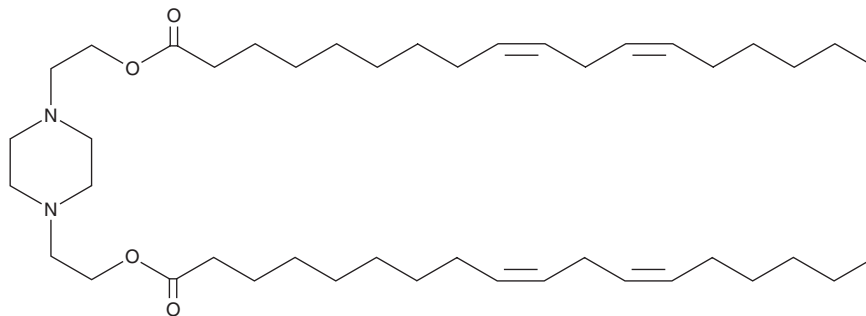
PRODUCT INFORMATION



AA3-DLin

Item No. 37903

CAS Registry No.: 3027565-91-6
Formal Name: piperazine-1,4-diylbis(ethane-2,1-diyl) (9Z,9'Z,12Z,12'Z)-bis(octadeca-9,12-dienoate)
MF: C₄₄H₇₈N₂O₄
FW: 699.1
Purity: ≥95%
Supplied as: A 25 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Description

AA3-DLin is an ionizable cationic amino lipid ($pK_a = 5.8$) that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA.¹ LNPs containing AA3-DLin and encapsulating mRNA for the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) spike glycoprotein induce immunogenicity in a mice.

Reference

1. Li, Z., Zhang, X.-Q., Ho, W., *et al.* Enzyme-catalyzed one-step synthesis of ionizable cationic lipids for lipid nanoparticle-based mRNA COVID-19 vaccines. *ACS Nano* **16(11)**, 18936–18950 (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, 11/06/2025

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