

PRODUCT INFORMATION



OF-C4-Deg-Lin

Item No. 37856

CAS Registry No.: 1853203-01-6
Formal Name: 9,12-octadecadienoic acid, (9Z,12Z)-1,1',1'',1'''-[(3,6-dioxo-2,5-piperazinediyl) bis(4,1-butanediyl)nitridi-4,1-butanediyl]] ester

MF: C₁₀₀H₁₇₆N₄O₁₀

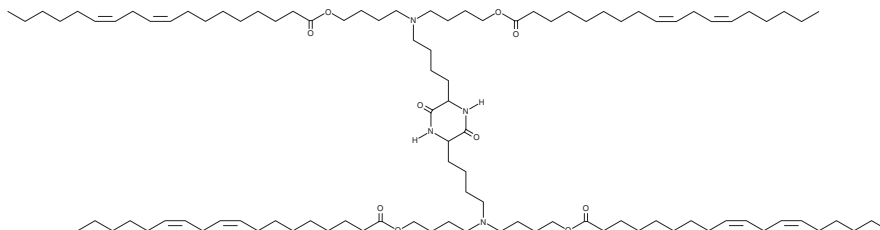
FW: 1,594.5

Purity: ≥90%

Supplied as: A 10 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

OF-C4-Deg-Lin is an ionizable cationic lipid.¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of siRNA and mRNA *in vitro* and *in vivo*. LNPs containing OF-C4-Deg-Lin and encapsulating an mRNA reporter selectively accumulate in mouse spleen after intravenous administration.

Reference

1. Fenton, O.S., Kauffman, K.J., McClellan, R.L., *et al.* Customizable lipid nanoparticle materials for the delivery of siRNAs and mRNAs. *Angew Chem. Int. Ed. Engl.* **57**(41), 13582-13586 (2018).

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