

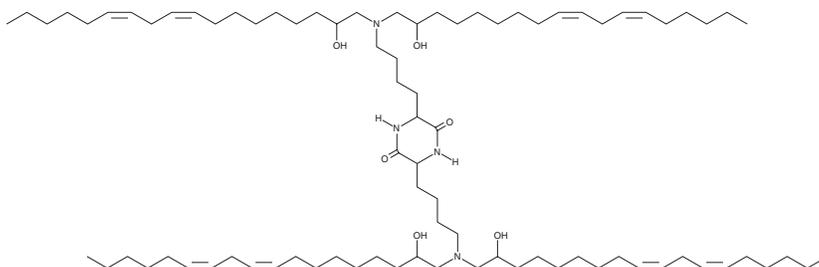
PRODUCT INFORMATION



OF-02

Item No. 37652

CAS Registry No.: 1883431-67-1
Formal Name: 3,6-bis[4-[bis[(9Z,12Z)-2-hydroxy-9,12-octadecadien-1-yl]amino]butyl]-2,5-piperazinedione
MF: C₈₄H₁₅₂N₄O₆
FW: 1,314.1
Purity: ≥90% (mixture of isomers)
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

OF-02 is an ionizable cationic amino lipid that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA.¹ LNPs containing OF-02 localize to the liver after intravenous administration in mice. Intravenous administration of LNPs containing OF-02 and encapsulating human erythropoietin (EPO) mRNA increases serum EPO levels in mice.

Reference

1. Fenton, O.S., Kauffman, K.J., McClellan, R.L., *et al.* Bioinspired alkenyl amino alcohol ionizable lipid materials for highly potent in vivo mRNA delivery. *Adv. Mater.* **28(15)**, 2939-2943 (2016).

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, 07/21/2025

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