

PRODUCT INFORMATION



RCB-4-8

Item No. 38803

CAS Registry No.: 2941228-91-5

Formal Name: bis((Z)-12-(((non-2-yn-1-yloxy)carbonyl)oxy)octadec-9-en-1-yl) 3,3'-((2-(diethylamino)ethyl)azanediyl)dipropionate

MF: $C_{68}H_{120}N_2O_{10}$

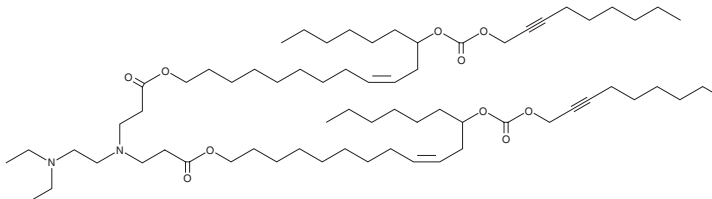
FW: 1,125.7

Purity: $\geq 95\%$

Supplied as: A 50 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

RCB-4-8 is an ionizable cationic lipid that has been used in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA.¹ RCB-4-8-containing LNPs improve lung transfection efficiency by approximately 100-fold over MC3-containing LNPs in mice.

Reference

1. Li, B., Manan, R.S., Liang, S.-Q., *et al.* Combinatorial design of nanoparticles for pulmonary mRNA delivery and genome editing. *Nat. Biotechnol.* Online ahead of print, (2023).

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, 11/12/2025

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