

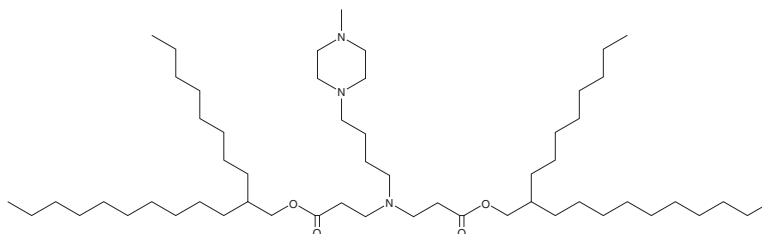
PRODUCT INFORMATION



Lipid C24

Item No. 37122

CAS Registry No.: 2767561-52-2
Formal Name: bis(2-octyldodecyl)
3,3'-((4-(4-methylpiperazin-1-yl)butyl)
azanediyldipropionate
MF: $C_{55}H_{109}N_3O_4$
FW: 876.5
Purity: $\geq 95\%$
Supplied as: A 5 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 C24 is a cationic ionizable lipid.¹ It has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).

Reference

1. Buschmann, M.D., Paige, M., Alishetty, S., *et al.* Ionizable lipids and methods of manufacture and use thereof. George Mason Research Foundation, Inc. and The Trustees of The University of Pennsylvania. WO2022/081750A1 (2021).

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