

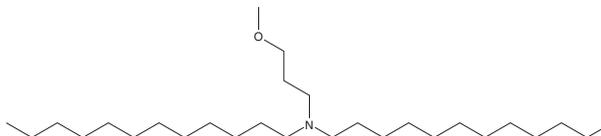
PRODUCT INFORMATION



AL-A12

Item No. 38001

CAS Registry No.: 3037065-56-5
Formal Name: N-dodecyl-N-(3-methoxypropyl)-1-dodecanamine
MF: C₂₈H₅₉NO
FW: 425.8
Purity: ≥95%
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

AL-A12 is an ionizable cationic amino lipid that has been used in the formation of lipid nanoparticles (LNPs).¹ LNPs containing AL-A12 and encapsulating minicircle DNA that encodes for GFP have been used to induce GFP expression in Huh7 cells. LNPs containing AL-A12 and encapsulating DNA that encodes for RFP have been used to induce RFP expression in endothelial cells and venous vasculature in a zebrafish embryo model.

Reference

1. Buck, J., Mueller, D., Mettal, U., *et al.* Improvement of DNA vector delivery of DOTAP lipoplexes by short-chain aminolipid. *ACS Omega* **5(38)**, 24724-24732 (2020).

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