

PRODUCT INFORMATION



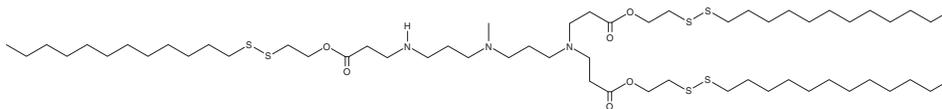
A1-28

Item No. 39590

CAS Registry No.: 3055572-07-8
Formal Name: bis(2-(dodecylsulfaneyl) ethyl) 3,3'-((4-methyl-11-oxo-12-oxa-15,16-dithia-4,8-diazaoctacosyl) azanediyl)dipropionate

Synonym: AMAP-O16B
MF: C₅₈H₁₁₅N₃O₆S₆
FW: 1,142.9
Purity: ≥95%

Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

A1-28 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. A1-28 is soluble in ethanol and chloroform.

Description

A1-28 is a disulfide bond-containing ionizable cationic lipid.¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of CRISPR complementary single-guide RNA (sgRNA) and Cas9 mRNA for genome editing *in vitro*.

Reference

1. Liu, J., Chang, J., Jiang, Y., *et al.* Fast and efficient CRISPR/Cas9 genome editing in vivo enabled by bio-reducible lipid and messenger RNA nanoparticles. *Adv. Mater.* **31(33)**, e1902575 (2019).

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, 01/26/2026

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