

PRODUCT INFORMATION



BAmP-O16B

Item No. 40473

Formal Name: bis(2-(dodecylsulfaneyl)ethyl)
3,3'-((3-(4-(3-((3-(2-(dodecylsulfaneyl)
ethoxy)-3-oxopropyl)amino)propyl)
piperazin-1-yl)propyl)azanediyl)
dipropionate

MF: C₆₁H₁₂₀N₄O₆S₆

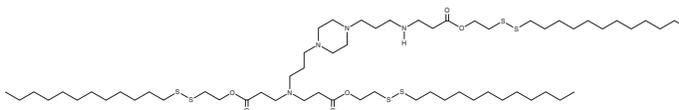
FW: 1,198.0

Purity: ≥95%

Supplied as: A 10 mg/ml solution in methyl acetate

Storage: -20°C

Stability: ≥3 years



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

Laboratory Procedures

BAmP-O16B 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. BAmP-O16B is soluble in dichloromethane and chloroform.

Description

BAmP-O16B is an ionizable cationic amino lipid that has been used in the generation of lipid nanoparticles (LNPs).¹ LNPs containing BAmP-O16B and encapsulating NSA-inactivated RNase A reduce the viability of SiHa and HeLa cancer cells, but not non-cancerous NIH3T3 and HEK293T cells, upon intracellular reactivation of RNase A by glutathione (GSH).

Reference

1. Zheng, Q., Ma, T., and Wang, M. Unleashing the power of proenzyme delivery for targeted therapeutic applications using biodegradable lipid nanoparticles. *Acc. Chem. Res.* **57**(2), 208-221 (2024).

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, 03/25/2026

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