

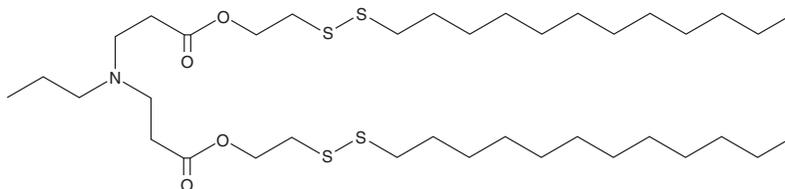
PRODUCT INFORMATION



PPPDA-O16B

Item No. 40517

Formal Name: bis(2-(dodecylsulfaneyl)ethyl)
3,3'-(propylazanediy)ldipropionate
MF: C₃₇H₇₃NO₄S₄
FW: 724.2
Purity: ≥98%
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

PPPDA-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. PPPDA-O16B is soluble in chloroform.

Description

PPPDA-O16B is a disulfide bond-containing ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for plasmid delivery *in vitro* and *in vivo*.¹ LNPs containing PPPDA-O16B and encapsulating plasmids encoding Ras/Rap1-specific endopeptidase (RRSP) increase RRSP levels and decrease Ras levels in, and reduce the viability of, HeLa cervical cancer cells. LNPs containing PPPDA-O16B and encapsulating plasmids encoding RRSP selectively localize to tumors over the heart, liver, spleen, lungs, and kidneys, as well as decrease tumor volume without reducing body weight, in an HCT116 colorectal cancer mouse xenograft model.

Reference

1. Li, W., Shao, L., Liu, J., et al. Intracellular delivery of bacterial effectors for cancer therapy using biodegradable lipid nanoparticles. *Biomater. Sci.* **11(9)**, 3172-3179 (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, 11/18/2025

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