

PRODUCT INFORMATION



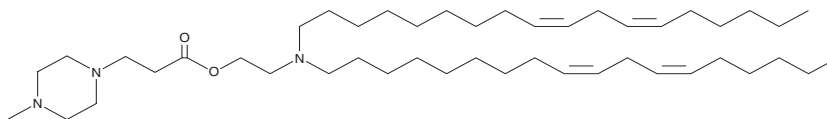
Lipid 10

Item No. 38705

CAS Registry No.: 2430034-02-7
Formal Name: 4-methyl-1-piperazinepropanoic acid,
2-[di-(9Z,12Z)-9,12-octadecadien-1-ylamino]ethyl ester

Synonym: EA-PIP
MF: C₄₆H₈₅N₃O₂
FW: 712.2

Purity: ≥95%
Supplied as: A 25 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.

Description

Lipid 10 is an ionizable cationic lipid ($pK_a = 6.2-6.5$).¹ Lipid nanoparticles (LNPs) containing lipid 10 and encapsulating siRNA for the gene encoding cytoskeleton-associated protein 5 (CKAP5) reduce its gene expression and protein levels *in vitro* and selectively reduce cell viability in a variety of genetically unstable cancer cell lines.² Administration of LNPs containing lipid 10 and encapsulating CKAP5 siRNA reduce tumor growth and increase survival in an NCI/ADR-RES human ovarian cancer mouse xenograft model.

References

1. Ramishetti, S., Hazan-Halevy, I., Chatterjee, S., *et al.* A combinatorial library of lipid nanoparticles for RNA delivery to leukocytes. *Adv. Mater.* **32(12)**, e1906128 (2020).
2. Chatterjee, S., Naidu, G.S., Hazan-Halevy, I., *et al.* Therapeutic gene silencing of CKAP5 leads to lethality in genetically unstable cancer cells. *Sci. Adv.* **9(14)**, eade4800 (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, 11/11/2025

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