

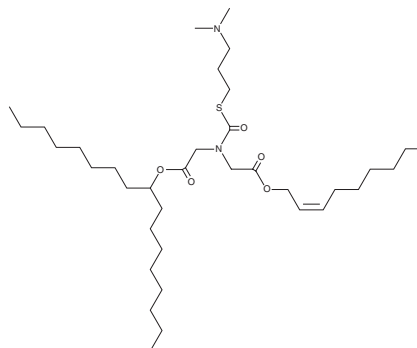
PRODUCT INFORMATION



Lipid 2,2(8,8) 4C CH₃

Item No. 37910

CAS Registry No.: 2230647-30-8
Formal Name: heptadecan-9-yl (Z)-N-(((4-(dimethylamino)butyl)thio)carbonyl)-N-(2-(non-2-en-1-yloxy)-2-oxoethyl)glycinate
Synonym: ATX-0114
MF: C₃₇H₇₀N₂O₅S
FW: 655.0
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.

Laboratory Procedures

Lipid 2,2(8,8) 4C CH₃ is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Lipid 2,2(8,8) 4C CH₃ is soluble in methanol and DMSO.

Description

Lipid 2,2(8,8) 4C CH₃ is an ionizable cationic lipid (pK_a = 6.69).¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of siRNA *in vivo*. LNPs containing lipid 2,2(8,8) 4C CH₃ and encapsulating siRNA targeting Factor VII decrease plasma Factor VII protein levels by 90% in mice.

Reference

1. Payne, J.E., Chivukula, P., Tanis, S.P., *et al.* Ionizable cationic lipid for RNA delivery. **WO2018/119163A1** (2018).

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, 11/06/2025

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