

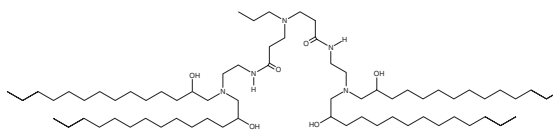
PRODUCT INFORMATION



C3-K2-E14

Item No. 39323

CAS Registry No.: 2933215-86-0
Formal Name: 3,3'-(propylazanediyl)bis(N-(2-(bis(2-hydroxytetradecyl)amino)ethyl)propanamide)
MF: $C_{69}H_{141}N_5O_6$
FW: 1,136.9
Purity: $\geq 98\%$
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥ 1 year



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

Laboratory Procedures

C3-K2-E14 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. Solvents such as chloroform purged with an inert gas can be used.

Description

C3-K2-E14 is an ionizable cationic lipid ($pK_a = 5.5$).¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA or siRNA *in vitro* and *in vivo*. LNPs containing C3-K2-E14 and encapsulating siRNA targeting the gene encoding colony-stimulating factor 1 (Csf1) reduce the percentage of circulating Ly6C^{hi} monocytes and increase the percentage of circulating Ly6C^{int} monocytes in mice.

Reference

1. Manning, A.M., Tilstra, G., Khan, A.B., *et al.* Ionizable lipid with supramolecular chemistry features for RNA delivery *in vivo*. *Small* **19**(41), e2302917 (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, 11/13/2025

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