

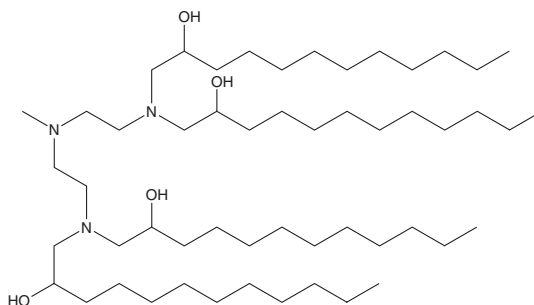
PRODUCT INFORMATION



C12-113

Item No. 39335

CAS Registry No.: 1220890-27-6
Formal Name: 1,1',1'',1'''-[(methylimino)bis(2,1-ethanediyl)nitri]o]tetrakis-2-dodecanol
MF: $C_{53}H_{111}N_3O_4$
FW: 854.5
Purity: $\geq 95\%$
Supplied as: A 25 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

C12-113 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 DMSO, purged with an inert gas can be used. The solubility of C12-113 in DMSO is approximately 20 mg/ml.

Description

C12-113 is an ionizable cationic lipidoid.¹ It has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA encoding severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) spike glycoprotein in mice.

Reference

1. Han, X., Alameh, M.-G., Butowska, K., et al. Adjuvant lipidoid-substituted lipid nanoparticles augment the immunogenicity of SARS-CoV-2 mRNA vaccines. *Nat. Nanotechnol.* (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, 03/26/2026

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