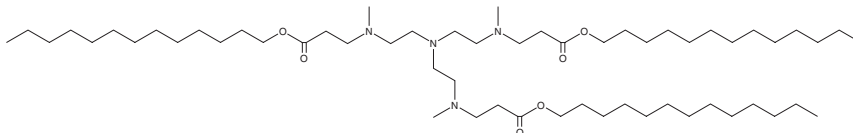


PRODUCT INFORMATION



304O₁₃
Item No. 38584

CAS Registry No.: 1566559-80-5
Formal Name: N,N',N''-(nitrilotri-2,1-ethanediyl) tris[N-methyl-β-alanine, 1,1',1''-tridecyl ester]
MF: C₅₇H₁₁₄N₄O₆
FW: 951.5
Purity: ≥95%
Supplied as: A 10 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.

Description

304O₁₃ is an ionizable cationic lipidoid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of siRNA *in vivo*.¹ LNPs containing 304O₁₃ and encapsulating siRNA targeting Factor VII mRNA decrease serum Factor VII levels in mice.

Reference

1. Whitehead, K.A., Dorkin, J.R., Vegas, A.J., *et al.* Degradable lipid nanoparticles with predictable *in vivo* siRNA delivery activity. *Nat. Commun.* **5**, 4277 (2014).

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, 03/16/2026

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