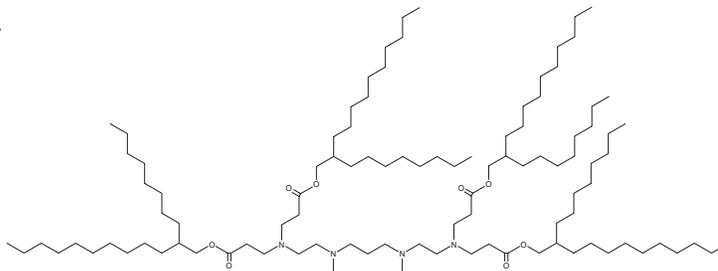


PRODUCT INFORMATION



503O_{8,12}
Item No. 44137

CAS Registry No.: 3067564-98-8
Formal Name: bis(2-octyldodecyl) 7,11-dimethyl-4,14-bis(3-((2-octyldodecyl)oxy)-3-oxopropyl)-4,7,11,14-tetraazaheptadecanedioate
MF: C₁₀₁H₂₀₀N₄O₈
FW: 1,598.7
Purity: ≥90%
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

503O_{8,12} 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. 503O_{8,12} is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

503O_{8,12} is an ionizable cationic lipidoid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ LNPs containing 503O_{8,12} and encapsulating an mRNA reporter traffic to the spleen in mice.

Reference

1. Petersen, D.M.S., Weiss, R.M., Hajj, K.A., *et al.* Branched-tail lipid nanoparticles for intravenous mRNA delivery to lung immune, endothelial, and alveolar cells in mice. *Adv. Healthc. Mater.* **13(22)**, e2400225 (2024).

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, 09/03/2025

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