

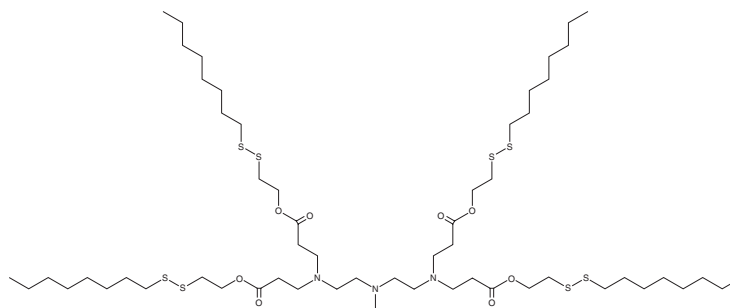
PRODUCT INFORMATION



113-O12B

Item No. 37671

CAS Registry No.: 2803699-72-9
Formal Name: tetrakis(2-(octyldisulfaneyl)ethyl) 3,3',3'',3'''-(((methylazanediy))bis(ethane-2,1-diyl))bis(azanetriyl))tetrapropionate
MF: C₅₇H₁₁₁N₃O₈S₈
FW: 1,223.0
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

113-O12B is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. 113-O12B is soluble in the organic solvent ethanol at a concentration of approximately 10 mg/ml.

Description

113-O12B is a disulfide bond-containing ionizable cationic lipidoid.¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA. Subcutaneous administration of LNPs containing 113-O12B and encapsulating an mRNA reporter accumulate specifically in mouse lymph nodes. LNPs containing 113-O12B and encapsulating mRNA encoding ovalbumin reduce tumor growth and increase the number of tumoral CD8⁺ T cells, M1-like macrophages, and activated dendritic cells in an ovalbumin-expressing B16/F10 murine melanoma model.

Reference

1. Chen, J., Ye, Z., Huang, C., *et al.* Lipid nanoparticle-mediated lymph node-targeting delivery of mRNA cancer vaccine elicits robust CD8⁺ T cell response. *Proc. Natl. Acad. Sci. USA* **119**(34), e2207841119 (2022).

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, 10/31/2023

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