

PRODUCT INFORMATION



VL422

Item No. 39874

CAS Registry No.: 2765519-47-7
Formal Name: octanedioic acid,

1,1'-[2-[[[3-(diethylamino)propoxy] carbonyl]oxy]methyl]-1,3-propanediyl] 8,8'-bis(1-octylnonyl) ester

MF: C₆₂H₁₁₇NO₁₁

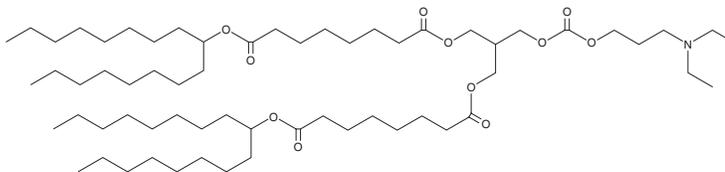
FW: 1,052.6

Purity: ≥95%

Supplied as: A 10 mg/ml 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

VL422 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. VL422 is soluble in chloroform.

Description

VL422 is an ionizable cationic lipid.¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of CRISPR complementary single-guide RNA (sgRNA) and *Cas9* mRNA for gene editing *in vitro* and *in vivo*. LNPs containing VL422 and encapsulating *Cas9* mRNA and sgRNA targeting the gene encoding angiopoietin-related protein 3 (ANGPTL3), a protein whose loss-of-function decreases LDL, HDL, and cholesterol plasma levels, induce a deletion in a premature stop codon in ANGPTL3 in the livers of cynomolgus monkeys.

Reference

1. Rajeev, K.G., Biswas, S., Kasiewicz, L.N., *et al.* Lipid formulations for gene editing. *Verve Therapeutics, Inc.* US2021/050511 (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, 01/26/2026

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