

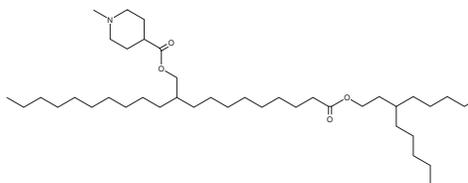
PRODUCT INFORMATION



L202

Item No. 37841

CAS Registry No.: 2170488-92-1
Formal Name: 1-methyl-4-piperidinecarboxylic acid,
2-[9-oxo-9-[(3-pentyl-octyl)oxy]nonyl]
dodecyl ester
MF: C₄₁H₇₉NO₄
FW: 650.1
Purity: ≥95%
Supplied as: A 5 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

L202 is an ionizable cationic lipid ($pK_a = 6.04$) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA.¹ Intramuscular administration of LNPs containing L202 and encapsulating nucleoside-modified mRNA encoding pre-fusion stabilized severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) spike glycoprotein, also known as surface glycoprotein, increase plasma levels of IgG antibodies targeting SARS-CoV-2 spike glycoprotein in mice and cynomolgus monkeys.

Reference

1. Suzuki, Y., Miyazaki, T., Muto, H., *et al.* Design and lyophilization of lipid nanoparticles for mRNA vaccine and its robust immune response in mice and nonhuman primates. *Mol. Ther. Nucleic Acids* **30**, 226-240 (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, 11/06/2025

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