

PRODUCT INFORMATION



LipidLaunch™ Ionizable Lipid LNP Screening Set Item No. 42832

The LipidLaunch™ Ionizable Lipid LNP Screening Set includes 12 ionizable cationic and cationic lipids to screen for use in lipid nanoparticle (LNP) formulations. The lipids are formulated at 10 mg/ml in ethanol for ease of addition to lipid mixes for LNP formulation. Lipids selected for this screening set include FDA-approved excipients, as well as some of the most commonly cited lipids used in nucleic acid-cargo LNPs. LNPs containing these lipids in combination with a sterol lipid, a phospholipid, and a PEGylated lipid have been used to encapsulate mRNA, plasmid DNA, siRNA, and other nucleic acid cargo for intracellular delivery. For more information, please see Cayman's LNP guide: Lipid Nanoparticle Formulation: Basic Concepts & Preparation Procedures. The composition of this lipid set may vary somewhat depending upon our inventory. For an accurate list of products in the current batch, please contact techserv@caymanchem.com.

Contents of LipidLaunch™ Ionizable Lipid LNP Screening Set

Item No.	Item Name	Size
33474	SM-102	5 mg
34337	ALC-0315	5 mg
34364	DLin-MC3-DMA	5 mg
36699	C12-200	5 mg
36700	cKK-E12	5 mg
37278	LP-01	5 mg
37667	Lipid A9	5 mg
15110	1,2-Dioleoyl-3-trimethylammoniumpropane (chloride)	5 mg
38320	Lipid CL1	5 mg
34363	DLin-KC2-DMA	5 mg
34372	Lipid 5	5 mg
35337	Lipid 29	5 mg

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, 01/13/2025

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