

PRODUCT INFORMATION



LipidLaunch™ Ionizable Lipid LNP Lung Screening Set Item No. 43435

The LipidLaunch™ Ionizable Lipid LNP Lung Screening Set includes 10 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 mixtures for LNP formulation. This screening set includes lipids reported to show delivery to the lung when 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. This screening set can be used in conjunction with LipidLaunch™ Helper Lipid Set (Item No. 42831) 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 Lung Screening Set

Item No.	Item Name	Size
34364	DLin-MC3-DMA	5 mg
37278	LP-01	5 mg
36699	C12-200	5 mg
15110	1,2-Dioleoyl-3-trimethylammoniumpropane (chloride)	50 mg
39284	244cis	5 mg
40420	IR-117-17	5 mg
42922	FO-32	5 mg
38155	4A3-SC8	5 mg
40493	Lipid 88	5 mg
46050	Lipid 392	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, 06/19/2026

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