

PRODUCT INFORMATION



LipidLaunch™ T Cell LNP Screening Set Item No. 43932

The LipidLaunch™ T Cell LNP Screening Set includes ten ionizable cationic lipids to screen for use in lipid nanoparticle (LNP) formulations for T cell transfection. 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 some of the most commonly cited ionizable lipids used in nucleic acid-cargo LNPs for transfection of activated T cells. LNPs containing these ionizable cationic 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 to T cells. For more information, please see Cayman's LNP guide: Lipid Nanoparticle Formulation: Basic Concepts & Preparation Procedures.

Contents of LipidLaunch™ T Cell LNP Screening Set

Item No.	Item Name	Size
33474	SM-102	5 mg
34364	DLin-MC3-DMA	5 mg
36699	C12-200	5 mg
36700	cKK-E12	5 mg
37667	Lipid A9	5 mg
38942	C14-4	5 mg
40805	C14-O2	5 mg
43084	E10i-494	5 mg
39783	Lipid C2	5 mg
40208	9322-O16B	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, 08/29/2025

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