

PRODUCT INFORMATION



LipidLaunch™ Helper Lipid Set Item No. 42831

The LipidLaunch™ Helper Lipid Set includes cholesterol, 1,2-dimyristoyl-*rac*-3-PEG(2000) (DMG-PEG(2000)), and 1,2-distearoyl-*sn*-glycero-3-PC (1,2-DSPC) for use in lipid nanoparticle (LNP) formulations. Cholesterol and 1,2-DSPC are formulated at 25 mM and DMG-PEG(2000) at 10 mg/ml in ethanol for ease of addition to LNP formulations. These lipids can be combined with cationic or ionizable cationic lipids to make a lipid mixture capable of encapsulating nucleic acids. Please note that warming to 37°C for up to 30 minutes may be required to ensure complete dissolution of these lipids in solvent. LNPs containing these lipids in combination with a cationic or ionizable cationic 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.

Contents of LipidLaunch™ Helper Lipid Set (Small)

Item No.	Item Name	Size
39189	1,2-Distearoyl- <i>sn</i> -glycero-3-PC (solution)	5 ml
39088	Cholesterol (solution)	5 ml
42746	DMG-mPEG(2000) (solution)	1 ml

Contents of LipidLaunch™ Helper Lipid Set (Large)

Item No	Item Name	Size
39189	1,2-Distearoyl- <i>sn</i> -glycero-3-PC (solution)	25 ml
39088	Cholesterol (solution)	50 ml
42746	DMG-mPEG(2000) (solution)	10 ml

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, 10/24/2025

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