

PRODUCT INFORMATION



LipidLaunch™ Phospholipid LNP Screening Set Item No. 43840

The LipidLaunch™ Phospholipid LNP Screening Set includes 12 phospholipids and one sphingolipid to screen for use in lipid nanoparticle (LNP) and liposomal formulations. Each lipid in the set is supplied in a vial containing 10 mg of lipid. Lipids selected for this screening set are some of the most cited phospholipids/sphingolipids used in nucleic acid-cargo LNPs and liposomal formulations. LNPs containing these lipids in combination with a cationic and/or ionizable cationic lipid, sterol lipid, and PEGylated lipid have been used to encapsulate mRNA, plasmid DNA, siRNA, and other nucleic acid cargo for intracellular delivery. In addition, liposomes consisting primarily of the phospholipids included in the set are commonly used for *in vitro* or *in vivo* delivery of lipophilic or hydrophobic compounds. For more information, please see Cayman's LNP guide: Lipid Nanoparticle Formulation: Basic Concepts & Preparation Procedures.

Contents of LipidLaunch™ Phospholipid LNP Screening Set

Item No.	Item Name	Size
15100	1,2-Distearoyl- <i>sn</i> -glycero-3-PC	10 mg
15095	1,2-Distearoyl- <i>sn</i> -glycero-3-PE	10 mg
15098	1,2-Dioleoyl- <i>sn</i> -glycero-3-PC	10 mg
15091	1,2-Dioleoyl- <i>sn</i> -glycero-3-PE	10 mg
15102	1-Palmitoyl-2-Oleoyl- <i>sn</i> -glycero-3-PC	10 mg
20067	1-Palmitoyl-2-Oleoyl- <i>sn</i> -glycero-3-PE	10 mg
10009473	1,2-Dipalmitoyl- <i>sn</i> -glycero-3-PC	10 mg
15092	1,2-Dipalmitoyl- <i>sn</i> -glycero-3-PE	10 mg
15097	1,2-Dimyristoyl- <i>sn</i> -glycero-3-PC	10 mg
15090	1,2-Dimyristoyl- <i>sn</i> -glycero-3-PE	10 mg
38152	1-Stearoyl-2-Oleoyl- <i>sn</i> -glycero-3-PC	10 mg
38153	1-Stearoyl-2-Oleoyl- <i>sn</i> -glycero-3-PE	10 mg
24345	Sphingomyelin (chicken egg)	10 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, 05/20/2025

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