

PRODUCT INFORMATION



LipidLaunch™ Cholesterol Analog LNP Screening Set Item No. 43889

The LipidLaunch™ Cholesterol Analog LNP Screening Set includes two types of cholesterol, five cholesterol analogs, and six bile acids to screen for use in lipid nanoparticle (LNP) formulations. Each lipid in the set is supplied in a vial containing 10 mg of lipid for ease of addition to lipid mixes for LNP formulation. Lipids selected for this screening set are some of the most cited cholesterol analogs and bile acids used in nucleic acid-cargo LNPs. LNPs containing these lipids in combination with a cationic and/or ionizable cationic lipid, phospholipid, and 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.

Contents of LipidLaunch™ Cholesterol Analog LNP Screening Set

Item No.	Item Name	Size
11031	6-ECDCA	10 mg
9003100	Cholesterol	10 mg
11686	Betulinic Acid	10 mg
11726	Oleanolic Acid	10 mg
39448	Cholesterol (plant)	10 mg
15121	Ursodeoxycholic Acid (sodium salt)	10 mg
34623	Norcholic Acid	10 mg
18079	Stigmasterol	10 mg
20098	7 α -hydroxy Cholesterol	10 mg
20250	Cholic Acid	10 mg
20253	Lithocholic Acid	10 mg
37386	24 α -ethyl Cholesterol	10 mg
10072	Ursolic Acid	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, 05/27/2025

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