

PRODUCT INFORMATION

LipidLaunch™ LNP Formulation Buffer (50 mM Acetate, pH 4.5)

Item No. 400813

Overview and Properties

Storage: -20°C (as supplied)
Stability: ≥2 years
Supplied in: 50 mM acetate, pH 4.5

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

LipidLaunch™ LNP Formulation Buffer (50 mM Acetate, pH 4.5) is an acidic aqueous buffer designed to optimally encapsulate nucleic acid cargo within lipid nanoparticles (LNPs). It consists of an RNase-free (by lateral flow assay) solution of 50 mM sodium acetate at pH 4.5. mRNA, or other nucleic acid cargo, can be diluted directly in LipidLaunch™ LNP Formulation Buffer (50 mM Acetate, pH 4.5) for subsequent use with a selected lipid mixture. For more information on methods and complementary products, please see Cayman's LNP guide: Lipid Nanoparticle Formulation: Basic Concepts & Preparation Procedures or Cayman's LNP resource center.

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, 07/31/2025

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