

PRODUCT INFORMATION



LipidLaunch™ ALC-0315 LNP (Luciferase)

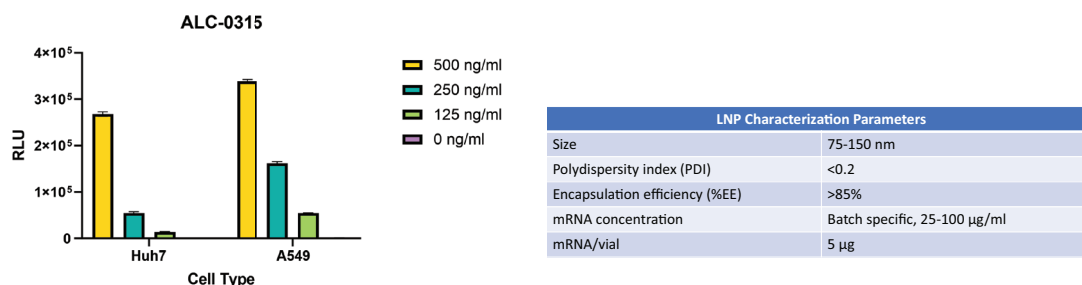
Item No. 40104

Overview and Properties

Storage: -80°C (as supplied)
Stability: ≥6 months
Supplied in: TBS, pH 7.5, with 10% sucrose

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

Images



Transfection of cells with luciferase-encapsulating ALC-0315 particles. Huh7 hepatocytes (left) and A549 lung epithelial cells (right) were incubated with Cayman's LipidLaunch™ ALC-0315 LNP (Luciferase) (Item No. 40104) at indicated RNA concentrations for 24 hours in cell culture media with 10% FBS. Luciferase expression (RLU) is shown.

Description

LipidLaunch™ ALC-0315 LNP (Luciferase) is a solution containing lipid nanoparticles (LNPs) composed of the ionizable cationic amino lipid ALC-0315 (Item No. 34337), cholesterol (Item Nos. 9003100 | 39088), the phospholipid 1,2-distearoyl-*sn*-glycero-3-PC (1,2-DSPC; Item Nos. 15100 | 39189), and the lipid excipient ALC-0159 (Item No. 34336) at a molar ratio of 46.3:42.7:9.4:1.6 and encapsulating mRNA encoding firefly luciferase protein. It is intended for proof-of-concept experiments to determine whether ALC-0315-based LNPs can effectively lead to the expression of a protein of interest in a target cell type, either *in vitro* or *in vivo*. Firefly luciferase is an enzyme that catalyzes the oxidation of ATP-dependent D-luciferin, resulting in chemiluminescence at a wavelength of approximately 560 nm.

Suggested *in vitro* use: Thaw LNPs on ice with occasional gentle swirling (*do not vortex*). Using a gentle pipetting technique, dilute 1:100-1:500 in complete cell culture media (with serum) and add to subconfluent cells in a luminescence-compatible tissue culture plate. Use a luciferase reporter assay substrate and read luminescence on a plate reader. Optimal conditions are highly dependent on cell type.

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, 04/16/2024

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