

PRODUCT INFORMATION



LipidLaunch™ SM-102 LNP (human cGAS)

Item No. 44930

Overview and Properties

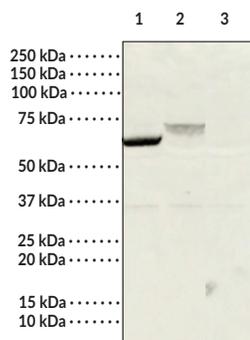
Storage: -80°C (as supplied)

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



Lane 1: 20 µg cGAS cell lysate (human)
Lane 2: 20 µg cGAS cell lysate (naked mole-rat)
Lane 3: 20 µg non-transfected cell lysate control

A549 cells were transfected with 250 ng/ml of LipidLaunch™ SM-102 LNP (human cGAS). Western blot analysis was performed on 20 µg of cell lysates using Cayman's His-Tag Recombinant Monoclonal Antibody - HRP Conjugated (Item No. 35511).

LNP Characterization Parameters	
Size	75-150 nm
Polydispersity index (PDI)	<0.25
Encapsulation efficiency (%EE)	>85%
mRNA concentration	Batch specific, 25-100 µg/ml
mRNA/vial	5 µg

Description

LipidLaunch™ SM-102 LNP (human cGAS) is a solution containing lipid nanoparticles (LNPs) composed of the ionizable cationic aminolipid SM-102 (Item No. 33474), cholesterol (Item No. 9003100), 1,2-distearoyl-*sn*-glycero-3-PC (1,2-DSPC; Item No. 15100), and the lipid excipient DMG-PEG(2000) (Item No. 33945) at a molar ratio of 50:38.5:10:1.5 and encapsulating mRNA encoding His-tagged human cyclic GMP-AMP (cGAMP) synthase (cGAS), a nucleotidyltransferase involved in the innate immune response.^{1,2} It is intended for the overexpression of His-tagged human wild-type cGAS *in vitro* and can be used in a variety of model cell lines.

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. Expression of His-tagged human cGAS may be detected between six and 48 hours after transfection. Optimal conditions are highly dependent on cell type.

References

1. Sun, L., Wu, J., Du, F., *et al.* Cyclic GMP-AMP synthase is a cytosolic DNA sensor that activates the type I interferon pathway. *Science* **339**(6121), 786-791 (2013).
2. Wu, J., Sun, L., Chen, X., *et al.* Cyclic GMP-AMP is an endogenous second messenger in innate immune signaling by cytosolic DNA. *Science* **339**(6121), 826-830 (2013).

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, 02/10/2026

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