

PRODUCT INFORMATION



BCP-NC2-C12

Item No. 42940

Formal Name: N¹,N³-bis(2-(didodecylamino)ethyl)bicyclo[1.1.1]pentane-1,3-dicarboxamide

MF: C₅₉H₁₁₆N₄O₂

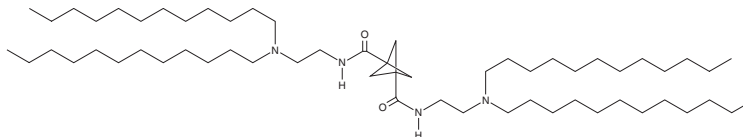
FW: 913.6

Purity: ≥98%

Supplied as: A 25 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Description

BCP-NC2-C12 is an ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ LNPs containing BCP-NC2-C12 accumulate primarily in the liver and spleen of mice. LNPs containing BCP-NC2-C12 and encapsulating mRNA encoding the Cas9 nuclease and small-guide RNA (sgRNA) targeting the gene encoding proprotein convertase subtilisin kexin type 9 (Pcsk9) induce insertion and deletion (indel) formation and reduce serum levels of Pcsk9 in mice.

Reference

1. Wu, S., Yang, Y., Lian, X., *et al.* Isosteric 3D bicyclo[1.1.1]pentane (BCP) core-based lipids for mRNA delivery and CRISPR/Cas gene editing. *J. Am. Chem. Soc.* **146**(50), 34733-34742 (2024).

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, 11/24/2025

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