

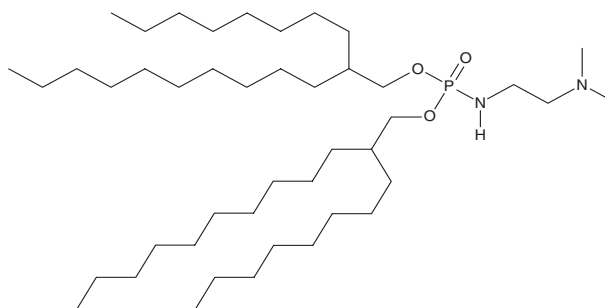
PRODUCT INFORMATION



BiP-20

Item No. 45618

Formal Name: bis(2-octyldodecyl)
(2-(dimethylamino)ethyl)
phosphoramidate
MF: C₄₄H₉₃N₂O₃P
FW: 729.2
Purity: ≥95%
Supplied as: A 50 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Description

BiP-20 is an ionizable cationic lipid that has been used in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA and single-guide RNA (sgRNA) *in vivo*.¹ LNPs containing BiP-20 are targeted to the liver. BiP-20-containing LNPs encapsulating Cas9 mRNA and transthyretin (TTR) sgRNA decrease serum levels of TTR in mice to a greater extent than LNPs containing LP-01 (Item No. 37278).

References

1. Jozić, A., Le Roux, C., Kim, J., *et al.* *In vivo* endosomal escape assay identifies mechanisms for efficient hepatic LNP delivery. *Nat. Biotechnol.* (2026).

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, 06/25/2026

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