

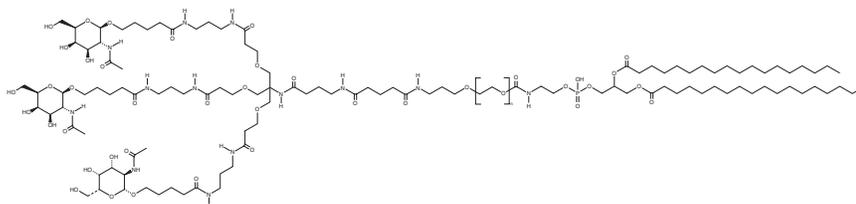
PRODUCT INFORMATION



DSPE-PEG(2000)-tri-GalNAc

Item No. 44551

Formal Name: 3-(((39-(((2R,3R,4R,5R,6R)-3-acetamido-4,5-dihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)-24,24-bis((3-((5-(((2R,3R,4R,5R,6R)-3-acetamido-4,5-dihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)pentanamido)propyl)amino)-3-oxopropoxy)methyl)-4,13,17,22,29,35-hexaoxo-5,8,26-trioxa-3,12,18,23,30,34-hexaazanonatriacontyl)oxy)(hydroxy)phosphoryl)oxy)propane-1,2-diyl distearate



MF: $(C_{2}H_{4}O)_{n}C_{115}H_{210}N_{13}O_{40}P$

Supplied as: A semi-solid

Storage: -20°C

Stability: ≥4 years

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

Laboratory Procedures

DSPE-PEG(2000)-tri-GalNAc is supplied as a solid. A stock solution may be made by dissolving the DSPE-PEG(2000)-tri-GalNAc in the solvent of choice, which should be purged with an inert gas. DSPE-PEG(2000)-tri-GalNAc is soluble in ethanol, chloroform, and DMSO. It is also soluble in water.

Description

DSPE-PEG(2000)-tri-GalNAc is a multivalent N-acetylgalactosamine-conjugated lipid composed of an O-deacetylated derivative of LNP GalNAc ligand 1 (Item No. 38952), a GABA linker, a PEG(2000) spacer, and 1,2-distearoyl-*rac*-glycero-3-phosphoethanolamine. Multivalent GalNAc-conjugated lipids have been used in the generation of lipid nanoparticles (LNPs) for targeted liver delivery *via* the asialoglycoprotein receptor (ASGPR).¹

Reference

1. Kasiewicz, L.N., Biswas, S., Beach, A., *et al.* GalNAc-Lipid nanoparticles enable non-LDLR dependent hepatic delivery of a CRISPR base editing therapy. *Nat. Commun.* **14**(1), 2776 (2023).

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, 10/13/2025

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