

PRODUCT INFORMATION



DMG-PEG(2000)

Item No. 33945

CAS Registry No.: 160743-62-4
Formal Name: α -[(2R)-2,3-bis[(1-oxotetradecyl)oxy]propyl]- ω -methoxy-poly(oxy-1,2-ethanediyl)

Synonyms: 1,2-Dimyristoyl-*rac*-glycero-3-methoxypolyethylene glycol-2000

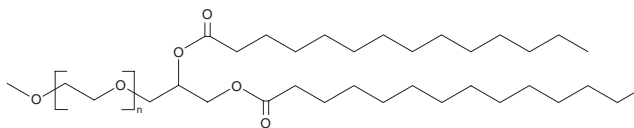
MF: $(C_2H_4O)_n C_{32}H_{62}O_5$

Purity: $\geq 90\%$

Supplied as: A crystalline 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

DMG-PEG(2000) is supplied as a crystalline solid. A stock solution may be made by dissolving the DMG-PEG(2000) in the solvent of choice, which should be purged with an inert gas. DMG-PEG(2000) is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of DMG-PEG(2000) in these solvents is approximately 1 mg/ml. DMG-PEG(2000) is also slightly soluble in DMSO.

Description

DMG-PEG(2000) is a lipid excipient that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).¹ Administration of fumarylacetoacetate hydrolase (FAH) mRNA in DMG-PEG(2000)-containing LNPs increases survival in a mouse model of hereditary tyrosinemia type 1 (HT-1). Formulations containing DMG-PEG(2000) have been used in the development of LNPs for the delivery of mRNA-based vaccines.

Reference

1. Cheng, Q., Wei, T., Jia, Y., *et al.* Dendrimer-based lipid nanoparticles deliver therapeutic FAH mRNA to normalize liver function and extend survival in a mouse model of hepatorenal tyrosinemia type I. *Adv. Mater.* **30(52)**, e1805308 (2018).

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, 10/13/2022

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