

PRODUCT INFORMATION



DMG-PEG(2000) (solution)

Item No. 42746

CAS Registry No.: 160743-62-4

Formal Name: α -[2,3-bis[(1-oxotetradecyl)oxy]propyl]- ω -methoxy-poly(oxy-1,2-ethanediyl)

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

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

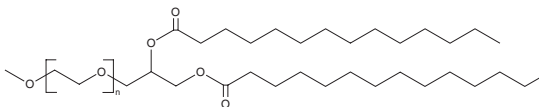
Purity: $\geq 90\%$

Supplied as: A 10 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Product may crash out of solution when cold. To restore the solution, gently warm and agitate the vial.



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

Description

DMG-PEG(2000) is a lipid excipient that has been used in combination with other lipids in the formation of lipid nanoparticles.¹ Administration of fumarylacetoacetate hydrolase (FAH) mRNA in DMG-PEG(2000)-containing lipid nanoparticles 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 lipid nanoparticles 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, 07/13/2026

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