

PRODUCT INFORMATION

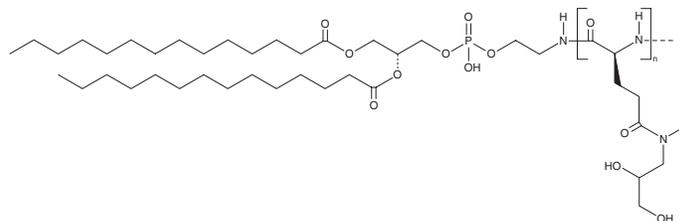


DMPE-PGA-diol40

Item No. 44172

Formal Name: (2R)-3-(((2-((2S)-5-((2,3-dihydroxypropyl)amino)-2-(methylamino)-5-oxopentanamido)ethoxy)(hydroxy)phosphoryl)oxy)propane-1,2-diyl ditetradecanoate

Synonyms: 1,2-Ditetradecanoyl-*sn*-glycero-3-Phosphoethanolamine-PGA-diol40, 1,2-Dimyristoyl-*sn*-glycero-3-Phosphoethanolamine-PGA-diol40, 1,2-Dimyristoyl-*sn*-glycero-3-Phosphoethanolamine-poly[L-glutamic acid-(3-aminopropane-1,2-diol)]₄₀, 1,2-DMPE-PGA-diol40, PE(14:0/14:0)-PGA-diol₄₀, PtdEtn-(1,2-Dimyristoyl)-PGA-diol40



MF: $(C_8H_{14}N_2O_4)_n C_{33}H_{65}NO_8P$

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

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

Laboratory Procedures

DMPE-PGA-diol40 is supplied as a solid. A stock solution may be made by dissolving the DMPE-PGA-diol40 in the solvent of choice, which should be purged with an inert gas. DMPE-PGA-diol40 is soluble in ethanol/water (80:20) at a concentration of approximately 25 mg/ml.

Description

DMPE-PGA-diol40 is a polymer and a poly[L-glutamic acid-3-aminopropane-1,2-diol]-conjugated form of 1,2-dimyristoyl-*sn*-glycero-3-PE (1,2-DMPE; Item No. 15090). It can be used as a shielding lipid in the generation of lipid nanoparticles (LNPs). This product was made using Curapath-patented technology.

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