

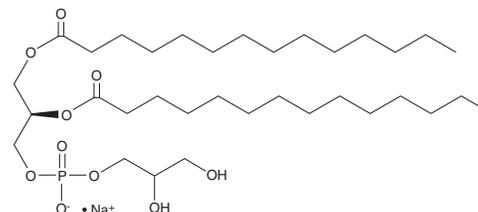
# PRODUCT INFORMATION



## 1,2-Dimyristoyl-*sn*-glycero-3-PG (sodium salt)

Item No. 15085

**CAS Registry No.:** 200880-40-6  
**Formal Name:** 1,2-dimyristoyl-*sn*-glycero-3-phosphoglycerol, monosodium salt  
**Synonyms:** 1,2-Dimyristoyl-*sn*-glycero-3-phospho-(1'-*rac*-glycerol), 1,2-DMPG  
**MF:** C<sub>34</sub>H<sub>66</sub>O<sub>10</sub>P • Na  
**FW:** 688.9  
**Purity:** ≥98%  
**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

1,2-Dimyristoyl-*sn*-glycero-3-PG (1,2-DMPG) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-DMPG (sodium salt) in the solvent of choice, which should be purged with an inert gas. 1,2-DMPG (sodium salt) is soluble in the organic solvent chloroform at a concentration of approximately 2 mg/ml.

### Description

1,2-DMPG is a phospholipid containing the saturated long-chain (14:0) myristic acid inserted at the *sn*-1 and *sn*-2 positions. It can be used in the generation of micelles, liposomes, and other types of artificial membranes.<sup>1-2</sup>

### References

1. Van Dijck, P.W., Ververgaert, P.H., Verkleji, A.J., *et al.* Influence of Ca<sup>2+</sup> and Mg<sup>2+</sup> on the thermotropic behaviour and permeability properties of liposomes prepared from dimyristoyl phosphatidylglycerol and mixtures of dimyristoyl phosphatidylglycerol and dimyristoyl phosphatidylcholine. *Biochim. Biophys. Acta* **406(4)**, 465-478 (1975).
2. Rodrigues, C., Gameiro, P., Reis, S., *et al.* Derivative spectrophotometry as a tool for the determination of drug partition coefficients in water/dimyristoyl-L- $\alpha$ -phosphatidylglycerol (DMPG) liposomes. *Biophys. Chem.* **94(1-2)**, 97-106 (2001).

#### 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, 11/11/2022

#### 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