

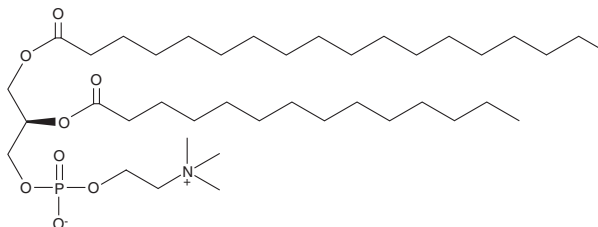
# PRODUCT INFORMATION



## 1-Stearoyl-2-Myristoyl-*sn*-glycero-3-PC

Item No. 15103

**CAS Registry No.:** 20664-02-2  
**Formal Name:** 1-stearoyl-2-myristoyl-*sn*-glycero-3-phosphatidylcholine  
**Synonyms:** 1,2-SMPC, 1-Stearoyl-2-Myristoyl-*sn*-glycero-3-Phosphocholine  
**MF:** C<sub>40</sub>H<sub>80</sub>NO<sub>8</sub>P  
**FW:** 734.1  
**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-Stearoyl-2-myristoyl-*sn*-glycero-3-PC (1,2-SMPC) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-SMPC in the solvent of choice, which should be purged with an inert gas. 1,2-SMPC is soluble in the organic solvent ethanol at a concentration of approximately 25 mg/ml.

### Description

1,2-SMPC belongs to the class of asymmetric phospholipids, which have two chains of unequal length attached to the glycerol backbone. 1,2-SMPC is abundant in biological membranes where it can interdigitate with equal length (symmetric) acyl chain phospholipids in bilayer membranes to potentially minimize the hydrophobic mismatch between acyl chains.<sup>1</sup>

### Reference

1. Høyrup, P., Mouritsen, O.G., and Jørgensen, K. Phospholipase A<sub>2</sub> activity towards vesicles of DPPC and DMPC-DSPC containing small amounts of SMPC. *Biochim. Biophys. Acta* **1515(2)**, 133-143 (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, 11/01/2022

### 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