

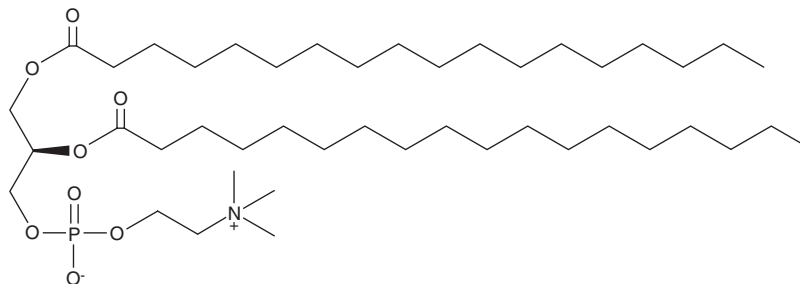
PRODUCT INFORMATION



1,2-Distearoyl-*sn*-glycero-3-PC

Item No. 15100

CAS Registry No.: 816-94-4
Formal Name: 1,2-distearoyl-*sn*-glycero-3-phosphatidylcholine
Synonyms: Coatsome MC 8080, 1,2-Distearoyl-*sn*-glycero-3-Phosphocholine, 1,2-DSPC
MF: C₄₄H₈₈NO₈P
FW: 790.2
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-Distearoyl-*sn*-glycero-3-PC (DSPC) is supplied as a crystalline solid. A stock solution may be made by dissolving the DSPC in the solvent of choice, which should be purged with an inert gas. DSPC is soluble in the organic solvent ethanol at a concentration of approximately 25 mg/ml.

Description

DSPC is a phospholipid containing the saturated long-chain (18:0) stearic acid inserted at the *sn*-1 and *sn*-2 positions. It is commonly used in the generation of micelles, liposomes, and other types of artificial membranes.^{1,2}

References

1. Hoyrup, P., Mouritsen, O.G., and Jorgensen, K. Phospholipase A₂ activity towards vesicles of DPPC and DMPC-DSPC containing small amounts of SMPC. *Biochim. Biophys. Acta.* **1515(2)**, 133-143 (2001).
2. Zhao, H., Joseph, J., Zhang, H., *et al.* Synthesis and biochemical applications of a solid cyclic nitron spin trap: A relatively superior trap for detecting superoxide anions and glutathyl radicals. *Free Radic. Biol. Med.* **31(5)**, 599-606 (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, 11/08/2022

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