

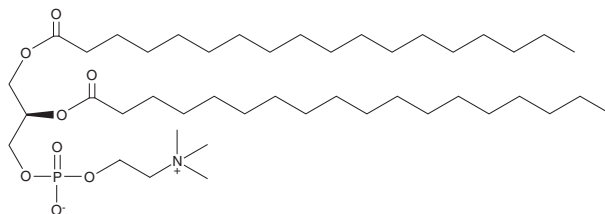
PRODUCT INFORMATION



1,2-Distearoyl-*sn*-glycero-3-PC (solution)

Item No. 39189

CAS Registry No.: 816-94-4
Formal Name: 1,2-distearoyl-*sn*-glycero-3-phosphatidylcholine
Synonyms: Coatsome MC 8080,
PC(18:0/18:0),
1,2-Distearoyl-*sn*-glycero-3-Phosphocholine,
1,2-DSPC,
18:0/18:0-PC
MF: C₄₄H₈₈NO₈P
FW: 790.2
Purity: ≥98%
Supplied as: A 25 mM solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

1,2-Distearoyl-*sn*-glycero-3-PC is a phospholipid containing stearic acid (Item No. 10011298) at the *sn*-1 and *sn*-2 positions. It has commonly been used in the generation of micelles, liposomes, and other types of artificial membranes.^{1,2}

References

1. Høyrup, P., Mouritsen, O.G., and Jørgensen, 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. Heberle, F.A. and Feigenson, G.W. Phase separation in lipid membranes. *Cold Spring Harb. Perspect. Biol.* **3**(4), 1-13 (2011).

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, 01/14/2025

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