

PRODUCT INFORMATION



Cardiolipin (18:0/18:0/18:0/18:0) (sodium salt)

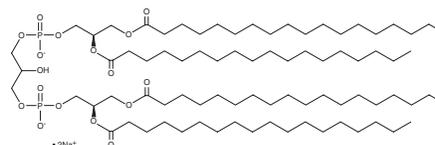
Item No. 39283

CAS Registry No.: 64069-64-3
Formal Name: octadecanoic acid, [R-(R*,R*)]-5,8,11-trihydroxy-5,11-dioxido-4,6,10,12-tetraoxa-5,11-diphosphapentadecane-1,2,14,15-tetrayl ester, disodium salt

Synonyms: 18:0 Cardiolipin, 72:0-CL, 4S-sn-CL, Diphosphatidylglycerol (18:0), Tetrastearoyl Cardiolipin

MF: C₈₁H₁₅₆O₁₇P₂ • 2Na
FW: 1,510.0

Purity: ≥98%
Supplied as: A 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

Cardiolipin (18:0/18:0/18:0/18:0) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the cardiolipin (18:0/18:0/18:0/18:0) (sodium salt) in the solvent of choice, which should be purged with an inert gas. Cardiolipin (18:0/18:0/18:0/18:0) (sodium salt) is soluble in methanol.

Description

Cardiolipin (18:0/18:0/18:0/18:0) is an anionic phospholipid containing the long-chain saturated fatty acid stearic acid (Item No. 10011298). It has been used to promote hydroxyapatite crystallization and supersaturation in a calcium solution.¹

Reference

1. Dalas, E., Ioannou, P.V., and Koutsoukos, P.G. Effect of fatty acyl and cation content of cardiolipins on in vitro calcification. *Langmuir* **5(1)**, 157-160 (1988).

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, 02/08/2024

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