

PRODUCT INFORMATION



Cardiolipin (16:0/18:1/16:0/18:1) (sodium salt)

Item No. 39282

CAS Registry No.: 2260669-35-8

Formal Name: 9-octadecenoic acid (9Z)-1,1'-
[[1R,13R)-4,7,10-trihydroxy-4,10-
dioxido-1,13-bis[[[(1-oxohexadecyl)
oxy]methyl]-3,5,9,11-tetraoxa-4,10-
diphosphatridecane-1,13-diyl] ester,
disodium salt

Synonyms: 16:0-18:1 Cardiolipin, 68:2-CL,
CL(1'-[16:0/18:1],3'-[16:0/18:1]),
Diphosphatidylglycerol (16:0/18:1)

MF: $C_{77}H_{144}O_{17}P_2 \cdot 2Na$

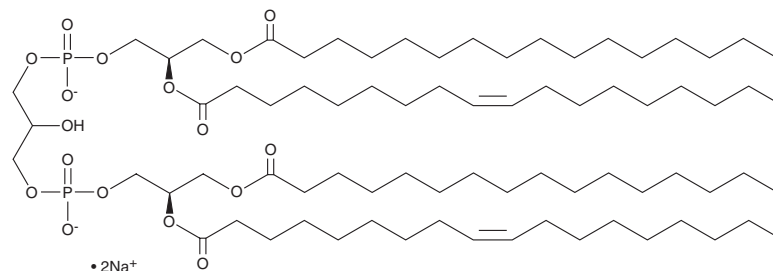
FW: 1,449.9

Purity: $\geq 95\%$

Supplied as: A 10 mg/ml solution in chloroform

Storage: $-20^{\circ}C$

Stability: ≥ 2 years



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

Description

Cardiolipin (16:0/18:1/16:0/18:1) is an anionic phospholipid containing the long-chain fatty acid palmitic acid (Item No. 10006627) and the monounsaturated fatty acid oleic acid (Item Nos. 90260 | 24659). It has been used in the synthesis of lipid nanodiscs for use in native mass spectrometry applications.¹

Reference

1. Kostelic, M.M., Zak, C.K., Jayasekera, H.S., *et al.* Assembly of model membrane nanodiscs for native mass spectrometry. *Anal. Chem.* **93**(14), 5972-5979 (2021).

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, 03/25/2026'

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