

PRODUCT INFORMATION



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

Item No. 39281

CAS Registry No.: 2260669-34-7

Formal Name: 9-hexadecenoic acid, 1,1',1'',1'''-[(2R,14R)-5,8,11-trihydroxy-5,11-dioxido-4,6,10,12-tetraoxa-5,11-diophosphapentadecane-1,2,14,15-tetrayl] ester, disodium salt

Synonyms: 16:1 Cardiolipin, 64:4-CL, Diphosphatidylglycerol (16:1), Tetrapalmitoleoyl Cardiolipin

MF: $C_{73}H_{132}O_{17}P_2 \cdot 2Na$

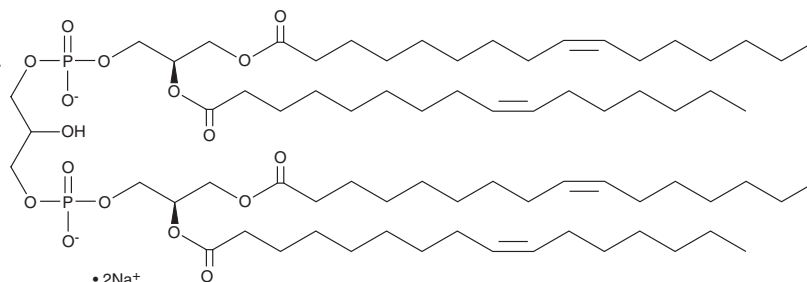
FW: 1,389.8

Purity: $\geq 95\%$

Supplied as: A 5 mg/ml solution in chloroform

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

Stability: ≥ 4 years



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

Description

Cardiolipin (16:1/16:1/16:1/16:1) is an anionic phospholipid containing the long-chain monounsaturated fatty acid (MUFA) palmitoleic acid (Item Nos. 10009871 | 21911). It increases the activation of, and cytokine production by, DN32.D3 type 1 natural killer T (NKT) cells when incubated with L-M(TK-) mouse fibroblasts expressing the lipid antigen-presenting molecule CD1d (L-CD1d cells) prior to their co-culture with the DN32.D3 NKT cells.¹

Reference

1. Lee, S., Sun, W., and Webb, T.J. Sphingosine kinase blockade leads to increased natural killer T cell responses to mantle cell lymphoma. *Cells* **9**(4), 1030 (2020).

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, 05/08/2026

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