

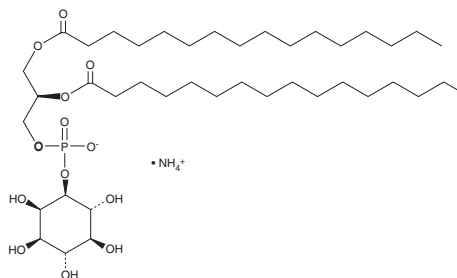
PRODUCT INFORMATION



PtdIns-(1,2-dipalmitoyl) (ammonium salt)

Item No. 10007710

CAS Registry No.: 34290-57-8
Formal Name: 1-(1,2R-dihexadecanoylphosphatidyl) inositol, monoammonium salt
Synonyms: DPPI, Phosphatidylinositol C-16, PI(16:0/16:0), 16:0/16:0 PI
MF: $C_{41}H_{78}O_{13}P \cdot NH_4$
FW: 828.1
Purity: $\geq 98\%$
Supplied as: A lyophilized powder
Storage: $-20^\circ C$
Stability: ≥ 5 years



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

Laboratory Procedures

PtdIns-(1,2-dipalmitoyl) (ammonium salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the PtdIns-(1,2-dipalmitoyl) (ammonium salt) in an organic solvent such as chloroform:methanol:water (4:3:1). The solubility of PtdIns-(1,2-dipalmitoyl) (ammonium salt) in this solvent is at least 1 mg/ml.

PtdIns-(1,2-dipalmitoyl) (ammonium salt) is not soluble in aqueous buffers, but can be dispersed in aqueous media by sonication or cosonication with detergents.

Description

PtdIns-(1,2-dipalmitoyl) is a derivative of phosphatidylinositol (PtdIns) that contains C16:0 fatty acyl chains at the *sn*-1 and *sn*-2 positions.

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, 07/11/2023

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