

PRODUCT INFORMATION



PtdIns-(1-palmitoyl-d₉, 2-palmitoyl) (sodium salt)

Item No. 28156

Formal Name: (R)-3-((hexadecanoyl-13,13,14,14,15,15,16,16,16-d₉)oxy)-2-(palmitoyloxy)propyl ((1S,2R,3R,4S,5S,6R)-2,3,4,5,6-pentahydroxycyclohexyl) phosphate, monosodium salt

Synonyms: DPPI-d₉, PI(16:0-d₉/16:0), 16:0-d₉/16:0 PI

MF: C₄₁H₆₉D₉O₁₃P • Na

FW: 842.1

Chemical Purity: ≥95% (PtdIns-(1,2-dipalmitoyl))

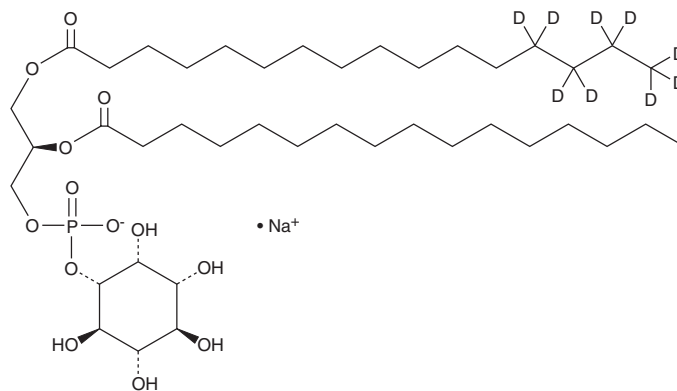
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

PtdIns-(1-palmitoyl-d₉, 2-palmitoyl) (sodium salt) is intended for use as an internal standard for the quantification of PtdIns-(1,2-dipalmitoyl) (Item No. 10007710) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

PtdIns-(1-palmitoyl-d₉, 2-palmitoyl) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the PtdIns-(1-palmitoyl-d₉, 2-palmitoyl) (sodium salt) in the solvent of choice, which should be purged with an inert gas. PtdIns-(1-palmitoyl-d₉, 2-palmitoyl) (sodium salt) is soluble in methanol. The solubility of PtdIns-(1-palmitoyl-d₉, 2-palmitoyl) (sodium salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PtdIns-(1-palmitoyl-d₉, 2-palmitoyl) is intended for use as an internal standard for the quantification of PtdIns-(1,2-dipalmitoyl) (Item No. 10007710) by GC- or LC-MS. 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, 02/27/2023

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