

L- α - Phosphatidylinositol- 3,4,5-P3 . 7Na

PKC substrate

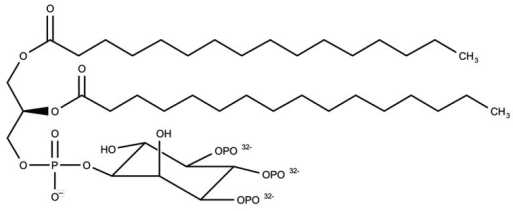
Activates Ca²⁺-insensitive PKC isotypes δ , ϵ and η . Binds to GRP1 *via* a PH domain.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PtdIns(3,4,5)P3, (1,2-dipalmitoyl)-, 1,2-Dipalmitoyl-sn-glycero-3-phospho-1-D-myo-inositol-3,4,5-trisphosphate . 7Na
Couple Target	PKC
Couple Type	Substrate
Formula	$C_{41}H_{75}O_{22}P_4 \cdot 7Na$
Formulation	Lyophilized.
MW	1204.8
Purity	≥98% (TLC)
Solubility	Soluble in water (10mg/ml) or in chloroform/methanol/water (1:1:0.3) (>10mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com