

L- α - Phosphatidylinositol- 3,4-P2

Activator of PKC and Akt

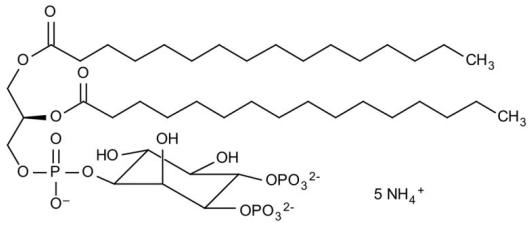
Activates Ca²⁺-insensitive PKC isotypes δ , ϵ and η . Activates Akt/PKB by direct interaction with the Akt PH domain.

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Hygroscopic.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
Couple Target	Akt, PKC
Couple Type	Activator
Formula	$C_{41}H_{76}O_{19}P_3 \cdot 5NH_4$
MW	966.0 . 90.2
Purity	≥98% (NMR)
Solubility	Soluble in chloroform/methanol/water, 50:50:10 (10mg/ml); not water soluble but resuspends, with vigorous vortexing, to ~0.5 mg/ml in water. See references for method of incorporation into mixed lipid vesicles by sonication, and for dissolving in a detergent-containing buffer.

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-be@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