

D-myo-Inositol 4,5-bisphosphate, L- α -Phosphatidyl-(1,2-dipalmitoyl)-

The presence of 4,5-bisphosphate group on the inositol ring has been found critical for binding specifically to PH domains that are commonly present in signalling proteins, also found in cytoskeleton.

Citations: 6

[View Online »](#)

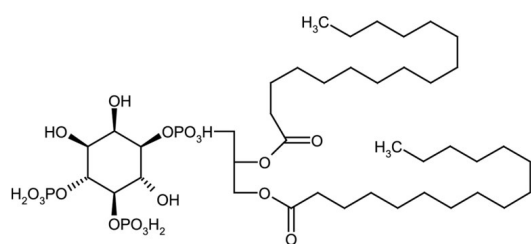
Ordering Information

[Order Online »](#)

ALX-307-039-C100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture. Hygroscopic.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PtdIns(4,5)P ₂ , (1,2-dipalmitoyl)-, L-α-Phosphatidyl-(1,2-dipalmitoyl)-D-myo-inositol 4,5-bisphosphate
Appearance	White to off-white powder.
Formula	C ₄₁ H ₈₁ O ₁₉ P ₃
Identity	Identity determined by 1H-NMR and 31P-NMR and MS.
MW	971.0
Purity	≥98% (HPLC)
Solubility	Soluble in water, DMSO or methanol.



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