

Phosphatidylethanol

Ultra-pure

In the presence of ethanol, phospholipase D converts phosphatidylcholine to phosphatidylethanol. This conversion is the basis of a sensitive and selective assay for PLD activity in intact cells. Phosphatidylethanol is a useful chromatographic standard for this assay.

Citations: 4

[View Online »](#)

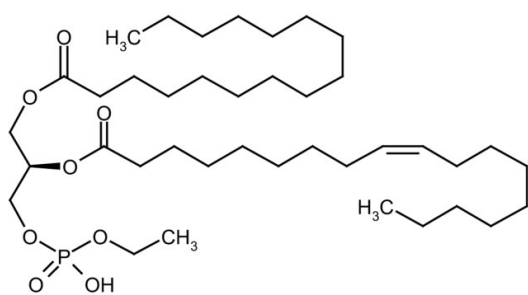
Ordering Information

[Order Online »](#)

BML-ST400-0010	10mg
BML-ST400-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Protect from oxygen. Store under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-Palmitoyl-2-oleoyl-sn-glycero-3-phosphoethanol, PEth-16:0/18:1
Appearance	Viscous oil or waxy solid.
Formula	$C_{39}H_{75}O_8P$
MW	703
Purity	≥98% (TLC)
Solubility	Soluble in chloroform.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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