

Oleyloxyethyl phosphorylcholine

Phospholipase inhibitor

A potent and specific substrate analog PLA_2 inhibitor ($\text{IC}_{50}=6.2\text{ }\mu\text{M}$ -porcine pancreatic PLA_2). It inhibits human recombinant sPLA_2 ($\text{IC}_{50}=13.7\text{ }\mu\text{M}$) selectively over cPLA_2 . At $1\text{ }\mu\text{M}$ it completely blocked the myogenic response in perfused dog renal arcuate arteries.

Citations: 4

[View Online »](#)

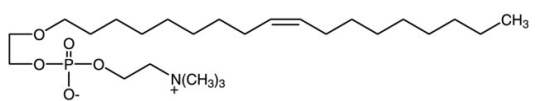
Ordering Information

[Order Online »](#)

BML-EI118-0010	10mg
BML-EI118-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Colorless oil or waxy solid.

CAS 96720-06-8

Couple Target Phospholipase

Couple Type Inhibitor

Formula $C_{25}H_{52}NO_5P$

MW 477.7

Purity ≥98% (HPLC)

Solubility Soluble in 100% ethanol.

Source Synthetic.

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)