

Thioetheramide-phosphorylcholine

Phospholipase inhibitor

A potent substrate analog inhibitor of snake venom PLA₂ (IC₅₀=2 µM). It inhibits human recombinant sPLA₂ (IC₅₀=0.9 µM) selectively over cPLA₂.

Citations: 2

[View Online »](#)

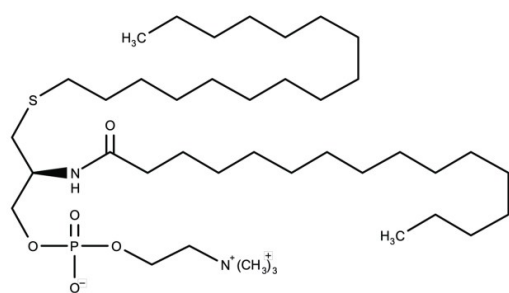
Ordering Information

[Order Online »](#)

BML-ST325-0005	5mg
BML-ST325-0025	25mg

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.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-Palmitylthio-2-palmitoylamido-1,2-dideoxy-sn-glycero-3-phosphorylcholine
Appearance	White solid.
CAS	116457-99-9
Couple Target	Phospholipase, PLA
Couple Type	Inhibitor
Formula	$C_{40}H_{83}N_2O_5PS$
MW	735.6
Purity	≥98% (TLC)
Solubility	Soluble in DMSO, 100% ethanol or methanol.

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-
fr@enzolifesciences.com](mailto:info-fr@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)