

PI 3K-SH2-OMT

PI3 kinase inhibitor

This unique SH2 domain ligand is derived from the sequence flanking Tyr⁷⁵¹ of the PDGF receptor. It inhibits the PI 3-kinase p85 C-terminal SH2 domain (IC_{50} =14.2 μ M). It contains the nonhydrolyzable phosphotyrosine mimetic L-O-(2-malonyl)tyrosine (OMT) and is provided as the trimethyl ester for greater cell permeability.

Ordering Information

[Order Online »](#)

BML-ST425-0500	0.5mg
-----------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow solid.
Formulation	Lyophilized.
MW	922.1
Purity	≥98% (TLC)
Sequence	Ac-Asp-Tyr(2-malonyl)-Val-Pro-Met-Leu-NH ₂ trimethyl ester
Solubility	Soluble in DMSO (0.5 mg/ml) or ethanol (0.3 mg/ml).



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)