

Wortmannin

PI3 kinase inhibitor

Potent and selective inhibitor of phosphatidyl-inositol 3-kinase. Active in purified preparations and cytosolic fractions ($IC_{50} = 5nM$) and is highly cell permeable. Inhibits fMLP induced PIP_3 and superoxide anion production ($IC_{50} = 50nM$) in guinea pig neutrophils and blocks the metabolic effects of insulin in isolated rat adipocytes without affecting the insulin receptor tyrosine kinase activity. Covalently binds to PI 3-kinase and is selective, inhibiting other kinases such as PI 4-kinase and myosin light chain kinase at concentrations 100-fold higher than that required for inhibition of PI 3-kinase. The half-life of wortmannin in culture media is quite short (8-13 minutes), care should be taken to account for this decomposition when planning experiments. Markedly potentiates the LPS-induced nitric oxide (NO) production from macrophages. Induces *in vivo* Alzheimer-like hyperphosphorylation in tau.

Citations: 73

[View Online »](#)

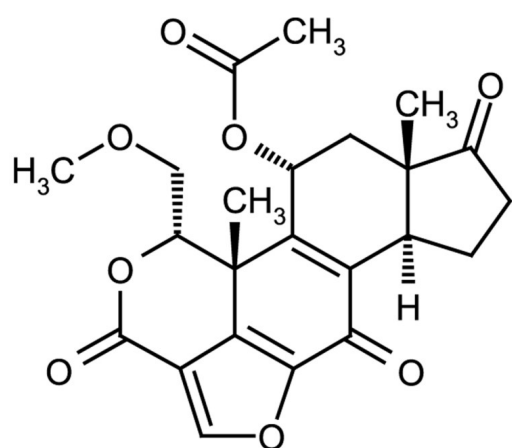
Ordering Information

[Order Online »](#)

BML-ST415-0001	1mg
BML-ST415-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Make solutions fresh immediately before use. DO NOT STORE SOLUTIONS.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	KY 12420
Appearance	White to off-white solid.
CAS	19545-26-7
Couple Target	PI3 kinase
Couple Type	Inhibitor
Formula	$C_{23}H_{24}O_8$
MI	14: 10053
MW	428.4
Purity	≥98% (HPLC, TLC)
RTECS	CB9641000
Solubility	Soluble in DMSO (50mg/ml), methanol (5mg/ml) or 100% ethanol (25mg/ml).
Source	Isolated from <i>Penicillium wortmannii</i>

Technical Info / Product Notes Replacement for **ADI-HPK-114**.



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)