

# 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 »](#)

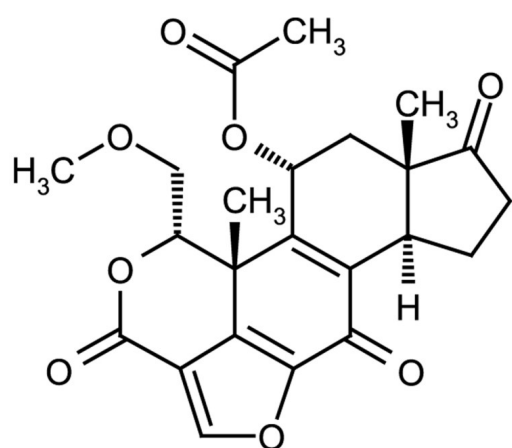
## Ordering Information

[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	<b>14:</b> 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**.



Last modified: May 29, 2024

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