

# PP1

## Src kinase inhibitor

A highly potent and selective inhibitor of Src family tyrosine kinases,  $IC_{50}$  =5 nM for p56<sup>lck</sup>; p59<sup>fyn</sup>T=6 nM; p60<sup>src</sup>=170 nM. Inhibition of Src family kinases by PP1 has been shown to: prevent  $\alpha$ -thrombin-stimulated Akt phosphorylation, delay PAF-mediated STAT3 phosphorylation, block  $Zn^{2+}$ -induced Ras activation, block adhesion-dependent tyrosine phosphorylation of p105<sup>HEF1</sup> <sup>5</sup>, and inhibit HLA-DR-induced tyrosine phosphorylation without affecting the activation of ERK1/2.

Citations: 23

[View Online »](#)

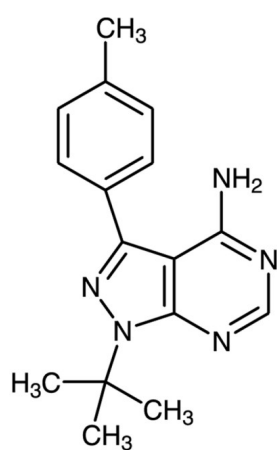
## Ordering Information

[Order Online »](#)

BML-EI275-0001	1mg
BML-EI275-0005	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	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Tan solid.
CAS	172889-26-8
Couple Target	Src kinase
Couple Type	Inhibitor
Formula	$C_{16}H_{19}N_5$
MW	281.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml), slightly soluble in 100% ethanol.



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)