PP1

Src kinase inhibitor

A highly potent and selective inhibitor of Src family tyrosine kinases, IC $_{50}$ =5 nM for p56 $^{\rm lck}$; p59 $^{\rm fynT}$ =6 nM; p60 $^{\rm src}$ =170 nM. Inhibition of Src family kinases by PP1 has been shown to: prevent a-thrombin-stimulated Akt phosphorylation, delay PAF-mediated STAT3 phosphorylation, block Zn²+-induced Ras activation, block adhesion-dependent tyrosine phosphorylation of p105 $^{\rm HEF1~5}$, and inhibit HLA-DR-induced tyrosine phosphorylation without affecting the activation of ERK1/2.

Citations: 23

View Online »

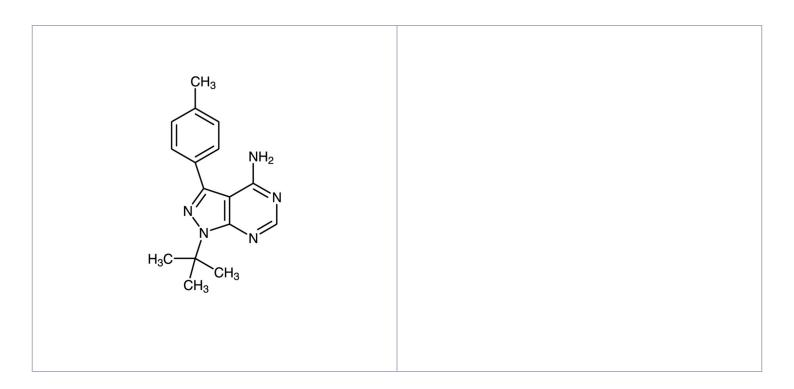
Ordering Information

Order Online »

BML-El275-0001	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)

Soluble in DMSO (25mg/ml), slightly soluble in 100%

ethanol.

Last modified: May 29, 2024

