

PP2

Src kinase inhibitor

A potent and selective inhibitor of the Src family tyrosine kinases. Inhibits Lck ($IC_{50}=4$ nM), Fyn ($IC_{50}=5$ nM), and Hck ($IC_{50}=5$ nM). It displays weaker inhibition of EGFR ($IC_{50}=480$ nM) and no inhibition of ZAP-70 (IC_{50} 100 μ M) and JAK2 (IC_{50} 50 μ M). PP2 is a useful tool for probing the involvement of Src family tyrosine kinases in cellular signaling.

Citations: 13

[View Online »](#)

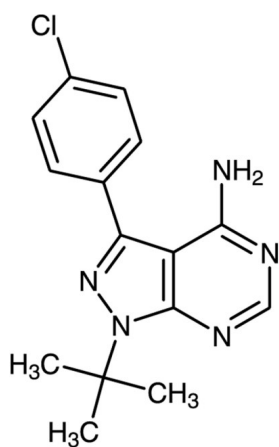
Ordering Information

[Order Online »](#)

BML-EI297-0001	1mg
BML-EI297-0005	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

Alternative Name	AG1879, 4-Amino-5-(4-chlorophenyl)-7-(t-butyl)pyrazolo[3,4-d]pyrimidine
Appearance	Off-white solid.
CAS	172889-27-9
Couple Target	Src kinase
Couple Type	Inhibitor
Formula	$C_{15}H_{16}ClN_5$
MW	301.8
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/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)