

Src I1

Inhibitor of Src kinase and p56lck tyrosine kinase

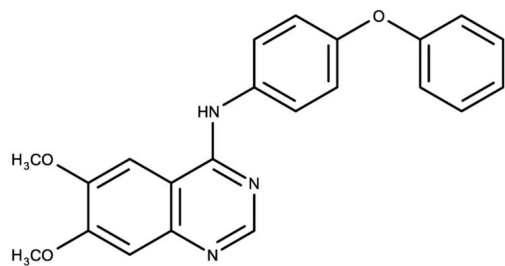
Potent, dual site (ATP domain and substrate-binding domain) Src and Lck kinase inhibitor (IC_{50} =44 and 88nM respectively). VEGFR2 IC_{50} =0.32 μ M; c-fms IC_{50} =30 μ M.

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	6,7-Dimethoxy-N-(4-phenoxyphenyl)-4-quinazolinamine
Appearance	Off-white solid.
CAS	179248-59-0
Couple Target	p56lck tyrosine kinase, Src kinase
Couple Type	Inhibitor
Formula	$C_{22}H_{19}N_3O_3$
MW	373.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (5mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com