

AG-1296

PDGF receptor tyrosine kinase inhibitor

A potent inhibitor of PDGF receptor tyrosine kinase ($IC_{50}=1\text{ }\mu\text{M}$). Also inhibits FGF receptor tyrosine kinase and c-kit. Induces apoptosis in a small-cell lung cancer cell line (H526).

Citations: 3

[View Online »](#)

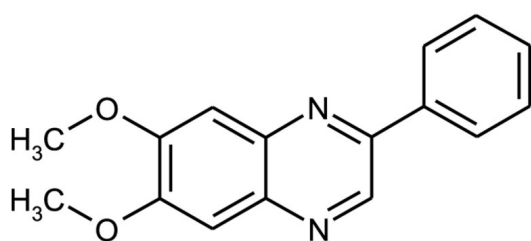
Ordering Information

[Order Online »](#)

BML-EI303-0005	5mg
BML-EI303-0025	25mg

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.
Handling	Store tightly sealed in the dark. Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tyrphostin AG 1296, 6,7-Dimethoxy-3-phenylquinoxaline
Appearance	Off-white to brownish solid.
CAS	146535-11-7
Couple Target	c-Kit, EGFR, PDGF receptor
Couple Type	Inhibitor
Formula	$C_{16}H_{14}N_2O_2$
MW	266.3
Purity	≥98% (TLC)
Solubility	Soluble in DMSO or dichloromethane.



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-
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)