

Tyrphostin 9

PDGF receptor inhibitor

Tyrphostin 9 is a selective inhibitor of the PDGF receptor tyrosine kinase ($IC_{50}=1.2\text{ }\mu\text{M}$). It is also a potent (10 nM) uncoupler of oxidative phosphorylation.

Citations: 10

[View Online »](#)

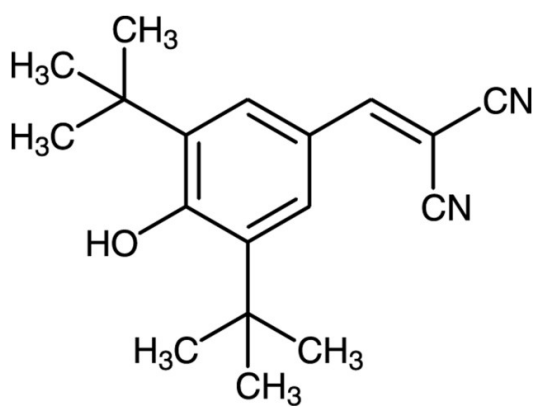
Ordering Information

[Order Online »](#)

BML-EI215-0050	50mg
BML-EI215-0250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Store tightly sealed in the dark. Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	RG-50872, Malonaben, SF 6847
Appearance	Yellow solid.
CAS	10537-47-0
Couple Target	PDGF receptor
Couple Type	Inhibitor
Formula	$C_{18}H_{22}N_2O$
MW	282.4
Purity	≥99% (HPLC)
RTECS	OO3737000
Solubility	Soluble in DMSO (>25mg/ml) or ethanol (20mg/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-fr@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