

Tyrphostin 47 (RG-50864) (AG-213)

EGFR inhibitor

A potent inhibitor of EGF receptor kinase activity (IC_{50} 's: EGFR = 2.4 μ M, PDGFR = 3.5 μ M, p210^{bcr-abl} = 5.9 μ M, insulin receptor = 640 μ M).

Citations: 5

[View Online »](#)

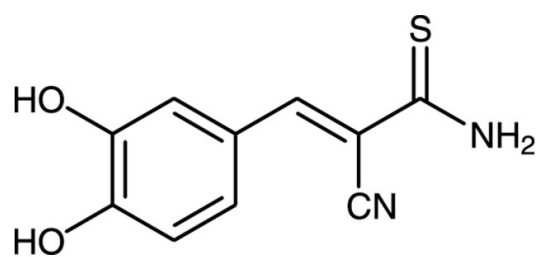
Ordering Information

[Order Online »](#)

BML-EI188-0005	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. Product is subject to oxidation. Make solutions fresh daily.
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

Appearance	Orange solid.
CAS	122520-86-9
Couple Target	EGFR
Couple Type	Inhibitor
Formula	$C_{10}H_8N_2O_2S$
MW	220.2
Purity	≥99% (TLC)
Solubility	Soluble at 20 mg/ml in 100% ethanol or DMSO.

Last modified: May 29, 2024



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)