

Tyrphostin 51

EGFR inhibitor

A potent inhibitor of EGF receptor kinase activity ($IC_{50}=0.8\text{ }\mu\text{M}$). Inhibition is mixed competitive with ATP and substrate.

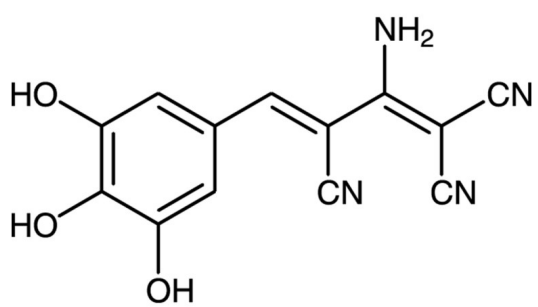
Ordering Information

[Order Online »](#)

BML-EI189-0005	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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Brown solid.
CAS	122520-90-5
Couple Target	EGFR
Couple Type	Inhibitor
Formula	$C_{13}H_8N_4O_3$
MW	268.2
Purity	≥99%
Solubility	Slightly soluble in water.



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