

SU11652

VEGFR inhibitor

Potent, ATP-competitive and cell permeable inhibitor of tyrosine kinases (VEGF-R2 IC_{50} =27nM, PDGF-R β IC_{50} =3nM). Also inhibits wild-type and all mutant forms of Kit in mast cells (IC_{50} 's=10 to 100nM) resulting in cell cycle arrest and apoptosis. Su11652 was also shown to induce autophagy in MCF-7 cells.

Citations: 3

[View Online »](#)

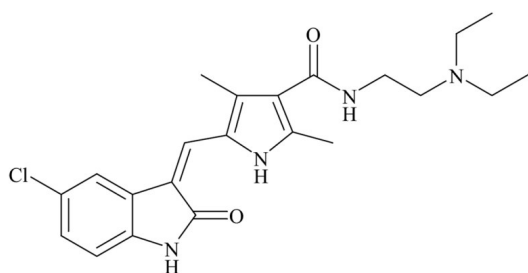
Ordering Information

[Order Online »](#)

BML-EI408-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months after receipt when stored at -20°C.
Handling	Solutions can be stored at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-[5-Chloro-2-oxo-1,2-dihydroindol-(3Z)-ylidenemethyl]-2,4-dimethyl-1H-pyrrole-3-carboxylic acid (2-diethylaminoethyl)amide
Appearance	Orange solid.
CAS	326914-10-7
Couple Target	VEGFR
Couple Type	Inhibitor
Formula	$C_{22}H_{27}ClN_4O_2$
MW	414.9
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>25 mg/ml).



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