

SU-5402

VEGFR2, FGFR1, and PDGFR β inhibitor

SU-5402 is a potent and selective inhibitor of multiple receptor tyrosine kinases, including vascular endothelial growth factor receptor 2 (VEGFR2), fibroblast growth factor receptor 1 (FGFR1), and platelet-derived growth factor receptor beta (PDGFRB). It has IC₅₀ values of 20 nM, 30 nM, and 510 nM for VEGFR2, FGFR1, and PDGFRB, respectively.

Key features and applications include:

- **Cancer Research:** SU-5402 is used to study the inhibition of angiogenesis and tumor growth by targeting VEGFR2, FGFR1, and PDGFRB pathways.
- **Stem Cell Research:** It is employed to maintain the pluripotency of embryonic stem cells and to study the differentiation processes.
- **Developmental Biology:** SU-5402 is utilized to investigate the roles of FGFR and VEGFR in embryonic development and tissue repair.

Relevant disease states include:

- **Cancer:** SU-5402 has shown antitumor properties and is studied for its potential to inhibit tumor growth in various cancers.
- **Neurodegenerative Diseases:** It is used in research to study its effects on neural stem cell differentiation, which can be relevant for conditions like Alzheimer's and Parkinson's diseases.
- **Developmental Disorders:** SU-5402 is studied for its role in inhibiting embryonic left-right determination, which can be relevant in understanding certain developmental disorders.

Ordering Information

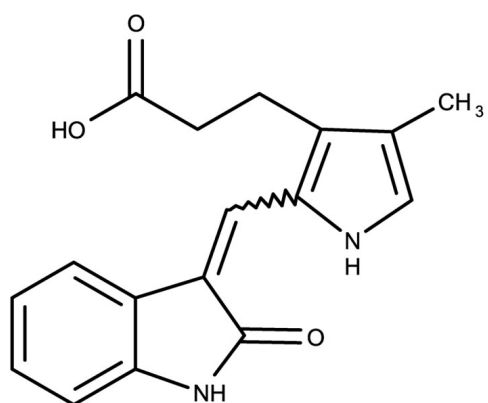
[Order Online »](#)

ENZ-CHM344-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20° for up to 3 months.
Handling	Provide appropriate exhaust ventilation at places where dust is formed. Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-[4-Methyl-2-(2-oxo-1,2-dihydro-indol-3-ylidenemethyl)-1H-pyrrol-3-yl]-propionic acid
Appearance	Orange solid.
CAS	215543-92-3
Couple Target	FGF receptor, PDGF receptor, VEGFR
Couple Type	Inhibitor
Formula	$C_{17}H_{16}N_2O_3$
Identity	Determined by NMR.
MW	296.32
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 10 mg/ml).



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)