

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

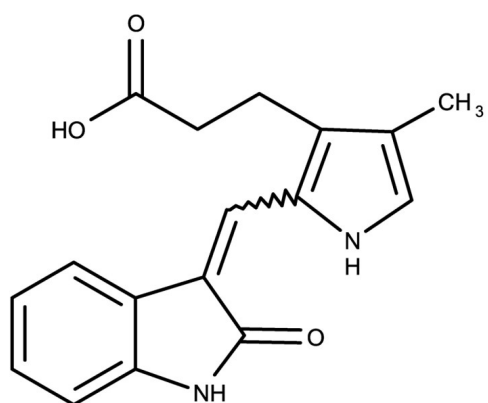
[Order Online »](#)

| | |
|-----------------|------|
| 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,
Last modified **March 27, 2025**
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