

MAZ51

VEGFR-3 inhibitor

Inhibits vascular endothelial cell growth factor-C (VEGF-C)-induced intracellular phosphorylation of VEGFR-3 (Flt-4) ($IC_{50}=1\mu M$).

Citations: 4

[View Online »](#)

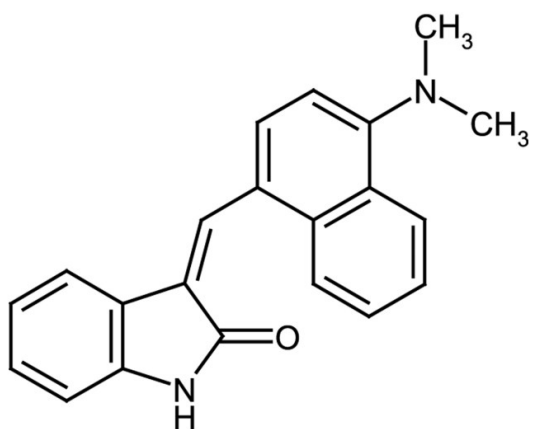
Ordering Information

[Order Online »](#)

ALX-270-342-M002	2mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-(4-Dimethylamino-naphthalen-1-yl-methylene)-1,3-dihydro-indol-2-one
Appearance	Orange crystals.
CAS	163655-37-6
Couple Target	VEGFR
Couple Type	Inhibitor
Formula	$C_{21}H_{18}N_2O$
Identity	Identity determined by MS, 1H- and 13C-NMR.
MW	314.4
Purity	≥98%
Solubility	Soluble in 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)