

ZM-323881

VEGFR inhibitor

Potent and selective inhibitor of the kinase activity of the human vascular endothelial growth factor receptor 2 (VEGFR2/KDR). Selectively inhibits VEGFR2 over VEGFR1 and other receptor tyrosine kinases with an IC_{50} = 2nM. Useful tool for exploring the biology of VEGFR receptors.

Citations: 3

[View Online »](#)

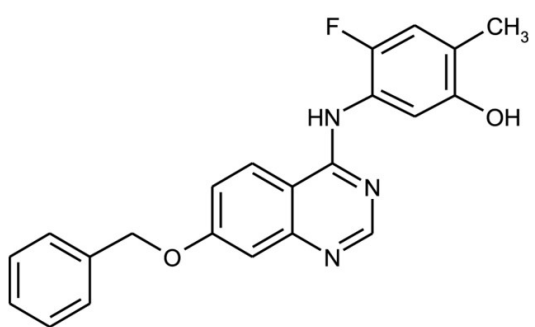
Ordering Information

[Order Online »](#)

BML-EI385-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 1 year as supplied. Stock solutions can be stored at -20°C for up to 3 months.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-((7-Benzyloxyquinazolin-4-yl)amino)-4-fluoro-2-methyl phenol
Appearance	White to off-white solid.
CAS	193001-14-8
Couple Target	VEGFR
Couple Type	Inhibitor
Formula	$C_{22}H_{18}FN_3O_2$
MW	375.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (3 mg/ml).
Source	Synthetic.

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)