

ZM-306416

VEGFR inhibitor

VEGF receptor tyrosine kinase inhibitor. Inhibits KDR ($IC_{50}=100\text{nM}$) and Flt ($IC_{50}=2\mu\text{M}$) tyrosine kinases.

Citations: 1

[View Online »](#)

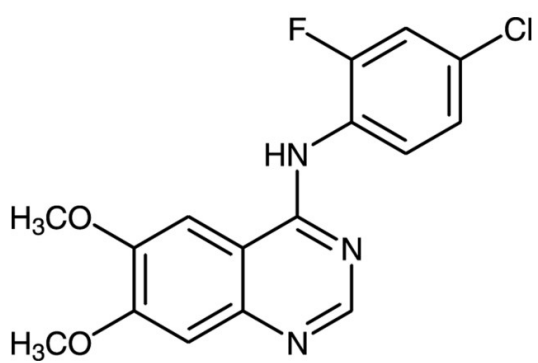
Ordering Information

[Order Online »](#)

BML-EI353-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white solid.
CAS	690206-97-4
Couple Target	VEGFR
Couple Type	Inhibitor
Formula	$C_{16}H_{13}ClFN_3O_2$
MW	333.7
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (3 mg/ml) or ethanol (1 mg/ml).

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-
be@enzolifesciences.com](mailto:info-be@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)