

Cabozantinib

Potent inhibitor of VEGFR-2, MET, Kit and FLT3-ITD

Cabozantinib is a very potent inhibitor, with an IC_{50} of less than 10 nM for VEGFR2, MET, Kit and FLT3-ITD.

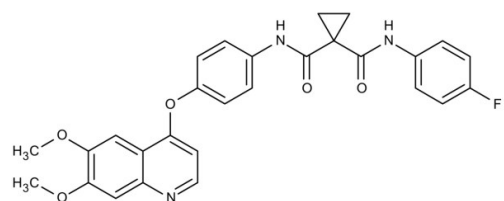
Ordering Information

[Order Online »](#)

ENZ-CHM182-0010	10mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	XL184, BMS-907351, N-(4-((6,7-Dimethoxyquinolin-4-yl)oxy)phenyl)-N-(4-fluorophenyl)cyclopropane-1,1-dicarboxamide
Appearance	Off-white to white powder.
CAS	849217-68-1
Couple Target	c-Kit, Flt, MET, VEGFR
Couple Type	Inhibitor
Formula	$C_{28}H_{24}FN_3O_5$
Identity	Determined by NMR.
MW	501.5
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (100mg/ml) or ethanol (1mg/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-
be@enzolifesciences.com](mailto:info-be@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)