

Motesanib

Cell-permeable inhibitor of VEGF receptors, PDGFR, and Kit receptor.

Motesanib potently inhibits angiogenesis and induces regression in tumor xenografts.

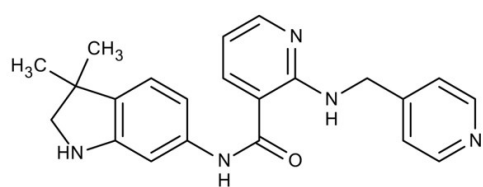
Ordering Information

[Order Online »](#)

ENZ-CHM147-0001	1mg
ENZ-CHM147-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AMG-706
Appearance	White powder solid.
CAS	453562-69-1
Couple Target	c-Kit, PDGF receptor, VEGFR
Couple Type	Inhibitor
Formula	$C_{22}H_{23}N_5O$
MW	373.5
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO or ethanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
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-fr@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