

Imatinib . mesylate

Tyrosine kinase inhibitor

Tyrosine kinase inhibitor. Highly specific for BCR-ABL, the enzyme associated with chronic myelogenous leukemia (CML) and certain forms of acute lymphoblastic leukemia (ALL). Also shown to inhibit the transmembrane receptor kit and platelet-derived growth factor (PDGF) receptors.

Citations: 24

[View Online »](#)

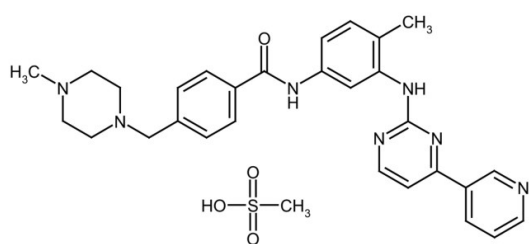
Ordering Information

[Order Online »](#)

ALX-270-492-M025	25mg
ALX-270-492-M500	500mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name STI-571, CGP-57148B

Appearance White to pale yellow crystalline powder.

CAS 220127-57-1

Couple Target Abl, PDGF receptor

Couple Type Inhibitor

Formula $C_{29}H_{31}N_7O \cdot CH_3SO_3H$

MI 14: 4902

MW 493.6 . 96.1

Purity ≥98% (HPLC)

Solubility Soluble in water (200mg/ml) or DMSO (100mg/ml) or DMF (10mg/ml) or PBS (2mg/ml). Sparingly soluble in ethanol (0.2mg/ml). Storing aqueous solutions for more than one day is not recommended.

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)