

# Dasatinib

## Broad-spectrum tyrosine kinase inhibitor

Dasatinib is a potent, orally active small-molecule inhibitor of multiple tyrosine kinases, including Bcr-Abl, Src family kinases, c-Kit, PDGFR, and EphA2. It inhibits Src, Abl, Lck, and c-Kit with an  $IC_{50}$  of 0.8 nM, 1 nM, 0.4 nM, and 79 nM, respectively. Originally developed as a second-generation treatment for chronic myeloid leukemia (CML), it is widely used in cancer research and cell signaling studies due to its broad kinase inhibition profile.

Key features and applications include:

- **Multi-Targeted Inhibition:** Blocks both receptor and non-receptor tyrosine kinases, including drug-resistant Bcr-Abl mutants.
- **Cancer Research:** Used to study leukemia, solid tumors, and kinase-driven malignancies.
- **Cell Signaling Studies:** Investigates pathways involving Src, Abl, and PDGFR signaling.
- **Immunology:** Modulates T-cell activation and natural killer (NK) cell function.
- **Fibrosis Models:** Explored for its anti-fibrotic effects in liver and lung disease models.
- **Autophagy & Apoptosis:** Induces programmed cell death and autophagy in various cancer cell lines.

Relevant disease states include:

- **Chronic Myeloid Leukemia (CML):** FDA-approved for treatment of CML, especially in cases resistant to imatinib.
- **Acute Lymphoblastic Leukemia (ALL):** Effective in Philadelphia chromosome-positive (Ph+) ALL.
- **Solid Tumors:** Investigated in breast, prostate, and pancreatic cancers for its anti-proliferative effects.
- **Fibrotic Diseases:** Studied for its potential to reduce fibrosis by targeting PDGFR and Src pathways.

- **Neurodegenerative Disorders:** Explored for neuroprotective effects via modulation of kinase signaling.

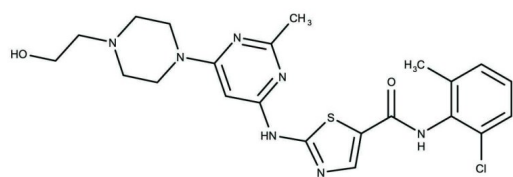
Ordering Information

Order Online »

ENZ-CHM333-0025	25mg
-----------------	------

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	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	BMS-354825, N-(2-Chloro-6-methylphenyl)-2-[[6-[4-(2-hydroxyethyl)-1-piperazinyl]-2-methyl-4-pyrimidinyl]amino]-5-thiazolecarboxamide
Appearance	White or off-white solid.
CAS	302962-49-8
Couple Target	Abl, c-Kit, p56lck tyrosine kinase, Src kinase
Couple Type	Inhibitor
Formula	$C_{22}H_{26}ClN_7O_2S$
Identity	Determined by NMR.
MW	488.01
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (200 mg/mL).



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)