

Lestaurtinib

FLT3/ TrkA inhibitor

Potent and selective FLT3 inhibitor. Inhibits RET and RET phosphorylation in medullary thyroid carcinoma cells. Suppresses JAK2/STAT5 signaling and the proliferation of primary erythroid cells from patients with myeloproliferative disorders. Potent Trk inhibitor. Cell permeable.

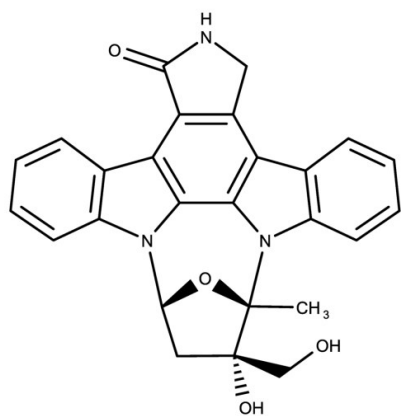
Ordering Information

[Order Online »](#)

ENZ-CHM467-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name KT5555, CEP-701

Appearance White or off-white solid.

CAS 111358-88-4

Couple Target Flt, Trk

Couple Type Inhibitor

Formula $C_{26}H_{21}N_3O_4$

Identity Determined by NMR.

MW 439.46

Purity ≥98% (TLC)

Solubility Soluble in DMSO (100 mg/ml) or ethanol (20 mg/ml).

Last modified: February 27, 2025



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)