

(R)-BI-2536

Potent and selective inhibitor of mammalian Plk1

Polo-like kinase 1 (Plk1) is a regulator of the cell cycle that has been implicated in the pathology of many cancers. BI-2536 is a potent and selective small-molecule inhibitor of mammalian Plk1. It has inhibitory activity at subnanomolar concentrations and inhibits tumor growth in multiple tumor lines with IC_{50} values below 1 μ M. BI-2536 also shows an >1,000-fold selectivity for PIK1 versus a large panel of other kinases. BI-2536 also demonstrated low K_d values against sister kinases Plk2 and Plk3. Preclinical studies in human cancer cell lines have shown that BI-2536 disrupts spindle assembly, resulting in mitotic arrest and inducing apoptosis. BI-2536 is currently in clinical trials against several types of solid tumor cancers.

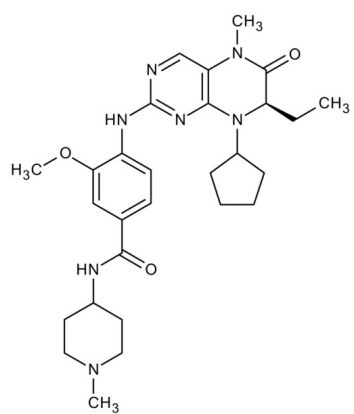
Ordering Information

[Order Online »](#)

ENZ-CHM160-0005	5mg
ENZ-CHM160-0010	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

Appearance	Off-white to yellow powder.
CAS	755038-02-9
Couple Target	PIK
Couple Type	Inhibitor
Formula	$C_{28}H_{39}N_7O_3$
Identity	Determined by NMR.
MW	521.7
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (21mg/ml) or ethanol (100mg/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-
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)