

mTOR inhibitor WYE-132

Highly potent, ATP-competitive, and specific mTOR kinase inhibitor. Antiproliferative.

A highly potent, ATP-competitive, and specific mTOR kinase inhibitor. Shows an anti-proliferative activity against a variety of tumor cell lines. WYE-132 causes inhibition of protein synthesis and cell size, induction of apoptosis, and cell cycle progression.

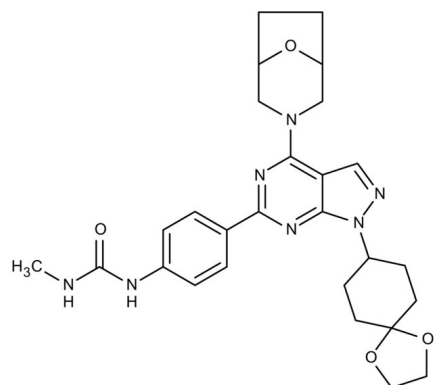
Ordering Information

[Order Online »](#)

ENZ-CHM140-0001	1mg
ENZ-CHM140-0005	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	1-{4-[1-(1,4-Dioxo-spiro[4.5]dec-8-yl)-4-(8-oxa-3-aza-bicyclo[3.2.1]oct-3-yl)-1H-pyrazolo[3,4-d]pyrimidin-6-yl]-phenyl}-3-methyl-urea, WYE-125132
Appearance	White powder solid.
CAS	1144068-46-1
Couple Target	mTOR
Couple Type	Inhibitor
Formula	$C_{27}H_{33}N_7O_4$
MW	519.6
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO.

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)