

mTOR inhibitor WYE-28

Cell-permeable, potent, selective, and ATP-competitive mTOR inhibitor.

Cell-permeable. A potent, selective, and ATP-competitive mTOR (mammalian target of rapamycin) inhibitor ($IC_{50} = 0.08$ nM). WYE-28 displays 75-fold selectivity over PI3K α ($IC_{50} = 6$ nM).

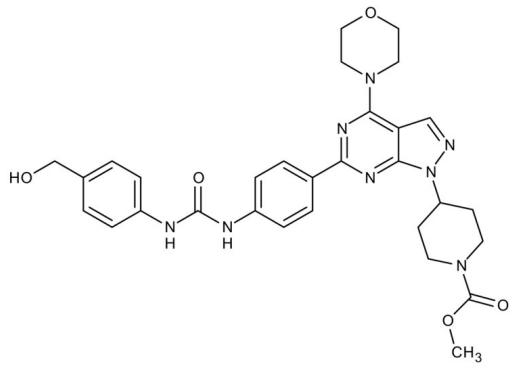
Ordering Information

[Order Online »](#)

ENZ-CHM137-0001	1mg
-----------------	-----

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	4-(6-{4-[3-(4-Hydroxymethyl-phenyl)-ureido]-phenyl}-4-morpholin-4-yl-pyrazolo[3,4-d]pyrimidin-1-yl)-piperidine-1-carboxylic acid methyl ester
Appearance	White powder solid.
CAS	1062172-60-4
Couple Target	mTOR
Couple Type	Inhibitor
Formula	$C_{30}H_{34}N_8O_5$
MW	586.6
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com