

Momelotinib

Selective inhibitor of JAK1 and JAK2 kinases including the JAK2 V617F mutant

Using ATP kinase assays, the IC_{50} values for CYT387 are as follows - JAK1 (11nM), JAK2, (18nM), JAK2 V617F (23nM) JAK3 (155nM), and TYK2 (17nM). In testing kinase arrays, CYT387 also had activity (defined as an $IC_{50} < 100nM$) detected against CDK2/cyclin A, MAPK8 (JNK1), PRKCN (PKD3), PRKD1 (PKC μ), ROCK2, and TBK1 kinase as well.

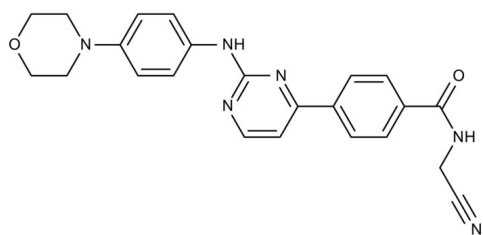
Ordering Information

[Order Online »](#)

ENZ-CHM197-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	CYT 387, N-(cyanomethyl)-4-[2-[[4-(4-morpholinyl)phenyl]amino]-4-pyrimidinyl]-benzamide
Appearance	Yellow to orange powder.
CAS	1056634-68-4
Couple Target	CDK, JAK, Jnk, PKD, ROCK, TANK-binding kinase, TYK
Couple Type	Inhibitor
Formula	$C_{23}H_{22}N_6O_2$
Identity	Determined by NMR.
MW	414.5
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (>20mM).

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)