

GSK-3 β inhibitor XVIII

Cell-permeable, potent and selective GSK-3 β inhibitor.

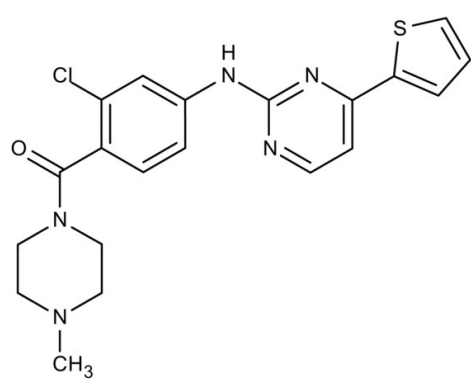
Ordering Information

[Order Online »](#)

ENZ-CHM134-0005	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	2-(Chloro-4-(4-thiophen-2-yl-pyrimidin-2-ylamino)-phenyl)-(4-methyl-pierazin-1-yl)-methanone
Appearance	White powder solid.
CAS	1139875-74-3
Couple Target	GSK
Couple Type	Inhibitor
Formula	$C_{20}H_{20}ClN_5OS$
MW	413.9
Purity	≥98% (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)