

GSK-3 inhibitor XIII

Potent inhibitor of GSK-3 ATP-binding site.

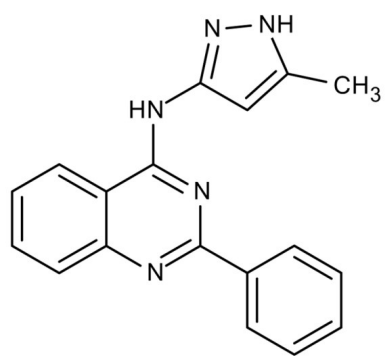
Ordering Information

[Order Online »](#)

| | |
|-----------------|-----|
| ENZ-CHM124-0001 | 1mg |
| ENZ-CHM124-0005 | 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 | (5-Methyl-1H-pyrazol-3-yl)-(2-phenylquinazolin-4-yl)amine |
| Appearance | White crystalline solid. |
| CAS | 404828-08-6 |
| Couple Target | GSK |
| Couple Type | Inhibitor |
| Formula | $C_{18}H_{15}N_5$ |
| MW | 301.4 |
| Purity | ≥99% (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