

# IWP-4

## Porcn inhibitor: Wnt/ $\beta$ -catenin pathway inhibitor

IWP-4 is an inhibitor of Wnt secretion and processing. It blocks Wnt-dependent signaling by inhibition of the O-palmitoyltransferase Porcn. Induces cardiomyocyte differentiation from human pluripotent stem cells. An important tool for probing the involvement of the Wnt pathway in physiological processes. Cell permeable.

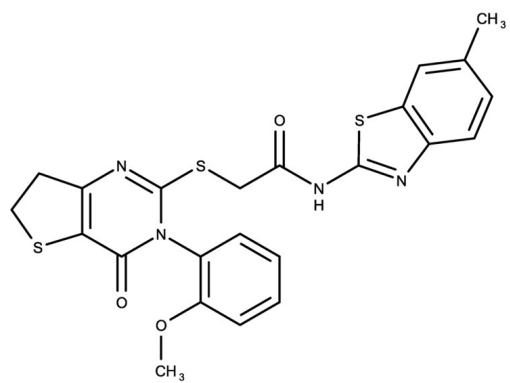
### Ordering Information

[Order Online »](#)

ENZ-CHM444-0025	25mg
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N-(6-Methyl-2-benzothiazolyl)-2-[[[3,4,6,7-tetrahydro-3-(2-methoxyphenyl)-4-oxothieno[3,2-d]pyrimidin-2-yl]thio]-acetamide

**Appearance** Pale yellow solid.

**CAS** 686772-17-8

**Couple Type** Inhibitor

**Formula**  $C_{23}H_{20}N_4O_3S_3$

**Identity** Determined by NMR.

**MW** 496.62

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO (5 mg/ml).

**Last modified: January 24, 2025**



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)