

# CCT036477

## Disrupts Wnt signaling

Inhibits Wnt signaling *via* blocking transcription at the  $\beta$ -catenin level independent of CBP or p300. Inhibited growth in several cancer cell lines (HT29, HCT116, SW480, and SNU475). CCT036447 was also active *in vivo* where it blocked development of zebrafish and *Xenopus* embryos and the expression of Wnt target genes.

Citations: 1

[View Online »](#)

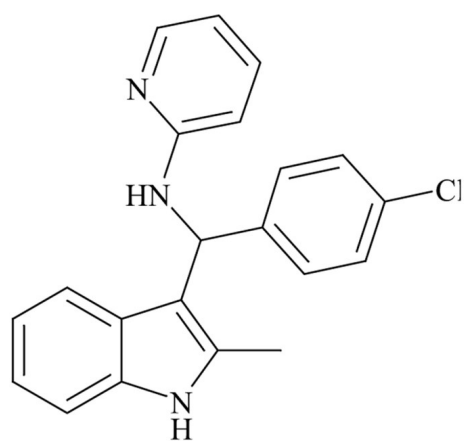
## Ordering Information

[Order Online »](#)

BML-WN114-0005	5mg
BML-WN114-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months after receipt when stored at -20°C. Solutions can be stored at -20°C for up to 3 months.
Handling	Solutions can be stored at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	N-((4-Chlorophenyl)(2-methyl-1H-indol-3-yl)methyl)pyridin-2-amine
Appearance	Yellow solid.
CAS	305372-78-5
Couple Target	Wnt
Couple Type	Inhibitor
Formula	$C_{21}H_{18}ClN_3$
MW	347.8
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml).



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)