

CCT036477

Disrupts Wnt signaling

Inhibits Wnt signaling *via* blocking transcription at the β -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 »](#)

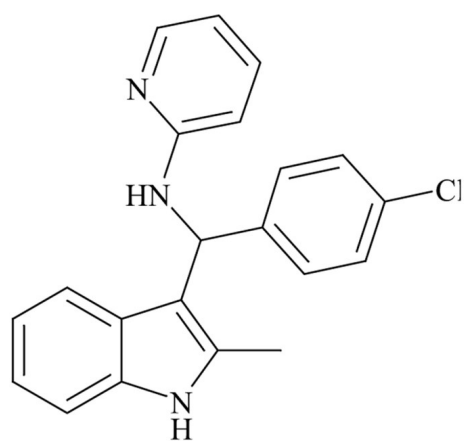
Ordering Information

[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).

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-fr@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