

Wnt-C59

PORCN inhibitor

Wnt-C59 is a highly potent and selective inhibitor of Porcupine (PORCN), a membrane-bound O-acyltransferase essential for Wnt protein palmitoylation and secretion. It has an IC₅₀ value of 74 pM in Wnt signaling reporter assays.

Key features and applications include:

- **Cancer Research:** Wnt-C59 is used to study the inhibition of Wnt signaling pathways, which are crucial in the development and progression of various cancers. It has been shown to block the growth of Wnt-driven mammary tumors in animal models.
- **Stem Cell Research:** This compound is employed to induce differentiation of human induced pluripotent stem cells (iPSCs) into cardiomyocytes and neural cells.
- **Developmental Biology:** Wnt-C59 is utilized to investigate the roles of Wnt signaling in embryonic development and tissue regeneration.

Relevant disease states include:

- **Cancer:** Wnt-C59 has shown potential in inhibiting the growth of various cancers by blocking Wnt signaling, which is often upregulated in tumors.
- **Neurodegenerative Diseases:** It is used in research to study its effects on neural differentiation, which can be relevant for conditions like Alzheimer's and Parkinson's diseases.
- **Developmental Disorders:** Wnt-C59 is studied for its role in inhibiting embryonic development processes, which can be relevant in understanding certain developmental disorders.

- GMP format available

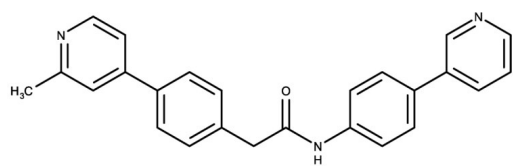
Ordering Information

[Order Online »](#)

ENZ-CHM345-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO or ethanol may be stored at -20°C for up to 1 month.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-(2-Methyl-4-pyridinyl)-N-[4-(3-pyridinyl)phenyl]benzeneacetamide, C59
Appearance	Tan solid.
CAS	1243243-89-1
Couple Target	Palmitoyltransferase
Couple Type	Inhibitor
Formula	$C_{25}H_{21}N_3O$
Identity	Determined by NMR.
MW	379.46
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 7 mg/ml) or in ethanol (up to 6 mg/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)