

PNU-74654

Catenin blocker

Anti-tumor agent that disrupts the canonical Wnt/ β -catenin pathway *via* blockage of the protein-protein interaction between β -catenin and TCF.

Citations: 3

[View Online »](#)

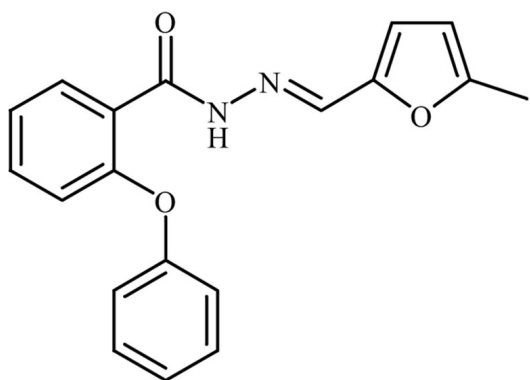
Ordering Information

[Order Online »](#)

BML-WN101-0005	5mg
BML-WN101-0025	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.
Handling	Solutions can be stored at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Phenoxybenzoic acid [(5-methyl-2-furanyl)methylene]hydrazide
Appearance	Off-white solid.
CAS	113906-27-7
Couple Target	Catenin
Couple Type	Blocker
Formula	$C_{19}H_{16}N_2O_3$
MW	320.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25 mg/ml) and in 100% ethanol (8 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)