

# BML-284

## Wnt activator

Wnt agonist that activates Wnt signaling in cellular assays and a *Xenopus* model. Induces  $\beta$ -catenin and TCF-dependent transcriptional activity in a dose-dependent manner ( $EC_{50}=0.7\mu\text{M}$ ). Wnt activation occurred in a GSK-3 $\beta$  independent manner.

Citations: 1

[View Online »](#)

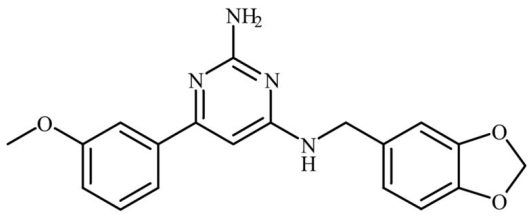
## Ordering Information

[Order Online »](#)

BML-WN105-0005	5mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	2-Amino-4-[3,4-(methylenedioxy)benzylamino]-6-(3-methoxyphenyl)pyrimidine
<b>Appearance</b>	Off-white solid.
<b>CAS</b>	853220-52-7
<b>Couple Target</b>	Wnt
<b>Couple Type</b>	Activator
<b>Formula</b>	$C_{19}H_{18}N_4O_3$
<b>MW</b>	350.4
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in DMSO (25 mg/ml) or in 100% ethanol (25 mg/ml).

Last modified: May 29, 2024



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-be@enzolifesciences.com](mailto:info-be@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)