

# IM-12

GSK-3 $\beta$  inhibitor (IC<sub>50</sub>=53nM (enzymatic), 3.8 $\mu$ M (cellular)). The compound significantly increases  $\beta$ -catenin levels, leads to TCF-induction and promotes neuronal differentiation in human neural progenitor cells *via* activation of canonical Wnt signaling.

Citations: 8

[View Online »](#)

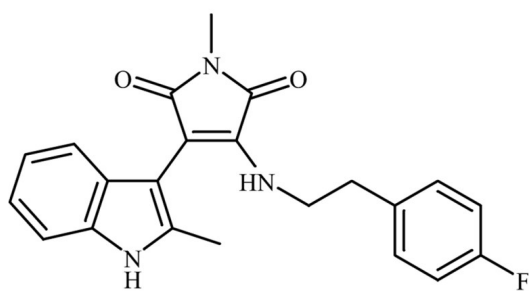
## Ordering Information

[Order Online »](#)

BML-WN102-0005	5mg
BML-WN102-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	3-(4-Fluorophenylethylamino)-1-methyl-4-(2-methyl-1H-indol-3-yl)-1H-pyrrole-2,5-dione
Appearance	Orange solid.
CAS	1129669-05-1
Couple Target	GSK
Couple Type	Inhibitor
Formula	$C_{22}H_{20}FN_3O_2$
MW	377.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25 mg/ml) or in 100% ethanol (5 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-  
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)