

# IQ-1

## Disrupts Wnt signaling

IQ-1 in conjunction with Wnt3a has been shown to be sufficient to maintain murine embryonic stem cell proliferation in the absence of serum. Binds to the PR72/PR130 subunit of protein phosphatase PP2A preventing  $\beta$ -catenin from switching co-activator usage from CBP to p300 leading to an increase in  $\beta$ -catenin/CBP-driven transcription at the expense of  $\beta$ -catenin/p300-driven transcription. Upregulates expression of Oct4 and Sox2 transcription factors.

Citations: 1

[View Online »](#)

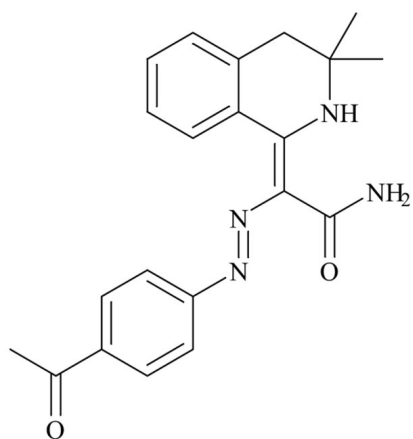
## Ordering Information

[Order Online »](#)

BML-WN115-0005	5mg
BML-WN115-0025	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.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2-(4-Acetylphenylazo)-2-(3,3-dimethyl-3,4-dihydro-2H-isoquinolin-1-ylidene)-acetamide
Appearance	Orange solid.
CAS	331001-62-8
Couple Target	Catenin
Formula	$C_{21}H_{22}N_4O_2$
MW	362.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or in 100% ethanol (10mg/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-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)