

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 β -catenin from switching co-activator usage from CBP to p300 leading to an increase in β -catenin/CBP-driven transcription at the expense of β -catenin/p300-driven transcription. Upregulates expression of Oct4 and Sox2 transcription factors.

Citations: 1

[View Online »](#)

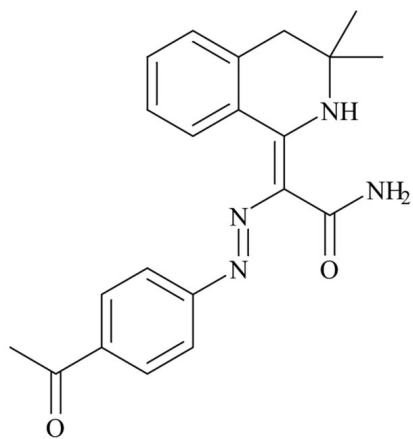
Ordering Information

[Order Online »](#)

BML-WN115-0005	5mg
----------------	-----

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).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com