

QS11

Wnt pathway activator

Synergizers with Wnt3a to activate the Wnt/ β -catenin pathway. Binds and inhibits GTPase activating protein of ADP-ribosylation factor 1 (ARFGAP1), thus activating ARF. It has also been suggested that a possible mechanism by which QS11 synergizes Wnt is through ARFGAP1 inhibition leading to increased β -catenin translocation via ARF activation.

Citations: 1

[View Online »](#)

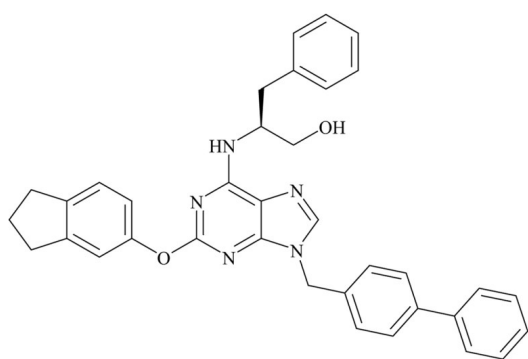
Ordering Information

[Order Online »](#)

BML-WN117-0005	5mg
BML-WN117-0025	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	(2S)-2-[2-(Indan-5-yloxy)-9-(1,1'-biphenyl-4-ylmethyl)-9H-purin-6-ylamino]-3-phenyl-propan-1-ol
Appearance	White solid.
CAS	944328-88-5
Couple Target	ADP-ribosylation factor, GTPase-activating protein
Couple Type	Activator, Inhibitor
Formula	$C_{35}H_{33}N_5O_2$
MW	567.7
Purity	≥98% (HPLC)
Solubility	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-
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)