

WAY-262611

DKK1 inhibitor

Wnt pathway agonist acting *via* inhibition of Dkk1 with no GSK-3 β inhibition. TCF-Luciferase EC₅₀=0.63 μ M. WAY262611 was able to enhance trabecular bone formation rate in the tibia of OVX rats following oral administration.

Citations: 2

[View Online »](#)

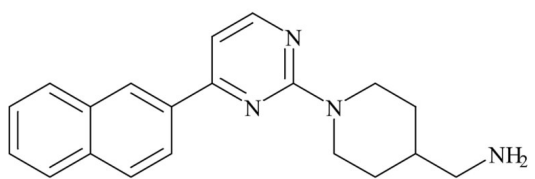
Ordering Information

[Order Online »](#)

BML-WN110-0005	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.
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	(1-(4-(Naphthalen-2-yl)pyrimidin-2-yl)piperidin-4-yl)methanamine
Appearance	Off-white solid.
CAS	123231-07-1
Couple Target	DKK
Couple Type	Inhibitor
Formula	$C_{20}H_{22}N_4$
MW	318.4
Purity	≥97% (TLC)
Solubility	Soluble in DMSO (30mg/ml) or in 100% ethanol (25mg/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