

NVP-TNKS656

TNKS2 inhibitor

NVP-TNKS657 is a potent and selective inhibitor of tankyrase enzymes, specifically targeting TNKS2 (IC₅₀ value of 6 nM). It exhibits over 300-fold selectivity against PARP1 and PARP2, making it directly applicable for research focused on the Wnt/ β -catenin signaling pathway.

Ordering Information

[Order Online »](#)

ENZ-CHM326-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

- GMP format available

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Powder: Stable at -20°C for 3 years or +4°C for 2 years; In solvent: Stable at -80°C for 2 years or -20°C for 1 year.
Handling	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Cyclopropylmethyl-2-[4-(4-methoxybenzoyl)piperidin-1-yl]-N-[(4-oxo-4,5,7,8-tetrahydro-4H-pyrano[4,3-d]pyrimidin-2-yl)methyl]acetamide, N-(Cyclopropylmethyl)-4-(4-methoxybenzoyl)-N-[(3,5,7,8-tetrahydro-4-oxo-4H-pyrano[4,3-d]pyrimidin-2-yl)methyl]-piperidineacetamide
Appearance	White to light yellow solid.
CAS	1419949-20-4
Couple Target	TNKS2
Couple Type	Inhibitor
Formula	$C_{27}H_{34}N_4O_5$
MW	494.58
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (35 mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-fr@enzolifesciences.com)
be@enzolifesciences.com

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
de@enzolifesciences.com

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

Last modified: November 4, 2024