

# NVP-TNKS656

## TNKS2 inhibitor

NVP-TNKS657 is a potent and selective inhibitor of tankyrase enzymes, specifically targeting TNKS2 (IC<sub>50</sub> value of 6 nM). It exhibits over 300-fold selectivity against PARP1 and PARP2, making it directly applicable for research focused on the Wnt/ $\beta$ -catenin signaling pathway.

### Ordering Information

[Order Online »](#)

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

### Manuals, SDS & CofA

[View Online »](#)

- GMP format available

## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	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
<b>Appearance</b>	White to light yellow solid.
<b>CAS</b>	1419949-20-4
<b>Couple Target</b>	TNKS2
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{27}H_{34}N_4O_5$
<b>MW</b>	494.58
<b>Purity</b>	≥99% (HPLC)
<b>Solubility</b>	Soluble in DMSO (35 mg/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)