# **PRODUCT** INFORMATION



NVP-TNKS656

Item No. 19930

| 1419949-20-4  |   |
|---|---|
| N-(cyclopropylmethyl)-4-(4-                                   | Q   |
| methoxybenzoyl)-N-[(3,5,7,8-tetrahydro-                       |   |
| 4-oxo-4H-pyrano[4,3-d]pyrimidin-2-yl)                         |   |
| methyl]-1-piperidineacetamide                                 |   |
| C <sub>27</sub> H <sub>34</sub> N <sub>4</sub> O <sub>5</sub> | н 🗸 🗸 🗸   |
| 494.6   | $\wedge$  |
| ≥98%  |   |
| λ <sub>max</sub> : 279 nm                                     |   |
| A crystalline solid   | $\checkmark \parallel \qquad \bigtriangledown$  |
| -20°C   | 0 V   |
| ≥4 years  |   |
|   | 1419949-20-4<br>N-(cyclopropylmethyl)-4-(4-<br>methoxybenzoyl)-N-[(3,5,7,8-tetrahydro-<br>4-oxo-4H-pyrano[4,3-d]pyrimidin-2-yl)<br>methyl]-1-piperidineacetamide<br>$C_{27}H_{34}N_4O_5$<br>494.6<br>≥98%<br>$\lambda_{max}$ : 279 nm<br>A crystalline solid<br>-20°C<br>≥4 years |

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

# Laboratory Procedures

NVP-TNKS656 is supplied as a crystalline solid. A stock solution may be made by dissolving the NVP-TNKS656 in the solvent of choice, which should be purged with an inert gas. NVP-TNKS656 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of NVP-TNKS656 in these solvents is approximately 33 mg/ml.

NVP-TNKS656 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NVP-TNKS656 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NVP-TNKS656 has a solubility of approximately 0.20 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

# Description

NVP-TNKS656 is an orally bioavailable inhibitor of tankyrase 2 (TNKS2;  $IC_{50} = 6 \text{ nM}$ ).<sup>1</sup> It is selective for TNKS2 over poly(ADP-ribose) polymerase (PARP) 1 and 2 ( $IC_{50}$ s = >19 and 32 nM, respectively). NVP-TNKS656 inhibits Wnt ligand-induced signaling with an  $IC_{50}$  value of 3.5 nM in an HEK293 cell reporter assay. In vivo, NVP-TNKS656 (350 mg/kg) reduces Axin 2 mRNA expression, a Wnt/β-catenin target gene, in MMTV-Wnt1 tumor bearing mice.

# Reference

1. Shultz, M.D., Cheung, A.K., Kirby, C.A., et al. Identification of NVP-TNKS656: The use of structureefficiency relationships to generate a highly potent, selective, and orally active tankyrase inhibitor. J. Med. Chem. 56(16), 6495-6511 (2013).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

## SAFETY DATA

SAFETY DATA This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

# WARRANTY AND LIMITATION OF REMEDY

uyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/30/2022

# CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD ANN ARBOR, MI 48108 · USA PHONE: [800] 364-9897 [734] 971-3335 FAX: [734] 971-3640 CUSTSERV@CAYMANCHEM.COM WWW.CAYMANCHEM.COM