

# PRODUCT INFORMATION

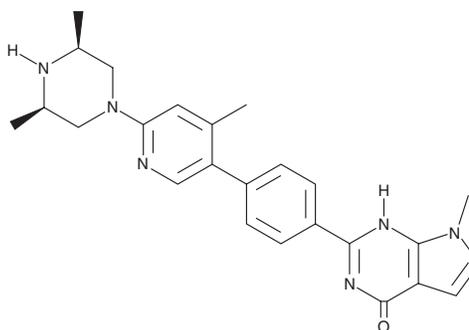


**AZ 6102**

Item No. 19898

**CAS Registry No.:** 1645286-75-4  
**Formal Name:** rel-2-[4-[6-[(3R,5S)-3,5-dimethyl-1-piperazinyl]-4-methyl-3-pyridinyl]phenyl]-3,7-dihydro-7-methyl-4H-pyrrolo[2,3-d]pyrimidin-4-one

**MF:** C<sub>25</sub>H<sub>28</sub>N<sub>6</sub>O  
**FW:** 428.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 327 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

## Laboratory Procedures

AZ 6102 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZ 6102 in the solvent of choice, which should be purged with an inert gas. AZ 6102 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of AZ 6102 in these solvents is approximately 10 and 15 mg/ml, respectively.

AZ 6102 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZ 6102 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AZ 6102 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

AZ 6102 is an inhibitor of tankyrases (TNKS) 1 and 2 with IC<sub>50</sub> values of 3 and 1 nM, respectively.<sup>1</sup> It demonstrates 100-fold selectivity for TNKS1/2 over PARP1, PARP2, and PARP6 (IC<sub>50</sub>s = 2, 0.5, and >3 μM, respectively).<sup>1</sup> By inhibiting TNKS1/2, AZ 6120 has been shown to inhibit Wnt signaling in DLD-1 cells with an IC<sub>50</sub> value of 5 nM.<sup>1</sup>

## Reference

1. Johannes, J. W., Almeida, L., Barlaam, B., *et al.* Pyrimidinone nicotinamide mimetics as selective tankyrase and Wnt pathways inhibitors suitable for *in vivo* pharmacology. *ACS Med. Chem. Lett.* **6**, 254-259 (2015).

### WARNING

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

### 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

Buyer 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, 11/04/2022

## 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