

# PRODUCT INFORMATION



## BMN 673

Item No. 19782

**CAS Registry No.:** 1207456-01-6  
**Formal Name:** (8S,9R)-5-fluoro-8-(4-fluorophenyl)-2,7,8,9-tetrahydro-9-(1-methyl-1H-1,2,4-triazol-5-yl)-3H-pyrido[4,3,2-de]phthalazin-3-one

**Synonym:** Talazoparib

**MF:** C<sub>19</sub>H<sub>14</sub>F<sub>2</sub>N<sub>6</sub>O

**FW:** 380.4

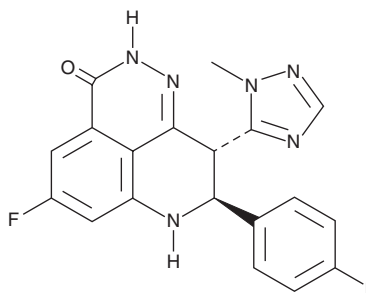
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 255, 309 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

BMN 673 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMN 673 in the solvent of choice, which should be purged with an inert gas. BMN 673 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BMN 673 in these solvents is approximately 0.25, 20, and 25 mg/ml, respectively.

BMN 673 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BMN 673 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BMN 673 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

BMN 673 is an orally available poly(ADP-ribose) polymerase (PARP) inhibitor with an IC<sub>50</sub> value of 0.57 nM.<sup>1</sup> It similarly binds to PARP1 and PARP2 with K<sub>i</sub> values of 1.2 and 0.85 nM, respectively, but has no effect on PARG.<sup>1</sup> BMN 673 selectively targets tumor cells with BRCA1, BRCA2, or PTEN gene mutations and elicits antitumor activity in xenografted tumors in mice.<sup>1</sup>

## Reference

1. Shen, Y., Rehman, F. L., Feng, Y., *et al.* BMN 673, a novel and highly potent PARP1/2 inhibitor for the treatment of human cancers with DNA repair deficiency. *Clin. Cancer. Res* **19(18)**, 5003-5015 (2013).

### 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, 09/29/2022

## 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