

# PRODUCT INFORMATION

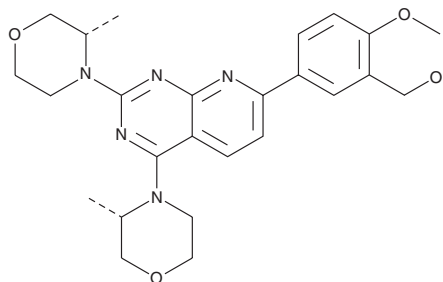


## AZD 8055

Item No. 16978

**CAS Registry No.:** 1009298-09-2  
**Formal Name:** 5-[2,4-bis[(3S)-3-methyl-4-morpholinyl]pyrido[2,3-d]pyrimidin-7-yl]-2-methoxybenzenemethanol

**Synonym:** CCG-168  
**MF:** C<sub>25</sub>H<sub>31</sub>N<sub>5</sub>O<sub>4</sub>  
**FW:** 465.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 211, 229, 281, 387 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

AZD 8055 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 8055 in the solvent of choice, which should be purged with an inert gas. AZD 8055 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AZD 8055 in these solvents is approximately 0.5, 1, and 10 mg/ml, respectively.

### Description

AZD 8055 is a potent, selective ATP-competitive inhibitor of mTOR, exhibiting an IC<sub>50</sub> value of 0.8 nM.<sup>1</sup> It is ~1,000-fold selective for mTOR over all PI3K isoforms and exhibits no activity against a panel of 260 kinases at concentrations up to 10 μM. AZD 8055 inhibits proliferation of A549 and H838 cells with IC<sub>50</sub> values of 53 and 20 nM, respectively. It inhibits growth of a variety of human tumor xenografts by at least 65% in mice at daily oral doses of 20 mg/kg.<sup>1</sup>

### Reference

1. Chresta, C.M., Davies, B.R., Hickson, I., *et al.* AZD8055 is a potent, selective, and orally bioavailable ATP-competitive mammalian target of rapamycin kinase inhibitor with *in vitro* and *in vivo* antitumor activity. *Cancer Res.* **70(1)**, 288-298 (2010).

**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, 09/30/2022

### 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