

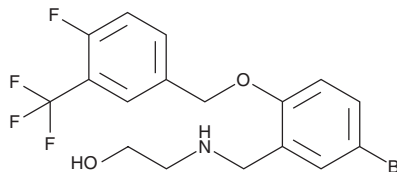
# PRODUCT INFORMATION



## AZ1 USP25/28 Inhibitor

Item No. 40335

**CAS Registry No.:** 2165322-94-9  
**Formal Name:** 2-[[[5-bromo-2-[[4-fluoro-3-(trifluoromethyl)phenyl]methoxy]phenyl]methyl]amino]-ethanol  
**Synonym:** AZ1 Ubiquitin-specific Protease 25/28 Inhibitor  
**MF:** C<sub>17</sub>H<sub>16</sub>BrF<sub>4</sub>NO<sub>2</sub>  
**FW:** 422.2  
**Purity:** ≥98%  
**Supplied as:** A 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

AZ1 USP25/28 Inhibitor is supplied as a solid. A stock solution may be made by dissolving the AZ1 USP25/28 inhibitor in the solvent of choice, which should be purged with an inert gas. AZ1 USP25/28 Inhibitor is soluble in acetonitrile and DMSO.

### Description

AZ1 USP25/28 Inhibitor is a dual inhibitor of ubiquitin-specific protease 25 (USP25) and USP28 (IC<sub>50</sub>s = 0.7 and 0.6 μM, respectively).<sup>1</sup> It is selective for USP25 and USP28 over USP2a (IC<sub>50</sub> = >100 μM), a panel of 22 other deubiquitinases at 10 μM, and a panel of eight caspases and cathepsins at 30 μM. AZ1 USP25/28 Inhibitor induces degradation of c-Myc in HCT116 cells in a concentration-dependent manner, an effect that can be reversed by the proteasome inhibitor MG132. It decreases the viability of several cancer cell lines (EC<sub>50</sub>s = 9.8-33.5 μM).

### Reference

1. Wrigley, J.D., Gavory, G., Simpson, I.A., *et al.* Identification and characterization of dual inhibitors of the USP25/28 deubiquitinating enzyme subfamily. *ACS Chem. Biol.* **12(12)**, 3113-3125 (2017).

#### 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, 03/08/2024

#### 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