

# PRODUCT INFORMATION

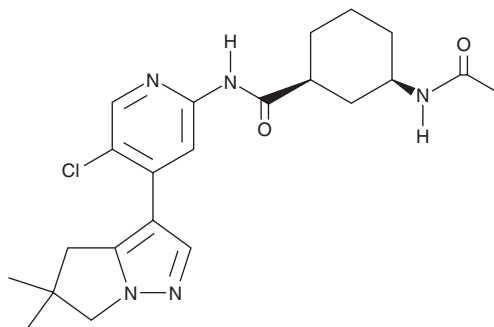


## AZD 4573

Item No. 33306

**CAS Registry No.:** 2057509-72-3  
**Formal Name:** (1S,3R)-3-(acetylamino)-N-[5-chloro-4-(5,6-dihydro-5,5-dimethyl-4H-pyrrolo[1,2-b]pyrazol-3-yl)-2-pyridinyl]-cyclohexanecarboxamide

**MF:** C<sub>22</sub>H<sub>28</sub>ClN<sub>5</sub>O<sub>2</sub>  
**FW:** 429.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 251 nm  
**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

AZD 4573 is supplied as a solid. A stock solution may be made by dissolving the AZD 4573 in the solvent of choice, which should be purged with an inert gas. AZD 4573 is soluble in organic solvents such as methanol and DMSO.

### Description

AZD 4573 is an inhibitor of cyclin-dependent kinase 9 (Cdk9; IC<sub>50</sub> = 14 nM).<sup>1</sup> It is selective for Cdk9 over Cdk1, Cdk7, Cdk4/6, and Cdk2 (IC<sub>50</sub>s = 370, 1,100, 1,100, and >10,000 nM, respectively). AZD 4573 (100 nM) induces apoptosis in MV4-11 acute myeloid leukemia (AML) cells. *In vivo*, AZD 4573 (15 mg/kg every other day), in combination with the Bcl-2 inhibitor venetoclax (ABT-199; Item No. 16233), reduces tumor volumes in SU-DHL-4 B cell lymphoma and OCI-AML-3 AML mouse xenograft models.

### Reference

1. Cidado, J., Boiko, S., Proia, T., *et al.* AZD4573 Is a highly selective CDK9 inhibitor that suppresses MCL-1 and induces apoptosis in hematologic cancer cells. *Clin. Cancer Res.* **26(4)**, 922-934 (2020).

#### 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, 09/25/2023

#### 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