

# PRODUCT INFORMATION

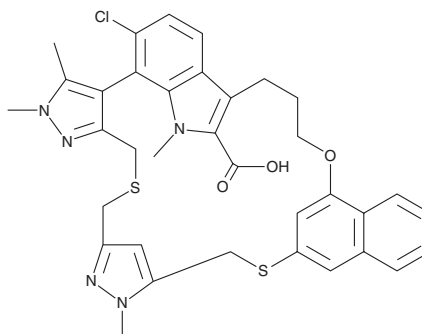


## AZD 5991

Item No. 28926

**CAS Registry No.:** 2143061-81-6  
**Formal Name:** (3aR)-5-chloro-2,11,12,24,27,29-hexahydro-2,3,24,33-tetramethyl-22H-9,4,8-(metheniminometheno)-14,20:26,23-dimetheno-10H,20H-pyrazolo[4,3-l][2,15,22,18,19]benzoxadithiadiazacyclohexacosine-32-carboxylic acid

**MF:** C<sub>35</sub>H<sub>34</sub>ClN<sub>5</sub>O<sub>3</sub>S<sub>2</sub>  
**FW:** 672.3  
**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

AZD 5991 is supplied as a solid. A stock solution may be made by dissolving the AZD 5991 in the solvent of choice, which should be purged with an inert gas. AZD 5991 is soluble in the organic solvent DMSO (sonicated) at a concentration of approximately 25 mg/ml.

### Description

AZD 5991 is an inhibitor of myeloid cell leukemia-1 (Mcl-1; IC<sub>50</sub> = 0.72 nM).<sup>1</sup> It is selective for Mcl-1 over Bcl-2, Bcl-xL, Bcl-W, and Bfl-1 (IC<sub>50</sub>s = 20, 36, 49, and 24 μM, respectively). AZD 5991 decreases the viability of MV4-11 acute myeloid leukemia (AML) and MOLP-8 multiple myeloma (MM) cells with EC<sub>50</sub> values of 24 and 33 nM, respectively. It disrupts formation of the Mcl-1-Bak complex and induces apoptosis in MOLP-8 cells when used at a concentration of 500 nM. AZD 5991 induces complete tumor regression in MV4-11 and MOLP-8 mouse xenograft models when administered at a single dose of 100 mg/kg.

### Reference

1. Tron, A.E., Belmonte, M.A., Adam, A., *et al.* Discovery of Mcl-1-specific inhibitor AZD5991 and preclinical activity in multiple myeloma and acute myeloid leukemia. *Nat. Commun.* **9**(1), 5341 (2018).

#### 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, 12/13/2022

#### 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