

PRODUCT INFORMATION

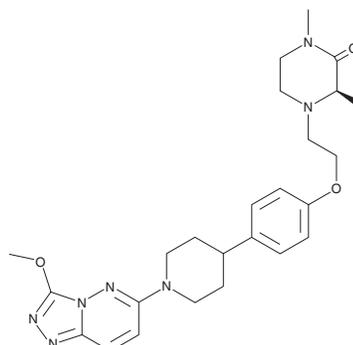


AZD 5153

Item No. 20864

CAS Registry No.: 1869912-39-9
Formal Name: (3R)-4-[2-[4-[1-(3-methoxy-1,2,4-triazolo[4,3-b]pyridazin-6-yl)-4-piperidinyl]phenoxy]ethyl]-1,3-dimethyl-2-piperazinone

MF: C₂₅H₃₃N₇O₃
FW: 479.6
Purity: ≥95%
UV/Vis.: λ_{max}: 239, 283, 303 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 5153 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 5153 in the solvent of choice, which should be purged with an inert gas. AZD 5153 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AZD 5153 in these solvents is approximately 20 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AZD 5153 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of AZD 5153 in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 5153 is an orally available, bivalent inhibitor of the bromodomain and extraterminal (BET) protein BRD4 (IC₅₀ = 5 nM).¹ It can simultaneously bind two bromodomains in BRD4, which has been shown to increase antitumor activity in multiple xenograft models of acute myeloid leukemia, multiple myeloma, and diffuse large B cell lymphoma.¹

Reference

1. Rhyasen, G.W., Hattersley, M.M., Yao, Y., *et al.* AZD5153: A novel bivalent BET bromodomain inhibitor highly active against hematologic malignancies. *Mol. Cancer Ther.* **15(11)**, 2563-2574 (2016).

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, 10/17/2022

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