

PRODUCT INFORMATION



AZD 4320

Item No. 41978

CAS Registry No.: 1357576-48-7

Formal Name: 4-[4-[(R)-(4'-chloro[1,1'-biphenyl]-2-yl)hydroxymethyl]-1-piperidinyl]-N-[[4-[[[(1R)-3-[(2-hydroxyethyl)methylamino]-1-[(phenylthio)methyl]propyl]amino]-3-[(trifluoromethyl)sulfonyl]phenyl]sulfonyl]-benzamide

MF: $C_{45}H_{48}ClF_3N_4O_7S_3$

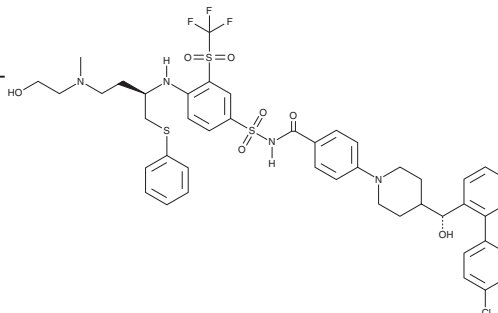
FW: 945.5

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

AZD 4320 is supplied as a solid. A stock solution may be made by dissolving the AZD 4320 in the solvent of choice, which should be purged with an inert gas. AZD 4320 is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

AZD 4320 is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 4320 is an inhibitor of the Bcl-2-Bim and Bcl-xL-Bim protein-protein interactions ($IC_{50}s = 3.9$ and 4.5 nM, respectively, in fluorescence polarization assays).¹ It induces apoptosis in MLL-AF9;Nras^{G12D} mouse acute myeloid leukemia (AML) cells overexpressing human Bcl-2 or Bcl-xL but not myeloid cell leukemia-1 (Mcl-1), Bfl-1, or Bcl-W when used at a concentration of $1 \mu M$. It also induces apoptosis in wild-type but not *Bax/Bak*^{-/-} SW48 cells. AZD 4320 (20 mg/kg) reduces tumor volume but induces transient thrombocytopenia in an RS4;11 acute lymphoblastic leukemia (ALL) mouse xenograft model.

Reference

1. Balachander, S.B., Criscione, S.W., Byth, K.F., *et al.* AZD4320, a dual inhibitor of Bcl-2 and Bcl-x_L, induces tumor regression in hematologic cancer models without dose-limiting thrombocytopenia. *Clin. Cancer Res.* **26**(24), 6535-6549 (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, 11/11/2024

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