

PRODUCT INFORMATION



WEHI-539 (hydrochloride)

Item No. 44330

CAS Registry No.: 2070018-33-4

Formal Name: 5-[3-[4-(aminomethyl)phenoxy]propyl]-2-[8-[2-(2-benzothiazolyl)hydrazinylidene]-5,6,7,8-tetrahydro-2-naphthalenyl]-4-thiazolecarboxylic acid, monohydrochloride

MF: $C_{31}H_{29}N_5O_3S_2 \cdot HCl$

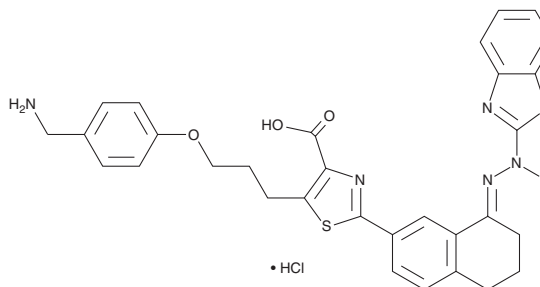
FW: 620.2

Purity: $\geq 95\%$

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

WEHI-539 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the WEHI-539 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. WEHI-539 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

WEHI-539 (hydrochloride) is slightly soluble 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

WEHI-539 is an inhibitor of Bcl-xL ($IC_{50} = 1.1$ nM).¹ It selectively binds to Bcl-xL ($K_d = 1.4$ nM) over Bcl-2, myeloid cell leukemia-1 (Mcl-1), Bcl-W, and Bfl-1, also known as Bcl-2-related protein A1 (K_d s = >750, >550, >550, and >550 nM, respectively). It selectively induces apoptosis in Mcl-1-deficient mouse embryonic fibroblasts (MEFs) over wild-type MEFs, Bcl-xL- or Bcl-2-deficient MEFs, and Mcl-1-overexpressing MEFs. WEHI-539 (1 μM) also induces apoptosis in SW480 colorectal cancer cells.²

References

1. Lessene, G., Czabotar, P.E., Sleebs, B.E., *et al.* Structure-guided design of a selective BCL-XL inhibitor. *ACS Med. Chem. Lett.* **5**(6), 662-667 (2014).
2. Scherr, A.L., Mock, A., Gdynia, G., *et al.* Identification of BCL-XL as highly active survival factor and promising therapeutic target in colorectal cancer. *Cell Death Dis.* **11**(10), 875 (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, 10/13/2025

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