

PRODUCT INFORMATION

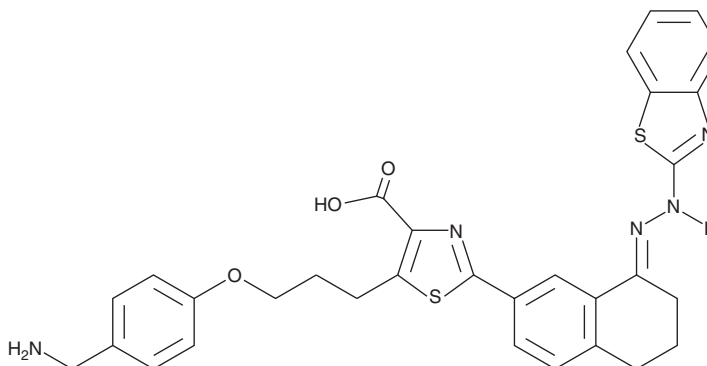


WEHI-539

Item No. 21478

CAS Registry No.: 1431866-33-9
Formal Name: 5-[3-[4-(aminomethyl)phenoxy]propyl]-2-[8-[2-(2-benzothiazolyl)hydrazinylidene]-5,6,7,8-tetrahydro-2-naphthalenyl]-4-thiazolecarboxylic acid

MF: C₃₁H₂₉N₅O₃S₂
FW: 583.7
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 252, 306, 335 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

WEHI-539 is supplied as a crystalline solid. A stock solution may be made by dissolving the WEHI-539 in the solvent of choice, which should be purged with an inert gas. WEHI-539 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of WEHI-539 in these solvents is approximately 1 and 15 mg/ml, respectively.

WEHI-539 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WEHI-539 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. WEHI-539 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

WEHI-539 is an inhibitor of Bcl-xL (IC₅₀ = 1.1 nM) that demonstrates more than 400-fold selectivity for Bcl-xL over Bcl-2, Mcl-1, Bcl-W, or A1.¹ It has been shown to induce apoptosis specifically in Mcl-1 deficient cells and does not affect Bcl-2 overexpressing cells.¹ WEHI-539 can be used as a tool for distinguishing the roles of Bcl-xL in both normal and cancer cells from those of the related prosurvival proteins such as Bcl-2 and Mcl-1.¹

Reference

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).

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, 11/09/2022

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