

PRODUCT INFORMATION



A-385358

Item No. 43417

CAS Registry No.: 406228-55-5

Formal Name: N-[[4-[[[(1R)-3-(dimethylamino)-1-[(phenylthio)methyl]propyl]amino]-3-nitrophenyl]sulfonyl]-4-(4,4-dimethyl-1-piperidiny)]-benzamide

MF: C₃₂H₄₁N₅O₅S₂

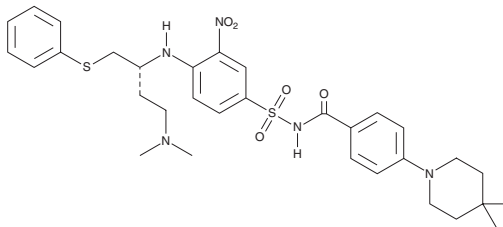
FW: 639.8

Purity: ≥95%

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

A-385358 is supplied as a solid. A stock solution may be made by dissolving the A-385358 in the solvent of choice, which should be purged with an inert gas. A-385358 is soluble (≥10 mg/ml) in acetonitrile and DMSO.

Description

A-385358 is an inhibitor of Bcl-xL ($K_i = 0.8$ nM).¹ It is selective for Bcl-xL over Bcl-2 ($K_i = 67$ nM). A-385358 decreases the proliferation of NCI H146 small cell lung carcinoma, MOLT-4 T lymphoblast, and CCRF CEM lymphoblastic leukemia cells in serum-free media (IC_{50} s = 0.35, 0.74, and 0.21 μ M, respectively). *In vivo*, A-385358 (100 mg/kg per day) in combination with paclitaxel (Item No. 10461) reduces tumor growth in an A549 lung cancer mouse xenograft model.

Reference

1. Shoemaker, A.R., Oleksijew, A., Bauch, J., *et al.* A small-molecule inhibitor of Bcl-XL potentiates the activity of cytotoxic drugs *in vitro* and *in vivo*. *Cancer Res.* **66(17)**, 8731-8739 (2006).

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, 05/16/2025

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