

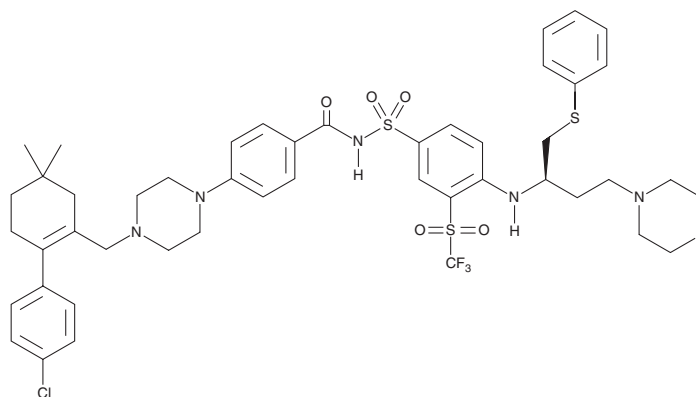
PRODUCT INFORMATION



ABT-263

Item No. 11500

CAS Registry No.: 923564-51-6
Formal Name: 4-[4-[[2-(4-chlorophenyl)-5,5-dimethyl-1-cyclohexen-1-yl]methyl]-1-piperazinyl]-N-[[4-[[[(1R)-3-(4-morpholinyl)-1-[(phenylthio)methyl]propyl]amino]-3-[(trifluoromethyl)sulfonyl]phenyl]sulfonyl]-benzamide
Synonyms: A-855071, Navitoclax
MF: C₄₇H₅₅ClF₃N₅O₆S₃
FW: 974.6
Purity: ≥98%
UV/Vis.: λ_{max}: 279, 320 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

ABT-263 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABT-263 in the solvent of choice, which should be purged with an inert gas. ABT-263 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ABT-263 in these solvents is approximately 0.5, 25, and 30, respectively.

ABT-263 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ABT-263 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ABT-263 has a solubility of approximately 0.3 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

ABT-263 is an inhibitor of the Bcl-2 family proteins Bcl-2, Bcl-xL, and Bcl-W (K_i s = <1, <0.5, and <1 nM, respectively).¹ It is selective for Bcl-2, Bcl-xL, and Bcl-W over Mcl-1 and A1 (K_i s = 550 and 354 nM, respectively). ABT-263 is cytotoxic in a panel of small cell lung cancer (SCLC), leukemia, and lymphoma cell lines with a mean EC_{50} value of 1 μ M. It induces apoptosis in influenza A-infected retinal pigment epithelial cells (EC_{50} = 0.08 μ M).² ABT-263 (100 mg/kg) induces tumor regression in NCI H146 SCLC and RS4;11 acute lymphocytic leukemia (ALL) mouse xenograft models.^{1,3}

References

1. Tse, C., Shoemaker, A.R., Adickes, J., *et al.* ABT-263: A potent and orally bioavailable Bcl-2 family inhibitor. *Cancer Res.* **68(9)**, 3421-3428 (2008).
2. Bulanova, D., Ianevski, A., Bugai, A., *et al.* Antiviral properties of chemical inhibitors of cellular anti-apoptotic Bcl-2 proteins. *Viruses* **9(10)**, E271 (2017).
3. Shoemaker, A.R., Mitten, M.J., Adickes, S., *et al.* Activity of the Bcl-2 family inhibitor ABT-263 in a panel of small cell lung cancer xenograft models. *Clin. Cancer Res.* **14(11)**, 3268-3277 (2008).

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, 11/18/2022

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