

ABT-263

Bcl-2 family inhibitor

ABT-263, also known as Navitoclax, is a potent and orally bioavailable inhibitor of the Bcl-2 family proteins, specifically targeting Bcl-2, Bcl-xL, and Bcl-w with K_i values of ≤ 0.5 nM, ≤ 1 nM, and ≤ 1 nM, respectively.

Key features and applications include:

- **High Selectivity:** Selectively inhibits Bcl-2, Bcl-xL, and Bcl-w over other Bcl-2 family proteins.
- **Potency:** Demonstrates strong inhibitory effects with low K_i values.
- **Induce Apoptosis:** Promotes apoptosis in various cancer cell lines, including small cell lung cancer (SCLC), leukemia, and lymphoma.
- **Enhance Chemotherapy:** Synergizes with chemotherapeutic agents to enhance their efficacy.
- **Reduce Tumor Growth:** Effective in mouse xenograft models of SCLC and acute lymphocytic leukemia (ALL).

Relevant disease states include:

- **Cancer:** Active in models of SCLC, leukemia, lymphoma, and other cancers.
- **Viral Infections:** Induces apoptosis in influenza A-infected retinal pigment epithelial cells.
- **Inflammation:** Reduces inflammation and viral-induced senescence in SARS-CoV-2 infection.

Ordering Information

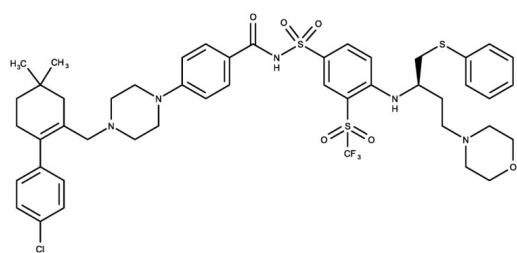
[Order Online »](#)

ENZ-CHM362-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20° for up to 3 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative 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, Navitoclax
Appearance	White solid.
CAS	923564-51-6
Couple Target	Bcl-2, Bcl-w, Bcl-X
Couple Type	Inhibitor
Formula	$C_{47}H_{55}ClF_3N_5O_6S_3$
Identity	Determined by NMR.
MW	974.61
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 25 mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: March 27, 2025