

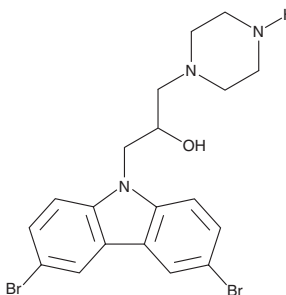
# PRODUCT INFORMATION



## Bax Channel Blocker

Item No. 37538

**CAS Registry No.:** 335165-68-9  
**Formal Name:** 3,6-dibromo- $\alpha$ -(1-piperazinylmethyl)-9H-carbazole-9-ethanol  
**Synonym:** BAI1  
**MF:** C<sub>19</sub>H<sub>21</sub>Br<sub>2</sub>N<sub>3</sub>O  
**FW:** 467.2  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Bax channel blocker is supplied as a solid. A stock solution may be made by dissolving the Bax channel blocker in the solvent of choice, which should be purged with an inert gas. Bax channel blocker is soluble in organic solvents such as DMSO, and dimethyl formamide. The solubility of Bax channel blocker in these solvents is approximately 10 and 5 mg/ml, respectively. Bax channel blocker is slightly soluble in ethanol.

### Description

Bax channel blocker is an inhibitor of Bax.<sup>1</sup> It binds to Bax in a cell-free assay ( $K_d = 15 \mu\text{M}$ ) and inhibits Bax-mediated membrane permeabilization induced by the membrane-targeted death ligand tBID in a liposomal release assay ( $IC_{50} = 3.3 \mu\text{M}$ ). Bax channel blocker inhibits apoptosis induced by TNF- $\alpha$ - and the protein synthesis inhibitor cycloheximide (Item No. 14126) in mouse embryonic fibroblasts (MEFs;  $IC_{50} = 1.8 \mu\text{M}$ ). It also inhibits cell death induced by the Bax activator BTS1 (Item No. 37305) in OCI-AML-3 cells.

### Reference

1. Garner, T.P., Amgalan, D., Reyna, D.E., *et al.* Small-molecule allosteric inhibitors of BAX. *Nat. Chem. Biol.* **15**(4), 322-330 (2019).

#### 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, 10/30/2023

#### 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