

# PRODUCT INFORMATION

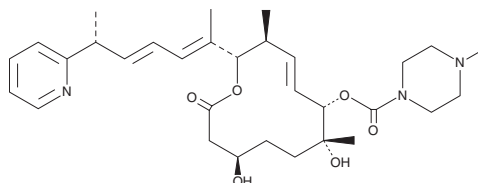


## H3B-8800

Item No. 42106

**CAS Registry No.:** 1825302-42-8  
**Formal Name:** 4-methyl-1-piperazinecarboxylic acid, (2S,3S,4E,6S,7R,10R)-7,10-dihydroxy-3,7-dimethyl-2-[(1E,3E,5R)-1-methyl-5-(2-pyridinyl)-1,3-hexadien-1-yl]-12-oxooxacyclododec-4-en-6-yl ester

**MF:** C<sub>31</sub>H<sub>45</sub>N<sub>3</sub>O<sub>6</sub>  
**FW:** 555.7  
**Purity:** ≥98%  
**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

H3B-8800 is supplied as a solid. A stock solution may be made by dissolving the H3B-8800 in the solvent of choice, which should be purged with an inert gas. H3B-8800 is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

### Description

H3B-8800 is an inhibitor of the splicing factor 3B (SF3B) complex.<sup>1</sup> It binds to SF3B complexes containing wild-type SF3B (IC<sub>50</sub> = 1.4 nM) or various SF3B mutations (IC<sub>50</sub>s = 0.3-1.8 nM) in cell-free assays. It inhibits splicing of canonical and aberrant pre-mRNA junctions induced by SF3B subunit 1 (SF3B1) in HEK293T nuclear extracts overexpressing SF3B1 with a lysine-to-glutamate mutation at residue 700 (SF3B1<sup>K700E</sup>). H3B-8800 selectively inhibits the growth of, and induces apoptosis in, Panc 05.04 pancreatic cancer cells expressing mutant SF3B over Panc 10.05, HAPF-II, and CFPAC-1 pancreatic cancer cells expressing wild-type SF3B in a concentration-dependent manner. It reduces tumor growth in a K562 chronic myeloid leukemia (CML) mouse xenograft model using K562 cells expressing wild-type SF3B1 when administered at a dose of 8 mg/kg or K562 cells expressing SF3B1<sup>K700E</sup> when administered at doses of 2, 4, or 8 mg/kg.

### Reference

1. Seiler, M., Yoshimi, A., Darman, R., *et al.* H3B-8800, an orally available small-molecule splicing modulator, induces lethality in spliceosome-mutant cancers. *Nat. Med.* **24**(4), 497-504 (2018).

#### 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, 11/11/2024

#### 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