

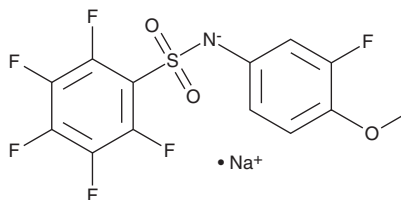
# PRODUCT INFORMATION



## Batabulin (sodium salt)

Item No. 33487

**CAS Registry No.:** 195533-98-3  
**Formal Name:** 2,3,4,5,6-pentafluoro-N-(3-fluoro-4-methoxyphenyl)-benzenesulfonamide, monosodium salt  
**Synonym:** T138067  
**MF:** C<sub>13</sub>H<sub>6</sub>F<sub>6</sub>NO<sub>3</sub>S • Na  
**FW:** 393.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 235 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

Batabulin (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the batabulin (sodium salt) in the solvent of choice, which should be purged with an inert gas. Batabulin (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of batabulin (sodium salt) in these solvents is approximately 30 mg/ml.

Batabulin (sodium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, batabulin (sodium salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Batabulin (sodium salt) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Batabulin is an inhibitor of tubulin polymerization.<sup>1</sup> It covalently binds to cysteine 239 (Cys<sup>239</sup>) of the β1, β2, and β4 tubulin isotypes and inhibits polymerization of bovine brain tubulin (IC<sub>50</sub> = 2 μM) in cell-free assays. Batabulin (30, 100, and 300 nM) induces cytoskeletal collapse and cell cycle arrest at the G<sub>2</sub>/M phase in MCF-7 breast cancer cells. It is cytotoxic against a panel of drug-sensitive and -resistant cancer cell lines (IC<sub>50</sub>s = 11-165 nM). Batabulin (40 mg/kg) reduces tumor volume in CCRF CEM lymphoblastic leukemia and CEM/VBL100 multidrug-resistant lymphoblastic leukemia mouse xenograft models.

### Reference

1. Shan, B., Medina, J.C., Santha, E., *et al.* Selective, covalent modification of β-tubulin residue Cys-239 by T138067, an antitumor agent with *in vivo* efficacy against multidrug-resistant tumors. *Proc. Natl. Acad. Sci. USA* **96**(10), 5686-5691 (1999).

#### 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