

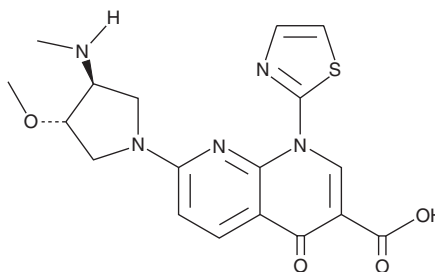
# PRODUCT INFORMATION



## Vosaroxin

Item No. 30614

**CAS Registry No.:** 175414-77-4  
**Formal Name:** 1,4-dihydro-7-[(3S,4S)-3-methoxy-4-(methylamino)-1-pyrrolidinyl]-4-oxo-1-(2-thiazolyl)-1,8-naphthyridine-3-carboxylic acid  
**Synonyms:** SNS-595, Voreloxin  
**MF:** C<sub>18</sub>H<sub>19</sub>N<sub>5</sub>O<sub>4</sub>S  
**FW:** 401.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 273, 345 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Vosaroxin is supplied as a crystalline solid. A stock solution may be made by dissolving the vosaroxin in the solvent of choice, which should be purged with an inert gas. Vosaroxin is slightly soluble in methanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of vosaroxin can be prepared by directly dissolving the crystalline solid in aqueous buffers. Vosaroxin is slightly soluble in PBS, pH 7.2. We do not recommend storing the aqueous solution for more than one day.

### Description

Vosaroxin is an inhibitor of topoisomerase II.<sup>1</sup> It induces relaxation of supercoiled DNA with an IC<sub>50</sub> value of 3.2 µg/ml in a cell-free assay. It induces DNA double-strand breaks in CCRF-CEM cells and cell cycle arrest at the G<sub>2</sub> phase in A549 cells in a concentration-dependent manner.<sup>2</sup> Vosaroxin inhibits proliferation in a panel of 15 cancer cell lines, including cancer cells resistant to doxorubicin (Item No. 15007), etoposide (Item No. 12092), or cisplatin (Item No. 13119), with IC<sub>50</sub> values ranging from 0.04 to 1.15 µM.<sup>3</sup> It inhibits tumor growth in multidrug-resistant SBC-3/ETP or PC-14/CDDP mouse xenograft models when administered at a dose of 20 mg/kg.

### References

1. Tsuzuki, Y., Tomita, K., Shibamori, K.-i., *et al.* Synthesis and structure-activity relationships of novel 7-substituted 1,4-dihydro-4-oxo-1-(2-thiazolyl)-1,8-naphthyridine-3-carboxylic acids as antitumor agents. Part 2. *J. Med. Chem.* **47**(8), 2079-2109 (2004).
2. Hawtin, R.E., Stockett, D.E., Byl, J.A.W., *et al.* Voreloxin is an anticancer quinolone derivative that intercalates DNA and poisons topoisomerase II. *PLoS One.* **5**(4), e10186 (2010).
3. Hoch, U., Lynch, J., Sato, Y., *et al.* Voreloxin, formerly SNS-595, has potent activity against a broad panel of cancer cell lines and in vivo tumor models. *Cancer Chemother. Pharmacol.* **64**(1), 53-65 (2009).

#### 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, 07/30/2020

#### 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