

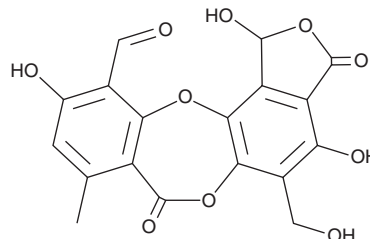
# PRODUCT INFORMATION



## Salazinic Acid

Item No. 29885

**CAS Registry No.:** 521-39-1  
**Formal Name:** 1,3-dihydro-1,4,10-trihydroxy-5-(hydroxymethyl)-8-methyl-3,7-dioxo-7H-isobenzofuro[4,5-b][1,4]benzodioxepin-11-carboxaldehyde  
**Synonym:** NSC 87509  
**MF:** C<sub>18</sub>H<sub>12</sub>O<sub>10</sub>  
**FW:** 388.3  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Usnea* sp.



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

### Laboratory Procedures

Salazinic acid is supplied as a solid. A stock solution may be made by dissolving the salazinic acid in the solvent of choice, which should be purged with an inert gas. Salazinic acid is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Salazinic acid is a depsidone lichen metabolite that has been found in *P. sulcata*.<sup>1</sup> It is active against *B. cereus*, *B. subtilis*, *S. aureus*, *P. aeruginosa*, *S. typhimurium*, *C. albicans*, and *A. niger* *in vitro* (MICs = 3.9-30.8 mM). Salazinic acid is cytotoxic to MM98, A431, and HaCaT cells in crystal violet (EC<sub>50</sub>s = 159, 2,870, and 48 μM, respectively) and neutral red uptake assays (EC<sub>50</sub>s = 1,925, 1,913, and 907 μM, respectively).<sup>2</sup> It increases the wound closure rate in scratch-wounded HaCaT monolayers and increases HaCaT cell migration in a transwell assay when used at a concentration of 30 μM.

### References

1. Candan, M., Yilmaz, M., Tay, T., *et al.* Antimicrobial activity of extracts of the lichen *Parmelia sulcata* and its salazinic acid constituent. *Z. Naturforsch. C J. Biosci.* **62(7-8)**, 619-621 (2007).
2. Burlando, B., Ranzato, E., Volante, A., *et al.* Antiproliferative effects on tumour cells and promotion of keratinocyte wound healing by different lichen compounds. *Planta. Med.* **75(6)**, 607-613 (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, 08/16/2023

#### 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