

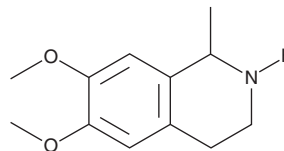
# PRODUCT INFORMATION



## (±)-Salsolidine

Item No. 32953

**CAS Registry No.:** 5784-74-7  
**Formal Name:** 1,2,3,4-tetrahydro-6,7-dimethoxy-1-methyl-isoquinoline  
**Synonym:** N-Norcarnegine  
**MF:** C<sub>12</sub>H<sub>17</sub>NO<sub>2</sub>  
**FW:** 207.3  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

(±)-Salsolidine is supplied as a crystalline solid. A stock solution may be made by dissolving the (±)-salsolidine in the solvent of choice, which should be purged with an inert gas. (±)-Salsolidine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (±)-salsolidine in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in 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 (±)-salsolidine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of (±)-salsolidine in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

(±)-Salsolidine is an alkaloid that has been found in *H. articulatum*.<sup>1</sup> It inhibits catechol-O-methyltransferase (COMT; K<sub>i</sub> = 190 μM).<sup>2</sup>

### References

1. El-Shazly, A. and Wink, M. Tetrahydroisoquinoline and β-carboline alkaloids from *Haloxylon articulatum* (Cav.) Bunge (Chenopodiaceae). *Z. Naturforsch C. J. Biosci.* **58(7-8)**, 477-480 (2003).
2. Sanft, K. and Thomas, H. Competitive inhibition of catechol-O-methyltransferase by the tetrahydroisoquinoline alkaloids salsolidine and 1-carboxysalsoline. *Z. Naturforsch C. J. Biosci.* **44(1-2)**, 173-176 (1989).

#### 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, 07/25/2024

#### 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