

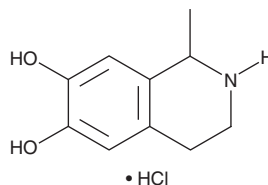
# PRODUCT INFORMATION



## (±)-Salsolinol (hydrochloride)

Item No. 15597

**CAS Registry No.:** 79923-51-6  
**Formal Name:** 1,2,3,4-tetrahydro-1-methyl-6,7-isoquinolinediol, monohydrochloride  
**MF:** C<sub>10</sub>H<sub>13</sub>NO<sub>2</sub> • HCl  
**FW:** 215.7  
**Purity:** ≥99%  
**UV/Vis.:** λ<sub>max</sub>: 288 nm  
**Supplied as:** A solid  
**Storage:** 4°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

(±)-Salsolinol (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the (±)-salsolinol (hydrochloride) in the solvent of choice. (±)-Salsolinol (hydrochloride) is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

(±)-Salsolinol is a metabolite of ethanol produced by the condensation of dopamine with acetaldehyde in the brain. It is thought to contribute to some aspects of alcohol's addictive properties in part through its ability to stimulate dopaminergic neurons in the posterior ventral tegmental area of the brain.<sup>1</sup> At 0.1 μM, (±)-salsolinol has been shown to: a) depolarize dopaminergic neurons, which increases their firing rate; b) activate μ-opioid receptors on GABAergic neurons, which disinhibits dopaminergic neurons; and c) activate dopamine1-like receptors, which enhances pre-synaptic glutamate transmission to dopaminergic neurons.<sup>1</sup>

### Reference

1. Xie, G., Krnjevic, K., and Ye, J.H. Salsolinol modulation of dopamine neurons. *Front. Behav. Neurosci.* 7(52), (2013).

#### 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, 12/12/2022

#### 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