

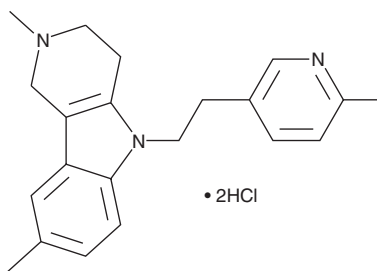
# PRODUCT INFORMATION



## Dimebolin (hydrochloride)

Item No. 10011349

**CAS Registry No.:** 97657-92-6  
**Formal Name:** 2,3,4,5-tetrahydro-2,8-dimethyl-5-[2-(6-methyl-3-pyridinyl)ethyl]-1H-pyrido[4,3-b]indole, dihydrochloride  
**Synonyms:** Dimebon, Latrepirdine  
**MF:** C<sub>21</sub>H<sub>25</sub>N<sub>3</sub> • 2HCl  
**FW:** 392.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 270 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

Dimebolin (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the dimebolin (hydrochloride) in the solvent of choice. Dimebolin (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of dimebolin (hydrochloride) in ethanol, DMSO, and DMF is approximately 3, 10, and 1 mg/ml, respectively.

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 dimebolin (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of dimebolin (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Dimebolin is a neuroprotective agent.<sup>1-3</sup> It binds to histamine H<sub>1</sub> and H<sub>2</sub> receptors (IC<sub>50</sub>s = 3.8 and 287 nM, respectively), as well as α<sub>1A</sub>-, α<sub>1B</sub>-, α<sub>1D</sub>-, and α<sub>2A</sub>-adrenergic receptors (ARs), imidazoline I<sub>2</sub> receptors, and the serotonin (5-HT) receptor subtypes 5-HT<sub>2A</sub>, 5-HT<sub>2C</sub>, 5-HT<sub>6</sub>, and 5-HT<sub>7</sub> (K<sub>s</sub> = 8-313 nM).<sup>4</sup> It also inhibits NMDA-evoked currents and voltage-gated calcium channels in mouse primary striatal neurons (IC<sub>50</sub>s = 10 and 50 μM, respectively).<sup>1</sup> Dimebolin (50 μM) inhibits glutamate-induced apoptosis in primary striatal neurons derived from the YAC128 transgenic mouse model of Huntington's disease. It inhibits cell death induced by amyloid-β (25-35) (Item No. 24155) in cerebellar granule cells when used at a concentration of 25 μM.<sup>2</sup> Dimebolin (1 mg/kg) inhibits decreases in two-way active avoidance reaction (TWAA) acquisition in a rat model of Alzheimer's disease induced by 1-ethyl-1-(2-hydroxyethyl) aziridinium (AF64A).<sup>3</sup>

### References

1. Wu, J., Li, Q., and Bezprozvanny, I. *Molecular Neurodegeneration* **3**(15), (2008).
2. Lermontova, N.N., Redkozubov, A.E., Shevtsova, E.F., et al. *B. Exp. Biol. Chem.* **132**(5), 1079-1083 (2001).
3. Lermontova, N.N., Lukoyanov, N.V., Serkova, T.P., et al. *B. Exp. Biol. Chem.* **129**(6), 640-642 (2000).
4. Okun, I., Tkachenko, S.E., Khvat, A., et al. *Curr. Alzheimer Res.* **7**(2), 97-112 (2010).

#### 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, 04/04/2024

#### 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