

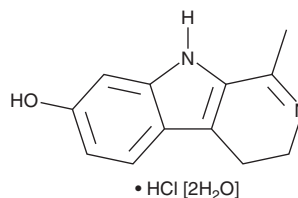
# PRODUCT INFORMATION



## Harmalol (hydrochloride hydrate)

Item No. 36090

**CAS Registry No.:** 6028-00-8  
**Formal Name:** 4,9-dihydro-1-methyl-3H-pyrido[3,4-b]indol-7-ol, monohydrochloride, dihydrate  
**Synonyms:** 11-hydroxy Harmalan, Harmidol  
**MF:**  $C_{12}H_{12}N_2O \cdot HCl [2H_2O]$   
**FW:** 272.7  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 216, 386 nm  
**Supplied as:** A solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years  
**Item Origin:** Plant/Unspecified sp.



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

### Laboratory Procedures

Harmalol (hydrochloride hydrate) is supplied as a solid. A stock solution may be made by dissolving the harmalol (hydrochloride hydrate) in the solvent of choice, which should be purged with an inert gas. Harmalol (hydrochloride hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of harmalol (hydrochloride hydrate) in these solvents is approximately 2, 1, and 10 mg/ml, respectively.

### Description

Harmalol is a  $\beta$ -carboline alkaloid and an active metabolite of harmaline that has been found in *P. harmala* and has diverse biological activities.<sup>1-5</sup> It is an inhibitor of dual-specificity tyrosine phosphorylation-regulated kinase 1A (DYRK1A;  $IC_{50} = 0.63 \mu M$ ) and monoamine oxidase A (MAO-A;  $IC_{50} = 0.66 \mu M$ ).<sup>3</sup> It is selective for DYRK1A over Cdk1, Cdk5, CK1 $\alpha$ 1, Clk4, DYRK2, Pim-1, and GSK3 $\beta$  but also inhibits DYRK1B and Clk1 at 10  $\mu M$ . It inhibits proliferation of H4 human glioblastoma cells ( $IC_{50}$ s = 23.7  $\mu M$ ). Harmalol (0.5-12.5  $\mu M$ ) reduces increases in the levels of the cytochrome P450 (CYP) isoform CYP1A1 induced by 2,3,7,8-tetrachlorodibenzo-*p*-dioxin (TCDD) in HepG2 cells and prevents TCDD-induced activation of the aryl hydrocarbon receptor (AhR) in guinea pig hepatic cytosolic extracts.<sup>4</sup> It reduces glutamate-induced cytotoxicity, cytochrome c release, caspase-3 activation, and the production of reactive oxygen species (ROS) in PC12 cells when used at a concentration of 25  $\mu M$ .<sup>5</sup> This product is also available as an analytical reference standard (Item No. 35145).

### References

1. Brierley, D.I. and Davidson, C. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **39**(2), 263-272 (2012).
2. Nikam, T.D., Nitnaware, K.M., and Ahire, M.L. *Natural products. Phytochemistry, botany and metabolism of alkaloids, phenolics and terpenes*. Ramawat, K.G. and Mérillon, J.-M., editors, Springer (2013).
3. Tarpley, M., Oladapo, H.O., Strepay, D., et al. *Eur. J. Pharm. Sci.* **162**, 105821 (2021).
4. El Gendy, M.A.M., Soshilov, A.A., Denison, M.S., et al. *Food Chem. Toxicol.* **50**(2), 353-362 (2012).
5. Han, E.S. and Lee, C.S. *Biomol. Ther. (Seoul)* **11**(2), 112-118 (2003).

#### 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, 12/19/2022

#### 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