

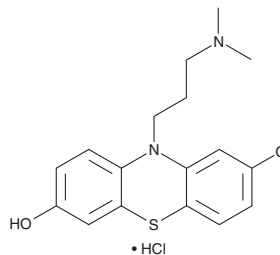
# PRODUCT INFORMATION



## 7-hydroxy Chlorpromazine (hydrochloride)

Item No. 33729

**CAS Registry No.:** 51938-11-5  
**Formal Name:** 8-chloro-10-[3-(dimethylamino)propyl]-10H-phenothiazin-3-ol, monohydrochloride  
**MF:** C<sub>17</sub>H<sub>19</sub>ClN<sub>2</sub>OS • HCl  
**FW:** 371.3  
**Purity:** ≥98%  
**Supplied as:** A 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

7-hydroxy Chlorpromazine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 7-hydroxy chlorpromazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 7-hydroxy Chlorpromazine (hydrochloride) is slightly soluble in methanol.

### Description

7-hydroxy Chlorpromazine is an active metabolite of the typical antipsychotic chlorpromazine (Item No. 16129).<sup>1</sup> *In vivo*, 7-hydroxy chlorpromazine (5 mg/kg) increases plasma prolactin levels in male rats. It also reverses D-amphetamine-induced depression of the dopaminergic neuronal firing rate in the rat substantia nigra in rats when administered at a dose of 1.6 mg/kg.<sup>2</sup>

### References

1. Meltzer, H.Y., Fang, V.S., Simonovich, M., *et al.* Effect of metabolites of chlorpromazine on plasma prolactin levels in male rats. *Eur. J. Pharmacol.* **41(4)**, 431-436 (1977).
2. Bunney, B.S. and Aghajanian, G.K. A comparison of the effects of chlorpromazine, 7-hydroxychlorpromazine and chlorpromazine sulfoxide on the activity of central dopaminergic neurons. *Life Sci.* **15(2)**, 309-318 (1974).

#### 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, 10/24/2022

#### 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