

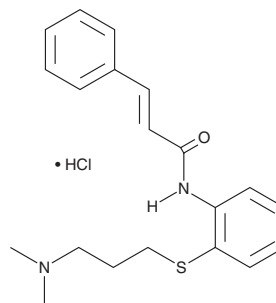
# PRODUCT INFORMATION



## Cinanserin (hydrochloride)

Item No. 30729

**CAS Registry No.:** 54-84-2  
**Formal Name:** N-[2-[[3-(dimethylamino)propyl]thio]phenyl]-3-phenyl-2-propenamamide, monohydrochloride  
**Synonym:** SQ 10,643  
**MF:** C<sub>20</sub>H<sub>24</sub>N<sub>2</sub>OS • HCl  
**FW:** 376.9  
**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

Cinanserin (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the cinanserin (hydrochloride) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Cinanserin is a serotonin (5-HT) receptor antagonist.<sup>1,2</sup> It inhibits 5-HT-induced effects on isolated rat uterus but not guinea pig ileum.<sup>1</sup> Cinanserin inhibits 5-HT-induced gross excitation, gastric mucosal erosion, and anaphylactoid edema in mice and rats. Cinanserin (12 and 36 mg/kg, i.p.) impairs acquisition of a running response in rats trained for food reward.<sup>2</sup> It inhibits severe acute respiratory syndrome coronavirus (SARS-CoV) 3C-like proteinase (3CL<sup>pro</sup>; IC<sub>50</sub> = 5 μM) and reduces SARS-CoV viral RNA levels in infected BHK-Rep-1 cells (IC<sub>50S</sub> = 19-34 μM).<sup>3</sup> It also inhibits viral replication of murine hepatitis virus (MHV) in CCL-9.1 murine liver epithelial cells (IC<sub>50</sub> = 31.25 μg/ml).<sup>4</sup>

### References

1. Rubin, B., Piala, J.J., Burke, J.C., *et al.* A new, potent, and specific serotonin inhibitor (SQ 10,643) 2'-(3-dimethylaminopropylthio) cinnamanilide hydrochloride: Antiserotonin activity on uterus and on gastrointestinal, vascular, and respiratory systems of animals. *Arch. Int. Pharmacodyn. Ther.* **152**, 132-143 (1964).
2. Rosen, A.J. and Cohen, M.E. The effects of cinanserin, a potent serotonin antagonist, on the acquisition of a running response in the rat. *Neuropharmacology* **12(6)**, 501-508 (1973).
3. Chen, L., Gui, C., Luo, X., *et al.* Cinanserin is an inhibitor of the 3C-like proteinase of severe acute respiratory syndrome coronavirus and strongly reduces virus replication in vitro. *J. Virol.* **79(11)**, 7095-7103 (2005).
4. Chiow, K.H., Phoon, M.C., Putti, T., *et al.* Evaluation of antiviral activities of *Houttuynia cordata* Thunb. extract, quercetin, quercetrin and cinanserin on murine coronavirus and dengue virus infection. *Asian Pac. J. Trop. Med.* **9(1)**, 1-7 (2016).

#### 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, 11/03/2022

#### 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