

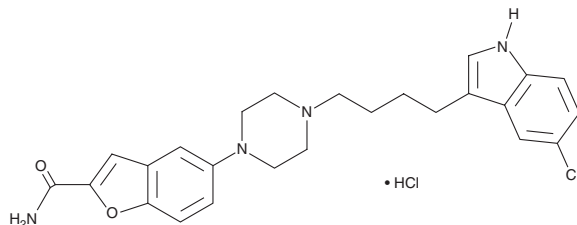
# PRODUCT INFORMATION



## Vilazodone (hydrochloride)

Item No. 21547

**CAS Registry No.:** 163521-08-2  
**Formal Name:** 5-[4-[4-(5-cyano-1H-indol-3-yl)butyl]-1-piperazinyl]-2-benzofurancarboxamide  
**Synonyms:** EMD 68843, SB 659746A  
**MF:** C<sub>26</sub>H<sub>27</sub>N<sub>5</sub>O<sub>2</sub> • HCl  
**FW:** 478.0  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 241, 271 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

Vilazodone (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the vilazodone (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Vilazodone (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of vilazodone (hydrochloride) in these solvents is approximately 30 and 5 mg/ml, respectively.

Vilazodone (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vilazodone (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Vilazodone (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Vilazodone is a selective serotonin reuptake inhibitor (SSRI) and a partial agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>1A</sub> (IC<sub>50</sub>s = 0.2 and 0.5 nM, respectively).<sup>1</sup> It increases extracellular 5-HT in the rat ventral hippocampus and frontal cortex when administered intraperitoneally at doses of 1 and 3 mg/kg. Vilazodone (1 mg/kg, i.p.) decreases immobility in the forced swim test in both rats and mice. Formulations containing vilazodone have been used in the treatment of depression.

### Reference

1. Page, M.E., Cryan, J.F., Sullivan, A., *et al.* Behavioral and neurochemical effects of 5-(4-[4-(5-Cyano-3-indolyl)-butyl]-butyl)-1-piperazinyl)-benzofuran-2-carboxamide (EMD 68843): A combined selective inhibitor of serotonin reuptake and 5-hydroxytryptamine<sub>1A</sub> receptor partial agonist. *J. Pharmacol. Exp. Ther.* **302**(3), 1220-1227 (2002).

#### 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, 10/04/2022

#### 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