

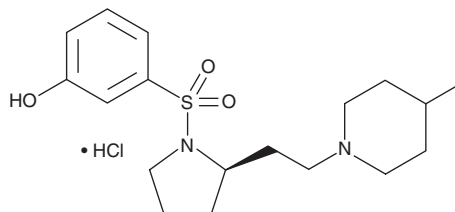
# PRODUCT INFORMATION



## SB-269970 (hydrochloride)

Item No. 17081

**CAS Registry No.:** 261901-57-9  
**Formal Name:** 3-[[[(2R)-2-[2-(4-methyl-1-piperidinyl)ethyl]-1-pyrrolidiny]sulfonyl]-phenol, monohydrochloride  
**MF:** C<sub>18</sub>H<sub>28</sub>N<sub>2</sub>O<sub>3</sub>S • HCl  
**FW:** 389.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 286 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

SB-269970 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SB-269970 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SB-269970 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SB-269970 (hydrochloride) in ethanol is approximately 5 mg/ml and approximately 50 mg/ml in DMSO and DMF.

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 SB-269970 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of SB-269970 (hydrochloride) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

The transduction of neurobehavioral effects by serotonin (5-hydroxy tryptamine; 5-HT) is mediated by at least seven major 5-HT receptor subtypes. SB-269970 is a potent 5-HT<sub>7A</sub> antagonist (pK<sub>i</sub> = 8.9) that demonstrates >50-fold binding selectivity over 5-HT<sub>5A</sub> and >250-fold selectivity over 5-HT<sub>1</sub>, 5-HT<sub>2</sub>, 5-HT<sub>4</sub>, 5-HT<sub>6</sub>, adrenergic α<sub>1</sub>, dopamine D<sub>2</sub>, and dopamine D<sub>3</sub> receptors.<sup>1</sup> It is also reported to block adrenergic α<sub>2</sub> receptors in guinea pig vas deferens.<sup>2</sup> SB-269970 has been used to target the 5-HT<sub>7</sub> receptor in the study of schizophrenia-like cognitive deficits.<sup>3</sup>

### References

1. Lovell, P.J., Bromidge, S.M., Dabbs, S., *et al.* A novel, potent, and selective 5-HT<sub>7</sub> antagonist: (R)-3-(2-(2-(4-Methylpiperidin-1-yl)ethyl)pyrrolidine-1-sulfonyl) phenol (SB-269970). *J. Med. Chem.* **43**(3), 342-345 (2000).
2. Foong, J.P.P. and Bornstein, J.C. 5-HT antagonists NAN-190 and SB 269970 block α<sub>2</sub>-adrenoceptors in the guinea pig. *NeuroReport* **20**(3), 325-330 (2009).
3. Nikiforuk, A., Kos, T., Fijal, K., *et al.* Effects of the selective 5-HT<sub>7</sub> receptor antagonist SB-269970 and amisulpride on ketamine-induced schizophrenia-like deficits in rats. *PLoS One* **8**(6), 1-12 (2013).

#### 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, 10/20/2022

#### 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