

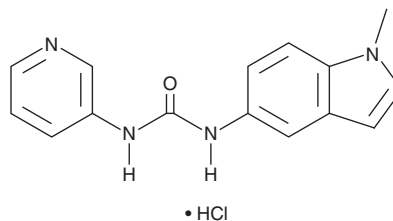
# PRODUCT INFORMATION



## SB-200646 (hydrochloride)

Item No. 30973

**CAS Registry No.:** 143797-62-0  
**Formal Name:** N-(1-methyl-1H-indol-5-yl)-N'-3-pyridinyl-urea, monohydrochloride  
**MF:** C<sub>15</sub>H<sub>14</sub>N<sub>4</sub>O • HCl  
**FW:** 302.8  
**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

SB-200646 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the SB-200646 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SB-200646 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

### Description

SB-200646 is a dual antagonist of the serotonin (5-HT) receptor subtypes 5-HT<sub>2C</sub> (K<sub>i</sub> = 125 nM in isolated rat cortex) and 5-HT<sub>2B</sub> (pA<sub>2</sub> = 7.5 in rat stomach fundus preparations).<sup>1</sup> It is selective for 5-HT<sub>2C</sub> and 5-HT<sub>2B</sub> over 5-HT<sub>1A</sub>, 5-HT<sub>2A</sub>, 5-HT<sub>1D</sub>, 5-HT<sub>3</sub>, benzodiazepine, GABA<sub>A</sub>, α<sub>1</sub>-, α<sub>2A</sub>-, α<sub>2B</sub>-, β<sub>1</sub>-, and β<sub>2</sub>-adrenergic, dopamine D<sub>1</sub> and D<sub>2</sub>, and histamine H<sub>1</sub> receptors (K<sub>i</sub>s = >10 μM for all). SB-200646 (20 and 40 mg/kg) reverses mCPP-induced hypophagia and decreases in social interaction time, indicating anxiolytic-like behavior, in rats.

### Reference

1. Kennett, G.A., Wood, M.D., Glen, A., *et al.* *In vivo* properties of SB 200646A, a 5-HT<sub>2C/2B</sub> receptor antagonist. *Br. J. Pharmacol.* **111(3)**, 797-802 (1994).

#### 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, 11/04/2022

#### 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