

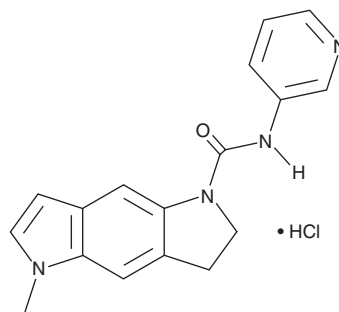
# PRODUCT INFORMATION



## SB-206553 (hydrochloride)

Item No. 34263

**CAS Registry No.:** 1197334-04-5  
**Formal Name:** 3,5-dihydro-5-methyl-N-3-pyridinyl-benzo[1,2-b:4,5-b']dipyrrole-1(2H)-carboxamide, monohydrochloride  
**MF:** C<sub>17</sub>H<sub>16</sub>N<sub>4</sub>O • HCl  
**FW:** 328.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 214, 236, 272, 325 nm  
**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-206553 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the SB-206553 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SB-206553 (hydrochloride) is soluble in the organic solvent DMSO.

### Description

SB-206553 is an antagonist of the serotonin (5-HT) receptor subtypes 5-HT<sub>2C</sub> (K<sub>i</sub> = 12.6 nM) and 5-HT<sub>2B</sub>.<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>1D</sub>, 5-HT<sub>1E</sub>, 5-HT<sub>1F</sub>, 5-HT<sub>2A</sub>, 5-HT<sub>3</sub>, 5-HT<sub>4</sub>, and 5-HT<sub>7</sub> receptors, as well as dopamine D<sub>2</sub>, D<sub>3</sub>, and D<sub>4</sub>, α<sub>1</sub>-adrenergic, histamine H<sub>1</sub>, and adenosine A<sub>1</sub> receptors (K<sub>i</sub>s = >1 μM for all). SB-206553 inhibits 5-HT-induced contraction of isolated rat stomach fundus, a tissue rich in 5-HT<sub>2B</sub> receptors, with an EC<sub>50</sub> value of 4.8 nM. It inhibits 5-HT-induced inositol phosphate release in HEK293 cells expressing human 5-HT<sub>2C</sub> receptors (EC<sub>50</sub> = 3.16 nM). SB-206553 (2 and 20 mg/kg) increases the time spent in active social interaction in a social interaction test and increases punished responding in the Geller-Seifter conflict test in rats, indicating anxiolytic-like activity. It also increases suppressed responding in a marmoset conflict model of anxiety when administered at doses of 15 and 20 mg/kg.

### Reference

1. Kennett, G.A., Wood, M.D., Bright, F., *et al.* *In vitro* and *in vivo* profile of SB 206553, a potent 5-HT<sub>2C</sub>/5-HT<sub>2B</sub> receptor antagonist with anxiolytic-like properties. *Br. J. Pharmacol.* **117**(3), 427-434 (1996).

#### 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, 12/02/2022

#### 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