

# PRODUCT INFORMATION



## SB-216641 (hydrochloride)

Item No. 32837

CAS Registry No.: 193611-67-5

Formal Name: N-[3-[2-(dimethylamino)ethoxy]-4-methoxyphenyl]-2'-methyl-4'-(5-methyl-1,2,4-oxadiazol-3-yl)-[1,1'-biphenyl]-4-carboxamide, monohydrochloride

MF: C<sub>28</sub>H<sub>30</sub>N<sub>4</sub>O<sub>4</sub> • HCl

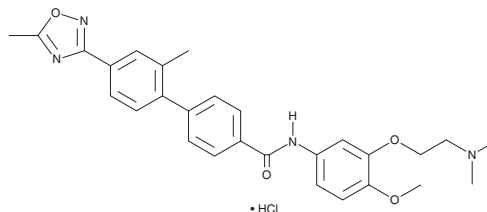
FW: 523.0

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-216641 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the SB-216641 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SB-216641 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

SB-216641 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT<sub>1B</sub>.<sup>1</sup> It binds to 5-HT<sub>1B</sub> receptors (K<sub>i</sub> = 1 nM) and is selective for 5-HT<sub>1B</sub> over 5-HT<sub>1A</sub>, 5-HT<sub>1D</sub>, 5-HT<sub>1E</sub>, and 5-HT<sub>1F</sub> receptors in radioligand binding assays (K<sub>i</sub>s = 500, 25, >1,000, and >1,000 nM, respectively). SB-216641 (1-1,000 nM) reduces 5-HT-induced inhibition of forskolin-induced cAMP production in CHO cells expressing human 5-HT<sub>1B</sub> receptors.

### Reference

1. Price, G.W., Burton, M.J., Collin, L.J., *et al.* SB-216641 and BRL-15572 – compounds to pharmacologically discriminate h5-HT1B and h5-HT1D receptors. *Naunyn Schmiedebergs Arch Pharmacol.* **356(3)**, 312-320 (1997).

#### 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