

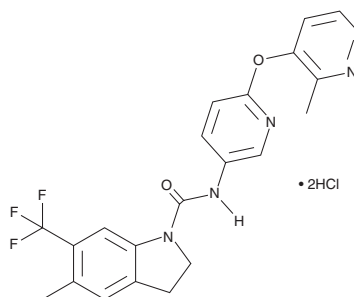
PRODUCT INFORMATION



SB-243213 (hydrochloride)

Item No. 36217

CAS Registry No.: 1780372-25-9
Formal Name: 2,3-dihydro-5-methyl-N-[6-[(2-methyl-3-pyridinyl)oxy]-3-pyridinyl]-6-(trifluoromethyl)-1H-indole-1-carboxamide, dihydrochloride
MF: $C_{22}H_{19}F_3N_4O_2 \cdot 2HCl$
FW: 501.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SB-243213 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the SB-243213 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SB-243213 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of SB-243213 (hydrochloride) in these solvents is approximately 100 and 50 mM, respectively.

Description

SB-243213 is an inverse agonist of the serotonin (5-HT) receptor subtype 5-HT_{2C}.¹ It is selective for 5-HT_{2C} over 5-HT_{2A} and 5-HT_{2B} receptors (K_i s = 0.001, 0.158, and 0.1 μM , respectively) and the cytochrome P450 (CYP) isoforms CYP1A2, CYP2C9, CYP2C19, CYP2D6, and CYP3A4 (IC_{50} s = >100, 23, >100, >100, and >100 μM , respectively). SB-243213 reduces the basal activity of 5-HT_{2C} in HEK293 cells expressing the human 5-HT_{2C} receptor (pK_B = 9.5). It reverses mCPP-induced hypolocomotion in rats (ID_{50} = 0.7 mg/kg). SB-243213 (0.2-5 mg/kg) has anxiolytic-like activity in the social interaction test in rats.

Reference

1. Bromidge, S.M., Dabbs, S., Davies, D.T., *et al.* Biarylcarbamoylindolines are novel and selective 5-HT_{2C} receptor inverse agonists: Identification of 5-methyl-1-[[2-[(2-methyl-3-pyridyl)oxy]-5-pyridyl]carbamoyl]-6-trifluoromethylindoline (SB-243213) as a potential antidepressant/anxiolytic agent. *J. Med. Chem.* **43**(6), 1123-1134 (2000).

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, 12/19/2022

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