

PRODUCT INFORMATION



S 32212 (hydrochloride)

Item No. 27639

CAS Registry No.: 847871-78-7
Formal Name: 1,2-dihydro-N-[4-methoxy-3-(4-methyl-1-piperazinyl)phenyl]-3H-benz[e]indole-3-carboxamide, monohydrochloride

MF: C₂₅H₂₈N₄O₂ • HCl

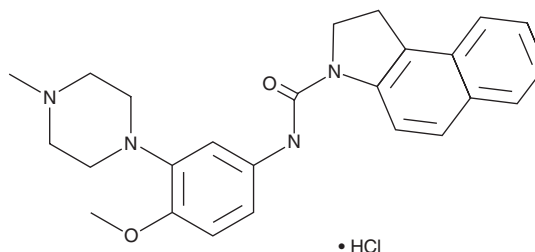
FW: 453.0

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

S 32212 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the S 32212 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. S 32212 (hydrochloride) is soluble in organic solvents such as DMSO. It is also soluble in water. The solubility of S 32212 (hydrochloride) in DMSO and water is approximately 25 mM. We do not recommend storing the aqueous solution for more than one day.

Description

S 32212 is an inverse agonist of the serotonin (5-HT) receptor subtypes 5-HT_{2CINI} and 5-HT_{2CVSV} (K_is = 6.6 and 8.9 nM, respectively).¹ It is also an antagonist of the 5-HT_{2A} receptor and the α_{2B}-adrenergic receptor (K_i = 5.8 nM for both). S 32212 is greater than 70-fold selective for these receptors in a panel of 80 receptors, enzymes, and ion channels. S 32212 reduces binding of GTPγS to Gα_q and decreases the activity of phospholipase C (PLC) in HEK293 cells expressing 5-HT_{2CINI} receptors (EC₅₀s = 38 and 18.6 nM, respectively) and in CHO cells expressing 5-HT_{2CVSV} receptors (EC₅₀s = 38 and 18.6 nM, respectively). S 32212 (2.5 mg/kg) decreases head twitching, penile erections, and drug discrimination induced by 5-HT receptor agonists in mice and rats. It reduces immobility time in the forced swim test and decreases marble burying in mice and rats when administered at doses of 10 and 40 mg/kg, indicating anti-depressant-like and anxiolytic activities.^{1,2}

References

1. Millan, M.J., Mannoury la Cour, C., Chanrion, B., *et al.* S32212, a novel serotonin type 2C receptor inverse agonist/α₂-adrenoceptor antagonist and potential antidepressant: I. A mechanistic characterization. *J. Pharmacol. Exp. Ther.* **340(3)**, 750-764 (2012).
2. Dekeyne, A., Brocco, M., Loiseau, F., *et al.* S32212, a novel serotonin type 2C receptor inverse agonist/α₂-adrenoceptor antagonist and potential antidepressant: II. A behavioral, neurochemical, and electrophysiological characterization. *J. Pharmacol. Exp. Ther.* **340(3)**, 765-780 (2012).

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, 12/22/2022

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