

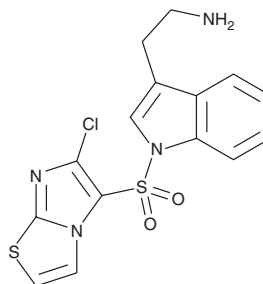
PRODUCT INFORMATION



WAY-181187

Item No. 23461

CAS Registry No.: 554403-49-5
Formal Name: 1-[(6-chloroimidazo[2,1-b]thiazol-5-yl)sulfonyl]-1H-indole-3-ethanamine
Synonym: SAX 187
MF: C₁₅H₁₃ClN₄O₂S₂
FW: 380.9
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

WAY-181187 is supplied as a solid. A stock solution may be made by dissolving the WAY-181187 in the solvent of choice, which should be purged with an inert gas. WAY-181187 is soluble (≥10 mg/ml) in DMSO.

Description

WAY-181187 is an agonist of the serotonin (5-HT) receptor subtype 5-HT₆.^{1,2} It increases intracellular cAMP levels in HeLa cells expressing 5-HT₆ (EC₅₀ = 6.5 nM).² WAY-181187 selectively binds to 5-HT₆ over 5-HT_{2B}, 5-HT_{2C}, and 5-HT₇ (K_is = 2,458, 124, and 679 nM, respectively). It is also selective for 5-HT₆ over a panel of 31 receptors and ion channels at 100 nM. It increases extracellular GABA levels in the rat frontal cortex when administered at a dose of 10 mg/kg. WAY-181187 (178 mg/kg) reduces schedule-induced polydipsia, an adjunctive model of compulsive disorder, in rats. It also reduces immobility in a modified forced swim test in rats when administered at a dose of 17 mg/kg, indicating antidepressant-like activity, and reduces the duration of defensive burying in rats at 10 and 17 mg/kg, indicating anxiolytic-like activity.³

References

1. Schechter, L.E., Lin, Q., Smith, D.L., *et al.* Neuropharmacological profile of novel and selective 5-HT₆ receptor agonists: WAY-181187 and WAY-208466. *Neuropsychopharmacology* **33**(6), 1323-1335 (2008).
2. Cole, D.C., Stock, J.R., Lennox, W.J., *et al.* Discovery of N1-(6-chloroimidazo[2,1-b][1,3]thiazole-5-sulfonyl)tryptamine as a potent, selective, and orally active 5-HT₆ receptor agonist. *J. Med. Chem.* **50**(23), 5535-5538 (2007).
3. Carr, G.V., Schechter, L.E., and Lucki, I. Antidepressant and anxiolytic effects of selective 5-HT₆ receptor agonists in rats. *Psychopharmacology (Berl)* **213**(2-3), 499-507 (2011).

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, 12/13/2024

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