

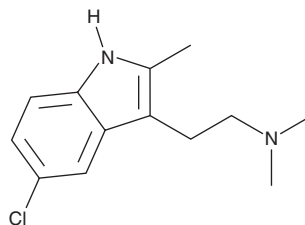
PRODUCT INFORMATION



ST1936

Item No. 39811

CAS Registry No.: 1210-81-7
Formal Name: 5-chloro-N,N,2-trimethyl-1H-indole-3-ethanamine
MF: C₁₃H₁₇ClN₂
FW: 236.7
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

ST1936 is supplied as a solid. A stock solution may be made by dissolving the ST1936 in the solvent of choice, which should be purged with an inert gas. ST1936 is soluble in the organic solvent acetonitrile.

Description

ST1936 is an agonist of the serotonin (5-HT) receptor subtype 5-HT₆.^{1,2} It selectively binds to 5-HT₆ (K_i = 31 nM) over 5-HT_{1A}, 5-HT_{1B}, 5-HT_{1D}, 5-HT_{2A}, 5-HT_{2B}, 5-HT_{2C}, 5-HT₃, 5-HT₄, 5-HT_{5A}, and the serotonin transporter (SERT; K_{iS} = >1,000 nM for all) but also binds to the 5-HT₇ receptor (K_i = 168 nM).¹ ST1936 induces cAMP accumulation in BHK fibroblasts (EC₅₀ = 16 nM).² It decreases the basal firing rate of dopaminergic neurons in the ventral tegmental area but not substantia nigra pars compacta in rats in a dose-dependent manner.³ ST1936 (10 and 30 mg/kg) prevents stress-induced decreases in vanilla sugar-sustained appetitive behavior, indicating anti-anhedonic-like activity, in rats.⁴ It induces self-administration in rats when administered at a dose of 1 mg/kg.⁵

References

1. Di Cesare, M.A., Minetti, P., Tarzia, G., *et al.* 5-halo-tryptamine derivatives used as ligands of the 5-HT₆ and/or 5-HT₇ serotonin receptors. **WO 03/000252 A1** (2003).
2. Mattsson, C., Sonesson, C., Sandahl, A., *et al.* 2-Alkyl-3-(1,2,3,6-tetrahydropyridin-4-yl)-1H-indoles as novel 5-HT₆ receptor agonists. *Bioorg. Med. Chem. Lett.* **15(19)**, 4230-4234 (2005).
3. Borsini, F., Bordi, F., Poggi, A., *et al.* Effects of ST1936, a selective serotonin-6 agonist, on electrical activity of putative mesencephalic dopaminergic neurons in the rat brain. *J. Psychopharmacol.* **29(7)**, 802-811 (2015).
4. Scheggi, S., Marchese, G., Borsini, F., *et al.* Effects of the 5-HT₆ receptor agonist ST 1936 on depression- and anhedonia-like experimental models. *Behav. Brain Res.* **224(1)**, 35-43 (2011).
5. Valentini, V., Piras, G., De Luca, M.A., *et al.* Evidence for a role of a dopamine/5-HT₆ receptor interaction in cocaine reinforcement. *Neuropharmacology* **65**, 58-64 (2013).

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, 11/15/2023

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