

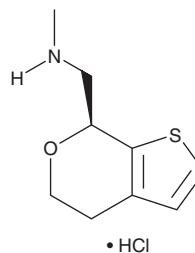
PRODUCT INFORMATION



SEP-363856 (hydrochloride)

Item No. 37399

CAS Registry No.: 1310422-41-3
Formal Name: 4,7-dihydro-N-methyl-5H-thieno[2,3-c]pyran-7-methanamine, monohydrochloride
Synonym: SEP-856
MF: C₉H₁₃NOS • HCl
FW: 219.7
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

SEP-363856 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the SEP-363856 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SEP-363856 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of SEP-363856 (hydrochloride) in these solvents is approximately 0.2 and 2 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of SEP-363856 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of SEP-363856 (hydrochloride) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

SEP-363856 is an agonist of trace amine-associated receptor 1 (TAAR1).¹ It increases intracellular cAMP levels in cells expressing TAAR1 (EC₅₀ = 0.14 μM). It is also an agonist of the serotonin (5-HT) receptor subtypes 5-HT_{1D} and 5-HT_{1A}, decreasing cAMP levels in cells expressing these receptors (EC₅₀s = 0.262 and 2.3 μM, respectively). SEP-363856 is selective for these receptors over the dopamine D_{2L} receptor (EC₅₀s = 10.44 and 8.02 μM in cAMP and arrestin recruitment assays, respectively), as well as α_{2A}- and α_{2B}-adrenergic, and 5-HT_{1B}, 5-HT_{2A}, 5-HT_{2C}, and 5-HT₇ receptors (EC₅₀s = 6.7->10 μM in functional assays). It reduces hyperactivity induced by phencyclidine (PCP) in mice when administered at 0.3, 1, and 3 mg/kg and reverses PCP-induced impairments in social interaction in rats. SEP-363856 (3, 10, and 30 mg/kg) increases prepulse inhibition of the acoustic startle response in mice. It also decreases the time spent immobile in the forced swim test in mice, indicating antidepressant-like activity.

Reference

1. Dedic, N., Jones, P.G., Hopkins, S.C., *et al.* SEP-363856, a novel psychotropic agent with a unique, non-D₂ receptor mechanism of action. *J. Pharmacol. Exp. Ther.* **371**(1), 1-14 (2019).

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, 04/14/2023

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