PRODUCT INFORMATION



Vortioxetine (hydrobromide)

Item No. 23694

| CAS Registry No.: | 960203-27-4 | Н |
|-------------------|---|-------------|
| Formal Name: | 1-[2-[(2,4-dimethylphenyl)thio]phenyl]- | |
| | piperazine, monohydrobromide | |
| MF: | $C_{18}H_{22}N_2S \bullet HBr$ | |
| FW: | 379.4 | N |
| Purity: | ≥95% | |
| UV/Vis.: | λ _{max} : 227 nm | S S |
| Supplied as: | A solid | |
| Storage: | 4°C | |
| Stability: | ≥4 years | ↔ • HBr ↔ ৲ |
| | | |

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

Laboratory Procedures

Vortioxetine (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the vortioxetine (hydrobromide) in the solvent of choice, which should be purged with an inert gas. Vortioxetine (hydrobromide) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of vortioxetine (hydrobromide) in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Vortioxetine (hydrobromide) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vortioxetine (hydrobromide) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Vortioxetine (hydrobromide) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Vortioxetine is a multimodal serotonergic agent that acts as a serotonin (5-HT; Item No. 14332) reuptake inhibitor (K_i = 1.6 nM) and binds to the 5-HT_{1A}, 5-HT_{1B}, 5-HT_{3A}, and 5-HT₇ receptors (K_is = 15, 33, 3.7, and 19 nM, respectively).¹ It also binds to the β_1 -adrenergic receptor (K_i = 46 nM). It acts as an antagonist at 5-HT_{3A} and 5-HT₇, a partial agonist at 5-HT_{1B}, and a full agonist at 5-HT_{1A} receptors. Vortioxetine is selective for the serotonin transporter (SERT) over the dopamine and norepinephrine transporters (IC₅₀s = 5.3, 890, and 140 nM, respectively). It has affinity for histamine H₂, melanocortin 4, β_2 -adrenergic, 5-HT_{2C}, and 5-HT₆ receptors with K_i values greater than 180 nM, but it is selective over the majority of targets in a panel of 75 ion channels, G protein-coupled receptors, enzymes, and transporters. Vortioxetine (5 or 10 mg/kg per day) increases 5-HT levels in vivo in rat brain and decreases occupancy at SERT. At acute and chronic doses of 5 mg/kg, but not 10 mg/kg, it increases the time mice spend in the center of the open field test, increases mobility in the forced swim test, and decreases the latency to feed in the novelty suppressed feeding test.² It also increases proliferation and survival of adult-born granule cells in the hippocampus. Formulations containing vortioxetine have been used in the treatment of major depressive disorder.

References

- 1. Bang-Andersen, B., Ruhland, T., Jørgensen, M. et al. Discovery of 1-[2-(2,4-dimethylphenylsulfanyl) phenyl]piperazine (Lu AA21004): A novel multimodal compound for the treatment of major depressive disorder. J. Med Chem. 54(9), 3206-3221 (2011).
- 2. Guilloux, J.P., Mendez-David, I., Pehrson, A. et al. Antidepressant and anxiolytic potential of the multimodal antidepressant vortioxetine (Lu AA21004) assessed by behavioural and neurogenesis outcomes in mice. Neuropharmacology 73, 147-159 (2013).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

subject to Cavman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information uyer agrees to purchase the material can be found on our website

Copyright Cayman Chemical Company, 01/27/2025

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