

# PRODUCT INFORMATION

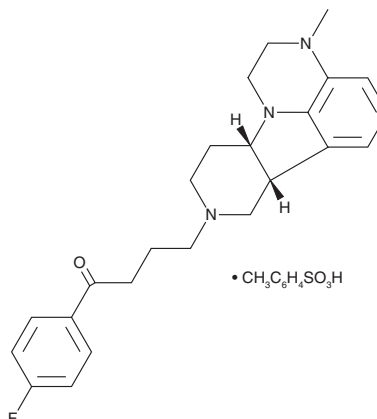


## Lumateperone (tosylate)

Item No. 34541

**CAS Registry No.:** 1187020-80-9  
**Formal Name:** 1-(4-fluorophenyl)-4-((6bR,10aS)-3-methyl-2,3,6b,9,10,10a-hexahydro-1H-pyrido[3',4':4,5]pyrrolo[1,2,3-de]quinoxalin-8(7H)-yl)butan-1-one 4-methylbenzenesulfonic acid

**Synonym:** ITI-007  
**MF:** C<sub>24</sub>H<sub>28</sub>FN<sub>3</sub>O • C<sub>7</sub>H<sub>8</sub>O<sub>3</sub>S  
**FW:** 565.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229 nm  
**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

Lumateperone (tosylate) is supplied as a solid. A stock solution may be made by dissolving the lumateperone (tosylate) in the solvent of choice, which should be purged with an inert gas. Lumateperone (tosylate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of lumateperone (tosylate) in these solvents is approximately 1, 10, and 25 mg/ml, respectively.

Lumateperone (tosylate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, lumateperone (tosylate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Lumateperone (tosylate) has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Lumateperone is an atypical antipsychotic.<sup>1</sup> It selectively binds to the serotonin (5-HT) receptor subtype 5-HT<sub>2A</sub> over 5-HT<sub>1A</sub> and 5-HT<sub>2C</sub> receptors (K<sub>i</sub>s = 0.54, 1,480, and 173 nM, respectively). Lumateperone also binds to dopamine D<sub>1</sub>, D<sub>2</sub>, and D<sub>4</sub> receptors (K<sub>i</sub>s = 52, 32, and 108 nM, respectively), α<sub>1A</sub>- and α<sub>1B</sub>-adrenergic receptors (K<sub>i</sub>s = 73 and 31 nM, respectively), and the 5-HT transporter (SERT; K<sub>i</sub> = 61 nM). It reduces head-twitch behavior induced by the 5-HT<sub>2A</sub> agonist quipazine in rats (ED<sub>50</sub> = 0.12 mg/kg). Formulations containing lumateperone have been used in the treatment of schizophrenia.

### Reference

1. Li, P., Zhang, Q., Robichaud, A.J., et al. Discovery of a tetracyclic quinoxaline derivative as a potent and orally active multifunctional drug candidate for the treatment of neuropsychiatric and neurological disorders. *J. Med. Chem.* **57**(6), 2670-2682 (2014).

**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, 10/04/2022

### 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