

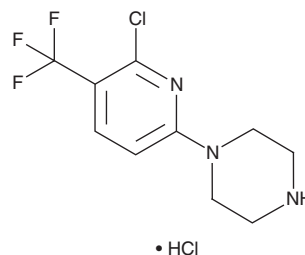
# PRODUCT INFORMATION



## ORG 12962 (hydrochloride)

Item No. 27675

**CAS Registry No.:** 210821-63-9  
**Formal Name:** 1-[6-chloro-5-(trifluoromethyl)-2-pyridinyl]-piperazine, monohydrochloride  
**Synonym:** Ro 44-6956  
**MF:** C<sub>10</sub>H<sub>11</sub>ClF<sub>3</sub>N<sub>3</sub> • HCl  
**FW:** 302.1  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

ORG 12962 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ORG 12962 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ORG 12962 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of ORG 12962 (hydrochloride) in DMSO and water is approximately 100 and 50 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

### Description

ORG 12962 is an agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>2C</sub> with an EC<sub>50</sub> value of 97.72 nM in a FLIPR-based intracellular calcium assay using CHO-K1 cells expressing the human receptor.<sup>1</sup> It is selective for 5-HT<sub>2C</sub> over 5-HT<sub>2A</sub> and 5-HT<sub>2B</sub> receptors (EC<sub>50</sub>s = 419.87 and 524.81 nM, respectively). ORG 12962 (1 and 3.2 mg/kg) has anti-aversive effects in a rat model of panic anxiety induced by dorsolateral periaqueductal gray matter (dPAG) stimulation.<sup>2</sup>

### References

1. Porter, R.H.P., Benwell, K.R., Lamb, H., *et al.* Functional characterization of agonists at recombinant human 5-HT<sub>2A</sub>, 5-HT<sub>2B</sub> and 5-HT<sub>2C</sub> receptors in CHO-K1 cells. *Br. J. Pharmacol.* **128**(1), 13-20 (1999).
2. Jenck, F., Moreau, J.-L., Berendsen, H.H.G., *et al.* Antiaversive effects of 5HT<sub>2C</sub> receptor agonists and fluoxetine in a model of panic-like anxiety in rats. *Eur. Neuropsychopharmacol.* **8**(3), 161-168 (1998).

#### 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, 02/10/2020

#### 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](http://WWW.CAYMANCHEM.COM)