

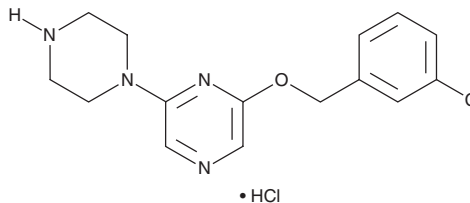
# PRODUCT INFORMATION



## CP 809,101 (hydrochloride)

Item No. 33283

**CAS Registry No.:** 1215721-40-6  
**Formal Name:** 2-[(3-chlorophenyl)methoxy]-6-(1-piperazinyl)-pyrazine, monohydrochloride  
**MF:** C<sub>15</sub>H<sub>17</sub>ClN<sub>4</sub>O • HCl  
**FW:** 341.2  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 248, 326 nm  
**Supplied as:** A crystalline 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

CP 809,101 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 809,101 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CP 809,101 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of CP 809,101 (hydrochloride) in DMSO and water is approximately 100 and 20 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

### Description

CP 809,101 is a potent agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>2C</sub>.<sup>1</sup> It is selective for human 5-HT<sub>2C</sub> and rat 5-HT<sub>2C</sub> over human 5-HT<sub>2A</sub>, 5-HT<sub>2B</sub>, and rat 5-HT<sub>2A</sub> receptors in a FLIPR membrane potential assay (EC<sub>50</sub>s = 0.11, 0.06, 153, 65.3, and 119 nM, respectively) as well as human dopamine D<sub>2</sub>, histamine H<sub>1</sub>, and α<sub>1</sub>- and α<sub>2</sub>-adrenergic receptors (K<sub>s</sub> = 872, 828, 217, and 956 nM, respectively). CP 809,101 inhibits the conditioned avoidance response, hyperactivity induced by PCP and D-amphetamine, and spontaneous locomotor activity in rats (ED<sub>50</sub>s = 4.8, 2.4, 2.9, and 2 mg/kg, respectively). It reduces prepulse inhibition deficits induced by apomorphine and improves novel object recognition in mice. CP 809,101 (0.01-1 μg/0.2 μl/site) reduces reinstatement of drug-seeking behavior in cocaine-primed rats when administered into the central amygdala but not the basolateral amygdala.<sup>2</sup> It also reduces reinstatement of food-seeking behavior in a rat model of dietary relapse.<sup>3</sup>

### References

1. Siuciak, J.A., Chapin, D.S., McCarthy, S.A., *et al.* CP-809,101, a selective 5-HT<sub>2C</sub> agonist, shows activity in animal models of antipsychotic activity. *Neuropharmacology* **52(2)**, 279-290 (2007).
2. Pockros-Burgess, L.A., Pentkowski, N.S., Der-Ghazarian, T., *et al.* Effects of the 5-HT<sub>2C</sub> receptor agonist CP809101 in the amygdala on reinstatement of cocaine-seeking behavior and anxiety-like behavior. *Int. J. Neuropsychopharmacol.* **17(11)**, 1751-1762 (2014).
3. Higgins, G.A., Silenieks, L.B., Altherr, E.B., *et al.* Lorcaserin and CP-809101 reduce motor impulsivity and reinstatement of food seeking behavior in male rats: Implications for understanding the anti-obesity property of 5-HT<sub>2C</sub> receptor agonists. *Psychopharmacology (Berl)* **233(14)**, 2841-2856 (2016).

#### 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, 11/17/2022

#### 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