

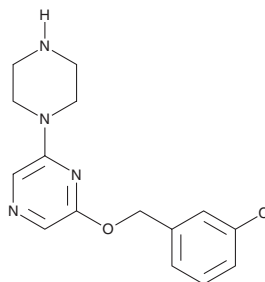
PRODUCT INFORMATION



CP 809,101

Item No. 24214

CAS Registry No.: 479683-64-2
Formal Name: 2-[(3-chlorophenyl)methoxy]-6-(1-piperazinyl)-pyrazine
MF: C₁₅H₁₇ClN₄O
FW: 304.8
Purity: ≥98%
UV/Vis.: λ_{max}: 249, 329 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 is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 809,101 in the solvent of choice. CP 809,101 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CP 809,101 in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

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 CP 809,101 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CP 809,101 in PBS, pH 7.2, is approximately 0.1 mg/ml. 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_{2C}.¹ It is selective for human 5-HT_{2C} and rat 5-HT_{2C} over human 5-HT_{2A}, 5-HT_{2B}, and rat 5-HT_{2A} receptors in a FLIPR membrane potential assay (EC₅₀s = 0.11, 0.06, 153, 65.3, and 119 nM, respectively) as well as human dopamine D₂, histamine H₁, and α₁- and α₂-adrenergic receptors (K_is = 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 (ED₅₀s = 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/side) reduces reinstatement of drug-seeking behavior in cocaine-primed rats when administered into the central amygdala but not the basolateral amygdala.² It also reduces reinstatement of food-seeking behavior in a rat model of dietary relapse.³

References

1. Xiang, Z., Pogozheva, I.D., Sorenson, N.B., *et al. Biochemistry* **46**(28), 8273-8287 (2007).
2. Hamberg, M. and Gotthammar, B. *Lipids* **8**(12), 733-744 (1973).
3. Kaminsky, Y.G., Marlatt, M.W., Smith, M.A., *et al. Exp. Neurol.* **221**(1), 26-37 (2010).

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, 10/18/2022

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