

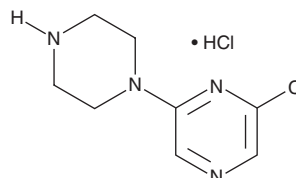
PRODUCT INFORMATION



MK-212 (hydrochloride)

Item No. 36089

CAS Registry No.: 61655-58-1
Formal Name: 2-chloro-6-(1-piperazinyl)-pyrazine,
monohydrochloride
MF: $C_8H_{11}ClN_4 \cdot HCl$
FW: 235.1
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 250, 329 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

MK-212 (hydrochloride) is supplied as a solid. Aqueous solutions of MK-212 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. MK-212 (hydrochloride) is slightly soluble in PBS (pH 7.2) and water. We do not recommend storing the aqueous solution for more than one day.

Description

MK-212 is an agonist of the serotonin (5-HT) receptor subtype 5-HT_{2C} ($EC_{50} = 0.028 \mu M$ in a calcium mobilization assay).¹ It is selective for 5-HT_{2C} over 5-HT_{2A} receptors ($EC_{50} = 0.42 \mu M$). MK-212 reduces food intake in fasted rats with an ED_{50} value of 5 mg/kg.² It reduces cocaine-primed reinstatement of cocaine-seeking behavior in rats when administered at doses of 0.32, 0.56, or 1 mg/kg.³

References

1. Jensen, A.A., Plath, N., Pedersen, M.H.F., *et al.* Design, synthesis, and pharmacological characterization of N- and O-substituted 5,6,7,8-tetrahydro-4H-isoxazolo[4,5-d]azepin-3-ol analogues: Novel 5-HT_{2A}/5-HT_{2C} receptor agonists with pro-cognitive properties. *J. Med. Chem.* **56**(3), 1211-1227 (2013).
2. J.C., H., Lawton, C.L., and Blundell, J.E. The 5-HT₂ receptor agonist MK-212 reduces food intake and increases resting but prevents the behavioural satiety sequence. *Pharmacol. Biochem. Behav.* **56**(1), 41-46 (1997).
3. Neisewander, J.L. and Acosta, J.I. Stimulation of 5-HT_{2C} receptors attenuates cue and cocaine-primed reinstatement of cocaine-seeking behavior in rats. *Behav. Pharmacol.* **18**(8), 791-800 (2007).

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, 12/19/2022

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