

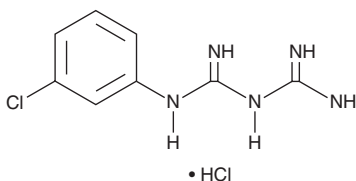
PRODUCT INFORMATION



1-(3-Chlorophenyl)biguanide (hydrochloride)

Item No. 34726

CAS Registry No.: 2113-05-5
Formal Name: N-(3-chlorophenyl)-imidodicyanimidic diamide, monohydrochloride
Synonyms: *m*-Chlorophenylbiguanide, *meta*-Chlorophenylbiguanide, *m*-CPBG, *meta*-CPBG
MF: C₈H₁₀ClN₅ • HCl
FW: 248.1
Purity: ≥95%
UV/Vis.: λ_{max}: 257 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

1-(3-Chlorophenyl)biguanide (*m*-CPBG) (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the *m*-CPBG in the solvent of choice, which should be purged with an inert gas. *m*-CPBG (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of *m*-CPBG (hydrochloride) in these solvents is approximately 5, 20, and 30 mg/ml, respectively.

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 *m*-CPBG (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of *m*-CPBG (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

m-CPBG is an agonist of the serotonin (5-HT) receptor subtype 5-HT₃.¹ It selectively binds 5-HT₃ over 5-HT_{1A} and 5-HT₂ receptors (K_s = 0.002, 10, and 10 μM, respectively) but also binds to high and low affinity sites on the dopamine transporter (DAT; IC₅₀s = 0.4 and 34.8 μM, respectively, in rat caudate putamen synaptosomal membranes).^{2,3} *m*-CPBG induces depolarization of isolated rat vagus nerve and stimulates inositol phosphate formation in rat frontocingulate cortical slices (EC₅₀s = 0.05 and 4.2 μM, respectively).^{1,4} It induces bradycardia, an effect that can be reversed by the 5-HT₃ receptor antagonist ondansetron, in anaesthetized cats (ED₅₀ = 20.3 nmol/kg).¹

References

1. Kilpatrick, G.J., Butler, A., Burrige, J., *et al.* 1-(*m*-chlorophenyl)-biguanide, a potent high affinity 5-HT₃ receptor agonist. *Eur. J. Pharmacol.* **182(1)**, 193-197 (1990).
2. Campbell, A.D., Womer, D.E., and Simon, J.R. The 5-HT₃ receptor agonist 1-(*m*-chlorophenyl)-biguanide interacts with the dopamine transporter in rat brain synaptosomes. *Eur. J. Pharmacol.* **290(2)**, 157-162 (1995).
3. Higgins, G.A., Joharchi, N., and Sellers, E.M. Behavioral effects of the 5-hydroxytryptamine₃ receptor agonists 1-phenylbiguanide and *m*-chlorophenylbiguanide in rats. *J. Pharmacol. Exp. Ther.* **264(3)**, 1440-1449 (1993).
4. Edwards, E., Hampton, E., Ashby, C.R., *et al.* 5-HT₃-like receptors in the rat medial prefrontal cortex: Further pharmacological characterization. *Brain Res.* **733(1)**, 21-30 (1996).

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, 10/17/2022

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