

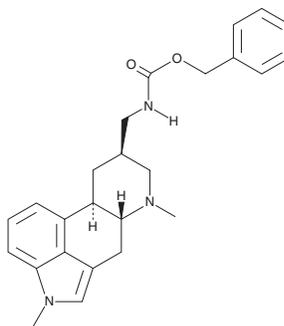
PRODUCT INFORMATION



Metergoline

Item No. 21336

CAS Registry No.: 17692-51-2
Formal Name: N-[[[(8β)-1,6-dimethylergolin-8-yl]methyl]-carbamic acid, phenylmethyl ester
MF: C₂₅H₂₉N₃O₂
FW: 403.5
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 289 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

Metergoline is supplied as a crystalline solid. A stock solution may be made by dissolving the metergoline in the solvent of choice. Metergoline is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of metergoline in these solvents is approximately 5, 25, and 30 mg/ml, respectively.

Metergoline is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, metergoline should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Metergoline has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Metergoline is a serotonin (5-HT) receptor antagonist that is active at 5-HT₁ (pK_i = 7.8 for 5-HT_{1B}), 5-HT₂, (pK_is= 8.64, 8.75, and 8.75 for 5-HT_{2A}, 5-HT_{2B}, and 5-HT_{2C}, respectively), and 5-HT₇ receptor subtypes.^{1,2} It blocks 5-HT₇ receptor-stimulated cAMP production (IC₅₀ = 321 nM) in HEK293 cells.³ Metergoline has been widely used to study serotonin receptor function, but it also dose-dependently inhibits voltage-dependent sodium and potassium channels transfected into *Xenopus* oocytes.^{4,5} Formulations containing metergoline have been studied for the treatment of hormone dysfunction, seasonal affective disorder, and anxiety.⁴

References

1. Millan, M.J., Newman-Tancredi, A., Lochon, S., *et al.* Specific labelling of serotonin 5-HT_{1B} receptors in rat frontal cortex with the novel, phenylpiperazine derivative, [³H]GR125,743. A pharmacological characterization. *Pharmacol. Biochem. Behav.* **71(4)**, 589-598 (2002).
2. Knight, A.R., Misra, A., Quirk, K., *et al.* Pharmacological characterisation of the agonist radioligand binding site of 5-HT_{2A}, 5-HT_{2B} and 5-HT_{2C} receptors. *Naunyn Schmiedeberg's Arch. Pharmacol.* **370(2)**, 114-123 (2004).
3. Knight, J.A., Smith, C., Toohey, N., *et al.* Pharmacological analysis of the novel, rapid, and potent inactivation of the human 5-Hydroxytryptamine₇ receptor by risperidone, 9-OH-Risperidone, and other inactivating antagonists. *Mol. Pharmacol.* **75(2)**, 374-380 (2009).
4. Yeom, H.D., and Lee, J.-H. Regulation of human Kv1.4 channel activity by the antidepressant metergoline. *Biol. Pharm. Bull.* **39(6)**, 1069-1072 (2016).
5. Lee, J.-H., Liu, J., Shin, M., *et al.* Metergoline inhibits the neuronal Nav1.2 voltage-dependent Na⁺ channels expressed in *Xenopus* oocytes. *Acta. Pharmacol. Sin.* **35(7)**, 862-868 (2014).

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, 12/22/2022

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