

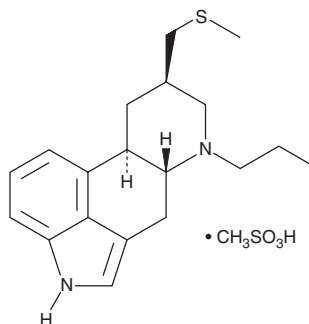
PRODUCT INFORMATION



Pergolide (mesylate)

Item No. 26085

CAS Registry No.: 66104-23-2
Formal Name: 8β-[(methylthio)methyl]-6-propyl-ergoline, monomethanesulfonate
Synonyms: LY127809, NSC 319773
MF: C₁₉H₂₆N₂S • CH₃SO₃H
FW: 410.6
Purity: ≥95%
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

Pergolide (mesylate) is supplied as a solid. A stock solution may be made by dissolving the pergolide (mesylate) in the solvent of choice, which should be purged with an inert gas. Pergolide (mesylate) is slightly soluble in methanol and DMSO.

Description

Pergolide is a potent dopamine D₁ and D₂ receptor agonist (K_is = 111 and 0.495 nM, respectively, for rat striatal receptors).¹ It depresses dopaminergic firing in paralyzed rats (ED₅₀ = <20 µg/kg), an effect that can be reversed by the dopamine D₂-selective antagonist spiperone or the dopamine D₁-selective antagonist SCH 23390 (Item No. 15631).² Pergolide (0.025 and 0.05 mg/kg) increases the volume threshold for inducing bladder contraction in a cynomolgus monkey model of Parkinson's disease induced by MPTP.³ It also reduces carrageenan-induced paw edema in adrenalectomized rats (ED₅₀ = 0.4 mg/kg).⁴

References

1. Miyagi, M., Itoh, F., Taya, F., *et al.* Dopamine receptor affinities *in vitro* and stereotypic activities *in vivo* of cabergoline in rats. *Biol. Pharm. Bull.* **19(9)**, 1210-1213 (1996).
2. Zhang, X.X., Jin, G.Z., and Wei, Y.F. Agonistic actions of pergolide on firing activity of dopamine neurons in substantia nigra compacta area. *Zhongguo Yao Li Xue Bao.* **16(5)**, 423-427 (1995).
3. Yoshimura, N., Mizuta, E., Yoshida, O., *et al.* Therapeutic effects of dopamine D₁/D₂ receptor agonists on detrusor hyperreflexia in 1-methyl-4-phenyl-1,2,3,6-tetrahydropyridine-lesioned parkinsonian cynomolgus monkeys. *J. Pharmacol. Exp. Ther.* **286(1)**, 228-233 (1998).
4. Bendele, A.M., Spaethe, S.M., Benslay, D.N., *et al.* Anti-inflammatory activity of pergolide, a dopamine receptor agonist. *J. Pharmacol. Exp. Ther.* **259(1)**, 169-175 (1991).

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