

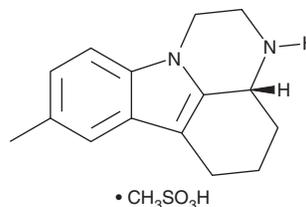
PRODUCT INFORMATION



(R)-(-)-Pirlindole (mesylate)

Item No. 43879

CAS Registry No.: 1821379-62-7
Formal Name: 2,3,3aR,4,5,6-hexahydro-8-methyl-1H-pyrazino[3,2,1-jk]carbazole, methanesulfonate
Synonym: (R)-Pirlindole
MF: C₁₅H₁₈N₂ • CH₃SO₃H
FW: 322.4
Purity: ≥98%
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

(R)-(-)-Pirlindole (mesylate) is supplied as a solid. A stock solution may be made by dissolving the (R)-(-)-pirlindole (mesylate) in the solvent of choice, which should be purged with an inert gas. (R)-(-)-Pirlindole (mesylate) is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

(R)-(-)-Pirlindole is an inhibitor of monoamine oxidase A (MAO-A; IC₅₀ = 0.43 μM).¹ It is selective for MAO-A over MAO-B. (R)-(-)-Pirlindole (50 mg/kg) decreases immobility time in the forced swim test in mice. It also inhibits reserpine-induced palpebral ptosis in mice when administered at a dose of 25 mg/kg.

Reference

1. Bruhwyler, J., Liégeois, J.F., Gérardy, J., *et al.* Comparative study of pirlindole, a selective RIMA, and its two enantiomers using biochemical and behavioural techniques. *Behav. Pharmacol.* **9(8)**, 731-737 (1998).

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, 06/04/2025

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