
Pargyline

(Catalog No. M55001)

Alternative Names

N-Methyl-N-(2-propynyl)benzylamine hydrochloride N-Methyl-N-propargylbenzylamine hydrochloride

Description

Pargyline is a mechanism-based inhibitor of LSD1 that effectively inhibits histone demethylation in vivo. Pargyline also selectively inhibits MAO-B.

CAS Number

306-07-0

Chemical Formula

C₁₁H₁₃N •HCl

Molecular Weight

195.69

Size

500 mg

Purity

>98%

Solubility

Soluble in water, DMSO, ethanol

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Pargyline	500 mg	M55001 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.