
Trans-2-Phenylcyclopropylamine hydrochloride

(Catalog No. M55000)

Alternative Names

5-[(2-hydroxy-5,8-dihydronaphthalen-1-yl)methyl]-6-phenyl-2-thioxo-2,3-dihy.

Alternative Names

2-PCPA

Description

trans-2-Phenylcyclopropylamine Tranylcyromine is an irreversible, mechanism-based inhibitor of LSD1 that effectively inhibits histone demethylation in vivo. Although not as selective, trans-2-Phenylcyclopropylamine Tranylcyromine also irreversibly inhibits MAO A and MAO B.

CAS Number

1986-47-6

Chemical Formula

C₉H₁₂ClN

Molecular Weight

169.7

Size

250 mg

Purity

>98%

Solubility

Water, DMSO

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Trans-2-Phenylcyclopropylamine hydrochloride	250 mg	M55000 -1

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