

Tranlycypromine (2-PCPA), hemisulfate salt

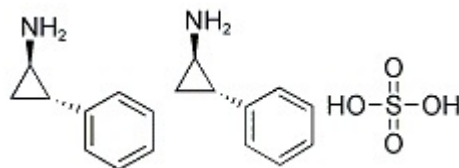
Catalog #: 27305

Lot #: 160429

Size: 100 mg

Structure:

CAS Registry #: 13492-01-8



Purity: >98%

Chemical Formula: C₁₈H₂₂N₂ · H₂SO₄

Molecular Weight: 364.46

Description: Tranlycypromine (2-PCPA) is an irreversible inhibitor of LSD1 histone demethylase and monoamine oxidases (MAO). The inhibitor targets monoamine oxidases A and B and has been used as a chemical scaffold to design new demethylase inhibitors. Tranlycypromine has been shown to increase Histone H3 lysine 4 (H3K4) methylation in P19 embryonal carcinoma cell.

Appearance: A white to off-white powder

Solubility: Soluble in water (25 mg/ml).

Biological Activity: Tranlycypromine inhibits LSD1 with an IC₅₀ value of 20.7 μM and a K_i value of 242.7 μM. It also irreversibly inhibits MAO A and MAO B with IC₅₀ values of 2.3 and 0.95 μM and K_i values of 101.9 and 16 μM, respectively.

Storage/Stability: Store solid form at +4°C. Hygroscopic. Solid form is stable at least 12 months from date of receipt, when stored as directed. Aqueous solutions are less stable; we recommend preparing and using stock solutions the same day.

Warnings: Hygroscopic. Protect from moisture. Protect from light.

Quality Control: The purity was determined by HPLC analysis.

Reference(s):

1. Schmidt, D.M.Z., and McCafferty, D.G. *Biochemistry* **46**:4408-4416 (2007).
2. Yang, M., *et al.* *Biochemistry* **46**:8058-65(2007).