

Tranlycypromine hemisulfate

Catalog No: 14046, 14047

Format: 250 mg, 50 mg

Chemical Properties:

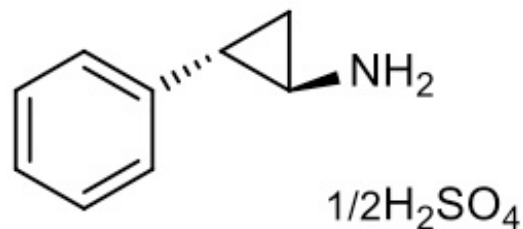
MW = 182.2

$C_9H_{11}N \cdot 1/2 H_2SO_4$

CAS 13492-01-8

Physical Properties: White solid

Names: Tranlycypromine hemisulfate, 2-PCPA, trans-2-phenylcyclopropanamine



Pharmacology: Irreversible inhibitor of lysine-specific demethylase 1 (LSD1/BHC110) (IC₅₀ less than 2 μM) (ref 1,2). Tranlycypromine treatment of P19 embryonal carcinoma cells resulted in global increase of H3K4 methylation along with transcriptional derepression of two BHC110 target genes, Egr1 and Oct4. When combined with CHIR99021, induces the generation of human-induced pluripotent stem cells in the absence of Sox2 (ref 3). Reduces proliferation, cell cycle progression and invasiveness of endometriotic stromal cells (ref 4). Upregulates CD86 expression in THP1 cells (ref 5). Cell permeable and active *in vivo*.

Solubilization: May be dissolved in water (25 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. NG Lee *et al. Bhem. Biol.* 2006, 13:563
2. DM Schmidt and DG McCafferty *Biochemistry* 2007, 46:4408
3. W Li *et al. Stem Cells* 2009, 27:2992
4. D Ding *et al. Fertil. Steril.* 2014, 101:740
5. JT Lynch *et al. Anal. Biochem.* 2013, 442:104

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of arrival.