

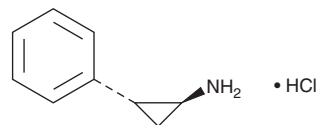
PRODUCT INFORMATION



Tranlycypromine (hydrochloride)

Item No. 10010494

CAS Registry No.: 1986-47-6
Formal Name: (1R,2S)-rel-2-phenyl-cyclopropanamine, monohydrochloride
Synonyms: 2-PCPA, *trans*-2-Phenylcyclopropylamine
MF: C₉H₁₁N • HCl
FW: 169.7
Purity: ≥98%
Supplied as: A crystalline 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

Tranlycypromine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the tranlycypromine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Tranlycypromine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of tranlycypromine (hydrochloride) in these solvents is approximately 5, 2.5, and 3 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of tranlycypromine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of tranlycypromine (hydrochloride) in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histones contain unstructured N-terminal residues that are the site of numerous post-translational modifications, involving acetylation, methylation, ubiquitination, and sumoylation to produce a specific gene regulatory outcome.¹ Complementary enzymes catalyze the addition and removal of these modifications as needed. The amine oxidase domain-containing enzyme lysine-specific demethylase 1 (LSD1) is a member of one of two classes of histone demethylases capable of demethylating lysine residues.¹ LSD1 shares similar catalytic sites with monoamine oxidases (MAO) A and B, the inhibition of which are used clinically to treat depression, anxiety, and Parkinson's disease.² Tranlycypromine is an irreversible, mechanism-based inhibitor of LSD1 with an IC₅₀ value of 20.7 μM and a K_i value of 242.7 μM that effectively inhibits histone demethylation *in vivo*.³ Although not as selective, tranlycypromine 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.³

References

1. Forneris, F., Binda, C., Vanoni, M.A., *et al.* Human histone demethylase LSD1 reads the histone code. *J. Biol. Chem.* **280(50)**, 41360-41365 (2005).
2. Lee, M.G., Wynder, C., Schidt, D.M., *et al.* Histone H3 lysine 4 demethylation is a target of nonselective antidepressive medications. *Chem. Biol.* **13(6)**, 563-567 (2006).
3. Schmidt, D.M.Z. and McCafferty, D.G. *trans*-2-Phenylcyclopropylamine is a mechanism-based inactivator of the histone demethylase LSD1. *Biochemistry* **46**, 4408-4416 (2007).

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, 10/25/2022

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