

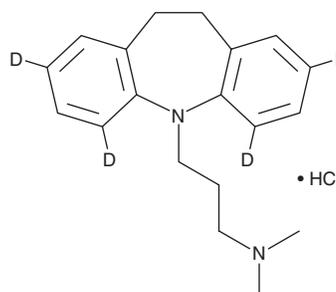
PRODUCT INFORMATION



Imipramine-d₄ (hydrochloride)

Item No. 43809

CAS Registry No.: 61361-33-9
Formal Name: 10,11-dihydro-N,N-dimethyl-5H-dibenz[b,f]azepine-2,4,6,8-d₄-5-propanamine, monohydrochloride
Synonym: Melipramine-d₄
MF: C₁₉H₂₀D₄N₂ • HCl
FW: 320.9
Chemical Purity: ≥98% (Imipramine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Supplied as: A 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

Imipramine-d₄ is intended for use as an internal standard for the quantification of imipramine (Item No. 15890) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Imipramine-d₄ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the imipramine-d₄ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Imipramine-d₄ (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of imipramine-d₄ (hydrochloride) in these solvents is approximately 25 mg/ml.

Description

Imipramine is a first generation tricyclic antidepressant.¹ It inhibits serotonin (5-HT) uptake by isolated human platelets by 54 and 79% when used at concentrations of 1 and 4 µg/ml, respectively, and inhibits norepinephrine uptake in rat brain by 63% when administered at a dose of 20 mg/kg.^{2,3} It also binds to histamine H₁, M₁ muscarinic acetylcholine, and α₁-adrenergic receptors (K_ds = 37, 46, and 32 nM, respectively).⁴ Imipramine (10 and 20 µM) prevents acid sphingomyelinase activation and subsequent ceramide release induced by infection with replication-deficient vesicular stomatitis virus pseudoviral particles (pp-VSV) presenting the severe acute respiratory coronavirus 2 (SARS-CoV-2) spike protein in Vero cells, an effect that can be overcome with exogenous application of C16 ceramide (d18:1/16:0) (Item No. 10681).⁵ Formulations containing imipramine have been used in the treatment of depression and as an adjunct in the treatment of childhood enuresis.

References

1. Spencer, P.S. *Br. J. Clin. Pharmacol.* **4(Suppl 2)**, 57S-68S (1977).
2. Yates, C.M., Todrick, A., and Tait, A.C. *J. Pharm. Pharmacol.* **16**, 460-463 (1964).
3. Glowinski, J. and Axelrod, J. *Nature* **204**, 1318-1319 (1964).
4. Cusack, B., Nelson, A., and Richelson, E. *Psychopharmacology (Berl.)* **114(4)**, 559-565 (1994).
5. Carpinheiro, A., Edwards, M.J., Hoffmann, M., et al. *Cell Rep. Med.* **1(8)**, 100142 (2020).

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, 06/16/2025

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