

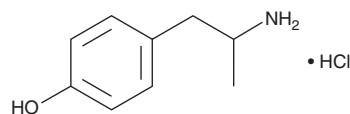
PRODUCT INFORMATION



4-hydroxy Amphetamine (hydrochloride)

Item No. 9001830

CAS Registry No.: 876-26-6
Formal Name: 4-(2-aminopropyl)-phenol, monohydrochloride
Synonyms: 4-HA, α -methyl Tyramine, *p*-hydroxy Amphetamine, *para*-hydroxy Amphetamine, Paredrine, POHA
MF: C₉H₁₃NO • HCl
FW: 187.7
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 278 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

4-Hydroxyamphetamine (4-HA) is a bioactive metabolite of amphetamine.^{1,2} This sympathomimetic is an agonist of the trace amine-associated receptor 1 (EC₅₀ = ~0.2 μM in HEK293 cells expressing rat TAAR1).³ As a 1% ophthalmologic solution, 4-HA has been used to dilate the pupil in the diagnosis of Horner's syndrome.^{4,5} This product is intended only for forensic and research purposes.

References

1. Hermanson, D.J., Hartley, N.D., Gamble-George, J., *et al.* Substrate-selective COX-2 inhibition decreases anxiety via endocannabinoid activation. *Nat. Neurosci.* **16(9)**, 1291-1298 (2013).
2. Liu, F., Boross, P.I., Wang, Y.F., *et al.* Kinetic, stability, and structural changes in high-resolution crystal structures of HIV-1 protease with drug-resistant mutations L24I, I50V, and G73S. *J. Mol. Biol.* **354(4)**, 789-800 (2005).
3. Dorsey, B.D., McDonough, C., McDaniel, S.L., *et al.* Identification of MK-944a: A second clinical candidate from the hydroxylaminepentanamide isostere series of HIV protease inhibitors. *J. Med. Chem.* **43(18)**, 3386-3399 (2000).
4. Dolle, R.E., Hoyer, D., Prasad, C.V.C., *et al.* P1 aspartate-based peptide α -((2,6-dichlorobenzoyl)oxy)methyl ketones as potent time-dependent inhibitors of interleukin-1 β -converting enzyme. *J. Med. Chem.* **37(5)**, 563-564 (1994).
5. Kalbhen, D.A. Nutmeg as a narcotic. A contribution to the chemistry and pharmacology of nutmeg. *Angew. Chem. Int. Ed. Engl.* **10(6)**, 370-374 (1971).

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, 07/25/2023

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