PRODUCT INFORMATION



N-hydroxy MDA (hydrochloride)

Item No. 13958

CAS Registry No.:	74341-83-6	
Formal Name:	N-hydroxy-a-methyl-1,3-benzodioxole-5-	
	ethanamine, monohydrochloride	
Synonyms:	MDOH,	н
	N-hydroxy-3,4-Methylenedioxyamphetamine	N
MF:	C ₁₀ H ₁₃ NO ₃ • HCl	
FW:	231.7	
Purity:	≥95%	
UV/Vis.:	λ _{max} : 236, 287 nm	• HCI
Supplied as:	A crystalline solid	
Storage:	-20°C	
Stability:	≥5 years	
Information represents the product energications. Batch energies analytical results are provided on each certificate of analysis		

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

Description

3,4-Methylenedioxyamphetamine (MDA) is psychedelic and entactogenic drug that is scheduled as a controlled substance in the United States.¹ N-hydroxy-MDA is an analog of MDA (Item No. 11554, hydrochloride form) which inhibits reuptake of noradrenaline (IC₅₀ = 87.8 μ M) more effectively than serotonin (IC₅₀ = 215 μ M).² Like MDA, N-hydroxy-MDA is regulated in the United States. This product is intended for forensic and research applications.

References

- 1. Baggott, M.J., Siegrist, J.D., Galloway, G.P., et al. Investigating the mechanisms of hallucinogen-induced visions using 3,4-methylenedioxyamphetamine (MDA): A randomized controlled trial in humans. PLoS One 5(12), 1-13 (2010).
- 2. Montgomery, T., Buon, C., Eibauer, S., et al. Comparative potencies of 3,4-methylenedioxymethamphetamine (MDMA) analogues as inhibitors of $[{}^{3}H]$ noradrenaline and $[{}^{3}H]$ 5-HT transport in mammalian cell lines. Br. J. Pharmacol. 152(7), 1121-1130 (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.

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