# **PRODUCT** INFORMATION



3,4-MDPA (hydrochloride)

Item No. 13972

CAS Registry No.:	74341-77-8		
Formal Name:	α-methyl-N-propyl-1,3-benzodioxole-5-		
	ethanamine, monohydrochloride		н
Synonym:	MDPR		Ï
MF:	$C_{13}H_{19}NO_2 \bullet HCI$		N N
FW:	257.8		Ť Ť Ì
Purity:	≥95%		
UV/Vis.:	λ <sub>max</sub> : 204, 236, 288 nm		
Supplied as:	A crystalline solid	• г	
Storage:	-20°C		
Stability:	≥5 years		

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

## Description

3,4-MDPA (hydrochloride) is a lesser known psychoactive substituted amphetamine of the phenethylamine class. Though little is known about the pharmacological properties, metabolism, and toxicity of 3,4-MDPA, it has been used as an internal standard for qualitative and quantitative analyses of hallucinogenic designer drugs in biological samples.<sup>1</sup>. This compound is intended for forensic and research purposes only.

## Reference

1. Pichini, S., Pujadas, M., Marchei, E., et al. Liquid chromatography-atmospheric pressure ionization electrospray mass spectrometry determination of "hallucinogenic designer drugs" in urine of consumers. J. Pharm. Biomed. Anal. 47(2), 335-342 (2008).

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