PRODUCT INFORMATION



5-methoxy EPT (hydrochloride)

Item No. 17298

CAS Registry No.:	2749303-04-4	
Formal Name:	N-ethyl-5-methoxy-N-propyl-1H-indole-	~
	3-ethanamine, monohydrochloride	
Synonym:	5-methoxy-N-ethyl-N-Propyltryptamine	
MF:	$C_{16}H_{24}N_2O \bullet HCI$	
FW:	296.8	
Purity:	≥98%	• HCl
Supplied as:	A neat solid	Н
Storage:	-20°C	
Stability:	≥1 year	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Description

5-methoxy EPT (hydrochloride) (Item No. 17298) is an analytical reference material that is classified as a tryptamine. It was detected in designer drugs purchased in an illegal market in Japan.¹ The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

Reference

1. Uchiyama, N., Miyazawa, N., Kawamura, M., et al. Analysis of newly distributed designer drugs detected in the products purchased in fiscal year 2008. Yakugaku Zasshi 130(2), 263-270 (2010).

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, 05/10/2022

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