

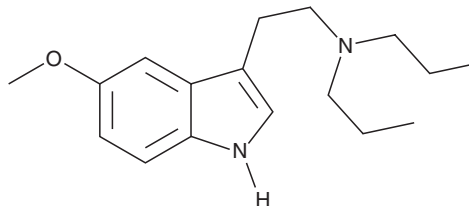
# PRODUCT INFORMATION



## 5-methoxy DPT

Item No. 17297

**CAS Registry No.:** 69496-75-9  
**Formal Name:** 5-methoxy-N,N-dipropyl-1H-indole-3-ethanamine  
**Synonyms:** 5-MeO DPT, 5-methoxy-N,N-Dipropyltryptamine  
**MF:** C<sub>17</sub>H<sub>26</sub>N<sub>2</sub>O  
**FW:** 274.4  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**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 DPT is a substituted form of DPT that is also an analog of 5-methoxy DiPT (Item No. 11865), a psychoactive monoamine re-uptake inhibitor.<sup>1,2</sup> The physiological and toxicological properties of this compound are not known. 5-methoxy DPT is regulated as a Schedule I compound in the United States. 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.

### References

1. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).
2. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11(5)**, 468-482 (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.

#### 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, 10/02/2020

#### 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