

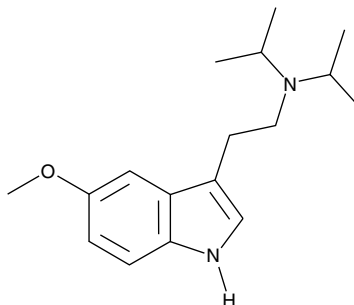
# Product Information



## 5-methoxy DiPT (exempt preparation)

Item No. 15697

**CAS Registry No.:** 4021-34-5  
**Formal Name:** 5-methoxy-N,N-bis(1-methylethyl)-1H-indole-3-ethanamine  
**Synonyms:** 5-methoxy-N,N-Diisopropyltryptamine, FOXY, 5-MeO DiPT  
**MF:** C<sub>17</sub>H<sub>26</sub>N<sub>2</sub>O  
**FW:** 274.4  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A solution in methanol  
**UV/Vis.:** λ<sub>max</sub>: 204, 224, 280 nm



### Laboratory Procedures

For long term storage, we suggest that 5-methoxy DiPT (exempt preparation) be stored as supplied at -20°C. It should be stable for at least two years.

5-methoxy DiPT (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 5-methoxy DiPT (exempt preparation) in these solvents is approximately 16, 3, and 5 mg/ml, respectively.

5-methoxy DiPT (exempt preparation) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 5-methoxy DiPT (exempt preparation) should be diluted with the aqueous buffer of choice. 5-methoxy DiPT (exempt preparation) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

5-methoxy DiPT is a tryptamine-type designer drug with pronounced psychoactive and physiological effects.<sup>1</sup> It potentially inhibits the re-uptake of monoamines (IC<sub>50</sub>s = 0.65, 2.2, and 8.2 μM for dopamine, serotonin, and norepinephrine, respectively) while not affecting their release from rat brain synaptosomes.<sup>2</sup> This product is intended for forensic applications.

### References

1. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11**, 468-482 (2010).
2. 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).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/15697](http://www.caymanchem.com/catalog/15697)

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 04/04/2014