

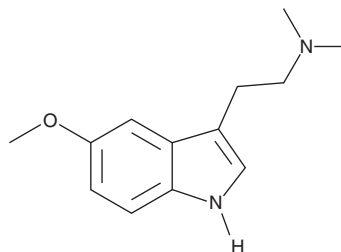
# PRODUCT INFORMATION



## 5-methoxy DMT

Item No. 11480

**CAS Registry No.:** 1019-45-0  
**Formal Name:** 5-methoxy-N,N-dimethyl-1H-indole-3-ethanamine  
**Synonyms:** 5-methoxy-N,N-Dimethyltryptamine, 5-MeO DMT  
**MF:** C<sub>13</sub>H<sub>18</sub>N<sub>2</sub>O  
**FW:** 218.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 278 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

5-methoxy DMT is a naturally occurring hallucinogenic indolealkylamine that potently activates serotonin (5-HT) receptors.<sup>1,2</sup> Although this compound is regulated in many countries, including the United States, it has been described as an 'emerging psychoactive substance' found in party pills and related mixtures.<sup>3</sup> As 5-methoxy DMT is inactivated by monoamine oxidases, inhibitors of monoamine oxidases are often combined with 5-methoxy DMT to prolong its activity.<sup>4,5</sup> The pharmacological and toxicological properties of this compound have been recently reviewed.<sup>6</sup> 5-methoxy DMT is regulated as a Schedule I compound in the United States. This product is intended for forensic applications.

### References

1. Glennon, R.A., Liebowitz, S.M., and Mack, E.C. Serotonin receptor binding affinities of several hallucinogenic phenylalkylamine and N,N-dimethyltryptamine analogues. *J. Med. Chem.* **21(8)**, 822-825 (1978).
2. Krebs-Thomson, K., Ruiz, E.M., Masten, V., *et al.* The roles of 5-HT<sub>1A</sub> and 5-HT<sub>2</sub> receptors in the effects of 5-MeO-DMT on locomotor activity and prepulse inhibition in rats. *Psychopharmacology (Berl)* **189(3)**, 319-329 (2006).
3. Bruno, R., Matthews, A.J., Dunn, M., *et al.* Emerging psychoactive substance use among regular ecstasy users in Australia. *Drug Alcohol Depend.* **124(1-2)**, 19-25 (2012).
4. Halberstadt, A.L., Buell, M.R., Masten, V.L., *et al.* Modification of the effects of 5-methoxy-N,N-dimethyltryptamine on exploratory behavior in rats by monoamine oxidase inhibitors. *Psychopharmacology (Berl)* **201(1)**, 55-66 (2008).
5. Shen, H.W., Wu, C., Jiang, X.L., *et al.* Effects of monoamine oxidase inhibitor and cytochrome P450 2D6 status on 5-methoxy-N,N-dimethyltryptamine metabolism and pharmacokinetics. *Biochem. Pharmacol.* **80(1)**, 122-128 (2010).
6. Shen, H.W., Jiang, X.L., Winter, J.C., *et al.* Psychedelic 5-methoxy-N,N-dimethyltryptamine: Metabolism, pharmacokinetics, drug interactions, and pharmacological actions. *Curr. Drug Metab.* **11(8)**, 659-666 (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, 07/26/2022

#### 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