

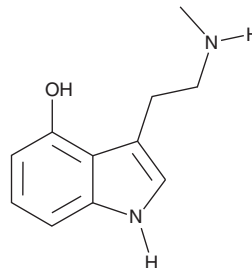
# PRODUCT INFORMATION



## Norpsilocin

Item No. 23586

**CAS Registry No.:** 28363-70-4  
**Formal Name:** 3-[2-(methylamino)ethyl]-1H-indol-4-ol  
**Synonyms:**  $\omega$ -N-methyl-4-Hydroxytryptamine,  
4-hydroxy-N-Methyltryptamine,  
4-hydroxy NMT  
**MF:** C<sub>11</sub>H<sub>14</sub>N<sub>2</sub>O  
**FW:** 190.2  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{\text{max}}$ : 224 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq 2$  years



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

### Description

Norpsilocin (Item No. 23586) is an analytical reference standard categorized as a tryptamine.<sup>1</sup> Norpsilocin has been found in *P. cubensis*, a mushroom with hallucinogenic properties. It is a potential metabolite of baeocystin (Item Nos. 37763 | 31583) based on the published metabolism of psilocybin (Item Nos. 9003134 | 14041).<sup>2</sup> At the time norpsilocin (Item No. 23586) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

### References

1. Lenz, C., Wick, J., and Hoffmeister, D. Identification of  $\omega$ -N-methyl-4-hydroxytryptamine (Norpsilocin) as a Psilocybe natural product. *J. Nat. Prod.* **80**(10), 2835-2838 (2017).
2. Dinis-Oliveira, R.J. Metabolism of psilocybin and psilocin: Clinical and forensic toxicological relevance. *Drug Metab. Rev.* **49**(1), 84-91 (2017).

#### 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, 07/11/2023

#### 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](http://WWW.CAYMANCHEM.COM)