

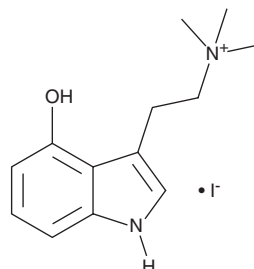
PRODUCT INFORMATION



4-hydroxy TMT (iodide)

Item No. 33670

CAS Registry No.: 2451455-99-3
Formal Name: 4-hydroxy-N,N,N-trimethyl-1H-indole-3-ethanaminium, monoiodide
Synonym: 4-hydroxy-N,N,N-Trimethyltryptamine
MF: C₁₃H₁₉N₂O • I
FW: 346.2
Purity: ≥98%
UV/Vis.: λ_{max}: 221 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

4-hydroxy TMT (iodide) (Item No. 33670) is an analytical reference standard categorized as a tryptamine.¹ 4-hydroxy TMT is an active metabolite of aeruginascin (Item Nos. 37762 | 31582). It does not induce the head-twitch response (HTR) in mice.² This product is intended for research and forensic applications.

References

1. Chadeayne, A.R., Pham, D.N.K., Reid, B.G., *et al.* Active metabolite of aeruginascin (4-hydroxy-N,N,N-trimethyltryptamine): Synthesis, structure, and serotonergic binding affinity. *ACS Omega* **5**(27), 16940-16943 (2020).
2. Glatfelter, G.C., Pottie, E., Partilla, J.S., *et al.* Structure-activity relationships for psilocybin, baeocystin, aeruginascin, and related analogues to produce pharmacological effects in mice. *ACS Pharmacol. Transl. Sci.* **5**(11), 1181-1196 (2022).

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, 02/14/2025

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