

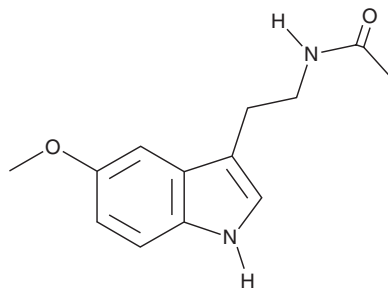
PRODUCT INFORMATION



Melatonin

Item No. 42524

CAS Registry No.: 73-31-4
Formal Name: N-[2-(5-methoxy-1H-indol-3-yl)ethyl]-acetamide
Synonyms: N-Acetyl-5-methoxytryptamine,
5-Methoxy-N-acetyltryptamine,
NSC 56423, NSC 113928
MF: C₁₃H₁₆N₂O₂
FW: 232.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

Melatonin (Item No. 42524) is an analytical reference standard categorized as a melatonin receptor agonist.¹ Melatonin has been found as an adulterant in products marketed as food supplements.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 14427).

References

1. Boutin, J.A., Audinot, V., Ferry, G., *et al.* Molecular tools to study melatonin pathways and actions. *Trends Pharmacol. Sci.* **26(8)**, 412-419 (2005).
2. Paiva, R., Correia, M., Delerue-Matos, C., *et al.* Adulteration of brain health (cognitive, mood, and sleep enhancement) food supplements by the addition of pharmaceutical drugs: A comprehensive review of analytical approaches and trends. *Foods* **13(6)**, 908 (2024).

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, 01/21/2024

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