

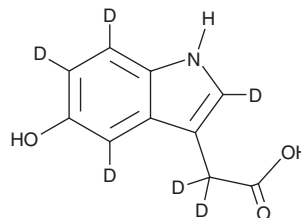
PRODUCT INFORMATION



5-hydroxy Indole-3-acetic Acid-d₆

Item No. 37199

CAS Registry No.: 2748469-82-9
Formal Name: 2-(5-hydroxy-1H-indol-3-yl-2,4,6,7-d₄)acetic-2,2-d₂ acid
Synonyms: 5-HIAA-d₆, 5-Hydroxyindoleacetic Acid-d₆, 5-hydroxy IAA-d₆
MF: C₁₀H₃D₆NO₃
FW: 197.2
Chemical Purity: ≥98% (5-hydroxy Indole-3-acetic Acid)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀
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.

Laboratory Procedures

5-hydroxy Indole-3-acetic acid-d₆ (5-HIAA-d₆) is intended for use as an internal standard for the quantification of 5-HIAA (Item No. 22889) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

5-HIAA-d₆ is supplied as a solid. A stock solution may be made by dissolving the 5-HIAA-d₆ in the solvent of choice, which should be purged with an inert gas. 5-HIAA-d₆ is soluble in organic solvents such as methanol, DMSO, and ethyl acetate.

Description

5-HIAA is a metabolite of the monoamine neurotransmitter serotonin (5-HT; Item No. 14332).¹ It is formed from serotonin *via* a 5-hydroxy indole-3-acetaldehyde intermediate by aldehyde dehydrogenase. 5-HIAA has been used as a biomarker for the detection of neuroendocrine tumors. Brain levels of 5-HIAA are decreased in patients with severe Alzheimer's disease.²

References

1. Corcuff, J.-B., Chardon, L., El Hajji Ridah, I., *et al.* Urinary sampling for 5HIAA and metanephrines determination: Revisiting the recommendations. *Endocr. Connect.* **6**(6), R87-R98 (2017).
2. Vermeiren, Y., Van Dam, D., Aerts, T., *et al.* Monoaminergic neurotransmitter alterations in postmortem brain regions of depressed and aggressive patients with Alzheimer's disease. *Neurobiol. Aging* **35**(12), 2691-2700 (2014).

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, 11/18/2022

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