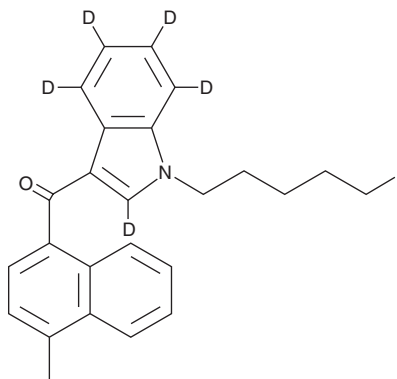


PRODUCT INFORMATION



MAM2201-d₅
Item No. 43606

CAS Registry No.: 2095544-38-8
Formal Name: [1-(5-fluoropentyl)-1H-indol-3-yl-2,4,5,6,7-d₅]
(4-methyl-1-naphthalenyl)-methanone
Synonyms: AM2201 4-methylnaphthyl analog-d₅,
JWH 122 N-(5-fluoropentyl) analog-d₅
MF: C₂₅H₁₉D₅FNO
FW: 378.5
Chemical Purity: ≥98% (MAM2201)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
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

MAM2201-d₅ (Item No. 43606) is intended for use as an internal standard for the quantification of MAM2201 (Item No. 9001219) 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).

MAM2201 is categorized as a synthetic cannabinoid.¹ MAM2201 has been found in Spice/K2-type herbal blends and has been associated with fatal overdose.^{2,3} Its use has been associated with impaired driving and motor vehicle collisions.⁴ MAM2201-d₅ is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Corli, G., Tirri, M., Bilel, S., *et al.* MAM-2201 acute administration impairs motor, sensorimotor, prepulse inhibition, and memory functions in mice: A comparison with its analogue AM-2201. *Psychopharmacology (Berl)* **240(7)**, 1435-1452 (2023).
2. Seely, K.A., Patton, A.L., Moran, C.L., *et al.* Forensic investigation of K2, Spice, and “bath salt” commercial preparations: A three-year study of new designer drug products containing synthetic cannabinoid, stimulant, and hallucinogenic compounds. *Forensic Sci. Int.* **233(1-3)**, 416-422 (2013).
3. Saito, T., Namera, A., Miura, N., *et al.* A fatal case of MAM-2201 poisoning. *Forensic toxicol.* **31**, 333-337 (2013).
4. Kaneko, S. Motor vehicle collisions caused by the ‘super-strength’ synthetic cannabinoids, MAM-2201, 5F-PB-22, 5F-AB-PINACA, 5F-AMB and 5F-ADB in Japan experienced from 2012 to 2014. *Forensic Toxicol.* **35(2)**, 244-251 (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, 06/05/2026

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