

PRODUCT INFORMATION



MAM2201

Item No. 9001219

CAS Registry No.: 1354631-24-5

Formal Name: [1-(5-fluoropentyl)-1H-indol-3-yl]
(4-methyl-1-naphthalenyl)-methanone

Synonyms: AM2201 4-methylnaphthyl analog,
JWH 122 N-(5-fluoropentyl) analog

MF: C₂₅H₂₄FNO

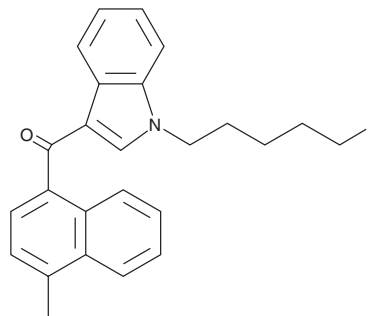
FW: 373.5

Purity: ≥98%

Supplied as: A neat solid

Storage: -20°C

Stability: ≥2 years



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

Description

MAM2201 (Item No. 9001219) is an analytical reference material 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 This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

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, 05/28/2025

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