

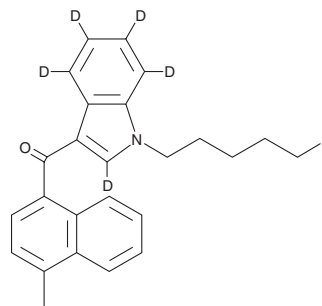
# PRODUCT INFORMATION



## MAM2201-d<sub>5</sub> (solution)

Item No. 18156

**CAS Registry No.:** 2095544-38-8  
**Formal Name:** [1-(5-fluoropentyl)-1H-indol-3-yl-d<sub>5</sub>]-  
(4-methyl-1-naphthalenyl)-methanone  
**Synonyms:** AM2201 4-methylnaphthyl analog,  
JWH 122 N-(5-fluoropentyl) analog  
**MF:** C<sub>25</sub>H<sub>19</sub>D<sub>5</sub>FNO  
**FW:** 378.5  
**Purity:** ≥98%  
**Supplied as:** A 100 µg/ml solution in methanol  
**Storage:** -20°C  
**Stability:** ≥3 years



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

### Description

MAM2201-d<sub>5</sub> is an analytical reference material intended for use as an internal standard for the quantification of MAM2201 (Item No. ISO00120) by GC- or LC-MS. AM2201 (Item No. 10707) is a potent synthetic cannabinoid (CB) which binds the CB<sub>1</sub> and CB<sub>2</sub> receptors with high affinity (K<sub>i</sub> = 1.0 and 2.6 nM, respectively).<sup>1</sup> MAM2201 is an analog of AM2201 that is methylated at the 4 position of the naphthyl group. The physiological and toxicological properties of this compound have not been delineated. This product is intended for research and forensic purposes.

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

### Reference

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832

#### 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, 08/05/2022

#### 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