

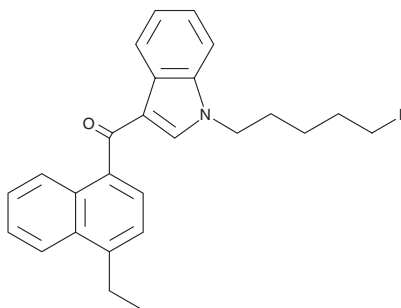
# PRODUCT INFORMATION



**EAM2201**

Item No. 11916

**CAS Registry No.:** 1364933-60-7  
**Formal Name:** (4-ethyl-1-naphthalenyl)  
[1-(5-fluoropentyl)-1H-indol-3-yl]-  
methanone  
**Synonym:** JWH 210 N-(5-fluoropentyl) analog  
**MF:** C<sub>26</sub>H<sub>26</sub>FNO  
**FW:** 387.5  
**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

M2201 (Item No. 10707) is a potent synthetic cannabinoid (CB) with K<sub>i</sub> values of 1.0 and 2.6 nM for the central CB<sub>1</sub> and peripheral CB<sub>2</sub> receptors, respectively.<sup>1</sup> It has been detected in herbal blends.<sup>2</sup> EAM2201 is an analog of AM2201 having an ethyl group added to the 4 position of the naphthyl group. The physiological properties of this compound are unknown. This product is intended for forensic and research applications.

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

## References

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832.
2. Simolka, K., Lindigkeit, R., Schiebel, H.-M., *et al.* Analysis of synthetic cannabinoids in "spice-like" herbal highs: Snapshot of the German market in summer 2011. *Anal. Bioanal. Chem.* [In press] 1-15 (2012).

### 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, 08/09/2019

## 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](http://WWW.CAYMANCHEM.COM)