

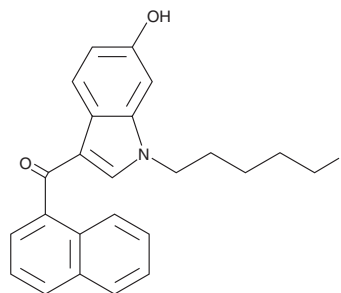
# PRODUCT INFORMATION



## AM2201 6-hydroxyindole metabolite

Item No. 11192

**CAS Registry No.:** 1427521-35-4  
**Formal Name:** (1-(5-fluoropentyl)-6-hydroxy-1H-indol-3-yl)(naphthalen-1-yl) methanone  
**MF:** C<sub>24</sub>H<sub>22</sub>FNO<sub>2</sub>  
**FW:** 375.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219, 294 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

AM2201 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> AM2201 6-hydroxyindole metabolite is an expected metabolite of AM2201 generated during phase I metabolism by cytochrome P450, suggesting that it would be detectable in blood and urine.<sup>2</sup> The biological properties of this compound are not known. This product is intended for forensic applications.

### References

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832.
2. Wintermeyer, A., Möller, I., Thevis, M., *et al.* *In vitro* phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**, 2141-53 (2010).

#### 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, 07/07/2023

#### 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)