

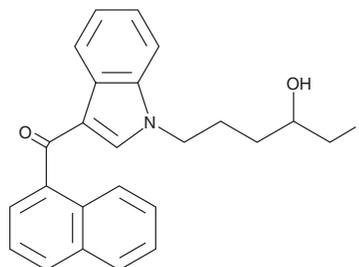
PRODUCT INFORMATION



AM2201 N-(4-hydroxypentyl) metabolite

Item No. 10203

CAS Registry No.: 1427521-34-3
Formal Name: (1-(5-fluoro-4-hydroxypentyl)-1H-indol-3-yl)(naphthalen-1-yl) methanone
MF: C₂₄H₂₂FNO₂
FW: 375.4
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 247, 315 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

AM2201 N-(4-hydroxypentyl) metabolite is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, dichloromethane, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of AM2201 N-(4-hydroxypentyl) metabolite in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO, dichloromethane, and DMF.

AM2201 N-(4-hydroxypentyl) metabolite is sparingly soluble in aqueous solutions. Therefore, further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

AM2201 is a potent synthetic cannabinoid (CB) with K_i values of 1.0 and 2.6 nM for the CB₁ and CB₂ receptors, respectively.¹ AM2201 N-(4-hydroxypentyl) metabolite is an expected urinary metabolite of AM2201, based on the known metabolism of similar compounds.² Its biological actions are unknown.

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, 11/20/2025

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