

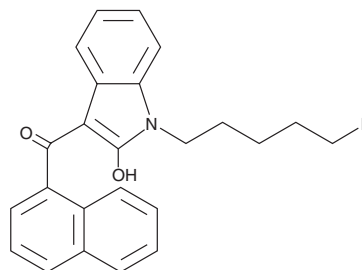
PRODUCT INFORMATION



AM2201 2-hydroxyindole metabolite

Item No. 11194

CAS Registry No.: 2749391-49-7
Formal Name: (1-(5-fluoropentyl)-2-hydroxy-1H-indol-3-yl)(naphthalen-1-yl) methanone
MF: C₂₄H₂₂FNO₂
FW: 375.4
Purity: ≥98%
Storage: -20°C
Stability: ≥1 year
Supplied as: A solution in acetonitrile
UV/Vis.: λ_{max}: 217, 265, 310 nm



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

Laboratory Procedures

AM2201 2-hydroxyindole metabolite is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of AM2201 2-hydroxyindole metabolite in these solvents is approximately 20 mg/ml.

AM2201 2-hydroxyindole metabolite is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of AM2201 2-hydroxyindole metabolite should be diluted with the aqueous buffer of choice. AM2201 2-hydroxyindole metabolite has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. 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 central CB₁ and peripheral CB₂ receptors, respectively.¹ AM2201 2-hydroxyindole metabolite is a potential monohydroxylated urinary metabolite of AM2201, based on the known metabolism of similar compounds.² Its biological actions are unknown. This product is intended for forensic and research applications.

References

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. **WO01/28557A1**, 1-25, PCT/US00/28832 (2001).
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, 05/23/2022

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