

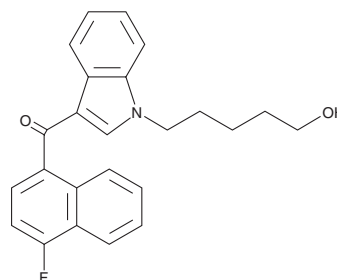
# PRODUCT INFORMATION



## JWH 412 N-(5-hydroxypentyl) metabolite

Item No. 17789

**CAS Registry No.:** 2748591-83-3  
**Formal Name:** (4-fluoronaphthalen-1-yl)(1-(5-hydroxypentyl)-1H-indol-3-yl)methanone  
**Synonym:** F2201 N-(5-hydroxypentyl) metabolite  
**MF:** C<sub>24</sub>H<sub>22</sub>FNO<sub>2</sub>  
**FW:** 375.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 317 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

JWH 412 (5-hydroxypentyl) metabolite (Item No. 17789) is an analytical reference standard that is structurally similar to known synthetic cannabinoids (CBs). It is a presumptive metabolite of JWH 412 (Item No. 11971), based on the metabolism of the structurally similar CBJWH 018 (Item No. 10900).<sup>1</sup> This product is intended for research and forensic applications.

### Reference

1. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200(1-3)**, 141-147 (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, 06/14/2023

#### 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