

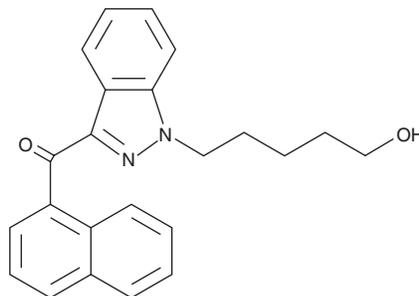
# PRODUCT INFORMATION



## THJ2201 N-(5-hydroxypentyl) metabolite

Item No. 9003736

**CAS Registry No.:** 1850409-16-3  
**Formal Name:** [1-(5-hydroxypentyl)-1H-indazol-3-yl]-1-naphthalenyl-methanone  
**Synonym:** THJ 018 N-(5-hydroxypentyl) metabolite  
**MF:** C<sub>23</sub>H<sub>22</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 358.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 220 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

THJ2201 N-(5-hydroxypentyl) metabolite (Item No. 9003736) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.<sup>1,2</sup> THJ2201 N-(5-hydroxypentyl) metabolite is a metabolite of THJ2201 (Item Nos. 14789 | 22860) and THJ 018 (Item No. 11962). This product is intended for research and forensic applications.

### References

1. Labutina, A.V. and Temerdashev, A.Z. Nontarget screening of the markers of synthetic cannabinoids in urine using HPLC-MS/MS. *J. Anal. Chem.* **70**, 1620-1628 (2015).
2. Diao, X., Wohlfarth, A., Pang, S., *et al.* High-resolution mass spectrometry for characterizing the metabolism of synthetic cannabinoid THJ-018 and its 5-fluoro analog THJ-2201 after incubation in human hepatocytes. *Clin. Chem.* **62(1)**, 157-169 (2016).

#### 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, 04/11/2025

#### 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