

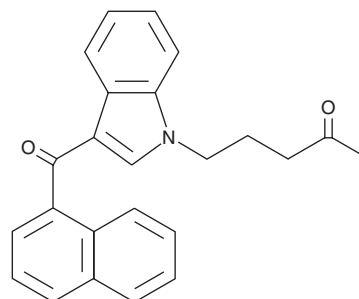
# PRODUCT INFORMATION



## JWH 018 N-(4-oxo-pentyl) metabolite

Item No. 13889

**CAS Registry No.:** 2748474-41-9  
**Formal Name:** 5-(3-(1-naphthoyl)-1H-indol-1-yl)pentan-2-one  
**Synonym:** JWH 018 4-keto metabolite  
**MF:** C<sub>24</sub>H<sub>21</sub>NO<sub>2</sub>  
**FW:** 355.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219, 246, 315 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

JWH 018 is a synthetic cannabinoid which has been identified in smoking mixtures.<sup>1,2</sup> JWH 018 N-(4-oxo-pentyl) metabolite is a potential metabolite of JWH 018 that is characterized by a carbonyl group at the 4 position of the alkyl chain. Its biological and toxicological properties are not known. This product is intended for forensic and research applications.

### References

1. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200**, 141-147 (2010).
2. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al.* Quantitative measurement of JWH-018 and JWH-073 metabolites excreted in human urine. *Anal. Chem.* **83(11)**, 4228-4236 (2011).

#### 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, 07/25/2023

#### 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