

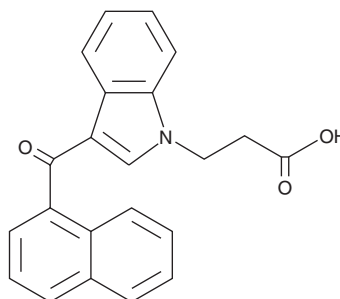
# PRODUCT INFORMATION



## JWH 018 N-propanoic acid metabolite

Item No. 14646

**CAS Registry No.:** 1451369-32-6  
**Formal Name:** 3-(1-naphthalenylcarbonyl)-1H-indole-1-propanoic acid  
**Synonyms:** AM2201 N-propanoic acid metabolite, JWH 073 N-propanoic acid metabolite  
**MF:** C<sub>22</sub>H<sub>17</sub>NO<sub>3</sub>  
**FW:** 343.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 218, 245, 313 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 (Item No. 10900) is a mildly selective agonist of peripheral cannabinoid (CB<sub>2</sub>) receptor, derived from the aminoalkylindole WIN 55,212-2. JWH 018 binds central cannabinoid (CB<sub>1</sub>) and CB<sub>2</sub> receptors with K<sub>i</sub> values of 9.0 and 2.94 nM, respectively.<sup>1</sup> JWH 018 is one of several synthetic CBs that have been included in smoking mixtures.<sup>2</sup> JWH 018 N-propanoic acid metabolite (Item No. 14646) is a major metabolite of JWH 018 that is detectable in urine.<sup>2</sup> The biological and toxicological properties of this compound have not been characterized. This analytical reference standard is intended for forensic and research applications only.

### References

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB<sub>1</sub> and CB<sub>2</sub> receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
2. Lovett, D.P., Yanes, E.G., Herbelin, T.W., *et al.* Structure elucidation and identification of a common metabolite for naphthoylindole-based synthetic cannabinoids using LC-TOF and comparison to a synthetic reference standard. *Forensic Sci. Int.* **226(1-3)**, 81-87 (2013).

#### 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, 06/21/2023

#### 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