

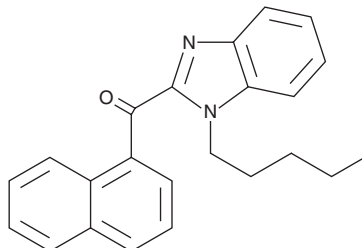
# PRODUCT INFORMATION



## JWH 018 benzimidazole analog

Item No. 15074

**CAS Registry No.:** 2316839-70-8  
**Formal Name:** naphthalen-1-yl(1-pentyl-1H-benzo[d]imidazol-2-yl)methanone  
**Synonym:** BIM-108  
**MF:** C<sub>23</sub>H<sub>22</sub>N<sub>2</sub>O  
**FW:** 342.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 219, 324 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 the peripheral cannabinoid (CB<sub>2</sub>) receptor with K<sub>i</sub> values of 9.0 and 2.94 nM for the central cannabinoid (CB<sub>1</sub>) and CB<sub>2</sub> receptors, respectively.<sup>1</sup> JWH 018 benzimidazole analog replaces the indole core of JWH 018 with a benzimidazole core. While the physiological and toxicological properties of this compound have not been tested, other benzimidazole derivatives demonstrate high selectivity for the CB<sub>2</sub> receptor.<sup>2</sup> This product is intended for forensic and research applications.

### 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. Pagé, D., Balaux, E., Boisvert, L., *et al.* Novel benzimidazole derivatives as selective CB<sub>2</sub> agonists. *Bioorg. Med. Chem. Lett.* **18(13)**, 3695-3700 (2008).

#### 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, 06/26/2023

#### 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