

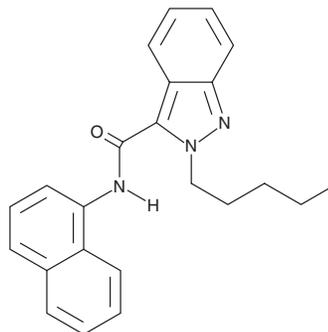
# PRODUCT INFORMATION



## NNEI 2'-indazole isomer

Item No. 17520

**CAS Registry No.:** 2365471-66-3  
**Formal Name:** N-1-naphthalenyl-2-pentyl-2H-indazole-3-carboxamide  
**Synonym:** MN-018 2'-indazole isomer  
**MF:** C<sub>23</sub>H<sub>23</sub>N<sub>3</sub>O  
**FW:** 357.5  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥6 years



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

### Description

NNEI (Item No. 15001) is an analog of JWH 018 (Item No. 10900), an aminoalkylindole-based synthetic cannabinoid that has been identified in smoking mixtures.<sup>1,2</sup> NNEI 2'-indazole isomer is an indazole analog of NNEI with the alkyl group at the two position. This compound is a close analog of MN-18 (Item No. 14817), which is also an indazole-based cannabinoid but has the alkyl group at the one position. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

### 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, 08/03/2022

#### 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