

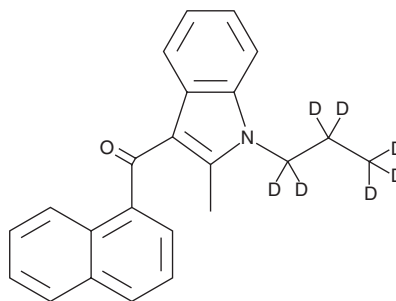
PRODUCT INFORMATION



JWH 015-d₇
Item No. 10660

CAS Registry No.: 1346601-16-8
Formal Name: (2-methyl-1-propyl-(1,1,2,2,3,3,3-d₇)-1H-indol-3-yl)-1-naphthalenyl-methanone
MF: C₂₃H₁₄D₇NO
FW: 334.5

Chemical Purity: ≥98% (JWH 015)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
UV/Vis.: λ_{max}: 219, 319 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 015-d₇ (Item No. 10660) is intended for use as an internal standard for the quantification of JWH 015 (Item No. 10009018) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

JWN 015 is categorized as a synthetic cannabinoid.¹ JWH 015 has been found in Spice/K2 herbal blends.² This product is intended for research and forensic applications.

References

1. Pertwee, R.G. Pharmacology of cannabinoid receptor ligands. *Curr. Med. Chem.* **6**(8), 635-664 (1999).
2. Seely, K.A., Patton, A.L., Moran, C.L., *et al.* Forensic investigation of K2, Spice, and "bath salt" commercial preparations: A three-year study of new designer drug products containing synthetic cannabinoid, stimulant, and hallucinogenic compounds. *Forensic Sci. Int.* **233**(1-3), 416-422 (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, 07/18/2023

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