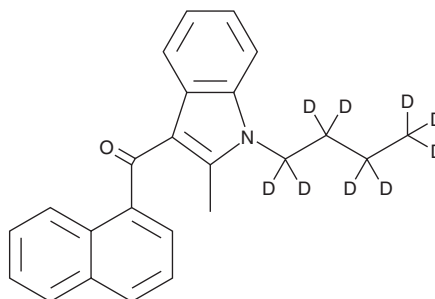


PRODUCT INFORMATION



JWH 016-d₉
Item No. 20850

CAS Registry No.: 2748469-65-8
Formal Name: (1-butyl-d₉-2-methyl-1H-indol-3-yl)-1-naphthalenyl-methanone
MF: C₂₄H₁₄D₉NO
FW: 350.5
Chemical Purity: ≥98% (JWH 016)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 218, 320 nm
Supplied as: A 1 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 016-d₉ (Item No. 20850) is intended for use as an internal standard for the quantification of JWH 016 (Item No. 10849) 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).

JWH 016 is categorized as a synthetic cannabinoid.^{1,2} It induces discriminative stimulus effects in rhesus monkeys comparable to those induced by Δ⁹-THC (Item Nos. ISO60157 | 12068).³ This product is intended for research and forensic 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 CB₁ and CB₂ receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
2. Wiley, J.L., Compton, D.R., Dai, D., *et al.* Structure-activity relationships of indole- and pyrrole-derived cannabinoids. *J. Pharmacol. Exp. Ther.* **285(3)**, 995-1004 (1998).
3. Wiley, J.L., Huffman, J.W., Balster, R.L., *et al.* Pharmacological specificity of the discriminative stimulus effects of Δ⁹-tetrahydrocannabinol in rhesus monkeys. *Drug Alcohol Depend.* **40(1)**, 81-86 (1995).

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, 11/12/2025

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