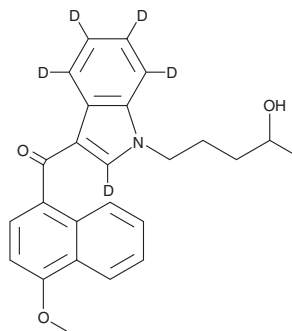


PRODUCT INFORMATION



JWH 081 N-(4-hydroxypentyl) metabolite-d₅ Item No. 14360

CAS Registry No.: 2747916-00-1
Formal Name: (1-(4-hydroxypentyl)-1H-indol-3-yl-2,4,5,6,7-d₅)
(4-methoxynaphthalen-1-yl)methanone
MF: C₂₅H₂₀D₅NO₃
FW: 392.5
Chemical Purity: ≥98% JWH 081 N-(4-hydroxypentyl) metabolite
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 212, 317 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 081 N-(4-hydroxypentyl) metabolite-d₅ (Item No. 14360) is intended for use as an internal standard for the quantification of JWH 081 N-(4-hydroxypentyl) metabolite (Item No. 14348) 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 081 N-(4-hydroxypentyl) metabolite is an expected product of the phase I metabolism of JWH 081 (Item No. 10579), based on studies on the metabolism of the related synthetic cannabinoids JWH 018 (Item No. 10900) and JWH 073 (Item No. 10904).^{1,2}

References

1. Wintermeyer, A., Möller, I., Thevis, M., *et al.* *In vitro* phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398**, 2141-53 (2010).
2. Chimalakonda, K.C., Moran, C.L., Kennedy, P.D., *et al.* Solid-phase extraction and quantitative measurement of omega and omega-1 metabolites of JWH-018 and JWH-073 in human urine. *Anal. Chem.* **83**(16), 6381-6388 (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, 06/14/2023

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