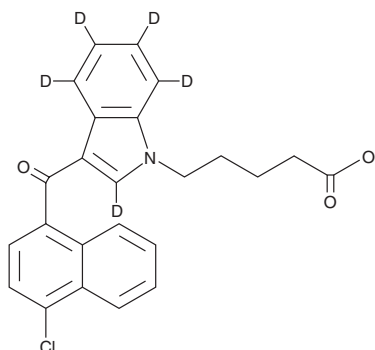


PRODUCT INFORMATION



JWH 398 N-pentanoic acid metabolite-d₅ Item No. 14372

CAS Registry No.: 2749394-76-9
Formal Name: 5-(3-(4-chloro-1-naphthoyl)-1H-indol-1-yl)-2,4,5,6,7-d₅)pentanoic acid
Synonym: JWH 398 N-(5-carboxypentyl) metabolite-d₅
MF: C₂₄H₁₅ClD₅NO₃
FW: 410.9
Chemical Purity: ≥98% (JWH 398 N-pentanoic acid metabolite)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 221, 313, 419 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

JWH 398 N-pentanoic acid metabolite-d₅ (Item No. 14372) is intended for use as an internal standard for the quantification of JWH 398 N-pentanoic acid metabolite (Item No. 10942) 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 398 (Item No. 13636) is a synthetic cannabinoid (CB) that acts as an agonist at both the central CB₁ receptor and the peripheral CB₂ receptor (K_is = 2.3 and 2.8 nM, respectively).¹ It has been detected as an additive of herbal mixtures.² JWH 398 N-pentanoic acid metabolite is an expected phase I metabolite of JWH 398. The biological and toxicological properties of this compound have not been characterized. This product is intended for forensic and research purposes.

References

- Huffman, J.W. Cannabimimetic indoles, pyrroles, and indenes: Structure-activity relationships and receptor interactions. *The cannabinoid receptors*. Reggio, P.H., editor, *Humana* (2009).
- Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-15 (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, 06/16/2023

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