

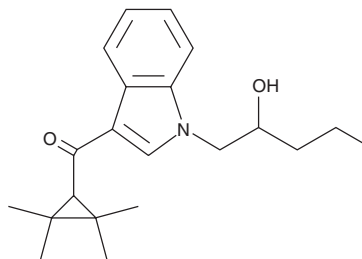
PRODUCT INFORMATION



UR-144 N-(2-hydroxypentyl) metabolite

Item No. 14470

CAS Registry No.: 2748288-94-8
Formal Name: (1-(2-hydroxypentyl)-1H-indol-3-yl)
(2,2,3,3-tetramethylcyclopropyl)methanone
MF: C₂₁H₂₉NO₂
FW: 327.5
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 256, 299 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

UR-144 N-(2-hydroxypentyl) isomer is a synthetic cannabinoid that differs from the parent compound, UR-144 (Item No. 11502), by the addition of a hydroxyl group to the alkyl chain. It is a potential metabolite of UR-144, based on the first pass metabolism of similar compounds by cytochrome P450 in the liver.^{1,2} The physiological and toxicological properties of this compound have not been studied. This product is intended for forensic and research applications.

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(5)**, 2141-53 (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, 06/19/2023

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