

PRODUCT INFORMATION

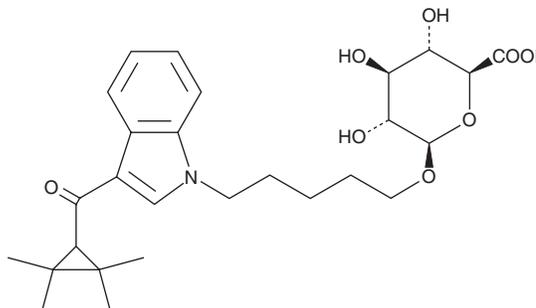


UR-144 N-(5-hydroxypentyl) β -D-Glucuronide

Item No. 11776

CAS Registry No.: 1630023-01-6
Formal Name: (2S,3S,4S,5R,6R)-3,4,5-trihydroxy-6-((5-(3-(2,2,3,3-tetramethylcyclopropane-1-carbonyl)-1H-indol-1-yl)pentyl)oxy)tetrahydro-2H-pyran-2-carboxylic acid

MF: C₂₇H₃₇NO₈
FW: 503.6
Purity: \geq 95%
UV/Vis.: λ_{max} : 215, 247, 305 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

UR-144 (Item No. 11502) is a potent synthetic cannabinoid (CB) which preferentially binds the peripheral CB₂ receptor ($K_i = 1.8$ nM) over the central CB₁ receptor ($K_i = 150$ nM).¹ UR-144 N-(5-hydroxypentyl) β -D-glucuronide is an expected major urinary metabolite of UR-144, based on the metabolism of similar cannabimimetics.^{2,3} The physiological and toxicological properties of this compound have not been tested. This product is intended for forensic and research applications.

References

1. Frost, J.M., Dart, M.J., Tietje, K.R., *et al.* Indol-3-ylcycloalkyl ketones: Effects of N1 substituted indole side chain variations on CB₂ cannabinoid receptor activity. *J. Med. Chem.* **53**, 295-315 (2010).
2. 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).
3. Chimalakonda, K.C., Moran, C.L., Kennedy, P.D., *et al.* Solid-phase extraction and quantitative measurement of ω and ω -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, 07/20/2023

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