

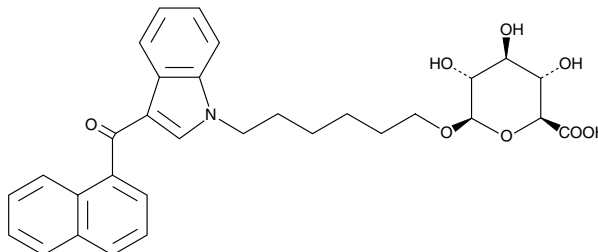
Product Information



JWH 019 N-(6-hydroxyhexyl) β -D-Glucuronide

Item No. 11466

CAS Registry No.: 1630023-00-5
Formal Name: (2S,3S,4S,5R,6R)-6-((6-(3-(1-naphthoyl)-1H-indol-1-yl)hexyloxy)-3,4,5-trihydroxytetrahydro-2H-pyran-2-carboxylic acid
MF: $C_{31}H_{33}NO_8$
FW: 547.6
Purity: $\geq 98\%$
Stability: ≥ 2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 218, 246, 314 nm



Laboratory Procedures

For long term storage, we suggest that JWH 019 N-(6-hydroxyhexyl) β -D-glucuronide be stored as supplied at -20°C . It should be stable for at least two years.

JWH 019 N-(6-hydroxyhexyl) β -D-glucuronide is supplied as a crystalline solid. A stock solution may be made by dissolving the JWH 019 N-(6-hydroxyhexyl) β -D-glucuronide in the solvent of choice. JWH 019 N-(6-hydroxyhexyl) β -D-glucuronide is soluble in DMSO. The solubility of JWH 019 N-(6-hydroxyhexyl) β -D-glucuronide in DMSO is approximately 10 mg/ml.

JWH 019 N-(6-hydroxyhexyl) β -D-glucuronide is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

JWH 019 (Item No. 13633) is a cannabimimetic indole that shows a high-affinity for both the central cannabinoid (CB_1) ($K_i = 9.8$ nM) and the peripheral cannabinoid (CB_2) ($K_i = 5.6$ nM) receptors.¹ JWH 019 N-(6-hydroxyhexyl) β -D-glucuronide is an expected major urinary metabolite of JWH 019, based on the metabolism of similar cannabimimetics.²⁻⁴ The physiological and toxicological properties of this compound have not been tested. This product is intended for forensic and research 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 CB_1 and CB_2 receptor binding. *Drug Alcohol Depend.* **60**, 133-140 (2000).
2. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200**, 141-147 (2010).
3. 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).
4. 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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/11466

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 11/20/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com