

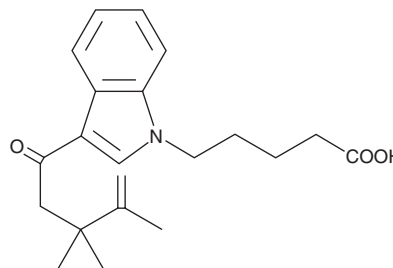
# PRODUCT INFORMATION



## UR-144 Degradant N-pentanoic acid metabolite

Item No. 9001453

**CAS Registry No.:** 1630022-97-7  
**Formal Name:** 3-(3,3,4-trimethyl-1-oxo-4-penten-1-yl)-1H-indole-1-pentanoic acid  
**MF:** C<sub>21</sub>H<sub>27</sub>NO<sub>3</sub>  
**FW:** 341.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 212, 246, 304 nm  
**Supplied as:** A solution in methanol  
**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 (Item No. 11502) is a potent synthetic cannabinoid (CB) that preferentially binds the peripheral CB<sub>2</sub> receptor (K<sub>i</sub> = 1.8 nM) over the central CB<sub>1</sub> receptor (K<sub>i</sub> = 150 nM).<sup>1</sup> UR-144 Degradant N-pentanoic acid metabolite (Item No. 9001453) is a potential phase I metabolite of a degradation product observed during GC-MS analysis of samples containing UR-144.<sup>2-4</sup> The biological activities of this compound have not been characterized. 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 CB<sub>2</sub> cannabinoid receptor activity. *J. Med. Chem.* **53(1)**, 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(5)**, 2141-2153 (2010).
3. 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).
4. Lovett, D.P., Yanes, E.G., Herbelin, T.W., *et al.* Structure elucidation and identification of a common metabolite for naphthoylindole-based synthetic cannabinoids using LC-TOF and comparison to a synthetic reference standard. *Forensic Sci. Int.* **226(1-3)**, 81-87 (2013).

#### 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, 07/11/2023

#### 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