

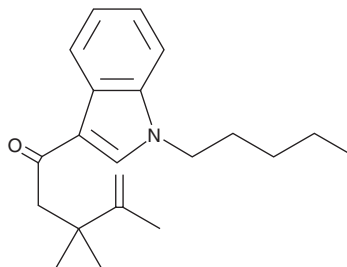
PRODUCT INFORMATION



UR-144 Degradant

Item No. 11928

CAS Registry No.: 1609273-88-2
Formal Name: 3,3,4-trimethyl-1-(1-pentyl-1H-indol-3-yl)pent-4-en-1-one
Synonyms: KM-X1 Degradant, UR-144
3,3,4-Trimethylpentenoyl isomer
MF: C₂₁H₂₉NO
FW: 311.5
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 247, 305 nm
Supplied as: A 10 mg/ml 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 synthetic cannabinoid (CB) analog of JWH 018 (Item No. 10900) in which the naphthalene ring is substituted with a tetramethylcyclopropyl group. This substitution results in a compound which is a potent agonist for the peripheral CB₂ receptor.¹ UR-144 Degradant is a common impurity observed during GC-MS analysis of samples containing UR-144. The opened ring of the degradant is presumed to be produced during heating of UR-144. This structure gives rise to a prominent fragment ion that is 15 amu greater than the base peak of UR-144. This pattern is consistent with McLafferty rearrangement of the degradant which does not occur with the parent compound.²

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. McLafferty, F.W. Mass spectrometric analysis. Molecular rearrangements. *Anal. Chem.* **31**(1), 82-87 (1959).

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, 11/21/2025

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