

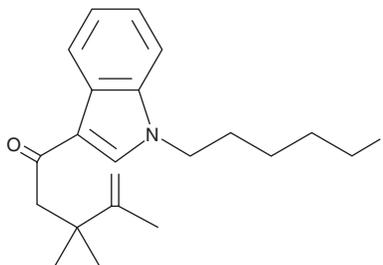
PRODUCT INFORMATION



XLR11 Degradant

Item No. 14055

CAS Registry No.: 1616469-09-0
Formal Name: 1-[1-(5-fluoropentyl)-1H-indol-3-yl]-3,3,4-trimethyl-4-penten-1-one
Synonyms: 5-FUR-144 Degradant, TCMP-2201 thermal isomer
MF: C₂₁H₂₈FNO
FW: 329.5
Purity: ≥98%
UV/Vis.: λ_{max}: 247, 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

XLR11 (Item No. 11565) is a synthetic cannabinoid (CB) which contains a tetramethylcyclopropyl group. In many synthetic CBs, this group serves to increase the affinity of the compound for the peripheral CB₂ receptor.¹ XLR11 degradant is a common impurity observed during GC-MS analysis of samples containing XLR11. The opened ring of the degradant is presumed to be produced during heating of XLR11. This structure gives rise to a prominent fragment ion that is 15 amu greater than the base peak of XLR11. 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, 07/25/2023

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