

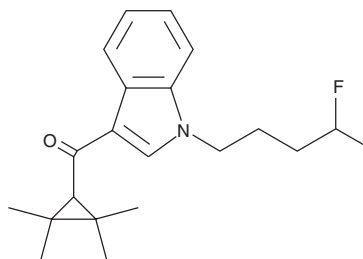
PRODUCT INFORMATION



XLR11 N-(4-fluoropentyl) isomer

Item No. 11769

CAS Registry No.: 2365471-50-5
Formal Name: (1-(4-fluoropentyl)-1H-indol-3-yl)
(2,2,3,3-tetramethylcyclopropyl)
methanone
MF: C₂₁H₂₈FNO
FW: 329.5
Purity: ≥95%
UV/Vis.: λ_{max}: 216, 246, 304 nm
Supplied as: A 1 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

XLR11 (Item No. 11565) is a synthetic cannabinoid (CB) featuring a tetramethylcyclopropyl group, which reportedly confers selectivity for the peripheral CB₂ receptor over the central CB₁ receptor.¹ XLR11 also has an N-(5-fluoropentyl) chain, which increases binding to both CB receptors.² XLR11 N-(4-fluoropentyl) isomer is an isomer of XLR11 which has the fluorine atom placed on the 4 position, rather than the terminal 5 position, of the alkyl group. The physiological and toxicological properties of this compound have not been evaluated. This product is intended for forensic and research purposes.

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. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832.

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, 11/20/2025

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