

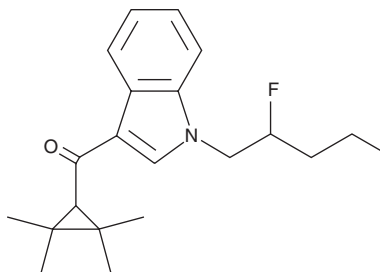
# PRODUCT INFORMATION



## XLR11 N-(2-fluoropentyl) isomer

Item No. 11767

**CAS Registry No.:** 1628690-25-4  
**Formal Name:** [1-(2-fluoropentyl)-1H-indol-3-yl]  
(2,2,3,3-tetramethylcyclopropyl)-methanone  
**MF:** C<sub>21</sub>H<sub>28</sub>FNO  
**FW:** 329.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 216, 244, 301 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) featuring a tetramethylcyclopropyl group, which reportedly confers selectivity for the peripheral CB<sub>2</sub> receptor over the central CB<sub>1</sub> receptor.<sup>1</sup> XLR11 also has an N-(5-fluoropentyl) chain, which increases binding to both CB receptors.<sup>2</sup> XLR11 N-(2-fluoropentyl) isomer differs structurally from XLR11 by having fluorine at the 2 position rather than the 5 position of the pentyl chain. The biological and toxicological properties of this compound have not been evaluated. 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**, 295-315 (2010).
2. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. *University of Connecticut*. **WO 01/28557 A1** (2001).

#### 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, 07/18/2023

#### 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](http://WWW.CAYMANCHEM.COM)