

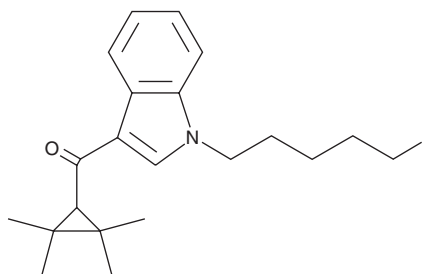
# PRODUCT INFORMATION



## XLR11 (exempt preparation)

Item No. 15033

**CAS Registry No.:** 1364933-54-9  
**Formal Name:** (1-(5-fluoropentyl)-1H-indol-3-yl)  
(2,2,3,3-tetramethylcyclopropyl)methanone  
**Synonyms:** 5-fluoro UR-144, 5-FUR-144  
**MF:** C<sub>21</sub>H<sub>28</sub>FNO  
**FW:** 329.5  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A solution in methanol  
**UV/Vis.:** λ<sub>max</sub>: 217, 246, 304 nm



### Description

XLR11 (exempt preparation) (Item No. 15033) is an analytical reference standard characterized as a synthetic cannabinoid.<sup>1</sup> XLR11 is regulated as a Schedule I compound in the United States. XLR11 (exempt preparation) (Item No. 15033) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

### Reference

1. Banister, S.D., Stuart, J., Kevin, R.C., *et al.* Effects of bioisosteric fluorine in synthetic cannabinoid designer drugs JWH-018, AM-2201, UR-144, XLR-11, PB-22, 5F-PB-22, APICA, and STS-135. *ACS Chem. Neurosci.* **6**(8), 1445-1458 (2015).

#### 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, 09/13/2015

#### 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