

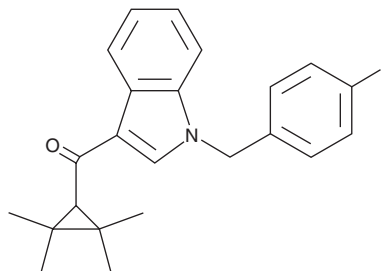
PRODUCT INFORMATION



FUB-144

Item No. 15537

CAS Registry No.: 2185863-15-2
Formal Name: [1-[(4-fluorophenyl)methyl]-1H-indol-3-yl]
(2,2,3,3-tetramethylcyclopropyl)-methanone
Synonym: FUB-UR-144
MF: C₂₃H₂₄FNO
FW: 349.5
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 245, 302 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 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 It also features an N-(5-fluoropentyl) chain, which increases binding to both CB receptors.² FUB-144 is a derivative of XLR11 in which the 5-fluoropentyl side chain is replaced by a 4-fluorobenzyl group, a functional group also found in the indazole-based synthetic cannabinoids AB-FUBINACA (Item No. 14292) and ADB-FUBINACA (Item No. 14039).^{3,4} The physiological and toxicological properties of this compound are not known. FUB-144 is regulated as a Schedule I compound in the United States. This product is intended for research and forensic 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 CB₂ cannabinoid receptor activity. *J. Med. Chem.* **53**(1), 295-315 (2010).
2. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. *University of Connecticut WO01/28557A1* (2001).
3. Uchiyama, N., Matsuda, S., Wakana, D., *et al.* New cannabimimetic indazole derivatives, N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-pentyl-1H-indazole-3-carboxamide (AB-PINACA) and N-(1-amino-3-methyl-1-oxobutan-2-yl)-1-(4-fluorobenzyl)-1H-indazole-3-carboxamide (AB-FUBINACA) identified as designer drugs in illegal products. *Forensic Toxicol.* **31**(1), 93-100 (2013).
4. Buchler, I.P., Hayes, M.J., Hedge, S.G., *et al.* Indazole derivatives. *Pfizer Inc. WO2009/106982A1* (2009).

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, 06/28/2023

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