

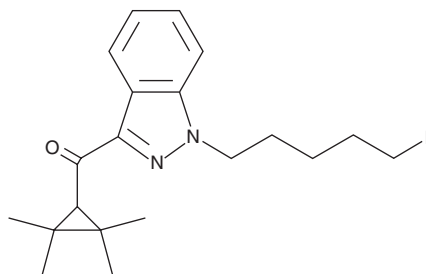
PRODUCT INFORMATION



FAB-144

Item No. 15155

CAS Registry No.: 2180935-79-7
Formal Name: [1-(5-fluoropentyl)-1H-indazol-3-yl]
(2,2,3,3-tetramethylcyclopropyl)-methanone
MF: C₂₀H₂₇FN₂O
FW: 330.4
Purity: ≥98%
UV/Vis.: λ_{max}: 241, 247, 307 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) with a tetramethylcyclopropyl group, which reportedly confers selectivity for the peripheral CB₂ receptor over the central CB₁ receptor.¹ FAB-144 is an analog of XLR11 that has an indazole, rather than an indole group. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

Reference

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).

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, 06/13/2023

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