

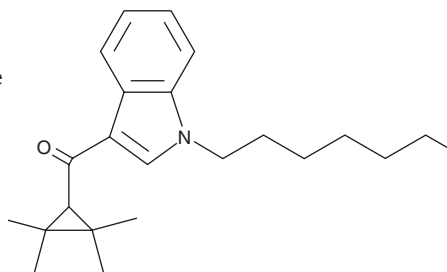
PRODUCT INFORMATION



UR-144 N-heptyl analog

Item No. 14231

CAS Registry No.: 1616469-06-7
Formal Name: (1-heptyl-1H-indol-3-yl)
(2,2,3,3-tetramethylcyclopropyl)-methanone
MF: C₂₃H₃₃NO
FW: 339.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 247, 304 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

UR-144 (Item No. 11502) is a potent synthetic cannabinoid (CB) which preferentially binds the peripheral CB₂ receptor (K_i = 1.8 nM) over the central CB₁ receptor (K_i = 150 nM).¹ UR-144 N-heptyl analog (Item No. 14231) differs from UR-144 by having two additional carbons in the alkyl group. The biological activities of this compound have not been characterized. 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, 06/15/2023

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