

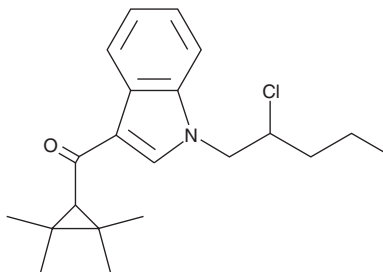
PRODUCT INFORMATION



UR-144 N-(2-chloropentyl) analog

Item No. 13115

CAS Registry No.: 2365471-32-3
Formal Name: (1-(2-chloropentyl)-1H-indol-3-yl)
(2,2,3,3-tetramethylcyclopropyl)
methanone
MF: C₂₁H₂₈ClNO
FW: 345.9
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 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

UR-144 N-(2-chloropentyl) analog is a synthetic cannabinoid which differs from the parent compound, UR-144 (Item No. 11502), by the addition of chloride to the 2 position of the pentyl chain. The physiological and toxicological properties of this compound have not been studied. This product is intended for forensic and research applications.

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, 07/25/2023

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