

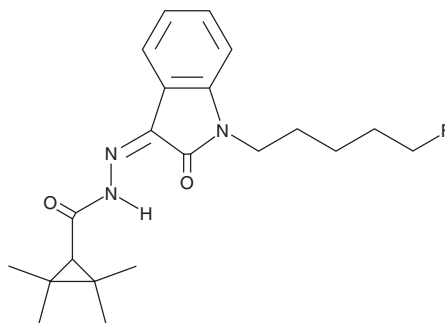
PRODUCT INFORMATION



Azidoindolene 1

Item No. 17083

CAS Registry No.: 1364933-69-6
Formal Name: 2,2,3,3-tetramethylcyclopropanecarboxylic acid (2Z)-2-[1-(5-fluoropentyl)-1,2-dihydro-2-oxo-3H-indol-3-ylidene]hydrazide
Synonym: F-TMCPRO-POXIZID
MF: 373.5
FW: C₂₁H₂₈FN₃O₂
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 250, 272, 324 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) and XLR11 (Item No. 11565) are potent synthetic cannabinoids (CBs) that have been identified as adulterants of herbal products.^{1,2} Azidoindolene 1 is structurally similar to UR-144, XLR11, and a number of additional synthetic CBs with a hydrazide linking the tetramethylcyclopropyl group to the aminoalkylindole group. The physiological and toxicological properties of this compound are not known. Typically, a tetramethylcyclopropyl group confers selectivity for the peripheral CB₂ receptor, and the addition of substituents at the N1-amine of the aminoalkylindole group is necessary for high affinity at either CB₁ or CB₂.³ This product is intended for forensic and research applications.

References

1. Uchiyama, N., Kawamura, M., Kikura-Hanajiri, R., *et al.* URB-754: A new class of designer drug and 12 synthetic cannabinoids detected in illegal products. *Forensic Sci. Int.* **227(1-3)**, 21-32 (2013).
2. Uchiyama, N., Matsuda, S., Kawamura, M., *et al.* Two new-type cannabimimetic quinolinyl carboxylates, QUPIC and QUCHIC, two new cannabimimetic carboxamide derivatives, ADB-FUBINACA and ADBICA, and five synthetic cannabinoids detected with a thiophene derivative a-PVT and an opioid receptor agonist AH-7921 identified in illegal products. *Forensic Toxicol.* **31(2)**, 223-240 (2013).
3. 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**, 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, 06/16/2023

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