

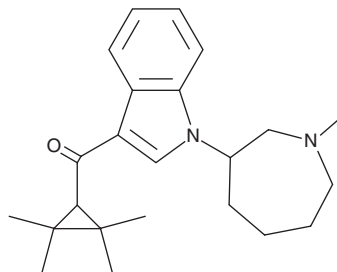
PRODUCT INFORMATION



AB-005 azepane isomer

Item No. 11877

CAS Registry No.: 1445751-74-5
Formal Name: [1-(hexahydro-1-methyl-1H-azepin-3-yl)-1H-indol-3-yl] (2,2,3,3-tetramethylcyclopropyl)-methanone
MF: C₂₃H₃₂N₂O
FW: 352.5
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

AB-005 (Item No. 11766) is a synthetic cannabinoid (CB) built on a 1-[(1-methylpiperidin-2-yl)methyl]-indole base that is characteristic of a series of potent CBs, including AM1220 (Item No. 9001055), AM1241 (Item No. 10010118), and AM1248 (Item No. 11282).¹ AB-005 azepane isomer is a variant of AB-005 in which the methylpiperidine group has been replaced with methylazepane. The physiological and toxicological properties of this compound have not been established. This product is intended for forensic and research applications.

Reference

1. Makriyannis, A. and Deng, H. Receptor selective cannabimimetic aminoalkylindoles. *Alix Yale & Ristas, LLP*. US2007024313134A1 (2007).

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, 07/24/2023

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