

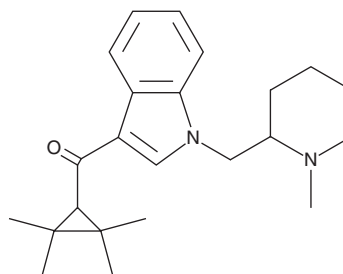
# PRODUCT INFORMATION



**AB-005**

Item No. 11766

**CAS Registry No.:** 895155-25-6  
**Formal Name:** [1-[(1-methyl-2-piperidiny)methyl]-1H-indol-3-yl](2,2,3,3-tetramethylcyclopropyl)-methanone  
**MF:** C<sub>23</sub>H<sub>32</sub>N<sub>2</sub>O  
**FW:** 352.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 246, 303 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 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).<sup>1</sup> The addition of a tetramethylcyclopropyl group has, in some cases, increased selectivity of synthetic CBs for the peripheral CB<sub>2</sub> receptor.<sup>2</sup> The physiological and toxicological properties of this compound have not been established. This product is intended for forensic and research applications.

## References

1. Makriyannis, A. and Deng, H. Receptor selective cannabimimetic aminoalkylindoles. *Alix Yale & Ristas, LLP. US2007024313134A1* (2007).
2. Poso, A. and Huffman, J.W. Targeting the cannabinoid CB<sub>2</sub> receptor: Modelling and structural determinants of CB<sub>2</sub> selective ligands. *Br. J. Pharmacol.* **153**(2), 335-346 (2008).

### 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, 07/18/2023

## 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](http://WWW.CAYMANCHEM.COM)