

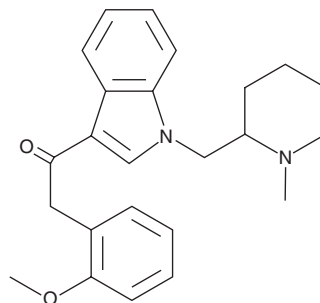
PRODUCT INFORMATION



Cannabipiperidiethanone

Item No. 11655

CAS Registry No.: 1345970-43-5
Formal Name: 2-(2-methoxyphenyl)-1-[1-[(1-methyl-2-piperidinyl)methyl]-1H-indol-3-yl]-ethanone
MF: C₂₄H₂₈N₂O₂
FW: 376.5
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 245, 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

Cannabipiperidiethanone (Item No. 11655) is a synthetic cannabinoid (CB) with structural similarities to JWH 250 (Item No. 13634) and AM2233 (Item No. 11008) that has been identified as a component of “herbal” blends illegally sold in Japan.¹ Cannabipiperidiethanone binds to the CB receptors CB₁ and CB₂ with IC₅₀ values of 591 and 968 nM, respectively.¹ Its reported affinities for CB₁ and CB₂ receptors are 2.3- and 9.4-fold lower than those of JWH 250, and 13- and 70-fold lower than those of (R)-(+)-WIN 55,212-2, respectively.¹ This product is intended for forensic and research applications.

Reference

1. Uchiyama, N., Kikura-Hanajiri, R., and Goda, Y. Identification of a novel cannabimimetic phenylacetylindole, cannabipiperidiethanone, as a designer drug in a herbal product and its affinity for cannabinoid CB₁ and CB₂ receptors. *Chem. Pharm. Bull. (Tokyo)* **59(9)**, 1203-1205 (2011).

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