

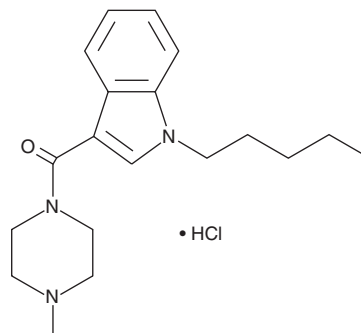
PRODUCT INFORMATION



Mepirapim (hydrochloride)

Item No. 15388

CAS Registry No.: 2365542-30-7
Formal Name: (4-methyl-1-piperazinyl)(1-pentyl-1H-indol-3-yl)-methanone, monohydrochloride
Synonym: JWH 018-4(methylpiperazine)
MF: C₁₉H₂₇N₃O • HCl
FW: 349.9
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 289 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

JWH 018 (Item No. 10900) is a potent synthetic cannabinoid (CB) that avidly binds both the central CB₁ receptor and the peripheral CB₂ receptor (K_is = 9.0 and 2.9 nM, respectively).¹ Structurally, JWH 018 is an aminoalkylindole with a naphthoyl group. Mepirapim is an analog of JWH 018 that has been identified in herbal blends in Japan.² It differs from JWH 018 by having a 4-methylpiperazine group in place of the naphthyl group. The physiological and toxicological properties of this compound are unknown. This product is intended for forensic and research applications.

References

1. Aung, M.M., Griffin, G., Huffman, J.W., *et al.* Influence of the N-1 alkyl chain length of cannabimimetic indoles upon CB₁ and CB₂ receptor binding. *Drug Alcohol Depend.* **60(2)**, 133-140 (2000).
2. Uchiyama, N., Shimokawa, Y., Matsuda, S., *et al.* Two new synthetic cannabinoids, AM-2201 benzimidazole analog (FUBIMINA) and (4-methylpiperazin-1-yl)(1-pentyl-1H-indol-3-yl)methanone (MEPIRAPIM), and three phenethylamine derivatives, 25H-NBOMe 3,4,5-trimethoxybenzyl analog, 25B-NBOMe, and 2C-N-NBOMe, identified in illegal products. *Forensic Toxicol.* (2013).

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, 06/27/2023

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