

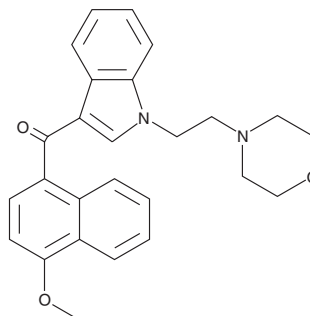
PRODUCT INFORMATION



JWH 198

Item No. 13981

CAS Registry No.: 166599-76-4
Formal Name: (4-methoxy-1-naphthalenyl)
[1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl]-methanone
MF: C₂₆H₂₆N₂O₃
FW: 414.5
Purity: ≥98%
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 198 is a synthetic cannabinoid (CB) which binds the central CB₁ receptor with high affinity (K_i = 10 nM).¹ This aminoalkylindole shares structural features with the antinociceptive CB₁ agonists pravadoline (Item No. 10006973) and WIN 55,212-2.¹ The physiological and toxicological properties of this compound have not been investigated. This product is intended for research and forensic applications.

Reference

1. Huffman, J.W. and Padgett, L.W. Recent developments in the medicinal chemistry of cannabimimetic indoles, pyrroles and indenenes. *Curr. Med. Chem.* **12**(12), 1395-1411 (2005).

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, 07/14/2023

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