

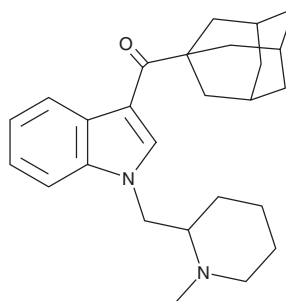
PRODUCT INFORMATION



AM1248

Item No. 11282

CAS Registry No.: 335160-66-2
Formal Name: [1-[(1-methyl-2-piperidinyl)methyl]-1H-indol-3-yl]tricyclo[3.3.1.1^{3,7}]dec-1-yl-methanone
MF: C₂₆H₃₄N₂O
FW: 390.6
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 242, 297 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

AM1248 is an adamantoylindole derivative with an N-methylpiperidin-2-ylmethyl substitution at the indole 1-position. It reportedly acts as a moderately potent agonist for both the central CB₁ and peripheral CB₂ cannabinoid receptors (K_is = 11.9 and 4.8 nM, respectively).¹ A related compound, 1-pentyl-3-(1-adamantoyl)indole, along with another synthetic cannabinoid, AM679, was identified as having been sold as a cannabinoid herbal blend known as “smart products” in Hungary in 2011.²

References

1. Makriyannis, A. and Deng, H. Receptor selective cannabimimetic aminoalkylindoles. US Patent 2007024313134A1 (2007), 1-16, 11/620,248.
2. Jankovics, P., Váradi, A., Tölgyesi, L., et al. Detection and identification of the new potential synthetic cannabinoids 1-pentyl-3-(2-iodobenzoyl)indole and 1-pentyl-3-(1-adamantoyl)indole in seized bulk powders in Hungary. *Forensic Sci. Int.* **214**, 27-32 (2012).

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