

PRODUCT INFORMATION



AM2232

Item No. 11503

CAS Registry No.: 335161-19-8

Formal Name: 3-(1-naphthalenylcarbonyl)-1H-Indole-1-pentanenitrile

MF: C₂₄H₂₀N₂O

FW: 352.4

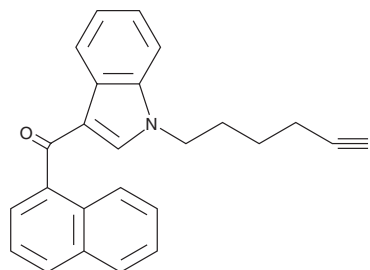
Purity: ≥95%

UV/Vis.: λ_{max}: 219, 311 nm

Supplied as: A 25 mg/ml solution in acetonitrile

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

AM2232 is a potent synthetic cannabinoid (CB) with K_i values of 0.28 and 1.48 nM for the CB₁ and CB₂ receptors, respectively.¹ The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

Reference

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832.

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, 11/21/2025

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