

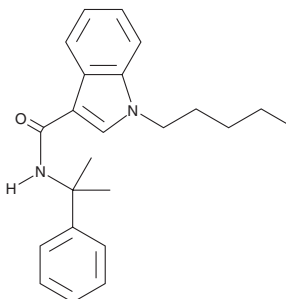
PRODUCT INFORMATION



CUMYL-PICA

Item No. 17211

CAS Registry No.: 1400742-32-6
Formal Name: N-(1-methyl-1-phenylethyl)-1-pentyl-1H-indole-3-carboxamide
MF: C₂₃H₂₈N₂O
FW: 348.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 290 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

AM2201 (Item No. 10707) is a potent synthetic cannabinoid (CB) developed on an indole base with K_i values of 1.0 and 2.6 nM for the central CB₁ and peripheral CB₂ receptors, respectively.¹ CUMYL-PICA is a derivative of AM2201 with a PINACA-like chain that lacks the alkyl-terminal fluorine atom and where the naphthoyl group has been modified with the addition of a benzyl group. The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

Reference

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. *University of Connecticut*. **WO01/28557A1** (2001).

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, 07/06/2023

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