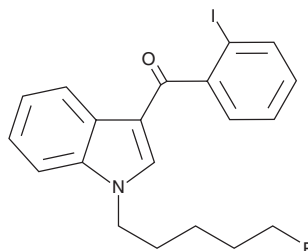


PRODUCT INFORMATION

AM694 (exempt preparation)

Item No. 15720

CAS Registry No.: 335161-03-0
 Formal Name: [1-(5-fluoropentyl)-1H-indol-3-yl]
 (2-iodophenyl)-methanone
 MF: C₂₀H₁₉FINO
 FW: 435.3
 Purity: ≥95%
 UV/Vis.: λ_{max}: 213, 251, 315 nm
 Supplied as: A solution in methanol
 Storage: -20°C
 Stability: ≥3 years



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

Description

AM694 is a potent synthetic cannabinoid (CB) with K_i values of 0.08 and 1.44 nM for the CB₁ and CB₂ receptors, respectively.¹ The physiological actions and metabolism of AM694 have not been characterized.

Reference

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. **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, 05/19/2025

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