

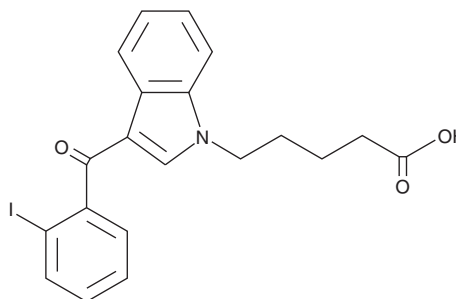
PRODUCT INFORMATION



AM694 N-pentanoic acid metabolite

Item No. 14359

CAS Registry No.: 1432900-96-3
Formal Name: 3-(2-iodobenzoyl)-1H-indole-1-pentanoic acid
MF: C₂₀H₁₈INO₃
FW: 447.3
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 251, 316 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

AM694 (Item No. 10567) is a potent synthetic cannabinoid (CB) with K_i values of 0.08 and 1.44 nM for the central CB₁ and peripheral CB₂ receptors, respectively.¹ AM694 N-pentanoic acid metabolite is a potential phase I metabolite of AM694, based on the known metabolism of similar compounds.² The physiological and toxicological properties of this compound have not been evaluated. This product is intended for forensic and research applications.

References

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. *University of Connecticut*. **WO01/28557A1** (2001).
2. Wintermeyer, A., Möller, I., Thevis, M., *et al.* In vitro phase I metabolism of the synthetic cannabimimetic JWH-018. *Anal. Bioanal. Chem.* **398(5)**, 2141-2153 (2010).

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, 06/16/2023

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