

PRODUCT INFORMATION

ADB-HEXINACA N-(6-hydroxyhexyl) metabolite

Item No. 35566

Formal Name: (S)-N-(1-amino-3,3-dimethyl-1-oxobutan-2-yl)-1-(6-hydroxyhexyl)-1H-indazole-3-carboxamide

MF: C₂₀H₃₀N₄O₃

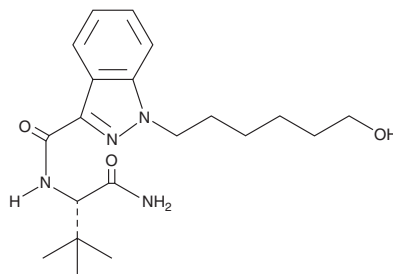
FW: 374.5

Purity: ≥98%

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

ADB-HEXINACA N-(6-hydroxyhexyl) metabolite (Item No. 35566) is an analytical reference standard that is categorized as a synthetic cannabinoid metabolite.¹ ADB-HEXINACA N-(6-hydroxyhexyl) metabolite is a metabolite of ABD-HEXINACA (Item No. 33820). This product is intended for research and forensic applications.

Reference

1. Baginski, S.R., Rautio, T., Nisbet, L.A., *et al.* The metabolic profile of the synthetic cannabinoid receptor agonist ADB-HEXINACA using human hepatocytes, LC-QTOF-MS and synthesized reference standards. *J. Anal. Toxicol.* (2023).

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, 09/27/2023

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