

PRODUCT INFORMATION

ADB-5'Br-PINACA N-(4-hydroxypentyl) metabolite

Item No. 40768

Formal Name: N-((S)-1-amino-3,3-dimethyl-1-oxobutan-2-yl)-5-bromo-1-(4-hydroxypentyl)-1H-indazole-3-carboxamide

MF: C₁₉H₂₇BrN₄O₃

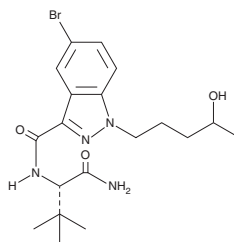
FW: 439.4

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Description

ADB-5'Br-PINACA N-(4-hydroxypentyl) metabolite (Item No. 40768) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. ADB-5'Br-PINACA N-(4-hydroxypentyl) metabolite is a potential metabolite of ADB-5'Br-PINACA (Item No. 37758) based on the published metabolism of ADB-PINACA (Item Nos. ISO60212 | ISO00150).¹ At the time ADB-5'Br-PINACA N-(4-hydroxypentyl) metabolite (Item No. 40768) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. Carlier, J., Diao, X., Scheidweiler, K.B., *et al.* Distinguishing intake of new synthetic cannabinoids ADB-PINACA and 5F-ADB-PINACA with human hepatocyte metabolites and high-resolution mass spectrometry. *Clin. Chem.* **63**(5), 1008-1021 (2017).

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, 01/13/2026

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