

PRODUCT INFORMATION

ADB-5'Br-BUTINACA 3,3-dimethyl butanoic acid metabolite

Item No. 37607

Formal Name: (S)-2-(5-bromo-1-butyl-1H-indazole-3-carboxamido)-3,3-dimethylbutanoic acid

Synonyms: ADB-5'Br-BINACA 3,3-dimethyl butanoic acid metabolite, ADB-BUT-5Br-INACA 3,3-dimethyl butanoic acid metabolite

MF: $C_{18}H_{24}BrN_3O_3$

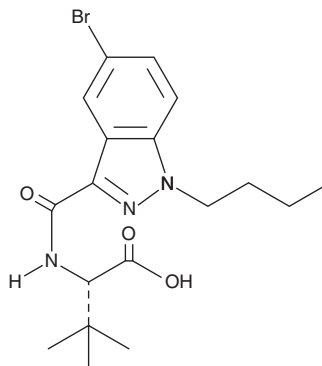
FW: 410.3

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 5 years



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

Description

ADB-5'Br-BUTINACA 3,3-dimethyl butanoic acid metabolite (Item No. 37607) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. ADB-5'Br-BUTINACA 3,3-dimethyl butanoic acid metabolite is a potential metabolite of ADB-5'Br-BUTINACA (Item No. 36770) based on the published metabolism of ADB-BUTINACA (Item No. 29350).¹ At the time ADB-5'Br-BUTINACA 3,3-dimethyl butanoic acid metabolite (Item No. 37607) 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. Kronstrand, R., Norman, C., Vikingsson, S., *et al.* The metabolism of the synthetic cannabinoids ADB-BUTINACA and ADB-4en-PINACA and their detection in forensic toxicology casework and infused papers seized in prisons. *Drug Test. Anal.* **14**(4), 634-652 (2021).

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, 07/07/2023

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