

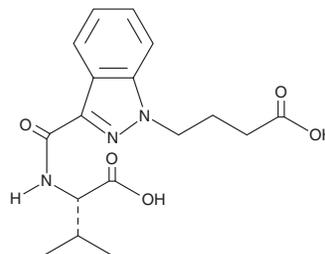
PRODUCT INFORMATION



4-cyano MMB-BUTINACA N-butanoic acid 3-methylbutanoic acid metabolite

Item No. 33836

Formal Name: (1-(3-carboxypropyl)-1H-indazole-3-carbonyl)-L-valine
Synonym: MMB-4CN-BUTINACA ester hydrolysis N-butanoic acid metabolite
MF: C₁₇H₂₁N₃O₅
FW: 347.4
Purity: ≥98%
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

4-cyano MMB-BUTINACA N-butanoic acid 3-methylbutanoic acid metabolite (Item No. 33836) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. 4-cyano MMB-BUTINACA N-butanoic acid 3-methylbutanoic acid metabolite is a potential metabolite of 4-cyano MMB-BUTINACA N-butanoic acid metabolite (Item No. 33835) and 4-cyano MMB-BUTINACA (Item No. 33334) based on the published metabolism of 5-fluoro AMB (Item No. 15489) and 4-cyano CUMYL-BUTINACA (Item No. 20194).^{1,2} At the time 4-cyano MMB-BUTINACA N-butanoic acid 3-methylbutanoic acid metabolite (Item No. 33836) 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.

References

1. Haschimi, B., Mogler, L., Halter, S., *et al.* Detection of the recently emerged synthetic cannabinoid 4F-MDMB-BINACA in "legal high" products and human urine specimens. *Drug Test Anal.* **11(9)**, 1377-1386 (2019).
2. Åstrand, A., Vikingsson, S., Lindstedt, D., *et al.* Metabolism study for CUMYL-4CN-BINACA in human hepatocytes and authentic urine specimens: Free cyanide is formed during the main metabolic pathway. *Drug Test Anal.* **10(8)**, 1270-1279 (2018).

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, 06/23/2023

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