

PRODUCT INFORMATION



4-cyano CUMYL-BUTINACA metabolite 10

Item No. 26399

Formal Name: 4-(3-((2-phenylpropan-2-yl)carbamoyl)-1H-indazol-1-yl)butanoic acid

Synonyms: CUMYL-4CN-BINACA N-butanoic Acid metabolite, CUMYL-CB-PINACA N-butanoic Acid metabolite, CUMYL-CYBINACA N-butanoic acid metabolite

MF: $C_{21}H_{23}N_3O_3$

FW: 365.4

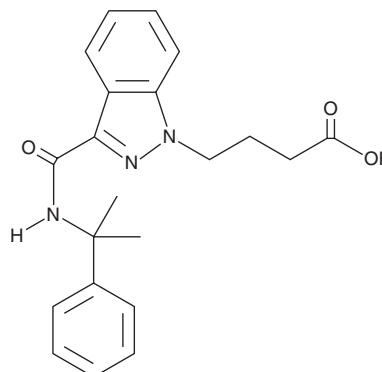
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 210, 303 nm

Supplied as: A solution in methyl acetate

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 CUMYL-BUTINACA metabolite 10 (Item No. 26399) is an analytical reference standard categorized as a synthetic cannabinoid metabolite.^{1,2} It is a metabolite of 4-cyano CUMYL-BUTINACA (Item No. 20194). This product is intended for research and forensic applications.

References

1. Staheli, S.N., Poetzsch, M., Veloso, V.P., *et al.* *In vitro* metabolism of the synthetic cannabinoids CUMYL-PINACA, 5F-CUMYL-PINACA, CUMYL-4CN-BINACA, 5F-CUMYL-P7AICA and CUMYL-4CN-B7AICA. *Drug Test Anal.* **10**(1), 148-157 (2018).
2. Öztürk, Y.E., Yeter, O., Öztürk, S., *et al.* Detection of metabolites of the new synthetic cannabinoid CUMYL-4CN-BINACA in authentic urine samples and human liver microsomes using high-resolution mass spectrometry. *Drug Test Anal.* **10**(3), 449-459 (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, 07/03/2023

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