

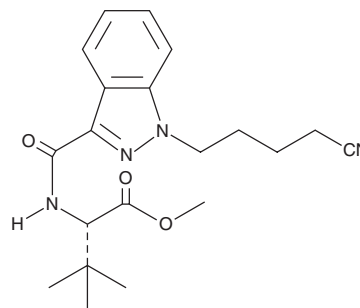
PRODUCT INFORMATION



4-cyano MDMB-BUTINACA

Item No. 26507

CAS Registry No.: 1185888-30-5
Formal Name: N-[[1-(4-cyanobutyl)-1H-indazol-3-yl]carbonyl]-3-methyl-L-valine, methyl ester
Synonym: 4CN-BINACA-ADB
MF: C₂₀H₂₆N₄O₃
FW: 370.5
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 302 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥5 years



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

Description

4-cyano MDMB-BUTINACA (Item No. 26507) is an analytical reference standard categorized as a synthetic cannabinoid.¹ This product is intended for research and forensic applications.

Reference

1. Buchler, I.P., Hayes, M.J., Hedge, S.G., et al. Indazole derivatives. *Pfizer Inc. WO 2009/106982 A1* (2009).

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