

PRODUCT INFORMATION



4-cyano CUMYL-BUT7AICA

Item No. 22818

CAS Registry No.: 2160555-53-1
Formal Name: 1-(4-cyanobutyl)-N-(1-methyl-1-phenylethyl)-1H-pyrrolo[2,3-b]pyridine-3-carboxamide

Synonyms: 4CN-CUMYL-B7AICA,
4CN-CUMYL-BUT7AICA,
CUMYL-4CN-B7AICA

MF: C₂₂H₂₄N₄O

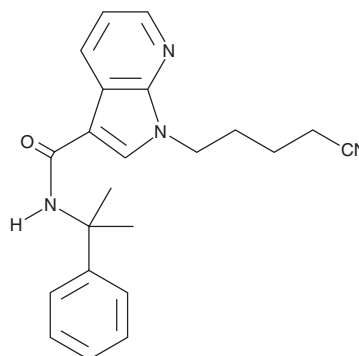
FW: 360.5

Purity: ≥98%

Supplied as: A neat solid

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 CUMYL-BUT7AICA (Item No. 22818) is an analytical reference material that is structurally similar to known synthetic cannabinoids. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

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, 08/28/2019

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