

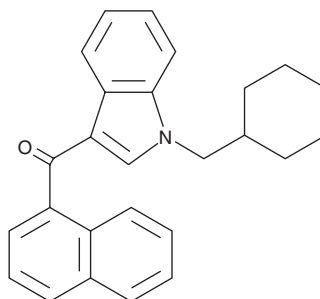
PRODUCT INFORMATION



NE-CHMIMO

Item No. 29301

CAS Registry No.: 1373876-11-9
Formal Name: [1-(cyclohexylmethyl)-1H-indol-3-yl]-1-naphthalenyl-methanone
Synonym: CHM-018
MF: C₂₆H₂₅NO
FW: 367.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 248, 316 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

NE-CHMIMO (Item No. 29301) is an analytical reference standard categorized as a synthetic cannabinoid.¹ It has been detected in human hair samples tested for verification of drug abstinence.² This product is intended for research and forensic applications.

References

1. Angerer, V., Bisel, P., Moosmann, B., *et al.* Separation and structural characterization of the new synthetic cannabinoid JWH-018 cyclohexyl methyl derivative NE-CHMIMO using flash chromatography, GC-MS, IR and NMR spectroscopy. *Forensic Sci. Int.* **266**, e93-e98 (2016).
2. Franz, F., Jechle, H., Angerer, V., *et al.* Synthetic cannabinoids in hair - pragmatic approach for method updates, compound prevalences and concentration ranges in authentic hair samples. *Anal. Chim. Acta.* **1006**, 61-73 (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, 07/28/2023

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