

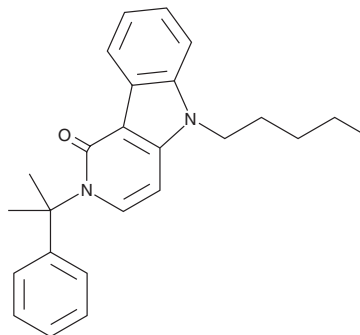
PRODUCT INFORMATION



CUMYL-PeGACLONE

Item No. 22696

CAS Registry No.: 2160555-55-3
Formal Name: 2,5-dihydro-2-(1-methyl-1-phenylethyl)-5-pentyl-1H-pyrido[4,3-b]indol-1-one
Synonym: SGT-151
MF: C₂₅H₂₈N₂O
FW: 372.5
Purity: ≥98%
UV/Vis.: λ_{max}: 252 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

CUMYL-PeGACLONE (Item No. 22696) is an analytical reference standard categorized as a synthetic cannabinoid.¹ It has been found in Spice-like herbal blends.² CUMYL-PeGACLONE is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Angerer, V., Mogler, L., Steitz, J.P., *et al.* Structural characterization and pharmacological evaluation of the new synthetic cannabinoid CUMYL-PEGACLONE. *Drug Test Anal.* (2017).
2. Ernst, L., Brandhorst, K., Papke, U., *et al.* Identification and quantification of synthetic cannabinoids in 'spicelike' herbal mixtures: Update of the German situation in early 2017. *Forensic Sci. Int.* **277**, 51-58 (2017).

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, 12/14/2023

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