

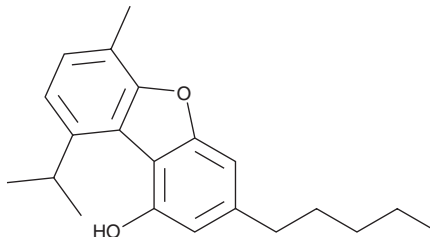
PRODUCT INFORMATION



Cannabifuran

Item No. 21108

CAS Registry No.: 56154-58-6
Formal Name: 6-methyl-9-(1-methylethyl)-3-pentyl-1-dibenzofuranol
Synonym: CBF
MF: C₂₁H₂₆O₂
FW: 310.4
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 232 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

Cannabifuran (Item No. 21108) is an analytical reference standard that is structurally similar to known phytocannabinoids. Cannabifuran has been found in *Cannabis* and *Cannabis* vape cartridges.^{1,2} This product is intended for research and forensic applications.

References

1. Lewis, M.M., Yang, Y., Wasilewski, E., *et al.* Chemical profiling of medical cannabis extracts. *ACS Omega* **2**(9), 6091-6103 (2017).
2. Guo, W., Vrdolijak, G., Liao, V.-C., *et al.* Major constituents of Cannabis vape oil liquid, vapor and aerosol in California vape oil cartridge samples. *Front. Chem.* **9**, 694905 (2021).

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, 02/14/2025

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