

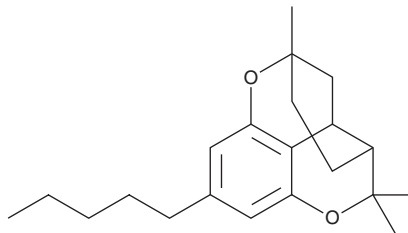
PRODUCT INFORMATION



Cannabicitran

Item No. 37475

CAS Registry No.: 31508-71-1
Formal Name: (6 α ,9 β ,10 α)-6a,7,8,9,10,10a-hexahydro-6,6,9-trimethyl-3-pentyl-1,9-epoxy-6H-dibenzo[b,d]pyran
Synonyms: CBT, CBT (citrin)
MF: C₂₁H₃₀O₂
FW: 314.5
Purity: \geq 98%
Supplied as: A solution in methanol
Storage: -20°C
Stability: \geq 5 year



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

Description

Cannabicitran (Item No. 37475) is an analytical reference standard that is structurally similar to known phytocannabinoids. Cannabicitran has been found in *Cannabis* and marijuana smoke condensate.^{1,2} This product is intended for research and forensic applications.

References

1. Bercht, C.A.L., Lousberg, R.J.J.C., Kupperts, F.J.E.M., *et al.* Cannabicitran: A new naturally occurring tetracyclic diether from Lebanese *Cannabis sativa*. *Phytochemistry* **13**(3), 619-621 (1974).
2. Novotný, M., Merli, F., Wiesler, D., *et al.* Fractionation and capillary gas chromatographic-mass spectrometric characterization of the neutral components in marijuana and tobacco smoke condensates. *J. Chromatogr.* **238**(1), 141-150 (1981).

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, 07/20/2023

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