

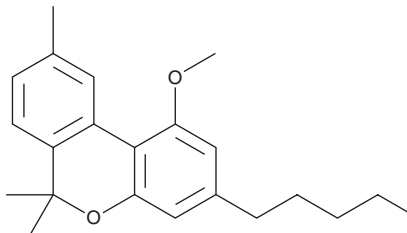
# PRODUCT INFORMATION



## Cannabinol methyl ether

Item No. 9002480

**CAS Registry No.:** 41935-92-6  
**Formal Name:** 1-methoxy-6,6,9-trimethyl-3-pentyl-6H-dibenzo[b,d]pyran  
**Synonym:** Cannabinol monomethyl ether, CBN monomethyl ether  
**MF:** C<sub>22</sub>H<sub>28</sub>O<sub>2</sub>  
**FW:** 324.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 218, 286 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Description

Cannabinol methyl ether (Item No. 9002480) is an analytical reference standard that is categorized as a phytocannabinoid. It can be isolated from *Cannabis* plants, derived from cannabinol (Item No. ISO60183), or synthesized.<sup>1-3</sup> The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

### References

1. Bercht, C.A., Lousberg, R.J., Küppers, F.J., *et al.* *Cannabis*. VII. Identification of cannabinol methyl ether from hashish. *J. Chromatogr.* **81(1)**, 163-166 (1973).
2. Lerner, M. Marihuana: Tetrahydrocannabinol and related compounds. *Science* **140(3563)**, 175-176 (1963).
3. Teske, J.A. and Deiters, A. A cyclotrimerization route to cannabinoids. *Org. Lett.* **10(11)**, 2195-2198 (2008).

#### 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, 07/25/2023

#### 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