

PRODUCT INFORMATION



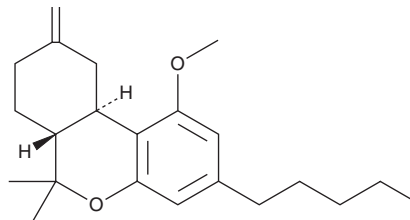
exo-THC methyl ether (exempt preparation)

Item No. 43654

CAS Registry No.: 27179-29-9

Formal Name: *trans*-6aR,7,8,9,10,10a-hexahydro-1-methoxy-6,6-dimethyl-9-methylene-3-pentyl-6H-dibenzo[b,d]pyran

Synonyms: (-)- Δ^7 -Tetrahydrocannabinol methyl ether,
 $\Delta^{9,11}$ -Tetrahydrocannabinol methyl ether,
 Δ^{11} -Tetrahydrocannabinol methyl ether,
(-)-*trans*- $\Delta^{9,11}$ -Tetrahydrocannabinol methyl ether,
exo-Tetrahydrocannabinol methyl ether,
(-)- Δ^7 -THC methyl ether,
(-)-*trans*- $\Delta^{9,11}$ -THC methyl ether



MF: $C_{22}H_{32}O_2$

FW: 328.5

Purity: $\geq 95\%$

Supplied as: A 1 mg/ml solution in acetonitrile

Storage: -20°C

Stability: ≥ 3 years

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

Description

exo-THC methyl ether (exempt preparation) (Item No. 43654) is an analytical reference standard that is structurally similar to known phytocannabinoids.¹ exo-THC methyl ether (exempt preparation) (Item No. 43654) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Magnuson, B.A., Roberts, A., and Nestmann, E.R. Critical review of the current literature on the safety of sucralose. *Food. Chem. Toxicol.* **106(Pt A)**, 324-355 (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, 04/25/2025

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