

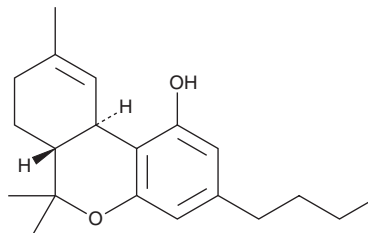
PRODUCT INFORMATION



Δ^9 -THCB

Item No. 29306

CAS Registry No.: 60008-00-6
Formal Name: *trans*-3-butyl-6aR,7,8,10a-tetrahydro-6,6,9-trimethyl-6H-dibenzo[b,d]pyran-1-ol
Synonyms: nor-THC, Δ^9 -Tetrahydrocannabinol-C4, Δ^9 -THC-butyl, Δ^9 -THC-C4, THCB
MF: C₂₀H₂₈O₂
FW: 300.4
Purity: ≥95%
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -80°C
Stability: ≥2 years



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

Description

Δ^9 -THCB (Item No. 29306) is an analytical reference standard categorized as a phytocannabinoid.¹⁻³ Δ^9 -THCB has been found in *Cannabis* samples seized by law enforcement. Δ^9 -THCB is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Smith, R.M. Identification of butyl cannabinoids in marijuana. *J. Forensic Sci.* **42(4)**, 610-618 (1997).
2. Harvey, D.J. Characterization of the butyl homologues of Δ^1 -tetrahydrocannabinol, cannabinol and cannabidiol in samples of cannabis by combined gas chromatography and mass spectrometry. *J. Pharm. Pharmacol.* **28(4)**, 280-285 (1976).
3. Linciano, P., Citti, C., Luongo, L., et al. Isolation of a high-affinity cannabinoid for the human CB₁ receptor from a medicinal *Cannabis sativa* variety: Δ^9 -Tetrahydrocannabinol, the butyl homologue of Δ^9 -tetrahydrocannabinol. *J. Nat. Prod.* **83(1)**, 88-98 (2019).

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, 01/05/2026

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