

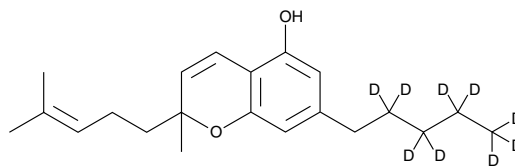
# Product Information



## (±)-Cannabichromene-d<sub>9</sub>

Item No. 16202

**Formal Name:** 2-methyl-2-(4-methyl-3-penten-1-yl)-7-pentyl-d<sub>9</sub>-2H-1-benzopyran-5-ol  
**Synonyms:** CBC-d<sub>9</sub>, Pentylcannabichromene-d<sub>9</sub>  
**MF:** C<sub>21</sub>H<sub>21</sub>D<sub>9</sub>O<sub>2</sub>  
**FW:** 323.5  
**Chemical Purity:** ≥98% (±)-Cannabichromene  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>9</sub>); ≤1% d<sub>0</sub>  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in methanol  
**UV/Vis.:** λ<sub>max</sub>: 229, 283 nm



### Laboratory Procedures

(±)-Cannabichromene-d<sub>9</sub> (CBC-d<sub>9</sub>) contains nine deuterium atoms on the pentyl chain. It is intended for use as an internal standard for the quantification of CBC (Item No. 9001880) by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that CBC-d<sub>9</sub> be stored as supplied at -20°C. It should be stable for at least one year.

CBC-d<sub>9</sub> is supplied as a solution in methanol. To change the solvent, simply evaporate the CBC-d<sub>9</sub> under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of CBC-d<sub>9</sub> in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

CBC-d<sub>9</sub> is used as an internal standard for the quantification of CBC by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

CBC is a non-psychoactive cannabinoid (CB) that has been shown to modulate the activity of proteins involved in nociceptive mechanisms through a non-CB mechanism of action.<sup>1</sup> It acts as an agonist of the transient receptor potential (TRP) channels of the vanilloid type-1 (EC<sub>50</sub> = 24.2 μM) and of the ankyrin type-1 (EC<sub>50</sub> = 90 nM) as well as inhibits proteins that facilitate anandamide inactivation (IC<sub>50</sub> = 12.3 μM).<sup>2</sup> In rodent models of pain, CBC demonstrates anti-inflammatory effects in the LPS-induced paw edema model at doses of 10-100 mg/kg and elicits antinociceptive responses in a tail flick test at 6 nM.<sup>1,3</sup>

### References

- DeLong, G.T., Wolf, C.E., Poklis, A., *et al.* Pharmacological evaluation of the natural constituent of *Cannabis sativa*, cannabichromene and its modulation by Δ<sup>9</sup>-tetrahydrocannabinol. *Drug Alcohol Depend.* **112**(1-2), 126-133 (2010).
- De Petrocellis, L., Ligresti, A., Moriello, A.S., *et al.* Effects of cannabinoids and cannabinoid-enriched *Cannabis* extracts on TRP channels and endocannabinoid metabolic enzymes. *Br. J. Pharmacol.* **163**(7), 1479-1494 (2011).
- Maione, S., Piscitelli, F., Gatta, L., *et al.* Non-psychoactive cannabinoids modulate the descending pathway of antinociception in anaesthetized rats through several mechanisms of action. *Br. J. Pharmacol.* **162**(3), 584-596 (2011).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/16202](http://www.caymanchem.com/catalog/16202)

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 02/19/2015