

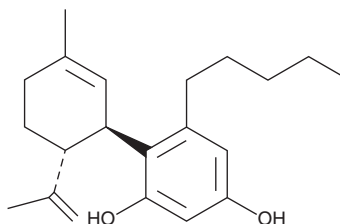
PRODUCT INFORMATION



Abnormal Cannabidiol

Item No. 10004259

CAS Registry No.: 22972-55-0
Formal Name: 4-[(1R,6R)-3-methyl-6-(1-methylethenyl)-2-cyclohexen-1-yl]-5-pentyl-1,3-benzenediol
Synonym: Abn-CBD
MF: C₂₁H₃₀O₂
FW: 314.5
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Abnormal cannabidiol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of abnormal cannabidiol in these solvents is at least 35 mg/ml.

Abnormal cannabidiol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of abnormal cannabidiol should be diluted with the aqueous buffer of choice. Abnormal cannabidiol has a solubility of 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Abnormal cannabidiol is a synthetic cannabidiol agonist of GPR55 (EC₅₀ = 2.5 μM in a GTPγS binding assay).¹ It is selective for GPR55 over cannabinoid (CB) receptor 1 (CB₁) and CB₂ (EC₅₀s = >30 μM for both in GTPγS binding assays). It is also a full agonist at GPR18, with an EC₅₀ value of 0.835 μM for inducing p44/42 MAPK phosphorylation in HEK293 cells expressing GPR18.³ Abnormal cannabidiol increases migration of BV-2 microglial cells (EC₅₀ = 0.6 μM).² It induces endothelium-dependent vasodilation via a CB₁-CB₂-NO-independent mechanism and inhibits vasoconstriction induced by endothelin 1 (ET-1; Item No. 24127) in retinal arterioles, an effect that can be blocked by the calcium-sensitive potassium channel blocker apamin (Item No. 17082).^{4,5} Abnormal cannabidiol also decreases blood pressure in anesthetized rats when administered intravenously at a dose of 20 μg/g or when administered directly into the rostral ventrolateral medulla at 0.4 and 0.8 μg.^{4,6}

References

1. Ryberg, E., Larsson, N., Sjögren, S., et al. *Br. J. Pharmacol.* **152**(7), 1092-1101 (2007).
2. McHugh, D., Page, J., Dunn, E., et al. *Br. J. Pharmacol.* **165**(8), 2414-2424 (2012).
3. Walter, L., Franklin, A., Witting, A., et al. *J. Neurosci.* **23**(4), 1398-1405 (2003).
4. Járai, Z., Wagner, J.A., Varga, K., et al. *Proc. Natl. Acad. Sci. U.S.A.* **96**(24), 14136-14141 (1999).
5. MacIntyre, J., Dong, A., Straiker, A., et al. *Eur. J. Pharmacol.* **735**(1), 105-114 (2014).
6. Penumarti, A., and Abdel-Rahman, A.A. *J. Pharmacol. Exp. Ther.* **349**(1), 29-38 (2014).

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, 03/28/2020

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