

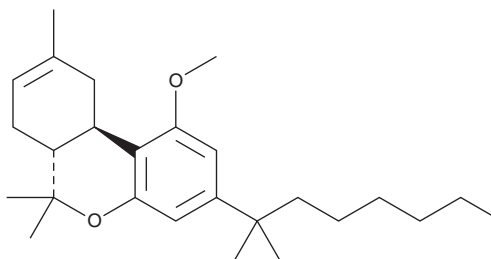
PRODUCT INFORMATION



L-759,633

Item No. 10009280

CAS Registry No.: 174627-50-0
Formal Name: 3-(1,1-dimethylheptyl)-6aR,7,10,10aR-tetrahydro-1-methoxy-6,6,9-trimethyl-6H-dibenzo[b,d]pyran
MF: C₂₆H₄₀O₂
FW: 384.6
Purity: ≥98%
Supplied as: A 10 mg/ml 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

L-759,633 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. Ethanol purged with an inert gas can be used. L-759,633 is soluble in ethanol at approximately 20 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of L-759,633 is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of L-759,633 in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-759,633 is a high-affinity peripheral cannabinoid receptor (CB₂)-selective receptor agonist with K_i values of 6.4 and 1,043 nM for CB₂ and central cannabinoid (CB₁) receptors, respectively. L-759,633 inhibits forskolin-stimulated cyclic AMP production in CHO cells transfected with CB₂ or CB₁ receptors with IC₅₀ values of 8.1 nM and 10 μM, respectively.¹

Reference

1. Ross, R.A., Brockie, H.C., Stevenson, L.A., *et al.* Agonist-inverse agonist characterization at CB₁ and CB₂ cannabinoid receptors of L759633, L759656 and AM630. *Br. J. Pharmacol.* **126**, 665-672 (1999).

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, 12/19/2025

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