

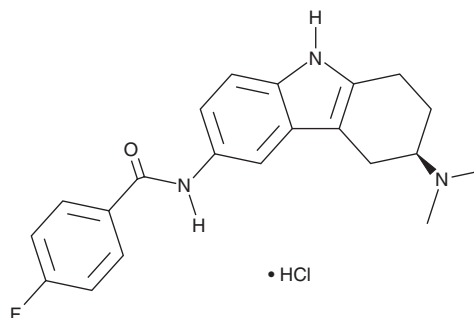
PRODUCT INFORMATION



LY344864 (hydrochloride)

Item No. 29496

CAS Registry No.: 1217756-94-9
Formal Name: N-[(3R)-3-(dimethylamino)-2,3,4,9-tetrahydro-1H-carbazol-6-yl]-4-fluoro-benzamide, monohydrochloride
MF: C₂₁H₂₂FN₃O • HCl
FW: 387.9
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 268 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

LY344864 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the LY344864 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. LY344864 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LY344864 (hydrochloride) in these solvents is approximately 10 and 3 mg/ml, respectively.

LY344864 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY344864 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LY344864 (hydrochloride) has a solubility of approximately 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

LY344864 is an agonist of the serotonin (5-HT) receptor subtype 5-HT_{1F} (K_i = 6 nM).¹ It is selective for 5-HT_{1F} over 5-HT_{1A-E}, 5-HT_{2A-C}, and 5-HT₇ receptors (K_is = 0.53-4.85 μM), as well as dopamine D₁ and D₂, GABA_A, histamine H₁, muscarinic, and α₁-, α₂-, and β-adrenergic receptors with K_i values ranging from 3.69 to greater than 100 μM. LY344864 inhibits forskolin-induced cAMP accumulation in L-M(TK-) cells expressing recombinant human 5-HT_{1F} receptors with an EC₅₀ value of 3 nM. It decreases electrically stimulated extravasation of plasma proteins in the dura mater in a rat trigeminal nerve model of migraine headache (ID₅₀s = 0.6 and 2.1 ng/kg for i.v. and oral administration, respectively).

Reference

1. Phebus, L.A., Johnson, K.W., Zgombick, J.M., *et al.* Characterization of LY344864 as a pharmacological tool to study 5-HT_{1F} receptors: Binding affinities, brain penetration and activity in the neurogenic dural inflammation model of migraine. *Life Sci.* **61**(21), 2117-2126 (1997).

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, 10/31/2022

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