

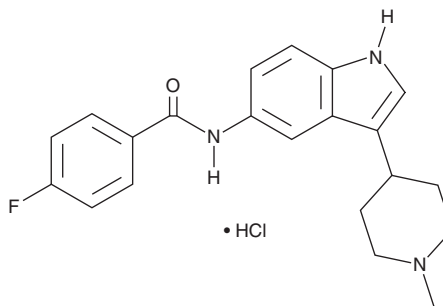
PRODUCT INFORMATION



LY334370 (hydrochloride)

Item No. 29481

CAS Registry No.: 199673-74-0
Formal Name: 4-fluoro-N-[3-(1-methyl-4-piperidiny)-1H-indol-5-yl]-benzamide, monohydrochloride
MF: C₂₁H₂₂FN₃O • HCl
FW: 387.9
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 267 nm
Supplied as: A crystalline 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

LY334370 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the LY334370 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. LY334370 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

LY334370 is an agonist of the serotonin (5-HT) receptor subtype 5-HT_{1F} (K_d = 0.446 nM).¹ It is selective for 5-HT_{1F} over other G protein-coupled 5-HT receptor subtypes with K_i values ranging from 16.4 to greater than 3,000 nM in radioligand binding assays. LY334370 inhibits forskolin-induced cAMP accumulation in mouse L-M(TK-) cell membranes expressing the recombinant human 5-HT_{1F} receptor (EC₅₀ = 1.51 nM).² It decreases electrically stimulated extravasation of plasma proteins in the dura mater in a guinea pig trigeminal nerve model of migraine headache.

References

1. Wainscott, D.B., Krushinski, J.H., Jr., Audia, J.E., *et al.* [3H]LY334370, a novel radioligand for the 5-HT_{1F} receptor. I. In vitro characterization of binding properties. *Naunyn Schmiedebergs Arch. Pharmacol.* **371(3)**, 169-177 (2005).
2. Johnson, K.W., Schaus, J.M., Durkin, M.M., *et al.* 5-HT_{1F} receptor agonists inhibit neurogenic dural inflammation in guinea pigs. *Neuroreport* **8(9-10)**, 2237-2240 (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, 10/21/2022

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