

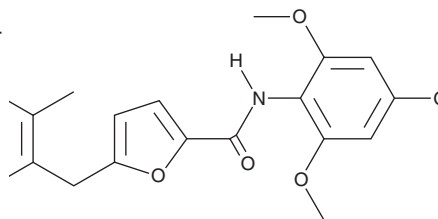
PRODUCT INFORMATION



AG-045572

Item No. 42244

CAS Registry No.: 263847-55-8
Formal Name: 5-[(5,6,7,8-tetrahydro-3,5,5,8,8-pentamethyl-2-naphthalenyl)methyl]-N-(2,4,6-trimethoxyphenyl)-2-furancarboxamide
MF: C₃₀H₃₇NO₅
FW: 491.6
Purity: ≥98%
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

AG-045572 is supplied as a solid. A stock solution may be made by dissolving the AG-045572 in the solvent of choice, which should be purged with an inert gas. AG-045572 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

AG-045572 is a nonpeptide antagonist of the gonadotropin-releasing hormone (GnRH) receptor ($K_s = 6, 2.2, \text{ and } 3.8 \text{ nM}$ for the human, mouse, and rat receptors, respectively).¹ It is selective for the GnRH receptor over a panel of 43 receptors, ion channels, and enzymes at 1 μM . AG-045572 prevents GnRH-induced increases in inositol phosphate production in HEK293 cells expressing the human GnRH receptor ($K_B = 25 \text{ nM}$). Oral or intravenous administration of AG-045572 decreases plasma luteinizing hormone (LH) levels in castrated rats. It inhibits GnRH-A-induced increases in plasma LH and testosterone levels in non-castrated rats when administered at a dose of 20 mg/kg.

Reference

1. Anderes, K.L., Luthin, D.R., Castillo, R., *et al.* Biological characterization of a novel, orally active small molecule gonadotropin-releasing hormone (GnRH) antagonist using castrated and intact rats. *J. Pharmacol. Exp. Ther.* **305**(2), 688-695 (2003).

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, 11/13/2024

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