

Human Recombinant H1 Histamine Receptor Stable Cell Line Cat. No. M00131

Version 05222014

1	Introduction	1
II	Background	1
Ш	Representative Data	2
IV	Thawing and Subculturing	2
V	References	3
	Limited Use License Agreement	4

I. INTRODUCTION

Catalog Number: M00131 Cell Line Name: HEK293/H1

Gene Synonyms: HRH1, H1-R, hisH1

Expressed Gene: Genbank Accession Number NM_ 000861, no expressed tags

Host Cell: HEK293

Quantity: Two vials of frozen cells (3×10⁶ per vial)

Stability: 16 passages

Application: Functional assay for H1 receptor

Freeze Medium: 45% culture medium, 45% FBS, 10% DMSO

Complete Growth Medium: DMEM, 10% FBS

Culture Medium: DMEM, 10% FBS, 200 µg/ml Zeocin

Mycoplasma Status: Negative

Storage: Liquid nitrogen immediately upon delivery

II. BACKGROUND

The H1 histamine receptor is expressed primarily in the lungs, vasculature, and brain. The H1 mediates the contraction of smooth muscles, neurotransmission in the central nervous system, the release of catecholamine from adrenal medulla, and increases in capillary permeability due to contraction of terminal venules. H1's role in inflammatory responses makes its antagonist suitable for treating allergies.



III. REPRESENTATIVE DATA

Concentration-dependent stimulation of intracellular calcium mobilization by Histamine in HEK293/H1 and HEK293 cells

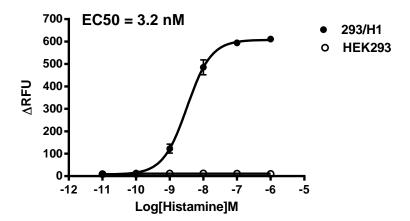


Figure Intracellular calcium response from HEK293 cells stably expressing human H1 histamine receptors and from untransfected control cells. Cells were loaded with Calcium-4 then stimulated with the indicated concentrations of histamine. Calcium responses were recorded on a FlexStation plate reader. Data represent the averge +/- standard deviation of triplicate determinations.

Note:

1. EC₅₀ value is calculated with four parameter logistic equation:

Y=Bottom + (Top-Bottom)/(1+10^((LogEC₅₀-X)*HillSlope))

X is the logarithm of concentration.

Y is RFU and starts at Bottom and goes to Top with a sigmoid shape.

2. Signal to background Ratio (S/B) = Top/Bottom

IV. THAWING AND SUBCULTURING

Thawing Protocol

- Remove the vial from liquid nitrogen tank and thaw cells quickly in a 37°C water-bath.
- 2. Just before the cells are completely thawed, decontaminate the outside of the vial with 70% ethanol and transfer the cells to a 15 ml centrifuge tube containing 9 ml of complete growth medium.
- 3. Pellet cells by centrifugation at 200 x g force for 5 min, and remove the medium.
- 4. Resuspend the cells in complete growth medium.
- 5. Transfer the cell suspension to a 10 cm dish with 10 ml of complete growth medium.
- 6. Grow the cells in incubator with 37°C, 5 %CO₂.
- 7. In the following day, replace the cells with fresh medium contains antibiotic.

Sub-culturing Protocol

Remove the culture medium from cells.



- 2. Wash cells with PBS (pH=7.4) to remove all traces of serum that contains trypsin inhibitor.
- 3. Add 2.0 ml of 0.05% (w/v) Trypsin- EDTA (GIBCO, Cat No. 25300) solution into 10 cm dish and observe the cells under an inverted microscope until cell layer is dispersed (usually within 3 to 5 minutes).
 Note: To avoid cells clumping, do not agitate the cells by hitting or shaking the dish while waiting for the cells detach.
 If cells are difficult to detach, please place the dish in 37°C incubator for ~2 min.
- 4. Add 6.0 to 8.0 ml of complete growth medium into dish and aspirate cells by gently pipetting.
- 5. Centrifuge the cells at 200 x g force for 5min, and remove the medium.
- 6. Resuspend the cells in culture medium and add the cells suspension to new culture dish.
- 7. Grow the cells in incubator with 37°C, 5 %CO₂.

V. REFERENCES

- 1. Sexton PM, Morfis M, Tilakaratne N, et.al (2006). Complexing receptor pharmacology: modulation of family B G protein-coupled receptor function by RAMPs. *Ann N Y Acad Sci.* 1070: 90–104.
- 2. Young A (2005). Amylin: Physiology and Pharmacology; Chapter 3: Receptor pharmacology. *Advances in Pharmacology* 52: 47–65.
- 3. *Calcitonin Receptors: Introduction. IUPHAR Database of Receptors and Ion Channels. International Union of Basic and Clinical Pharmacology.
- 4. McLatchie LM, Fraser NJ, Main MJ, et. al(1998). RAMPs regulate the transport and ligand specificity of the calcitonin-receptor-like receptor. *Nature* 393 (6683): 333–9.
- 5. Foord SM, Marshall FH (1999). RAMPs: accessory proteins for seven transmembrane domain receptors. *Trends Pharmacol. Sci.* 20 (5): 184–7.

GenScript USA Inc.

860 Centennial Ave. Piscataway, NJ 08854 Toll-Free: 1-877-436-7274

Tel: 1-732-885-9188, Fax: 1-732-210-0262

Email: product@genscript.com
Web: http://www.genscript.com

For Research Use Only.

- 860 Centennial Ave., Piscataway, NJ 08854, USA -



Limited Use License Agreement

This is a legal agreement between you (Licensee) and GenScript USA Inc. governing use of GenScript's stable cell line products and protocols provided to licensee. By purchasing and using the stable cell line, the buyer agrees to comply with the following terms and conditions of this label license and recognizes and agrees to such restrictions:

- The products are not transferable and will be used at the site where they were purchased. Transfer to another site
 owned by buyer will be permitted only upon written request by buyer followed by subsequent written approval by
 GenScript.
- 2) The purchaser cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party.
- 3) The products sold by GenScript are for laboratory and animal research purposes only. The products are not to be used on humans, for consumption, or for any unlawful uses.

GenScript USA Inc. will not assert against the buyer a claim of infringement of patents owned or controlled by GenScript USA Inc. and claiming this product based upon the manufacture, use or sale of a clinical diagnostic, therapeutic and vaccine, or prophylactic product developed in research by the buyer in which this product or its components has been employed, provided that neither this product nor any of its components was used in the manufacture of such product. For information on the use of this product for other purposes, contact Marketing Department, GenScript USA Inc., 860 Centennial Avenue, Piscataway, New Jersey 08840, U.S.A. Phone: 1-732-885-9188. Fax: 1-732-210-0262. Email: marketing@genscript.com.