

# Human Recombinant KiSS1-derived Peptide Receptor Stable Cell Line Cat. No. NA Version 06112020

I.	Introduction	
II.	Background1	
III.	Representative Data	
IV.	Thawing and Subculturing	
٧.	References	
VI.	Limited Use License Agreement	

#### I. INTRODUCTION

Catalog Number: M00190

Cell Line Name: CHO-K1/EDG2

Gene Synonyms: KISS1R; AXOR12; GPR54; HOT7T175

Expressed Gene: Genbank Accession Number NM\_032551; no expressed tags

Host Cell: CHO-K1

Culture Properties: Adherent

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

Stability: 16 passages

Application: Functional assay for KiSS1 receptor

Freeze Medium: 45% culture medium, 45% FBS (Cat. #10099-141, Gibco), 10% DMSO (Cat.

#D2650, Sigma)

Complete Growth Medium: Ham's F12 (Cat. #11765, Life Technologies), 10% FBS Culture Medium: Ham's F12, 10% FBS, 400 µg/ml G418 (Cat. #10131-035, Gibco)

Mycoplasma Status: Negative\*

Storage: Liquid nitrogen immediately upon delivery

#### II. BACKGROUND

The KiSS1-derived peptide receptor (also known as GPR54 or the Kisspeptin receptor) is a G<sub>q</sub>-coupled receptor which binds the peptide hormone kisspeptin (metastin). The KiSS1 gene inhibits metastatic activity of melanoma and other tumor cell lines, and clinical evidence supports a role for KiSS1 in inhibition of metastasis in human cancer. Kisspeptins and GPR54 also play a central role in hypothalamic regulation of puberty, by directly governing the release of gonadotropin-releasing hormone from the hypothalamus. In addition, mutations in GPR54 in mice and humans result in hypogonadotropic hypogonadism.

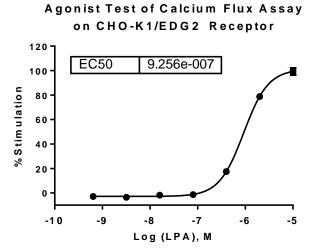


<sup>\*</sup> The mycoplasma test was performed with MycoAlert™ PLUS Mycoplasma Detection Kit of Lonza.

# III. REPRESENTATIVE DATA

Concentration-dependent stimulation of intracellular calcium mobilization by LPA in CHO-

# K1/EDG2



**Figure 1:** LPA-induced concentration-dependent stimulation of intracellular calcium mobilization in CHO-K1/EDG2 cells. The cells were loaded with Calcium-4 prior to stimulation with EDG2 receptor agonist, LPA. The intracellular calcium change was measured by FLIPR. The relative fluorescent units (RFU) were plotted against the log of the cumulative doses (5-fold dilution) of LPA (Mean  $\pm$  SD, n = 2). The EC50 of LPA on EDG2/CHO-K1 cells was 0.93  $\mu$ M. The S/B of LPA on EDG2 in CHO-K1 cells was 31.

### Note:

- EC<sub>50</sub> value is calculated with four parameter logistic equation:
   Y=Bottom + (Top-Bottom)/(1+10^((LogEC<sub>50</sub>-X)\*HillSlope))
   X is the logarithm of concentration. Y is the response 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**

- 1. Remove the vial from liquid nitrogen tank and thaw cells quickly in a 37°C water-bath.
- 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 for 5 min, and remove the medium.
- 4. Resuspend the cells with 1 ml complete growth medium.
- 5. Transfer the cell suspension to a 10 cm dish containing 10 ml complete growth medium.
- 6. Transfer the dish into an incubator of 37°C, 5% CO<sub>2</sub>.



7. Add antibiotic into the medium on the next day.

### **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 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 during incubation. 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 complete growth medium into dish and aspirate cells by gently pipetting.
- 5. Centrifuge the cells at 200 X g for 5min, and remove the medium.
- 6. Resuspend the cells in culture medium and transfer the cells to a new culture dish.
- 7. Transfer the dish into an incubator of 37°C, 5% CO<sub>2</sub>.

Subcultivation Ratio: 1:3 to 1:8

Medium Renewal: Every 2 to 3 days

#### V. REFERENCES

- 1. Lee DK, Nguyen T, O'Neill GP, et al., Discovery of a receptor related to the galanin receptors (1999). *FEBS Lett.* 446 (1): 103–7.
- 2. Kotani M, Detheux M, Vandenbogaerde A, et al. The metastasis suppressor gene KiSS-1 encodes kisspeptins, the natural ligands of the orphan G protein-coupled receptor GPR54 (2001). *J. Biol. Chem.* 276 (37): 34631–6.
- 3. Shahab M, Mastronardi C, Seminara SB, et al. (2005) Increased hypothalamic GPR54 signaling: a potential mechanism for initiation of puberty in primates. *Proceedings of the National Academy of Sciences of the United States of America*. 102 (6): 2129–34.
- 4. Colledge WH (2005) GPR54 and puberty. Trends Endocrinol. Metab. 15 (9): 448–53.

GenScript USA Inc.

860 Centennial Ave., Piscataway, NJ 08854 Tel: 732-885-9188, 732-885-9688

> Fax: 732-210-0262, 732-885-5878 Email: <u>info@genscript.com</u>

Web: <a href="http://www.genscript.com">http://www.genscript.com</a>
For Research Use Only.



# **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.