

Human Recombinant Glucagon Receptor Stable Cell Line

Technical Manual No. TM0587

Version 10132010

I	Introduction	1
II	Background	1
	Representative Data	2
	Thawing and Subculturing	
	References	
	Limited Use License Agreement	4

I. Introduction

Catalog Number: M00422 Cell Line Name: HEK293/GCGR/Gα15 Gene Synonyms: GCGR; GGR; MGC138246 Expressed Gene: Genbank Accession Number NM_000160; no expressed tags Host Cell: HEK293 Quantity: Two vials of frozen cells (3×10⁶ per vial) Stability: 16 passages Application: Functional assay for glucagon receptor Freeze Medium: 45% culture medium, 45% FBS, 10% DMSO Complete Growth Medium: DMEM, 10% FBS Culture Medium: DMEM, 10% FBS, 50 µg/ml Hygromycin B, 300 µg/ml G418 Mycoplasma Status: Negative

Storage: Liquid nitrogen immediately upon delivery

II. Background

Glucagon regulates blood glucose via control of hepatic glycogenolysis and gluconeogenesis and via regulation of insulin release from the β cell. Pharmacological administration of glucagon increases blood glucose in normal and diabetic subjects, and produces positive inotropic and chronotropic cardiovascular effects, relaxation of smooth muscle in the gastrointestinal tract and stimulation of growth hormone secretion. The actions of glucagon are mediated via a single adenylate cyclase-coupled glucagon receptor that also couples to the phospholipase C-inositol phosphate (PLC-IP) pathway leading to Ca²⁺ release from intracellular stores.

§: GenScript employs a PCR-based method to test the mycoplasma. The test covers 11 of the most common strains of mycoplasma, (covering approximately 95% of M. fermentans, M. hyorhinis, M. arginini, M. orale, M. salivarium, M. hominis, M. pulmonis, M. arthritidis, M. neurolyticum, M. hyopneumoniae and M. capricolum) and one species Ureaplasma (U. urealyticum), with sufficient sensitivity and specificity.

-1-							
860 Centennial Ave., Piscataway, NJ 08854, USA							
Tel: 1-732-885-9188	Fax: 1-732-210-0262	Email: info@genscript.com	Web: www.genscript.com				

.



III. Representative Data

Concentration-dependent stimulation of intracellular calcium mobilization by Glucagon in

HEK293/GCGR/Ga15 and HEK293 cells

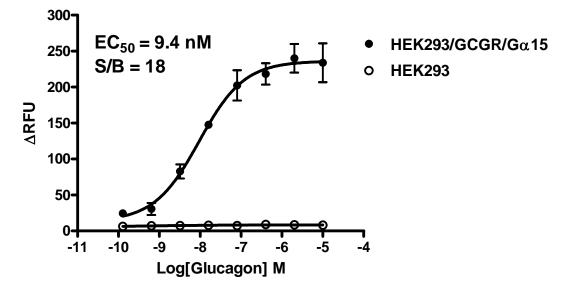


Figure 1. Glucagon-induced concentration-dependent stimulation of intracellular calcium mobilization in HEK293/GCGR/G α 15 and HEK293 cells. The cells were loaded with Calcium-4 prior to stimulation with a GCGR receptor agonist, Glucagon. The intracellular calcium change was measured by FlexStation. The relative fluorescent units (RFU) were plotted against the log of the cumulative doses (5-fold dilution) of Glucagon (Mean ± SD, n = 2). The EC₅₀ of Glucagon on GCGR co-expressing with G α 15 in HEK293 cells was 9.4 nM. The S/B of Glucagon on GCGR co-expressing with G α 15 in HEK293 cells was 18.

Notes:

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 the response

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

- 1. 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 discard the medium.
- 4. Resuspend the cells in complete growth medium.
- 5. Add 10 ml of the cell suspension in a 10 cm dish.
- 6. Add Hygromycin B and G418 to concentrations of 50 μg/ml and 300 μg/ml respectively the following day.

-2-							
860 Centennial Ave., Piscataway, NJ 08854, USA							
Tel: 1-732-885-9188	Fax: 1-732-210-0262	Email: info@genscript.com	Web: www.genscript.com				



Subculturing: Protocol

- 1. Remove and discard culture medium.
- 2. Wash cells with PBS (pH=7.4) to remove all traces of serum that contains trypsin inhibitor.
- Add 2.0 ml of 0.05% (w/v) Trypsin- EDTA (GIBCO, Cat No. 25300) solution to 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 clumping, do not agitate the cells by hitting or shaking the dish while waiting for the cells to detach. Cells that are difficult to detach may be placed at 37°C to facilitate dispersal.
- 4. Add 6.0 to 8.0 ml of complete growth medium and aspirate cells by gently pipetting, centrifuge the cells 200 x g force for 5min, and discard the medium.
- 5. Resuspend the cells in culture medium and add appropriate aliquots of the cell suspension to new culture vessels.
- 6. Incubate cultures at 37°C.

Subcultivation Ratio: 1:3 to 1:8 weekly. Medium Renewal: Every 2 to 3 days

V. References

- 1. Moore, M. C. and Cherrington, A. D., Autoregulation of hepatic glucose production(1998). *Eur. J. Endocrinol.*. 138: 240-248
- 2. Samols, E. and Marks, V., Interrelationship of glucagon, insulin, and glucose: The insulinogenic effect of glucagon(1966). *Diabetes*. 15: 855-865
- 3. Muhlhauser, I. et al., Incidence and management of severe hypoglycemia in 434 adults with insulin-dependent diabetes mellitus(1985). *Diabetes Care.* 8: 268-273
- 4. White, C. M., A review of potential cardiovascular uses of intravenous glucagon administration (1999). *J. Clin. Pharmacol.* 39: 442-447
- 5. Wakelam, M. J. and Houslay, M. D., Activation of two signal-transduction systems in hepatocytes by glucagon(1986). *Nature*. 323: 68-71

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: <u>http://www.genscript.com</u>

For Research Use Only.

-3-



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 or its components are used for research purposes only.

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 purposes other than research, contact Marketing Department, GenScript USA Inc., 860 Centennial Avenue, Piscataway, New Jersey 08854, U.S.A. Phone: 1-732-885-9188. Fax: 1-732-210-0262. Email: marketing@genscript.com.

Fax: 1-732-210-0262