

---

**Human Recombinant Parathyroid Hormone Receptor 1 Stable Cell Line**  
**Cat. No. M00315** **Version 07292020**

---

I. Introduction .....	1
II. Background .....	1
III. Representative Data .....	2
IV. Thawing and Subculturing .....	2
V. References .....	3
VI. Limited Use License Agreement.....	4

---

## **I. Introduction**

Catalog Number: M00315

Cell Line Name: CHO-K1/PTH1/Gα15

Gene Synonyms: MGC138426; MGC138452; PFE; PTHR; PTHR1

Expressed Gene: Genbank Accession Number NM\_000316; no expressed tags

Host Cell: CHO-K1/Gα15

Culture Properties: Adherent

Quantity: Two vials of frozen cells (>1×10<sup>6</sup> per vial)

Stability: 16 passages

Application: Functional assay for PTH1 receptor (Calcium flux assay, cAMP assay)

Freeze Medium: 45% culture medium, 45% FBS (Cat. #10099-141, Gibco), 10% DMSO (Cat. #D2650, Sigma)

Complete Growth Medium: Ham's F-12K (Kaighn's) (Cat. #21127, Life Technologies), 10% FBS

Culture Medium: Ham's F-12K (Kaighn's), 10% FBS, 100 µg/ml Hygromycin B (Cat. #10687010, Invitrogen), 400 µg/ml G418 (Cat. #10131-035, Gibco)

Mycoplasma Status: Negative\*

Storage: Liquid nitrogen immediately upon receipt

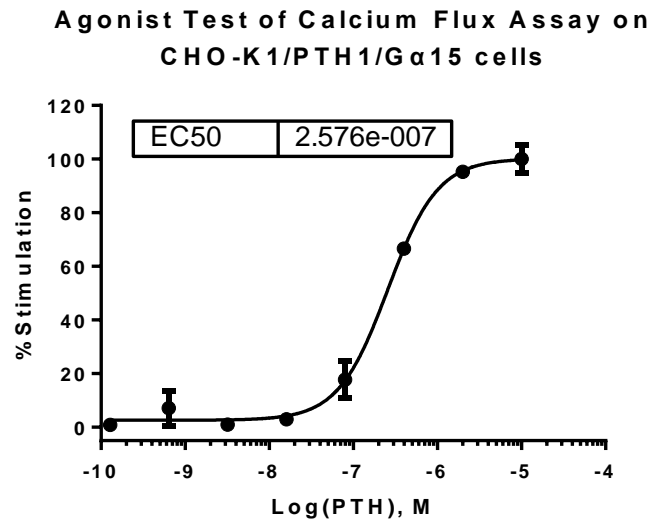
## **II. Background**

Parathyroid hormone (PTH)/PTH-related peptide (PTHrP) receptor (PTHR1) is a G protein coupled receptor which mediates the actions of both amino-terminal PTH and PTHrP fragments. The most abundant expression of PTHR1 is found in renal tubular cells and in osteoblasts, where the PTH/PTHrP receptor mediates the endocrine actions of PTH, and in prehypertrophic chondrocytes of the metaphyseal growth plate, where it mediates the autocrine/paracrine actions of PTHrP. Intact PTH (PTH 1-84) is compounded by a peptide of 84 amino acids (AA), the amino-terminal sequence, constituted by the first 34 AA (N-terminal structure), is necessary for its action.

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

### III. Representative Data

#### Concentration-dependent stimulation of intracellular calcium mobilization by PTH (1-34) in CHO-K1/PTH1/Gα15 cells



**Figure 1.** PTH (1-34)-induced concentration-dependent stimulation of intracellular calcium mobilization in CHO-K1/PTH/Gα15 cells. The cells were loaded with Calcium-4 prior to being stimulated with agonist PTH (1-34). The intracellular calcium change was measured by FLIPR<sup>TETRA</sup>. The relative fluorescent units (RFU) were plotted against the log of the cumulative doses of PTH (1-34) (Mean ± SD, n = 2). The EC<sub>50</sub> of PTH (1-34) on this cell was 0.26 μM.

#### Notes:

- EC<sub>50</sub> value is calculated with four parameter logistic equation:  

$$Y = \text{Bottom} + \frac{(\text{Top} - \text{Bottom})}{(1 + 10^{-(\text{LogEC}_{50} - X) \cdot \text{HillSlope}})}$$
 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.

### IV. Thawing and Subculturing

#### Thawing: Protocol

- Remove the vial from liquid nitrogen tank and thaw the 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.
- Pellet cells by centrifugation at 200 x g force for 5 min, and discard the medium.
- Resuspend the cells in complete growth medium.
- Add 10 ml of the cell suspension in a 10 cm dish.
- Add Hygromycin B and G418 to concentrations of 100 μg/ml and 400 μg/ml respectively the following day.

---

## 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.
3. Add 2.0 ml of 0.05% (w/v) Trypsin- EDTA (GIBCO, Cat No. 25300) solution to a 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

Medium Renewal: Every 2 to 3 days

## V. References

1. C Leroy, D Manen<sup>1</sup>, R Rizzoli<sup>1</sup>, M Lombès<sup>2</sup> and C Silve (2004) Functional importance of Myc-associated zinc finger protein for the human parathyroid hormone (PTH)/PTH-related peptide receptor-1 P2 promoter constitutive activity. *Journal of Molecular Endocrinology* 32, 99–113
2. C. de La Piedra, E. Fernández, M.<sup>a</sup> L. González Casaus and E. González Parra (2008) Differences in the function of parathyroid peptides. What are we measuring? *Nefrología*; 28 (2) 123-128

**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](mailto:product@genscript.com)  
Web: <http://www.genscript.com>  
**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:

- 1) 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 Ave. Piscataway, NJ 08854, U.S.A. Phone: 1-732-885-9188. Fax: 1-732-210-0262. Email: [marketing@genscript.com](mailto:marketing@genscript.com).