

---

**Human Recombinant PAR4 Proteinase-activated Receptor Stable Cell Line**  
**Cat. No. M00206** **Version 06232020**

---

|  |   |
|--|---|
| 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: M00206

Cell Line Name: CHO-K1/PAR4

Gene Synonyms: PAR4, PAR-4, F2RL3

Expressed Gene: Genbank Accession Number NM\_003950; no expressed tags

Host Cell: CHO-K1

Culture Properties: Adherent

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

Stability: More than 16 passages

Application: Functional assay for PAR4 receptor (Calcium flux assay)

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

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

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

Mycoplasma Status: Negative\*

Storage: Liquid nitrogen immediately upon receipt

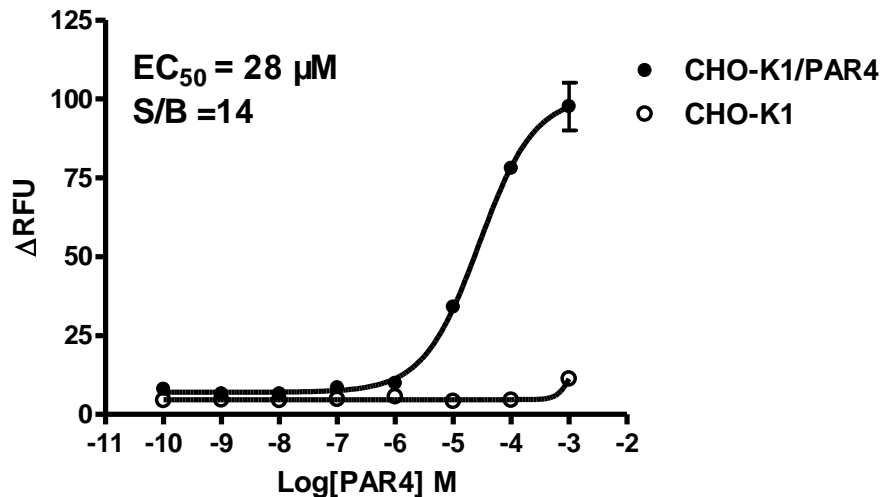
**II. BACKGROUND**

Protease-activated receptor (PAR)-4 is a member of a unique family of GPCRs. The protease-activated receptors (PARs) that are activated by proteolytic cleavage of the N-terminal domain of the receptor reveal a tethered ligand. The PAR family consists of 4 receptors; PAR1 and PAR3 are activated by thrombin, and PAR2 and PAR4 are activated by several serine proteases (Macfarlane et al., 2001). PAR4 is a recently identified low-affinity thrombin receptor that plays a pathophysiological role in many types of tissues including the lung. Mice lacking PAR4 are protected from mesenteric arteriole thrombosis, indicating that PAR4 is a potential target for treatment of thrombosis in humans.

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

### III. REPRESENTATIVE DATA

Concentration-dependent stimulation of intracellular calcium mobilization by PAR4 in CHO-K1/PAR4 cells



**Figure 1:** PAR4-induced concentration-dependent stimulation of intracellular calcium mobilization in CHO-K1/PAR4 cells. The cells were loaded with Calcium-4 prior to being stimulated with a PAR4 receptor agonist, PAR4. The intracellular calcium change was measured by FLIPR. The relative fluorescent units (RFU) were normalized and plotted against the log of the cumulative doses (10-fold dilution) of PAR4 (Mean  $\pm$  SD, n = 2). The EC<sub>50</sub> of PAR4 on CHO-K1/PAR4 cells was 28  $\mu$ M.

#### Note:

- 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 and starts at Bottom and goes to Top with a sigmoid shape.
- 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.
- 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 for 5 min, and remove the medium.
- Resuspend the cells with 1 ml complete growth medium.
- Transfer the cell suspension to a 10 cm dish containing 10 ml complete growth medium.
- Transfer the dish into an incubator of 37°C, 5% CO<sub>2</sub>.
- Add antibiotic into the medium on the next day.

### Sub-culturing Protocol

1. 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. Mao Y, Jin J, Daniel JL, Kunapuli SP (2009) Regulation of plasmin-induced protease-activated receptor 4 activation in platelets. *Platelets*. 2009 May; 20(3):191-8.
2. Nieman MT. (2008) Protease-activated receptor 4 uses anionic residues to interact with alpha-thrombin in the absence or presence of protease-activated receptor 1. *Biochemistry*. 2008 Dec 16; 47(50):13279-86
3. Ando S, Otani H, Yagi Y, Kawai K, Araki H, Nakamura T, Fukuhara S, Inagaki C. (2007) Protease-activated receptor 4-mediated Ca<sup>2+</sup> signaling in mouse lung alveolar epithelial cells. *Life Sci*. 2007 Aug 16; 81(10):794-802. Epub 2007 Aug 17

**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).