



## MMP9 Stably Expressing HEK293 Stable Cell Line

Catalog Number: EL-1003 (For Research Use Only)

### Introduction

Matrix metalloproteinase (MMP)-9, one of the most widely investigated MMPs, regulates pathological remodeling processes that involve inflammation and fibrosis in cardiovascular disease and cancer. It has been shown that MMP9 plays an important role in several aspects of central nervous system activity. MMP-9 expression and activity is a common endpoint measured. Little information is available on the underlying mechanisms of MMP-9 function. Therefore, MMP-9 overexpression cell line will be beneficial and highlights for unrevealing MMP-9 function. Signosis has developed MMP9 Stably Expressing HEK 293 Cell Line, which allows you to study the roles MMP9 playing on a wide variety of biological actions and functionalities.

### Provided Materials

One vial of  $2 \times 10^6$  cells, at passage 4, in Freezing Media.

**IMPORTANT:** store the frozen cells in liquid nitrogen until you are ready to thaw and propagate them.

### Handling cells upon arrival

It is strongly recommended that you propagate the cells by following instructions as soon as possible upon arrival \*\*.

**IMPORTANT:** An adequate number of frozen stocks must be made from early passages as cells will undergo genotypic changes. Genetic instability in transfected cells will result in a decreased responsiveness over time in normal cell culture conditions.

### Required Cell Culture Media

- **Complete Growth Media**  
In 450mL of DMEM, add 50mL FBS (10% final) and 5mL Penicillin/Streptomycin (1% final).
- **Freezing Media**  
Add 10% DMSO (final) to Complete Growth Media and sterile filter. Make fresh each time.

### Materials required but not provided

(can be substituted with a comparable third-party product)

Materials	Product number
Dulbecco's Modified Eagles Medium (DMEM)	Hyclone SH30243.01
Fetal Bovine Serum (FBS)	Fisherbrand P/N 03-600-511
Penicillin/Streptomycin	Hyclone P/N SV30010
Trypsin	Hyclone P/N SH30236.02
Phosphate-buffered saline (PBS)	Cellgro P/N 21-040-CV
DMSO	Sigma P/N D8418
96-well white plate	Greiner Bio-One P/N 655098

### Initial Culture Procedure

1. Quickly thaw cells in a 37 °C water bath with careful agitation. Remove from the bath as soon as the vial is thawed.
2. Transfer cells to a 15ml centrifuge tube containing 7ml of pre-warmed Complete Growth Media.
3. Centrifuge tube at 1200-1500 RPM for 5 minutes
4. Remove supernatant and resuspend cells with 1ml Complete Growth Media.
5. Transfer cells to a T75cm<sup>2</sup> tissue culture flask or 100 mm culture dish containing 8-12ml of Complete Growth Media.
6. Place the flask with cells in a humidified incubator at 37 °C with 5% CO<sub>2</sub>.

### Subculture Procedure

- A sub-cultivation ratio of 1:3 to 1:4 is recommended with media changes every 2 to 3 days.

**Preparing frozen stocks**

*This procedure is designed for 60mm<sup>2</sup>dish or T25cm<sup>2</sup> flask. Scale volumes accordingly to other vessels.*

1. When cells reach 1-1.5x10<sup>6</sup>/ml, freeze down cells.
2. Transfer cells to a 15ml conical centrifuge tube and centrifuge at 1200-1500 RPM for 5 minutes to collect the cells into a pellet.
3. Carefully aspirate the media and resuspend cells in 1ml freezing media and gently resuspend by pipetting up and down.
4. Transfer 1mL of cells into a cryogenic vial.
5. Place the cryogenic vial in a freezing container (e.g. Nalgene # 5100-0001) and store it at -80°C freezer overnight.
6. Transfer cells to liquid nitrogen for long-term storage.

---

\*\* Signosis products are warranted for 30 days from the date of shipment, and this warranty is valid only if the product is stored and handled according to this user manual or product information sheet and shipped directly by Signosis.