

GSK3B Stably Expressing Cell Line

HeLa Catalog number: EL-1001 HEK293 Catalog number: EL-1002 (For Research Use Only)

Introduction

GSK3β is a multifunctional kinase known to be a key regulator of a wide range of cellular functions. GSK3β is involved in modulating a variety of functions including cell signaling, growth metabolism, and various transcription factors regulating the survival or apoptosis. β-Catenin, a key component in Wnt signaling pathway, is phosphorylated by GSK3β. Inhibition of GSK3ß activity leads to stabilization and accumulation of \beta-catenin in the cytosol, which is shuttled into the nucleus and regulates gene expression. Dysregulation of GSK3β expression leads to many pathological conditions, including Alzheimer's disease, inflammation, diabetes, and cancer. Many studies have supported GSK3β playing a role as a tumor suppressor and tumor promoter. Signosis provided GSK-3β Stably Expressing HeLa or HEK293 Cell Line for GSK3β signaling pathway study.

Provided Materials

One vial of 2 x 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 results 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

- Quickly thaw cells in a 37 °C water bath with careful agitation. Remove from the bath as soon as the vial is thawed.
- 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.
- Transfer cells to a T75cm² tissue culture flask or 100 mm culture dish containing 8-12ml of Complete Growth Media.
- Place the flask with cells in a humidified incubator at 37 °C with 5% CO₂.

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 60mm²dish or T25cm² flask. Scale volumes accordingly to other vessels.

- 1. When cells reach $1-1.5 \times 10^6$ /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.
- 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.