

HLA-A:02:01 Stably expressing HeLa Cell Line

Catalog #: HLA-0009

ISO 13485

Introduction

HeLa cells are the oldest and most commonly used human cancer cell lines in biomedical research. It is an immortal cell line derived from cervical cancer cells taken from a lady called Henrietta Lacks in 1951, which was found to be remarkably durable and prolific. HeLa- HLA-A:02:01 stable cell line is transformed from the HeLa cell line and stably expresses the HLA-A:02:01 protein. The expression of HLA-A:02:01 has been validated by western blot.

Protein Accession Number: P04439

Provided Materials

One vial of 2×10^6 cells, at P4 in Freezing Media.**IMPORTANT:** store the frozen cells in liquid nitrogen until you are ready to thaw and propagate them.

Additional Materials Required

1. Dulbecco's Modified Eagle Medium (DMEM)
2. Fetal Bovine Serum (FBS)
3. Penicillin/Streptomycin
4. Trypsin
5. Phosphate-buffered saline (PBS)
6. DMSO
7. 96-well white plate

Handling Cells Upon Arrival

It is strongly recommended that you propagate the cells 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.

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 T75cm² 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% 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 60mm² dish or T25cm² flask. Scale volumes according to other vessels.

1. When cells reach 1-1.5x10⁶/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 and store it at -80 °C freezer overnight.
6. Transfer cells to liquid nitrogen for long-term storage.

This product is for research use only.