

Jurkat Cell Line

Catalog Number: PC-006 (For Research Use Only)

Jurkat cells (Clone E6-1 ATCC® TIB-152TM) are an immortalized line of T lymphocyte cells that are used to study acute T cell leukemia, T cell signaling, and the expression of various chemokine receptors susceptible to viral entry, particularly HIV. Jurkat cells are also useful in science because of their ability to produce interleukin 2. Their primary use, however, is to determine the mechanism of differential susceptibility of cancers to drugs and radiation.

Product

One vial of 2 x 10⁶ cells 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.

Required Cell Culture Media

• Complete Growth Media

In 450mL of RPMI, add 50mL FBS (10% final) and 5mL Penicillin/Streptomycin (1% final).

• 2x 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)

- RPMI -- Hyclone P/N SH30027.01
- 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*

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.
- 2. Wipe the vial with 70% ethanol and transfer cell suspension into a 15ml centrifuge tube containing 9.0 mL warm complete culture medium.
- 3. Spin at 200 g, for 5minutes at room temperature.
- Resuspend cell pellet with the appropriate volume of complete medium and transfer the cell suspension into a T25 culture flask.
- 5. Place the flask with cells in a humidified incubator at 37°C with 5% CO₂.
- **6.** The next day, replace media with fresh Complete Growth Medium.
- 7. When cells are 6-8x10⁵ cells/ml, split them 1:4 with fresh media. Add appropriate aliquots of the cell suspension to new culture vessels (T25 = 10 ml; T75 = 50 ml; T150 = 100 ml maximum volume). Grow cells to no more than 8 x 10⁵ cells/ml. Disperse clumps gently for counting.

Preparing frozen stocks

This procedure is designed for 100mm²dish or T75cm² flask. Scale volumes accordingly to other vessels.

- When cells reach 90-100% confluency, freeze down cells.
- 2. Transfer cells to a 15ml conical centrifuge tube and centrifuge at 200 g for 5 minutes to collect the cells into a pellet.
- Carefully aspirate the media and resuspend cells in 0.5mL complete growth media.
- Add 0.5mL of 2X freezing media and gently resuspend by pipetting up and down.
- 5. Transfer 1mL of cells into a cryogenic vial.
- Place the cryogenic vial in a freezing container (Nalgene # 5100-0001) and store it at -80°C freezer overnight.
- Transfer cells to liquid nitrogen for long-term storage.