

# NF-kB Luciferase Reporter Jurkat Stable Cell Line (For Research Use Only)

Catalog Number: SL-0050

#### Introduction

NF-κB is a critical regulator of inflammatory responses, proliferation, and differentiation of T-cells. The aberrant activation of NFkB can contribute to the development of autoimmunity, chronic inflammation, or lymphoid cancer. Jurkat cells are human T lymphocyte cells widely used to study T cell signaling. Signosis developed a stable Jurkat NFkB-luciferase reporter stable cell line, which can be used for easily monitoring the activation of NFkB activation in T cells through sensitive luciferase analysis. This cell line was established by transfection using a pTA-NFkB-luciferase reporter vector, along hygromycin expression vector followed hygromycin selection. The hygromycin resistant clones were subsequently screened for luciferase activity induced by TNFa treatment

# **Product description**

Signosis has developed NF-kB luciferase reporter Jurkat stable cell line by co-transfecting NF-kB luciferase reporter vector and hygromycin expression vector. The hygromycin-resistant clones were subsequently screened for TNF $\alpha$  luciferase activity. The cell line can be used as a reporter system for monitoring the activation of NF-kB triggered by stimuli treatments, gene overexpression, and gene knockdown.

# Materials provided

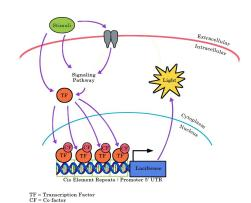
One vial of 4 x 10<sup>6</sup> cells, at passage 4, in Freezing Media. **IMPORTANT**: store the frozen cells in liquid nitrogen until you are ready to thaw and propagate them.

# Handling cells upon arrival



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

**IMPORTANT**: It is imperative that an adequate number of frozen stocks be made from early passages as cells may undergo genotypic changes. Possible genetic instability in transfected cells may results in a



decreased responsiveness over time in normal cell culture conditions.

# Required Cell Culture Media

# • Complete Growth Media

In 450mL of RPMI-1640, 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)

producty		
Materials	Product number	
RPMI-1640 Medium	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	Cellgro P/N 21-040-	
(PBS)	CV	
DMSO	Sigma P/N D8418	
96-well white plate	Greiner Bio-One P/N	
	655098	
Luciferase substrate	Signosis P/N LUC015	
Cell lysis buffer	Signosis P/N LS-001	
Hygromycin B	Toku-E P/N H010	

#### **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.
- 2. Transfer cells to 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 100mm<sup>2</sup> tissue culture dish (or T-75cm<sup>2</sup> flask) containing 10ml of Complete Growth Media.
- Place the dish with cells in a humidified incubator at 37 °C with 5% CO<sub>2</sub>.

#### Cell maintenance

- 1. After cells recovered and are growing well (after at least one passage), maintain and subculture the cells in Complete Growth Media with 100-200µg/ml hygromycin B.
- Pass the cells every 3 days by inoculating 5x10<sup>5</sup>/ml. Do not allow the cell concentration to exceed 3x10<sup>6</sup>/ ml.

# Preparing frozen stocks

This procedure is designed for 60mm<sup>2</sup>dish or T-25cm<sup>2</sup> flask. Scale volumes accordingly to other vessels.

- 1. When cells reach 2-3x10<sup>6</sup>cells/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 at a density of 5-7x10<sup>6</sup>/ml in freezing media and gently resuspend by pipetting up and down.
- 4. Aliquot 1ml of cells into cryogenic vials.
- Place the cryogenic vial in a freezing container (Nalgene # 5100-0001) and store it at -80°C freezer overnight.
- **6.** Transfer cells to liquid nitrogen for long-term storage.

#### Assay procedure

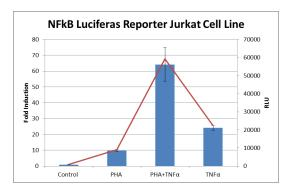
The following procedure should be followed as a guideline. You will need to optimize the assay conditions based on your experimental setup.

- 1. Centrifuge cells at 1000-1500 RPM for 5 minutes
- 2. Remove supernatant and resuspend NF-kB Jurkat cells at 1-1.2x10<sup>6</sup> cells/ml in pre-warmed RPMI medium+0.1%FBS.
- 3. Add 90µl of cell suspension (~100,000 cells) per well of a 96 well white plate.
- Add 10µl of 10x stock of inducers per well and 10µl PBS or endotoxin-free water as a negative control.

- Note: For PHA stimulation,  $1\mu$  of 10x stock PHA can be added (to get a final concentration of  $50\mu g/ml$ ) and incubate at  $37^{\circ}C$  in a  $CO_2$  incubator for 4 hrs. before adding inducers.
- 5. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for the appropriate time to produce maximal induction
- 6. Slowly discard 80μl of the media by using a pipette. *Note: Do not disturb the cells on the bottom of each well.* 
  - Optional Step: Centrifuge the plate at 1200 RPM for 1 minute to settle the cells onto the bottom of each well, then discard 80uL of the media.
- 7. Add 20μl of 2x lysis buffer to each well (To prepare 2x lysis buffer, add two volumes of 5x lysis buffer to three volumes of distilled water).
- **8.** Incubate cells in lysis buffer for 15-30 minutes at room temperature with gentle agitation.
- Transfer cell lysate solution from each well to 1.5 ml centrifuge tubes.
- 10. Centrifuge the tubes at full speed for 1 minute.
- 11. Carefully pipet 20µl of the supernatant back to the 96 well white plate.

Note: It is very important to centrifuge the cell lysate and test the supernatant only.

- **12.** Add 50μl of luciferase substrate to each well and gently pipette up and down.
- 13. Immediately read the plate in a luminometer.



0.1x106 Cells were plated on a 96-well plate and incubate in  $100\mu l$  media +0.1% FBS. Additions to wells were as follows: (Control ) No addition; (PHA) 50µg/ml incubated for 20 hours; (PHA+TNF $\alpha$ ) PHA50µg/ml incubated for 4 hours then TNF $\alpha$ 20ng/ml was added and continue incubating for 16 hours, (TNF $\alpha$ ) TNF $\alpha$ 20ng/ml incubated for 16 hours.

# Signosis Luciferase Reporter Stable Cell Lines

For a complete list of cell lines please visit our website at <a href="http://www.signosisinc.com/category/cell-based-assays">http://www.signosisinc.com/category/cell-based-assays</a>

Transcription Factor	Pathway	Cell Line	Cat #
NFkB	NFkB	Hela; human cervical cancer	SL0001
NFkB	NFkB	NIH/3T3; mouse fibroblast	SL0006
NFkB	NFkB	HEK293; human embryonic kidney	SL0012
NFkB	NFkB	MCF-7; human breast cancer	SL0013
NFkB	NFkB	A549; human lung cancer	SL0014
NFkB	NFkB	HepG2; human river cancer	SL0017
NFkB	NFkB	MEF; murine embryonic fibroblast	SL0033
NFAT	Calcium Signaling	Jurkat; human T lymphocytes	SL0032
NFAT	Calcium Signaling	Hela; human cervical cancer	SL0018
p53	p53	Hela; human cervical cancer	SL0011
p53	p53	RKO; human colon cancer	SL0007
SMAD	TGFbeta	HepG2; human river cancer	SL0016
SMAD	TGFbeta	NIH/3T3; mouse fibroblast	SL0030
NRF2	Antioxidant Response	MCF7; human breast cancer	SL0010
STAT1	JAK-STAT	Hela; human cervical cancer	SL0004
STAT3	JAK-STAT	Hela; human cervical cancer	SL0003
HIF	Hypoxia Response	NIH/3T3; mouse fibroblast	SL0005
HIF	Hypoxia Response	Hela; human cervical cancer	SL0023
HIF	Hypoxia Response	Neuro2a; mouse neuroblastoma	SL0027
ER	Estrogen Receptor Signaling	T47D; human breast cancer	SL0002
AR	Androgen Receptor Signaling	MDA-MB-453; human breast cancer	SL0008
GR	Glucocorticoid Receptor Signaling	MDA-MB-453; human breast cancer	SL0009
GR	Glucocorticoid Receptor Signaling	Hela; human cervical cancer	SL0021
AP-1	JNK, ERK, MAPK Signaling	Hela; human cervical cancer	SL0019
CREB	cAMP, PICA, CaMK Signaling	HEK293; human embryonic kidney	SL0020
CREB	cAMP, PICA, CaMK Signaling	NIH/3T3; mouse fibroblast	SL0031
CHOP	Unfolded Protein Response, ER stress	Mia-Paca2; human pancreatic cancer	SL0025
TCF/LEF	Wnt/b-catenin	HEK293; human embryonic kidney	SL0015
TCF/LEF	Wnt/b-catenin	Hela; human cervical cancer	SL0022
TCF/LEF	Wnt/b-catenin	CHO-KI; Chinese Hamster Ovary	SL0028
ELK	MAPK Signaling	HEK293; human embryonic kidney	SL0040
ELK	MAPK Signaling	Hela; human cervical cancer	SL0041
IRF	Immune Response Pathway	HEK293; human embryonic kidney	SL0035

<sup>\*\*</sup> 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.