

Datasheet for HEK293T/NFAT-Luc Cell Line

Catalog number: SL402

Product: HEK293T cell line stably expressing NFAT transcriptional response (TRE) gene.

Description: The NFAT transcriptional response (TRE) cell line assays the status of the

PKC/Ca++ signaling pathway through the activity of the transcription factor NFAT. The NFAT transcriptional TRE cell line contains a minimal promoter and tandem repeats of the NFAT transcriptional response element upstream

of a secreted Gaussia luciferase reporter gene.

Quantity: 1 vial of 2 x 10⁶ cells; frozen

Shipping conditions: Dry ice

Storage conditions: Liquid nitrogen vapor phase. Remove the item from the dry ice packaging

and check all items for damage and leakage. Place immediately into storage at

or below -140 °C, preferably into the liquid nitrogen vapor phase, until use.

Transgene integration:



TPS-NFAT-TRE-miniCMV-GLuc-SV40-Puro

Source of parental line:

HEK293T

Organism: *Homo sapiens,* human Tissue: kidney Cell type: Epithelial

Quality control: >95% viability before freezing. All cells were tested and found to be free

of mycoplasma, bacteria, viruses, and other toxins.

Safety instructions: To ensure safety, protective gloves, clothing, and a face mask should be worn

when handling frozen vials. Some leakage may occur into the vial during storage. The liquid nitrogen will be converted to gas upon thawing. Due to the nature of nitrogen gas, pressure may build within the vial and possibly result in the vial

exploding or losing its cap. This may cause flying debris.



Thawing procedure: The vial of cells should be thawed in a 37 °C water bath with gentle agitation.

For optimal performance, the vial should be thawed in under two minutes. Ensure that the cap of the vial did not loosen upon thawing, and re-tighten as needed. Spray the vial with 70% EtOH and wipe off. Repeat. Using aseptic technique, add the contents of the vial to 9 ml of complete growth medium (without selection). Centrifuge for 5 min. at 125 x g. Aspirate the medium, being careful not to disturb the pellet. Resuspend in 10 mL of complete growth medium, and place into a culture vessel of your choice. Only add selection to the medium after 24 hours in culture.

Culture conditions:

Complete Growth Medium

The base medium for this cell line is DMEM. For optimal growth and maintenance of selection, add the following components to the base medium: dialyzed fetal bovine serum to a final concentration of 10%.

Selection: Puromycin to a final concentration of 1 μg/mL

Culture temperature: 37 °C with 5% CO₂

Subculture: Replace culture medium with selection-free medium and incubate for up to 6 hours. Rinse the cells with PBS without cations, digest cells with 0.25% (w/v) Trypsin-EDTA (0.53 mM) solution and split at 1:3 to 1:10 ratio.

Cryopreservation: Freeze slowly in complete growth medium supplemented with 5% (v/v) DMSO.

Mycoplasma: Negative

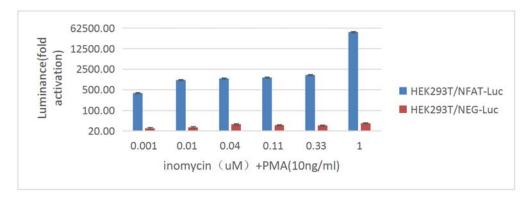
(MycoAllert Mycoplasma Detection Kit from Lonza)

Product QC:

Dose curve

Split cells from Master Cell Bank clone into 96 well plate, HEK293T/NEG-Luc cell line as negative control, cells were incubated for 24 hours, followed by 18 hours of treatment with PMA and ionomycin. PMA:10ng/ml, ionomycin(uM): 0, 0.01, 0.04, 0.11, 0.33, 1. A Gaussia Luciferase Assay was performed.





Citation of product: If use of this item results in a publication, please use this information: HEK293T/NFAT-Luc Cell Line (SL402, GeneCopoeia, Inc., Rockville, MD).

Limited Use License

A limited use license is granted to the Buyer of the Product. The Product shall be used by the Buyer for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Copyright ©2016 GeneCopoeia, Inc.

TRSCL-DS-062816

GeneCopoeia, Inc. 9620 Medical Center Drive, #101 Rockville, MD 20850 USA

Tel: 301-762-0888; Fax: 301-762-3888 Email: support@genecopoeia.com Web: www.genecopoeia.com