THP1-Lucia[™] ISG Cells

Interferon Regulatory Factor-Inducible Reporter Monocytes

Catalog code: thpl-isg https://www.invivogen.com/thp1-lucia-isg

> For research use only Version 23E29-MM

PRODUCT INFORMATION

Contents and Storage

• 3-7 x 10⁶ of THP1-Lucia[™] ISG cells in a cryovial or shipping flask. IMPORTANT: If cells provided in a cryovial are not frozen upon arrival, contact InvivoGen immediately.

• 1 ml of Zeocin[®] (100 mg/ml). Store at 4 °C or at -20 °C.*

• 1 ml Normocin[™] (50 mg/ml), a formulation of three antibiotics active against mycoplasmas, bacteria and fungi. Store at -20°C.* *The expiry date is specified on the product label.

Note: Data sheets for all components are available on our website.

• **1 tube of QUANTI-Luc™ 4 Reagent**, a Lucia luciferase detection reagent (sufficient to prepare 25 ml). Store at -20 °C. Avoid repeated freeze-thaw cycles. Note: This product is photosensitive and should be protected from light.

Handling Frozen Cells Upon Arrival

Cells must be thawed immediately upon receipt and grown according to handling procedures (as described on the next page) to ensure the best cell viability and proper assay performance.

Note: **Avoid freezing cells upon receipt** as it may result in irreversible damage to the cell line.

Disclaimer: We cannot guarantee cell viability if the cells are not thawed immediately upon receipt and grown according to handling procedures.

IMPORTANT: For cells that arrive in a shipping flask please refer to the enclosed 'cell recovery procedure'.

Quality Control

- Reporter activity has been validated using functional assays.
- The stability for 20 passages following thawing has been verified.
- THP1-Lucia[™] ISG cells are guaranteed mycoplasma-free.

Cell Line Stability

Cells will undergo genotypic changes resulting in reduced responsiveness over time in normal cell culture conditions. Genetic instability is a biological phenomenon that occurs in all stably transfected cells. Therefore, it is critical to prepare an adequate number of frozen stocks at early passages.

USE RESTRICTIONS

These cells are distributed for research purposes only.

This product is covered by a Limited Use License. By use of this product, the buyer agrees to the terms and conditions of all applicable Limited Use Label Licenses. For non-research use, such as screening, quality control or clinical development, contact info@invivogen.com.

PRODUCT DESCRIPTION

THP1-Lucia[™] ISG (interferon-stimulated genes) cells were derived from the human monocytic cell line THP-1, which represents a model of choice to study the activation and signaling of cytosolic DNA sensors (CDSs). Indeed, THP-1 cells have been shown to express all the CDSs identified so far1-3, with the exception of DAI4. THP1-Lucia™ ISG cells express the secreted luciferase Lucia reporter gene under the control of an IRF-inducible promoter. This composite promoter is comprised of five IFN-stimulated response elements (ISRE) fused to an ISG54 minimal promoter, which is unresponsive to activators of the NF-kB or AP-1 pathways. As a result, THP1-Lucia™ ISG cells allow the monitoring of the IRF pathway by determining the activity of Lucia luciferase. The levels of IRF induced Lucia luciferase in the cell culture supernatant are readily assessed with QUANTI-Luc[™] 4 Lucia/Gaussia, a Lucia and Gaussia luciferase detection reagent. THP1-Lucia[™] ISG cells are highly responsive to CDS ligands, such as transfected double-stranded nucleic acids

THP1-Lucia[™] ISG cells are resistant to Zeocin[®].

1. Zhang Z. *et al.*, 2011. The helicase DDX41 senses intracellular DNA mediated by the adaptor STING in dendritic cells. Nat Immunol.12(10):959-65. 2. Veeranki S. *et al.*, 2011. IFI16 protein mediates the anti-inflammatory actions of the type-I interferons through suppression of activation of caspase-1 by inflammasomes. PLoS One. 6(10):e27040. 3. Arakawa R. *et al.*, 2010. Characterization of LRRFIP1. Biochem Cell Biol. 88(6):899-906. 4. Lippmann J. *et al.*, 2010. IFNbeta responses induced by intracellular bacteria or cytosolic DNA in different human cells do not require ZBP1 (DLM-1/DAI). Cell Microbiol. 10(12):2579-88.

SAFETY CONSIDERATIONS

Biosafety Level 1

HANDLING PROCEDURES

Required Cell Culture Medium

• Growth Medium: RPMI 1640, 2 mM L-glutamine, 25 mM HEPES, 10% heat-inactivated fetal bovine serum (30 min at 56°C), 100 µg/ml Normocin[™], Pen-Strep (100 U/ml-100 µg/ml)

Initial culture of all THP-1 derived cells must be performed in growth medium containing 20% heat-inactivated FBS.

Note: The use of Normocin[™] together with Pen-Strep is required to keep the cells free of microbial contaminants. Contamination of this cell line may activate TLRs resulting in differentiation of the monocytes and activation of the reporter gene.

- Freezing Medium: 95% fetal bovine serum (FBS), 5% DMSO
- Test Medium: RPMI 1640, 2 mM L-glutamine, 25 mM HEPES, 10% heat-inactivated fetal bovine serum, Pen-Strep (100 U/ml-100 µg/ml)

Required Selective Antibiotic $Zeocin^{(\!R\!)}$

TECHNICAL SUPPORT InvivoGen USA (Toll-Free): 888-457-5873 InvivoGen USA (International): +1 (858) 457-5873 InvivoGen Europe: +33 (0) 5-62-71-69-39 InvivoGen Asia: +852 3622-3480 E-mail: info@invivogen.com



Any questions about our cell lines? Visit our FAQ page.



Initial Culture Procedure

The first propagation of cells should be for generating stocks for future use. This ensures the stability and performance of the cells for subsequent experiments.

1. Thaw the vial by gentle agitation in a 37° C water bath. To reduce the possibility of contamination, keep the O-ring and cap out of the water. Thawing should be rapid.

2. Remove the vial from the water bath as soon as the contents are thawed, and decontaminate by dipping in or spraying with 70% ethanol.

Note: All steps from this point should be carried out under strict aseptic conditions.

3. Transfer cells in a vial containing 15 ml of pre-warmed growth medium (with 20% heat-inactivated FBS). Do not add selective antibiotics until the cells have been passaged twice.

4. Centrifuge vial at 150 x g (RCF) for 10 minutes.

5. Remove supernatant containing the cryoprotective agent and resuspend cells with 1 ml of growth medium (with 20% heat-inactivated FBS). 6. Transfer the vial contents to a 25 cm² tissue culture flask containing 5 ml of growth medium (with 20% heat-inactivated FBS). 7. Place the culture at 37°C in 5% CO₂.

Frozen Stock Preparation

1. Resuspend cells at a density of $5-7 \times 10^{\circ}$ cells/ml in freezing medium prepared extemporaneously with cold FBS.

2. Aliquot 1 ml cells into cryogenic vials.

3. Place vials in a freezing container and store at -80°C overnight.

4. Transfer vials to liquid nitrogen for long term storage.

Note: If properly stored, cells should remain stable for years.

Cell Maintenance

1. After cells have recovered and are growing well (after at least one passage), maintain and subculture the cells in growth medium. To maintain selection pressure, add $100 \,\mu$ g/ml of Zeocin[®] to the growth medium every other passage.

2. Pass the cells every 3 days by inoculating 5 x 10 $^{\rm s}$ cells/ml. Do not allow the cell concentration to exceed 2 x 10 $^{\rm o}$ cells/ml.

Notes: To ensure the best results:

- Use THP1-Lucia[™] ISG cells with less than 20 passages.

- Handling of cells should be as short as possible to prevent any damage resulting from the prolonged stay at room temperature without 5% CO_2 .

PMA-INDUCED DIFFERENTIATION (OPTIONAL)

Following Phorbol 12-myristate 13-acetate (PMA) treatment, THP1-Lucia[™] ISG cells are more sensitive to IFN-inducers, such as transfected Poly(dA:dT), while the response to TLR ligands, such as Pam3CSK4, is diminished.

Day 1

1. Add 180 μl of THP1-Lucia $^{\rm m}$ ISG cell suspension per well of a 96-well plate (1 x 10⁵ cells/well).

2. Treat THP1-Lucia^M ISG cells with 20 µl of PMA (final concentration 20-50 ng/ml) for 3 hours at 37°C in 5% CO₂.

3. Wash cells gently with pre-warmed and add 200 μI pre-warmed growth medium.

Day 4

4. Wash cells with pre-warmed PBS and add 180 μl growth medium.

5. Add 20 µl of an IFN inducer, such as Poly(dA:dT)/LyoVec™.

6. Incubate overnight at 37°C in 5% CO₂.

DETECTION OF IFN INDUCTION Sample Preparation

1. Resuspend all powdered samples in endotoxin-free water to avoid activation of TLR4 of the THP1-Lucia[™] ISG cell line.

2. Warm the samples at 37°C before use.

Notes:

Avoid testing of pure samples soluble only in ethanol or DMSO: these solutions are toxic to the cell line and can result in false negative results.
Samples containing a phosphatase activity cannot be tested as they can result in false positive results (like serum not previously heat-inactivated).

IFN Induction

1. Centrifuge cells at 150 x g (RCF) for 10 minutes or 300 x g (RCF) for 5 minutes.

2. Remove supernatant and resuspend THP1-Lucia^M ISG cells at 5x10⁵ cells/ml in fresh, pre-warmed growth medium.

3. Add 20 µl of sample per well including Poly(dA:dT)/LyoVec[™] as the positive control and endotoxin free water as a negative control (use new tips for each well to avoid cross-contamination).

4. Add 180 μl of cell suspension (1x10 $^{\rm s}$ cells) per well of a flat-bottom 96-well plate.

5. Incubate the plate at 37°C in a CO₂ incubator for 18-24 h.

6. Prepare QUANTI-Luc[™] 4 Reagent working solution following the instructions on the pouch.

7. Set the luminometer with the following parameters: 50 μ l of injection, end-point measurement with a 4 second start time and 0.1 second reading time.

8. Pipet 10 µl of THP1-Lucia[™] ISG cell culture supernatant per well in a 96-well white (opaque) or black plate, or a luminometer tube.

9. Prime the injector with QUANTI-Luc[™] 4 Reagent working solution and proceed with the measurement.

RELATED PRODUCTS

Product	Description	Cat. Code
Normocin™	Antimicrobial agent	ant-nr-1
Pam3CSK4	TLR2 ligand	tlrl-pms
Poly(dA:dT)/LyoVec™	RIG-I & CDS ligand	tlrl-patc
PMA	Phorbol myrisate acetate	tlrl-pma
QUANTI-Luc™ 4 Lucia/Gaussia	Luminesence detection kit	rep-qlc4lg1
Zeocin®	Selection antibiotic	ant-zn-1



Any questions about our cell lines? Visit our FAQ page.



QUANTI-Luc[™] 4 Reagent

A coelenterazine-based luminescence assay reagent

https://www.invivogen.com/quanti-luc

For research use only Version 23C24-AK

PRODUCT INFORMATION

Contents

• 1 tube of QUANTI-Luc[™] 4 Reagent (20X)

One tube of QUANTI-Luc[™] 4 Reagent is sufficient for 5 x 96-well plates (25 ml standard Flash/end-point detection).

Note: This sample cannot be sold separately from the QUANTI-Luc™ 4 Lucia/Gaussia kit.

QUANTI-Luc[™] 4 Lucia/Gaussia comprises two liquid components:

• QUANTI-Luc[™] 4 Reagent 20X (coelenterazine substrate)

• QUANTI-Luc[™] 4 Stabilizer 25X (optimized Glow assay reagent) Find more information at <u>https://www.invivogen.com/quanti-luc</u>.

Storage and Stability

- Store QUANTI-Luc[™] 4 Reagent at -20°C for up to 12 months.

– After preparation, the working solution is stable for 48 hours at 4° C and 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

Note: This product is photosensitive and should be protected from light.

Quality Control

Each lot is thoroughly tested to ensure the absence of lot-to-lot variation.

- Physicochemical characterization (pH, appearance).

- Functional assays using recombinant Lucia protein or reporter cells.

DESCRIPTION

QUANTI-Luc[™] 4 Reagent is a component of the QUANTI-Luc[™] 4 Lucia/Gaussia kit. It contains the coelenterazine substrate for the detection of secreted Lucia or Gaussia activity in live-cell supernatants, and of intracellular Renilla after cell lysis. The light signal produced correlates to the amount of luciferase protein expressed. It is quantified using a luminometer and expressed as relative light units (RLUs).

METHODS

Preparation of QUANTI-Luc[™] 4 Reagent working solution (1X):

1. Dilute the total volume of the 20X tube (1.25 ml) of Reagent into 23.75 ml of sterile water to obtain 25 ml of working solution. 2. Vortex **very briefly** (a few seconds).

3. Use the working solution immediately or store until required for use. QUANTI-Luc[™] 4 Reagent working solution can be stored for 48 hours at 4°C or 1 month at -20°C.

Flash detection of luciferase activity from cell culture medium:

To obtain **end-point readings** using a luminometer **with an injector**. 1. Set the luminometer with the following parameters: $50 \ \mu$ l of injection, end-point measurement with a 4 second start time and 0.1 second reading time.

2. Pipet 10-20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.

3. Prime the injector with QUANTI-Luc[™] 4 Reagent 1X and proceed **immediately** with the measurement.

To obtain **end-point readings** using a luminometer **without injectors**.

- 1. Set the luminometer with a 0.1 second reading time.
- 2. Pipet 10-20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.
- 3. Add 50 µl of QUANTI-Luc[™] 4 Reagent 1X to each well or tube.
- 4. Gently tap the plate several times to mix (do not vortex).
- 5. Proceed **immediately** with the measurement.

RELATED PRODUCTS

Product	Cat. Code
QUANTI-Luc™ 4 Lucia/Gaussia Kit 500 tests 2 x 500 tests 5 x 500 tests	rep-qlc4lg1 rep-qlc4lg2 rep-qlc4lg5

TECHNICAL SUPPORT InvivoGen USA (Toll-Free): 888-457-5873 InvivoGen USA (International): +1 (858) 457-5873 InvivoGen Europe: +33 (0) 5-62-71-69-39 InvivoGen Asia: +852 3622-3480 E-mail: info@invivogen.com

