

# Jurkat-Lucia™ NFAT-CD16 Cells

NFAT-CD16 Lucia Luciferase Reporter T Lymphocytes

Catalog code: jknl-nfat-cd16

<https://www.invivogen.com/jurkat-lucia-nfat-adcc-adcp-cells>

For research use only

Version 22C10-NJ

## PRODUCT INFORMATION

### Contents and Storage

- 1 vial of Jurkat-Lucia™ NFAT-CD16 cells (3-7 x 10<sup>6</sup> cells)

**IMPORTANT:** If cells provided in a cryovial are not frozen upon arrival, contact InvivoGen immediately.

- 1 ml of Blasticidin (10 mg/ml). Store at 4°C or at -20°C.\*
- 1 ml of Zeocin® (100 mg/ml). Store at 4°C or at -20°C.\*
- 1 ml of Normocin™ (50 mg/ml), a formulation of three antibiotics active against mycoplasmas, bacteria and fungi. Store at -20°C.\*
- 1 pouch of QUANTI-Luc™. Store QUANTI-Luc™ pouch at -20°C for 12 months. Reconstituted QUANTI-Luc™ medium is stable for 1 week at 4°C or for 1 month at -20°C. Keep reconstituted QUANTI-Luc™ away from light.

\*The expiry date is specified on the product label.

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

### 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

- Human CD16A expression has been verified by flow-cytometry.
- Induction of antibody-dependent cellular cytotoxicity (ADCC) has been validated using InvivoGen's anti-hCD20-hIgG1 antibody and Raji-Null target cell line.
- The stability for 20 passages following thawing has been verified.
- Jurkat-Lucia™ NFAT-CD16 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.

1. Leibson P.J., 1997. Signal transduction during natural killer cell activation: inside the mind of a killer. *Immunity*. 6:655. 2. Bruhns P. et al., 2009. Specificity and affinity of human Fcγ receptors and their polymorphic variants for human IgG subclasses. *Blood*. 113(16):3716. 3. Nagelkerke S.Q. et al., 2019. Genetic variation in low-to-medium-affinity Fcγ receptors: functional consequences, disease associations, and opportunities for personalized medicine. *Front. Immunol.* 10:2237. 4. Shaw J-P. et al., 1998. Identification of a putative regulator of early T cell activation genes. *Science*. 241:202.

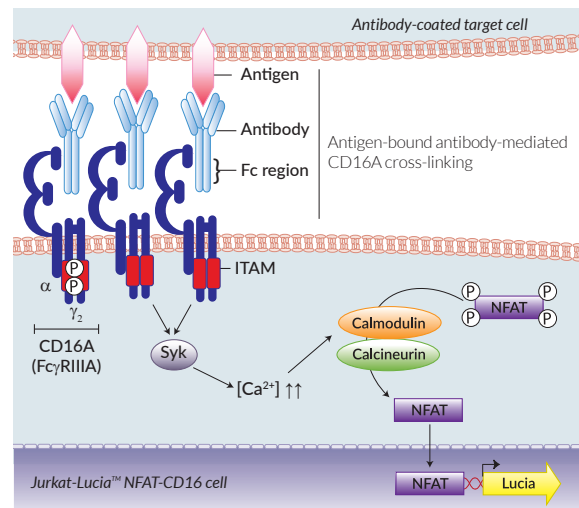
## INTRODUCTION

Antibody-dependent cellular cytotoxicity (ADCC) is an immune mechanism through which Fc receptor-bearing effector cells can recognize and kill antibody (Ab)-coated target cells expressing antigens on their surface. ADCC is triggered by the cross-linking between antigen-bound Abs and the Fc receptor CD16A (FcγRIIIA) at the surface of immune effector cells, such as Natural Killer cells. These interactions induce the increase of intracellular calcium concentrations and the translocation of the NFAT transcription factor to the nucleus, where it can bind to the promoter regions of ADCC relevant genes<sup>1</sup>.

CD16A features allelic polymorphisms among the human population, notably at position 158 in the mature protein (or position 161 in the full protein)<sup>2</sup>. The V158 allotype is reported to have higher affinity for monoclonal immunoglobulin G (IgGs) than the F158 allotype<sup>2,3</sup>.

## PRODUCT DESCRIPTION

Jurkat-Lucia™ NFAT-CD16 cells were engineered from the human T-lymphocyte Jurkat cell line. Jurkat cells naturally express a functional NFAT pathway<sup>4</sup>. Jurkat-Lucia™ NFAT-CD16 cells stably express the cell surface Fc receptor CD16A (FcγRIIIA; V158 allotype<sup>2,3</sup>) and the Lucia luciferase reporter gene under the control of an ISG54 minimal promoter fused to six NFAT response elements. These cells have been functionally tested with various target cells and specific monoclonal Ab isotype combinations. Antibodies displaying lower EC<sub>50</sub> have higher ADCC potency. These cells are resistant to Blasticidin and Zeocin®.



## 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](mailto:info@invivogen.com)



Any questions about our cell lines?

Visit our FAQ page.

 **InvivoGen**  
www.invivogen.com

## 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](mailto:info@invivogen.com).

## SAFETY CONSIDERATIONS

Biosafety Level 1

## HANDLING PROCEDURES

### Required Cell Culture Medium

- **Growth Medium:** IMDM, 2 mM L-glutamine, 25 mM HEPES, 10% heat-inactivated fetal bovine serum (FBS; 30 min at 56 °C), 100 µg/ml Normocin™, Pen-Strep (100 U/ml-100 µg/ml)
- **Freezing Medium:** 90% FBS, 10% DMSO
- **Test Medium:** IMDM, 2 mM L-glutamine, 25 mM HEPES, 10% heat-inactivated FBS, Pen-Strep (100 U/ml-100 µg/ml) **without Normocin, Blastcidin, and Zeocin®**

### Required Selective Antibiotics

- Blastcidin and Zeocin®

### 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 must 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 larger vial containing 15 ml of pre-warmed growth medium. **Do not add selective antibiotics until the cells have been passaged twice.**
4. Centrifuge cells at 150 x g (RCF) for 10 min.
5. Remove supernatant containing the cryoprotective agent and resuspend cells with 1 ml of growth medium.
6. Transfer the vial contents to a T-25 culture flask containing 5 ml of growth medium **without selective antibiotics.**
7. Place the culture at 37°C in 5% CO<sub>2</sub>.

### Frozen Stock Preparation

1. Resuspend cells at a density of 5-7 x 10<sup>6</sup> cells/ml in freezing medium freshly prepared 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. Jurkat-Lucia™ NFAT-CD16 cells grow in suspension.
2. After cells have recovered (after at least one passage), maintain and subculture the cells in growth medium. To maintain selection pressure, add 10 µg/ml of Blastcidin and 100 µg/ml of Zeocin® to the growth medium every other passage.
3. Pass the cells every 3 days by inoculating 2-5 x 10<sup>5</sup> cells/ml. Do not allow the cell concentration to exceed 2 x 10<sup>6</sup> cells/ml.

*Note: The average doubling time for the Jurkat-Lucia™ NFAT-CD16 cells is ~ 48 hours using the conditions described above.*

## Cell-Handling Recommendations

To ensure the best results, use Jurkat-Lucia™ NFAT CD16 cells with less than 20 passages.

## APPLICATION

Jurkat-Lucia™ NFAT-CD16 cells have been designed as effector reporter cells for InvivoGen's antibody-dependent cellular cytotoxicity (ADCC) assay using our expanding collection of Raji-derived target cells. For more information, visit <https://www.invivogen.com/raji-derived-target-cells>.

## ADCC REPORTER ASSAYS

Below is a protocol to perform an ADCC assay with Raji-Null cells which constitutively express human CD20 at the cell surface.

### Cell Preparation

Pass effector and target cells 2 days prior to the reporter assay.

1. Day -2: Resuspend Jurkat-Lucia™ NFAT-CD16 cells at 5 x 10<sup>5</sup> cells/ml, and Raji-Null cells at 4 x 10<sup>5</sup> cells/ml in pre-warmed test medium.
2. Incubate at 37°C in a CO<sub>2</sub> incubator for 48 h.
3. Day 0: Centrifuge Raji-Null cells at 300 x g (RCF) for 5 min.
4. Remove supernatant and resuspend at 1.1 x 10<sup>6</sup> cells/ml in pre-warmed test medium.

*Note: In steps 5 & 6, Jurkat-Lucia™ NFAT-CD16 cells should be prepared just prior to their addition to the antibody-coated target cells.*

5. Centrifuge Jurkat-Lucia™ NFAT-CD16 cells at 300 x g (RCF) for 5 min.
6. Remove supernatant and resuspend at 2.2 x 10<sup>6</sup> cells/ml in fresh, pre-warmed test medium.

*IMPORTANT: To ensure reproducible results, homogenize the cell suspensions.*

### ADCC Induction

Below is a protocol for end-point readings using a luminometer. This protocol can be adapted for use with kinetic measurements.

1. Add 20 µl of test anti-hCD20 mAb per well including a positive control (e.g. Anti-hCD20 IgG1) and a negative control (e.g. Anti-β-galactosidase IgG1).

*Note: We recommend to prepare 1:4 or 1:2 dilution series.*

2. Add 90 µl of Raji-Null cell suspension (~100,000 cells) per well of a flat-bottom 96-well plate.
3. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 1 h.
4. Add 90 µl of Jurkat-Lucia™ NFAT CD16 cell suspension (~200,000 cells) per well.
5. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 6 h.
6. Prepare QUANTI-Luc™ following the instructions on the data sheet.
7. Transfer 20 µl of co-incubated Raji-Null and Jurkat-Lucia™ NFAT CD16 cell supernatant into a 96-well white (opaque) or black plate, or a luminometer tube.
8. Add 50 µl of QUANTI-Luc™ per well.
9. Proceed **immediately** with the measurement.

## RELATED PRODUCTS

Product	Description	Cat. Code
Anti-β-Gal-hlgG1	Control antibody	bgal-mab1
Anti-hCD20-hlgG1	Anti-hCD20 antibody	hcd20-mab1
Blastcidin	Selection antibiotic	ant-bl-05
QUANTI-Luc™	Lucia detection medium	rep-qlc1
Raji-Null cells	ADCC target cell line	raji-null
Zeocin®	Selection antibiotic	ant-zn-05

## 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](mailto:info@invivogen.com)



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

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)

# QUANTI-Luc™

A coelenterazine-based luminescence assay reagent

Catalog code: rep-qlc1, rep-qlc2

<https://www.invivogen.com/quantiluc>

For research use only

Version 19A04-MM

## PRODUCT INFORMATION

### Contents

QUANTI-Luc™ is provided as packs of individually sealed pouches.

- rep-qlc1: 2 pouches of QUANTI-Luc™
- rep-qlc2: 5 pouches of QUANTI-Luc™

Each pouch contains everything needed to prepare 25 ml of reagent allowing the preparation of 500 wells of a 96-well plate.

### Storage and Stability

- Store QUANTI-Luc™ pouches at -20°C for 12 months.
- Reconstituted QUANTI-Luc™ is stable for 1 week at 4°C and for 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

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

## DESCRIPTION

QUANTI-Luc™ is an assay reagent containing all the components required to quantitatively measure the activity of Lucia luciferase and other coelenterazine-utilizing luciferases. QUANTI-Luc™ contains the coelenterazine substrate and stabilizing agents for the luciferase reaction. The light signal produced is quantified using a luminometer and expressed as relative light units (RLU). The signal produced correlates to the amount of luciferase protein expressed, indicating promoter activity in the reporter assay.

QUANTI-Luc™ is optimized for use with Lucia luciferase reporter cell lines. Lucia luciferase is a secreted coelenterazine luciferase encoded by a synthetic gene. As Lucia luciferase is secreted, it can be directly measured in the cell culture medium using bioluminescent assays.

InvivoGen provides a recombinant Lucia luciferase protein (see Related Products) which is a positive control for QUANTI-Luc™. A dilution series of the recombinant Lucia luciferase protein can also be used to determine the linear range of the assay.

## METHODS

### Preparation of QUANTI-Luc™

1. Pour the pouch contents into a 50 ml screw cap tube.
2. Add 25 ml of sterile water.
3. Swirl product gently until powder is completely dissolved.
4. Use QUANTI-Luc™ assay solution immediately or store until required for use. Reconstituted QUANTI-Luc™ can be stored for 1 week at 4°C and for 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

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

### 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 µl 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 the QUANTI-Luc™ assay solution 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™ assay solution 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	Catalog Code
QUANTI-Luc™ Gold (For standard and HTS assays)	rep-qlcg1
pSelect-zeo-Lucia™ (expression plasmid)	psetz-lucia
Recombinant Lucia luciferase protein	rec-lucia
<b>Reporter Cells</b>	
THP1-Dual™ (IRF-Lucia/NF-κB-SEAP) Cells	thpd-nfis
THP1-Lucia™ NF-κB Cells	thp1-nfkb

For a complete list of InvivoGen's Lucia luciferase Reporter Cell Lines visit <https://www.invivogen.com/lucia-reporter-cells>.

## 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](mailto:info@invivogen.com)

