

## Protocol

# IncuCyte® – iQue3 Workflow for the Immune Cell Killing Assay

This protocol provides a detailed workflow for the combination of the IncuCyte and iQue3 to facilitate measurement of immune cell mediated killing of adherent or non-adherent target tumor cells and T-cell activation (TCA). This assay workflow describes how the two platforms can be used both simultaneously and concurrently to provide a deeper understanding of immune cell killing (ICK). The workflow allows for (1) direct measurements of tumor cell death

with a no-wash, mix-and-read protocol using the IncuCyte ICK application and co-culture methodology; (2) quantification of morphological and spatial data using the IncuCyte Cell-by-Cell analysis software and (3) quantification of T-cell activation markers and cytokine release using the Intellicyt Human T Cell Activation Cell and Cytokine Profiling Kit (TCA kit). The ICK assay can be directly transferred between the two platforms with minimal sample manipulation.

### Required materials

- Target cells of interest: cells are required to be stably expressing the NuLight Green Lentivirus (Essen BioScience Cat. No. 4475)
- Poly-L-ornithine (Sigma Cat. No. P4957) for coating plate when using non-adherent target cells
- IncuCyte Annexin V Red Reagent (Essen BioScience Cat. No. 4641)
- Immune (effector) cells of interest
- Effector cell culture media
- Effector cell activator (e.g. CD3/CD28 Dynabeads®)
- Human T Cell Activation Cell and Cytokine Profiling Kit-TCA kit (Intellicyt Cat. No. 90560)
- QBeads® if temporal cytokine analysis desired (Intellicyt Assay Builder: IFN $\gamma$  option 1 and TNF $\alpha$  option 2).
- 96-well flat-bottom microplate (e.g. Corning Cat. No. 3595)
- 96-well v-bottom microplate (e.g. Costar Cat. No. 3363)
- Accutase (e.g. Gibco Cat. No. A1110501)

### Suggested materials

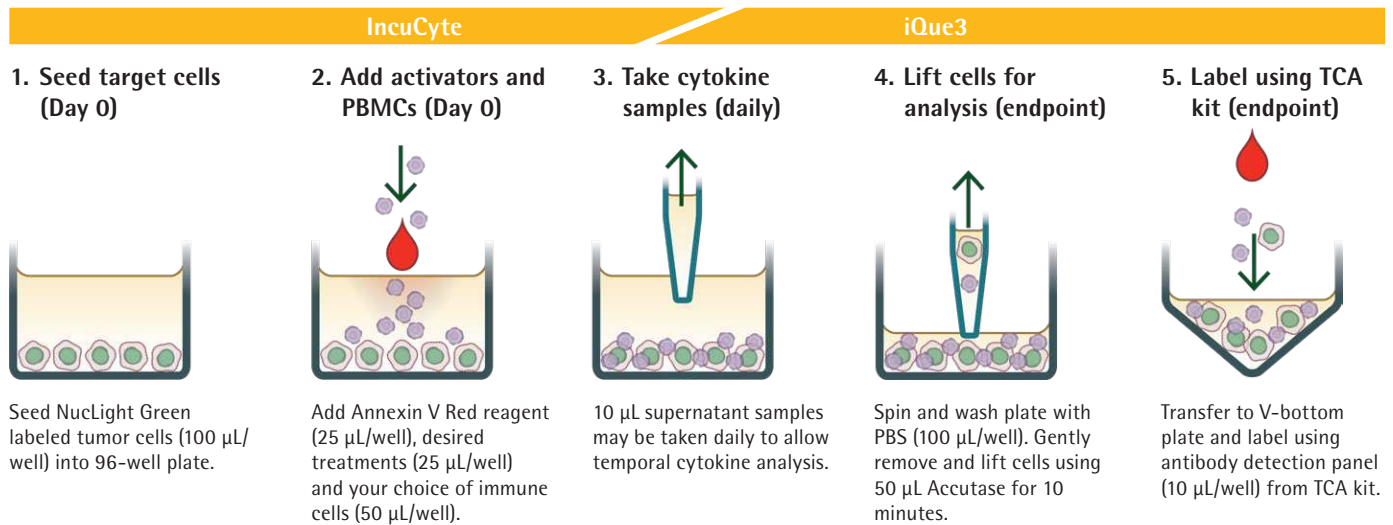
- Plate washer: a number of points in the protocol require a wash step and while plates can be 'flicked' in order to remove unwanted supernatant, a plate washer is highly recommended.

NOTE: There is no recommended plate washer and instrument settings must be optimized before use.

## General guidelines

- Following cell seeding, place plates at ambient temperature (15 minutes for adherent cell lines and 45 minutes for non-adherent cell lines) to ensure homogenous cell settling.
- It is strongly recommended that the following controls are included on the plate: (1) a non-activated PBMC control in co-culture with target cells (2) target cell in mono-culture (3) activated PBMC in mono-culture.
- Remove bubbles from all wells by gently squeezing a wash bottle (containing 70–100% ethanol with the inner straw removed) to blow vapor over the surface of each well.
- After placing the plate in the IncuCyte® Live-Cell Analysis System, allow the plate to warm to 37° C for 30 minutes prior to scanning.
- For the final cytokine sample the 10 µl volume must be taken prior to cell transfer and assayed in a separate plate to the antibody labeled cells.
- Cytokine standards require fresh cell media to ensure the reproducibility and reliability of your data.
- The optional Cell Proliferation and Encoder Dye (B/Green) from TCA kit cannot be included in this experiment due to the use of NucGreen target cells.

## Immune cell killing of adherent or non-adherent tumor cells protocol



### 1. Seed target cells

#### 1.1 Coat plate for non-adherent cells:

Coat a 96-well flat bottom plate with relevant coating matrix. It is recommended that wells are coated using 50 µL of 0.01% poly-L-ornithine solution. Coat plates for 1 hour at RT, remove solution from wells, allow plates to dry for 30 minutes prior to cell addition.

- 1.2 **Seed target cancer cells:** An appropriate density of cells must be seeded (100 µL/well). 2,000 to 5,000 cells/well for adherent cells (seeded in the morning) and 10,000 to 20,000 cells/well for non-adherent (seeded prior to assay) are reasonable starting points.

**NOTE:** NucLight Green expressing cells must be used to allow integration of IncuCyte and iQue3. NucLight Red cells are not compatible with the Human T Cell Activation Cell and Cytokine Profiling Kit (TCA kit).

### 2. Reagent and treatment preparation

- 2.1 Prepare the following reagents in medium:
- Test materials (e.g. T-cell stimulants, antibodies, cytokines; 25 µL/well, prepared 8x FAC).
  - Apoptosis detection reagent, IncuCyte Annexin V Red (Cat. No. 4641): solubilize Annexin V by adding 100 µL of complete medium or PBS. The reagent can then be diluted as required in complete medium containing at least 1 mM CaCl<sub>2</sub> (25 µL/well at 8x FAC, for a final dilution of 1:200).
- 2.2 Add all reagents to assay plate to achieve a volume of 150 µL/well.

### 3. Add immune cells

- 3.1 Prepare an effector cell suspension at an appropriate density. It is recommended that different target-to-effector ratios are tested (e.g., 1:3, 1:5).
- NOTE:** Cell Proliferation and Encoder Dye (B/Green) from the TCA kit cannot be included in this assay due to the presence of NucGreen Target cells.
- 3.2 Seed 50 µL/well to achieve a total assay volume of 200 µL. Allow plates to settle on level surface at RT for 30 minutes.
- 3.3 Place the assay plate into the IncuCyte Live-Cell Analysis System, allow to warm to 37° C for 30 minutes. Schedule 24 hour repeat scanning:
- Scan Type: Non-adherent Cell-by-Cell or Standard (4-images per well)
  - Objective: 20x
  - Channel selection: Phase + Green + Red (if including apoptosis reagent)
  - Scan interval: 3 hours

#### 4. Optional: Daily cytokine samples

- 4.1 Combine lyophilized standards from the QBeads kit and add 200 µL fresh culture media incubate at RT for 15 minutes without trituration.
- 4.2 Prepare an 11-point 1:3 serial dilution of solubilised standards. Plate 10 µL of each dilution into duplicate wells of a V-bottom plate, with concentration going from low to high (Rows A and B, 2-12). Place 10 µL of media only in 1st wells as a control.
- 4.3 10 µL of supernatant from each sample on the assay plate should be transferred to the remaining empty wells of the V-bottom plate, i.e. Rows C-H.  
**NOTE:** Ensure removal and replacement of plate from IncuCyte will not interrupt scheduled 3 hour scans.
- 4.4 Capture beads are diluted by 18 times volume with fresh media. Add 190 µL/well and incubate at RT in the dark for 1 hour.
- 4.5 Centrifuge at 300xg 5 minutes and aspirate supernatant. Re-suspend beads in residual liquid by strong shaking (3000 rpm 1 minute).
- 4.6 Add 10 µL/well Cytokine Detection, followed by a quick spin (300xg 5 seconds) and brief shake (2000 rpm 20 seconds). Leave in the dark at RT for 2 hours.
- 4.7 Add 100 µL/well wash buffer, spin at 300xg 5 minutes and aspirate.
- 4.8 Re-suspend beads in residual liquid and add 20 µL/well wash buffer.
- 4.9 Run plate on iQue3 with TCA template. Optimization of protocol may be necessary to achieve required results. A minimum sip time of 4 seconds is recommended.

#### 5. Endpoint T-cell subset and cytokine analysis

- 5.1 For endpoint cytokine analysis, remove 10 µL of supernatant from each sample well and follow section 4.
- 5.2 Centrifuge plate at 300xg for 5 minutes and remove supernatant gently using multichannel pipette, leaving at least 50 µL media in well.
- 5.3 Wash once with PBS and centrifuge at 300xg for 5 minutes to ensure all cells remain in the wells.
- 5.4 Gently remove PBS and add 50 µL/well Accutase for 10 minutes at 37° C followed by a 2 minute shake at 1400 rpm.
- 5.5 Add 100 µL media to quench Accutase, triturate sample and then transfer to a V-bottom plate.  
**NOTE:** Use a microscope to check that cells have been lifted and transferred from the cell culture plate, if cells still remain repeat steps 5.3-5.4.
- 5.6 Centrifuge plate at 300xg for 5 minutes, aspirate supernatant and re-suspend in 10 µL of fresh media.
- 5.7 Dilute Cell Viability Dye 1:500 into Antibody Detection Panel. Add 10 µL/well followed by quick spin (300xg 5 seconds), brief shake (2000 rpm 20 seconds). Leave at RT in the dark 1 hour.
- 5.8 Add 100 µL wash buffer and centrifuge at 300xg for 5 minutes.
- 5.9 Re-suspend cells in residual liquid and add 20 µL/well wash buffer.
- 5.10 Run plate on iQue3 with TCA kit template. Optimization may be necessary to achieve required results. A minimum sip time of 4 seconds is recommended.

#### 6. Analysis using ForeCyt

- 6.1 TCA kit comes with a templated analysis. This includes automated compensation that should not require changes for this assay.
- 6.2 A change in the gating should be included in order to exclude your target cells from the gates:
  - a. All events: select cell populations using SSC-H vs FSC-H.
  - b. Single cells: identify single cells using FSC-H vs FSC-A.
  - c. NucGreen exclusion: BL1-H vs FSC-H will identify two populations. Low green are effector cells and high green target cells.
  - d. Live/Dead cells: RL1-H vs SSC-H can then be used to exclude any dead cells.
  - e. Continuation of TCA kit gates: copy gates onto new population to allow exclusion of target cells.

A complete suite of cell health and T-cell biology applications is available to fit your experimental needs. Find more information at [www.sartorius.com/incucyte](http://www.sartorius.com/incucyte) and [www.sartorius.com/intellicyt](http://www.sartorius.com/intellicyt)

For additional product or technical information, please email us at [AskAScientist@sartorius.com](mailto:AskAScientist@sartorius.com)

Essen Bioscience,  
a Sartorius Company  
300 West Morgan Road  
Ann Arbor, Michigan, 48108  
USA

[www.sartorius.com/incucyte](http://www.sartorius.com/incucyte)  
[www.sartorius.com/intellicyt](http://www.sartorius.com/intellicyt)  
Email: [AskAScientist@sartorius.com](mailto:AskAScientist@sartorius.com)  
USA +1.734.769.1600  
UK +44.1707.358688  
Japan, Australia and other countries  
around the world: +65.6872.3966