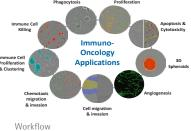
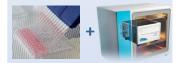
## Quantitative Live-Cell Imaging Assays for Immmunotherapy: Chemotaxis, Immune Cell Killing and Phagocytosis

G. Lovell, N. Bevan, C. Szybut, N. Rodinova, D. Appledorn, M. Tikhoneko, L. O'Clair, T. O'Callaghan, T.Dale, and D. Trezise Essen BioScience R&D, Welwyn Garden City, UK and Ann Arbor, MI, USA

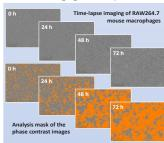


## Micro-titre Plate Models



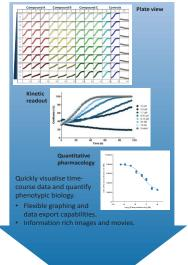
- Live cell protocols enable real time measures within your incubator - novel instrumentation, reagents & consumables.
- Simple mix-and-read methods no wash, no fix, no cell lifting.
- Analyse up to 6 plates in the IncuCyte<sup>®</sup> system.

Automated Imaging and Analysis



- Time-lapse images taken from every well and automatically analysed.
- Non-perturbing, non-invasive and direct measures of phenotypic cell biology.
- High definition phase contrast image processing and fluorescence object quantification.

## Real Time Quantification



Summary and Impact

- For the body to defend and fight against cancer, immune cells must recognise, engage, destroy and ultimately remove unwanted tumour cells. Understanding
- these processes and interactions at the cellular level is central to identifying and validating new drug targets and cellular therapy approaches.
- Essen BioScience offers a flexible range of phenotypic assays to explore all aspects of the immuno-oncology research area. All of these assays are based on noninvasive live-cell analysis of cells in 96-well micro-plates using an IncuCyte<sup>®</sup> ZOOM live cell analysis system.
- Here we describe a cluster of new assays for quantifying immune cell biology and interactions with tumour cells.

Immune Cell Killing

- Phase-contrast/fluorescence images from cells gathered over time are processed to measure phenotypes, such as immune cell proliferation, migration & death.
  Experiments are performed with low cell numbers using simple mix and read formats which are non-perturbing to the cell model.
  - Experiments are performed with low centralities using simple mix and read formats which are non-performing to the centrological model. Each of these approaches provide a full time-course of the biology; images and time-lapse movies provide credibility and valuable biological insight. Added throughput and automated image analysis enhances productivity.

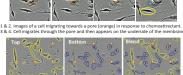


PRINCIPLE: Cells seeded onto the top surface of the IncuCyte™ ClearView membrane insert and migrate towards chemoattractant in the lower reservoir plate.

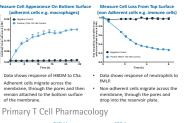
Cell movement monitored over time, imaging every 0.5-2 h. IncuCyte® 200M live cell analysis system images and quantifies the top and bottom surface of the membrane. Cell movement is detected as a loss of cell area from the top surface or an increase in cell area on the bottom surface.

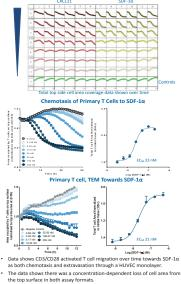
APPLICATIONS: Measure migration, invasion or trans-cellular migration (e.g. TEM) with both adherent and non-adherent cells.

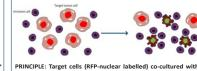
## Visualise and Quantify Images



Automated image processing separates cells located on the top surface (yellow) and bottom surface (blue) of the membrane, pores shown (orange). Images are processed ccquisition.



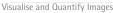


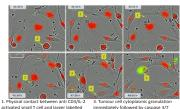


PRINCIPLE: Target cells (KPP-nuclear Tabelled) co-cultured with immune cells (T cells, NK, PBMC), with various activators and the IncuCyte® ZOOM Caspase 3/7 apoptosis reagent (green).

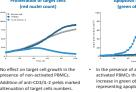
Cells monitored over time, imaging every 2 h. IncuCyte® ZOOM live cell analysis system images and quantifies phase and fluorescence images Target cell number is quantified as the number of red objects (nuclei), apoptosis by counting the green-labelled nuclei.

APPLICATIONS: Measure T-cell killing/ADCC in adherent or non adherent target cells.

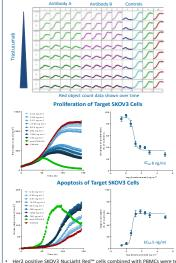




Wated shall i Cell and larger labelled
 minicusery nonwea or cospase or
 VO3 Nuccight Red" cell. T cell division.
 Iumour cells under attack from T cells.
 Proliferation of target cells
 Apoptosis of target cells



Antibody-Dependent Cell Cytotoxicity



 Her2 positive SKOV3 Nuclight Red<sup>™</sup> cells combined with PBMCs were tested in the presence of Trastuzumab to induce antibody-dependent cell-mediated cytotoxicity (ADCC).
 Concentration\_dependent decrease in proliferation and increase in apontosis

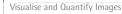
Concentration-dependent decrease in proliferation and increase in apoptosi
 No response was seen in Her2 negative cells (A549; lung carcinoma).

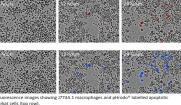


PRINCIPLE: pHrodo<sup>®</sup> labelled cells added to phagocytes (e.g. macrophages), phagocytosis is initiated following receptor activation and pHrodo<sup>®</sup> labelled cells are engulfed.

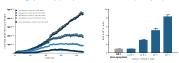
Cells monitored over time, imaging every 15-30 min. Little or no fluorescence is measured while target cells are in the extracellular environment (pH 7.4). Once in the acidic phagosome (pH 4.5-5.5) there is an increase in pHrodo<sup>®</sup> fluorescence.

APPLICATIONS: Measure phagocytosis, efferocytosis and ADCP of pHrodo® labelled target cells.



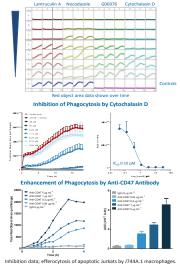


jurkat cells (top row). Masking of red fluorescent objects (blue) for analysis and enabling quantification (bottom row). Phagocytosis of anontotic pHrodo<sup>®</sup> labelled lurkat cells (efferocytosis)



 Addition of apoptotic cells to the J744A.1 macrophages results in an increase in fluorescent area as labelled cells transition into the acidic phagosome .
 Amplitude of fluorescent signal is proportional to the number of target cells added.
 The fluorescent signal is minimal with the addition of non-anontotic lurkat cells.

Modulation of Phagocytosis



Inhibition data; efferocytosis of apoptotic Jurkats by J744A.1 macrophages. Enhancement data; phagocytosis of CCRF-CEM cells in the presence of increasing concentrations of anti-CD47 by BMDM. Inclusion of anti-CD47 binds to the "don't eat me" signal on CCRF-CEM to promote phagocytosis by the macrophages.