

1. ABSTRACT

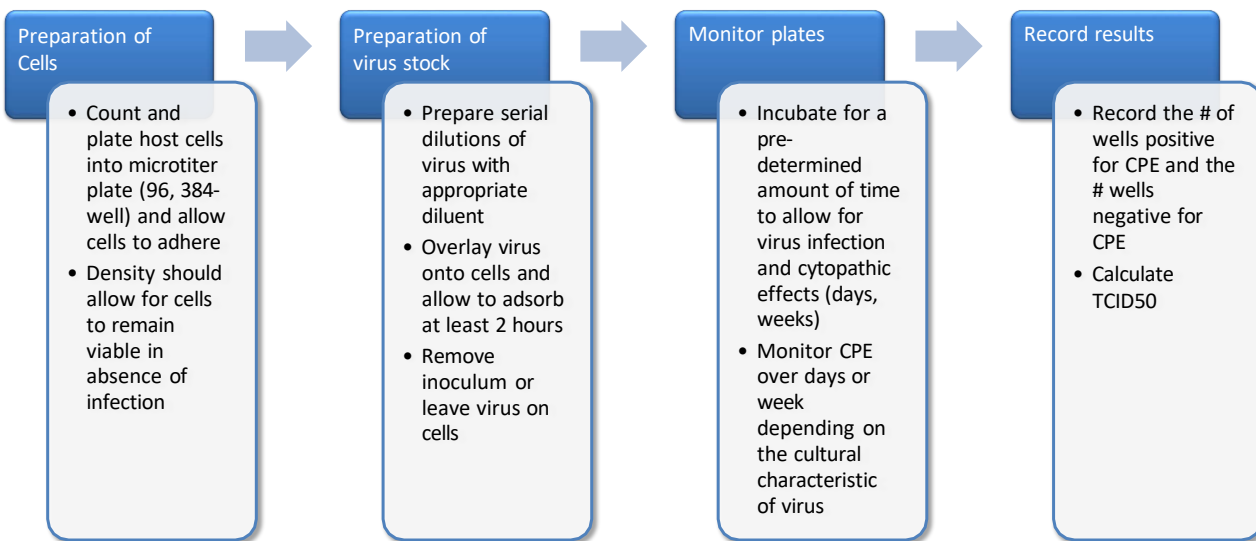
Qualitative assessment of pathogenic infections, as well as the protective effects of candidate therapeutics, often relies on the observation of specific pathological changes within the host cell. These phenotypic changes, such as cell rounding, swelling/shrinking, granularity, etc., are known as a Cytopathic Effect (CPE) and can be visualized via light microscopy. As the magnitude and localization of the CPE may vary considerably, careful examination of replicate samples at various titers is required for reliable, qualitative results. This subjective approach which is specific to the infectious agent as well as the host cell is tedious, time consuming and low throughput, requiring manual well-by-well examination by highly skilled personnel.

Revvity's Celigo imaging cytometer has been applied to provide automated, rapid assessment of viral infectivity in a range of plate formats. Using f-theta optics, Celigo provides high quality, whole well images using brightfield and/or fluorescent illumination. Automated segmentation and analysis provides quantitative output of CPE based on characteristic changes to the host cell monolayer.

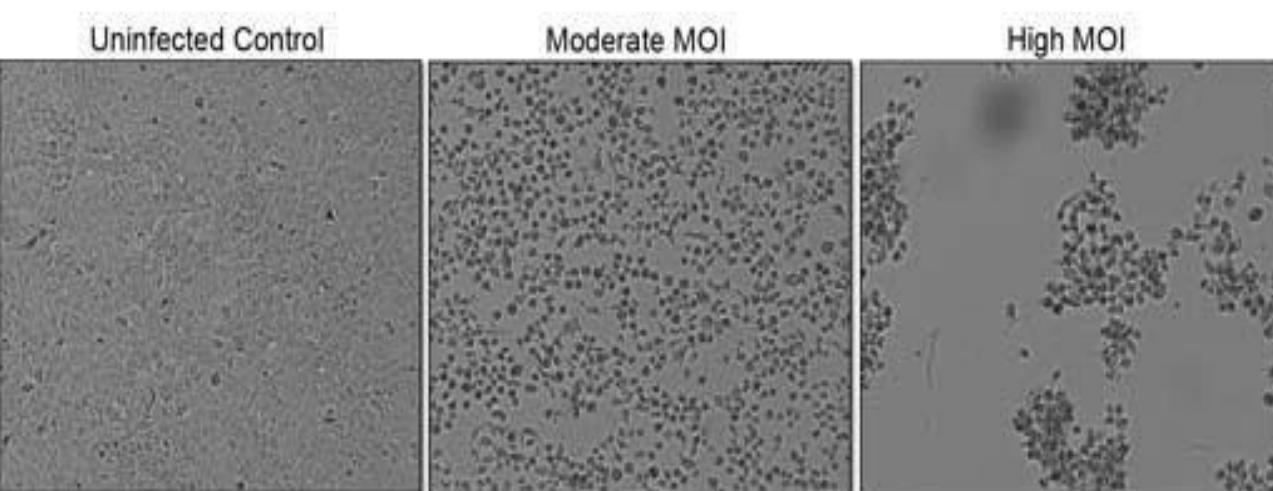
Celigo provides several key benefits;

- Objective segmentation and quantitative output of magnitude of infection
- Automated sample analysis reduces time, labor and variability
- High throughput and scaleable – less than 5 min for 96 or 384 well plate
- Capture high resolution, whole well images for documentation or manual assessment of CPE
- Supports related fluorescent based functional assays relevant to infection (e.g. expression analysis, apoptosis)

2. TCID₅₀ ASSAY PROTOCOL

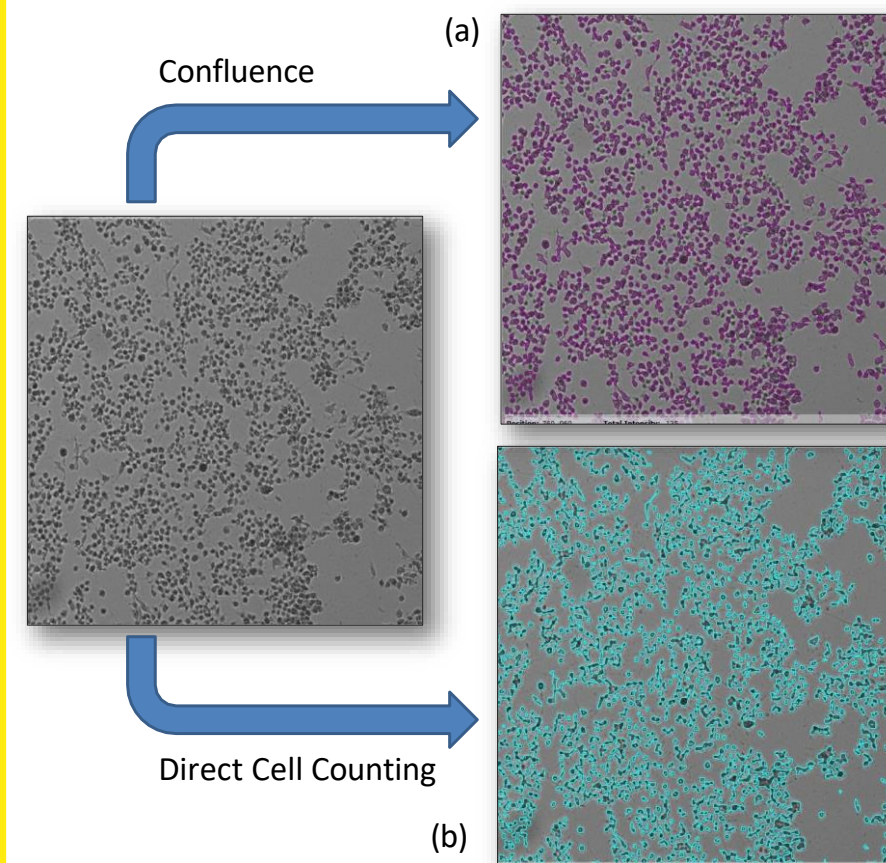


3. BRIGHT FIELD IMAGING



- BVDV (Bovine Viral Diarrhea Virus) Infection of BK6 (Bovine kidney) cells
- 96-well plates were scanned on the Celigo using the Cell Counting Application
- A dark focus was selected to better separate the healthy monolayer from the infected cells

4. DETERMINING BEST ANALYSIS METHOD TO SCORE CPE



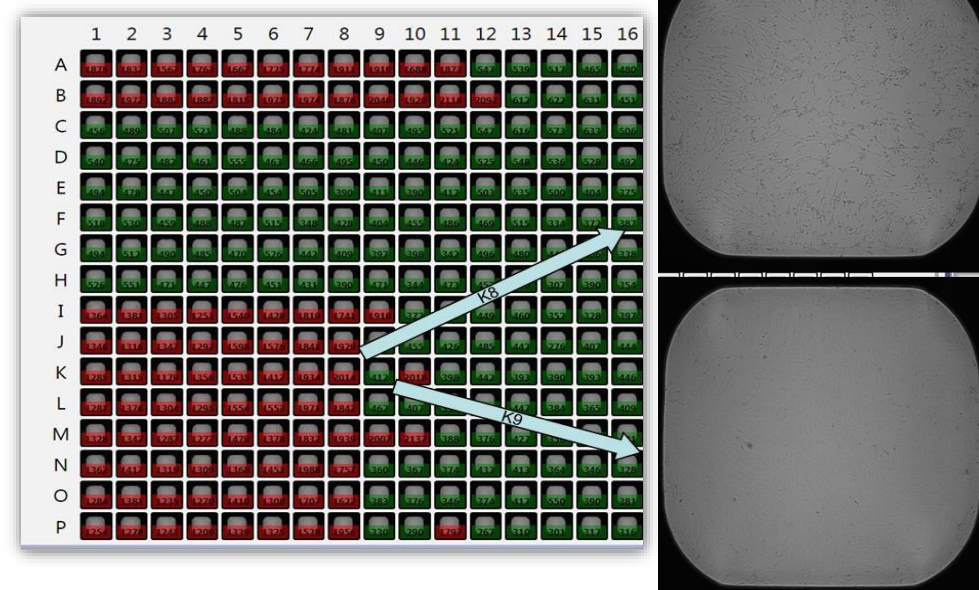
- Each cell/virus pair will have distinct morphological changes
- Analysis setting can be created and saved for each pair
- CPE was scored + based on the appearance of dark, rounded cells
- In this example, the reduction in confluence (a) and the increase in sick/dying cells (b) scored equally well

5. RESULTS TAB DISPLAYS PLATE AND WELL LEVEL

384 Well Plate (Corning 3712)
Plate set up: Rows A&B, Titrated virus control
Rows C-H, Negative control
Rows I-P, Virus inoculum (1) 5×10^{-1} → (16) 1×10^{-8}

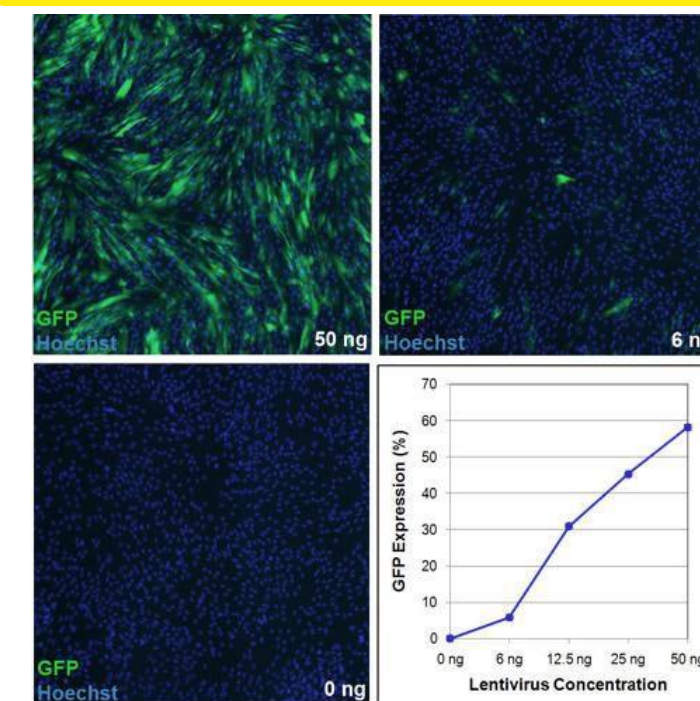
Plate Level view

Well Level view



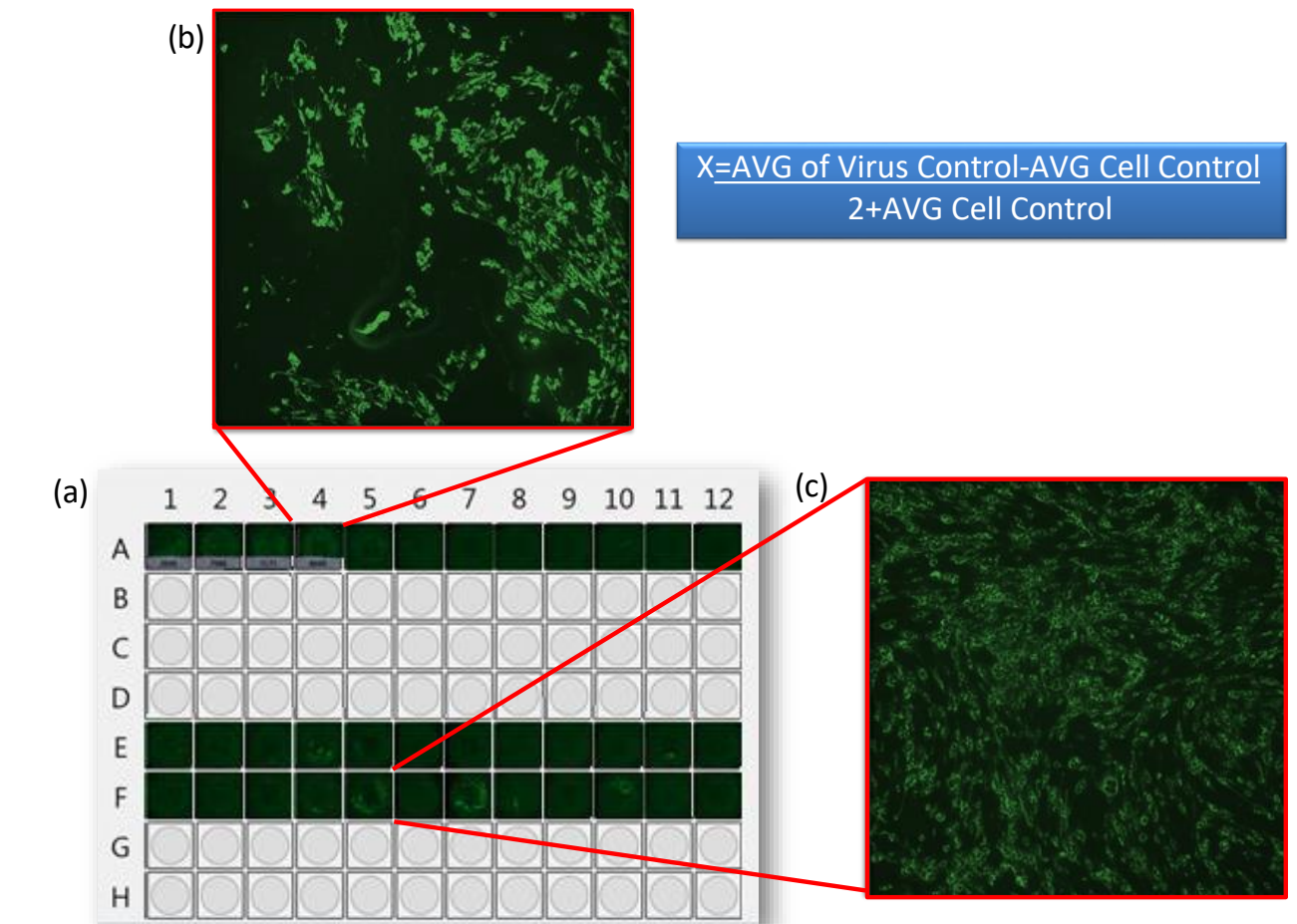
- Bovine Viral Diarrhea (BVD) infection of Bovine Turbinate cells (BT2)
- Threshold function indicates red wells (infected) vs. green wells (uninfected)
- Breakpoint of infection is clearly visible at plate level
- Well level data is exported as CSV
- Celigo results correlated 100% with assessment via light microscope

6. MOI BY GFP-INDICATOR VIRUS



- Known concentrations of viral stock (ng/ml) are used to infect primary fibroblasts. Infectious titre is calculated by measuring the % of infected cells/dose.
- Viability and overall cell health can also be determined by staining with propidium iodide for sick/dying cells (not shown)
- The Celigo can scan and report up to 4 separate channels at one time (Brightfield, Red, Green, Blue)

7. FLUORESCENT TITERS FROM IFA



- Celigo allows for multiple analyses/plate to be carried out in one experiment. (a) Result thumbnails following scan are displayed. Zoomed images of infection by BVD1 (b) and BVD2 (c) demonstrate the varied morphologies of infected cells as well as the localization of the viral antigen
- Antibodies directed against viral antigen are often used as indicators of infection. Unlike TCID₅₀ where a single event of infection is scored as a positive, these assays measure % infection. Intensity of FL signal may also be used to indicate strength of infection.

8. CALCULATION OF TCID₅₀

Definition: 50% tissue culture infectious dose of a virus

Spearman-Kärber formula:

$$M = xk + d [0.5 - (1/n)(r)]$$

xk = dose of the highest dilution.
 r = sum of the number of "n" responses.
 d = spacing between dilutions.
 n = wells per dilution.

Reed and Muench formula:

% infection immediately above 50% - 50%

% infected immediately above 50% - % infected immediately below 50%

- Determination of TCID₅₀ MOI and neutralization assays are easily calculated by exporting the well-level data as CSV into a number of online calculators

9. CONCLUSION

The Celigo[®] S Imaging cytometer is a benchtop *in situ* cellular analysis system that rapidly provides high integrity whole well images for routine brightfield and fluorescent cellular analysis with a simple workflow.

- Settings for each cell type and virus can be saved and re-used to assure objective results
- Additional information such as cell health, cell cycle status can be determined from the same plates using Revvity designed applications.
- Time course studies can be carried out by scanning plates over several days/weeks allowing for additional information than may be generated by carrying out end-point assays alone.
- User-friendly and intuitive software allows even those with little imaging experience to generate valuable data.

