

# Performance of a Label-Free Image-Based 2D Scratch Wound Healing Assay to Monitor Cell Migration and its Inhibition

## Author

Leonie Rieger and Brad Larson  
Agilent Technologies, Inc.

Ulf Rädler  
ibidi GmbH

## Introduction

Cell migration, the random movement of cells, plays a central role in physiological and pathological processes. Cell migration is an integral part of basic biological processes such as tissue formation, wound healing, and the immune response, as well as processes such as cancer invasion and metastasis.<sup>1,2</sup> The basic principal of cell movement is a cyclical 5-step procedure starting with cell polarization and protrusion of the leading edge. In the second step, the leading edge attaches to the substrate, followed by the proteolytic degradation of physical barriers, for example, tissue components and actomyosin contraction. Finally, the cell rear slides forward. The main modification of collective cell movement is the remaining intercellular cell-cell connections of the leading edge. Therefore, the principle steps are simultaneously coordinated.<sup>2</sup> An advanced knowledge and understanding of the mechanism of collective cell migration is crucial to develop novel therapeutic strategies to control, for example, cellular invasion, the key role in metastasis of cancer.

A powerful tool to investigate cell migration is the wound healing assay. After creating an artificial gap into a confluent cell monolayer, either in the form of a scratch or by removing a growth barrier, the assay follows cell migration from the edges over time until the establishment of new cell-cell contact, and the gap is closed again. The extent of wound healing and wound healing rate provide information about the migration characteristics, cell health, and ability to interact with other cells, in addition to how test-molecules can affect each of these parameters.

In this application note experiments were performed to track the ability of various cell types to move into a created gap, or heal the wound. In addition, inhibition of wound healing with Cytochalasin D is also shown. Culture-Inserts and 35 mm  $\mu$ -Dishes, provided by ibidi, were incorporated to perform the wound healing experiment process. Wound closure was monitored using an Agilent BioTek Lionheart FX automated microscope, and cellular analysis of the kinetic images captured was automatically performed using Agilent BioTek Gen5 microplate reader and imager software.

## Materials and methods

### Materials

**ibidi Culture-Insert:** The ibidi Culture-Insert was mainly developed to replace classical scratch assays. Silicon inserts, placed on a cell culture surface, provide two cell culture reservoirs. These two chambers are separated with a 500  $\mu\text{m}$  thick wall. The culture insert sticks to the bottom of a  $\mu\text{-Dish}$  or 24-well plate due to a specially designed surface. Due to the adhesive design, the cells are not able to grow under the wall. For the experiments performed, the Culture-Inserts 2-well in a  $\mu\text{-Dish}^{35\text{ mm, high}}$  (part number 81176) were used and provided by ibidi GmbH (Munich, Germany).

**Cells:** HT-1080 fibrosarcoma cells (part number CCL-121) and immortalized keratinocytes (part number CRL-2309) were purchased from ATCC (Manassas, VA). RFP expressing human neonatal dermal fibroblasts (part number cAP-0008RFP) were obtained from Angio-Proteomie (Boston, MA).

**Wound healing inhibitor:** The compound Cytochalasin D (part number 1233, R&D Systems; Minneapolis, MN) was used to inhibit cell migration due to its interfering with actin polymerization and potent disruption of actin filament function.<sup>3</sup>

### Agilent BioTek Lionheart FX automated microscope:

The Lionheart FX automated microscope with augmented microscopy provides digital microscopy with high magnification images, using up to 100x objectives. In addition, the microscopy module provides microscopy in fluorescence, brightfield, color brightfield, and phase contrast. Due to environmental control, including temperature, gas control, and a humidity chamber, the system is optimized for long-term kinetic assays and live-cell imaging. The instrument was used to monitor wound closure in phase contrast.

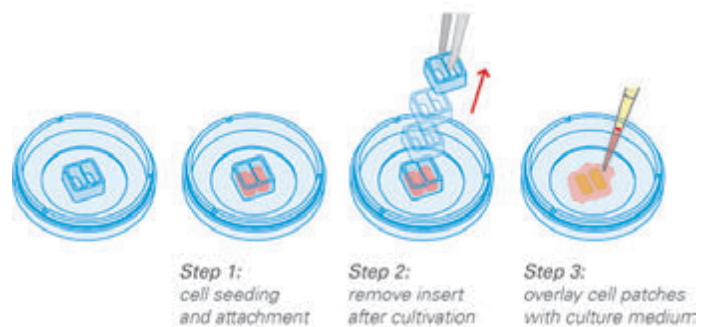
### Agilent BioTek Gen5 microplate reader and imager software:

For automated image capturing and analysis, the Gen5 3.0 microplate reader and imager software was used. Image acquisition is completely automated from sample translation, focusing, and exposure control. The included cell analysis tool calculates precise values in variable parameters, for example, object area and size, as well as the object intensity. Using the customized transformation tool, the calculated values can be converted into the preferred presentation method.

**MetaVi Labs cell analysis software:** To validate the data provided and calculated by Gen5, a second independent analysis tool was used. Images captured over the total incubation period were uploaded to their website. Pixel resolution values were also required by the software for calculation. After the images were analyzed, a detailed report was provided by the software including graphs showing different wound healing determinations.

### Methods

**Cell preparation:** HT-1080 fibrosarcoma cells, immortalized keratinocytes, and the neonatal dermal fibroblasts were cultured in a T75-Flask until they reached 80% confluency. Accutase was used to detach the cells from the flask. Each cell type was resuspended to a concentration of  $5.0 \times 10^5$  cells/mL and 70  $\mu\text{L}$  of cell suspension (35,000 cells per reservoir) were added to each chamber of the insert (Figure 1). The cells were allowed to attach to the coated bottom of the dish and to reach confluency overnight. The plastic inserts were removed with sterile forceps. After washing the cell patch with 2 mL of fresh culture medium, it was replaced with 2 mL of culture medium containing 10,000, 100, or 0 nM Cytochalasin D.



**Figure 1.** Performed experimental workflow for the wound healing assay in the ibidi Culture-Insert. The medium used in step 3 contained either 10,000, 100, or 0 nM Cytochalasin D.

**Imaging:** The dishes were placed in the  $\mu\text{-Dish}$  microscopy rack (ibidi, part number 80035) and fixed with the magnetic lid for additional stabilization. The rack was placed in the Agilent BioTek Lionheart FX, previously set to 37  $^{\circ}\text{C}$  and 5%  $\text{CO}_2$ . Imaging of the wound area was performed every 30 minutes over a 20-hour period using phase contrast and a 4x objective.

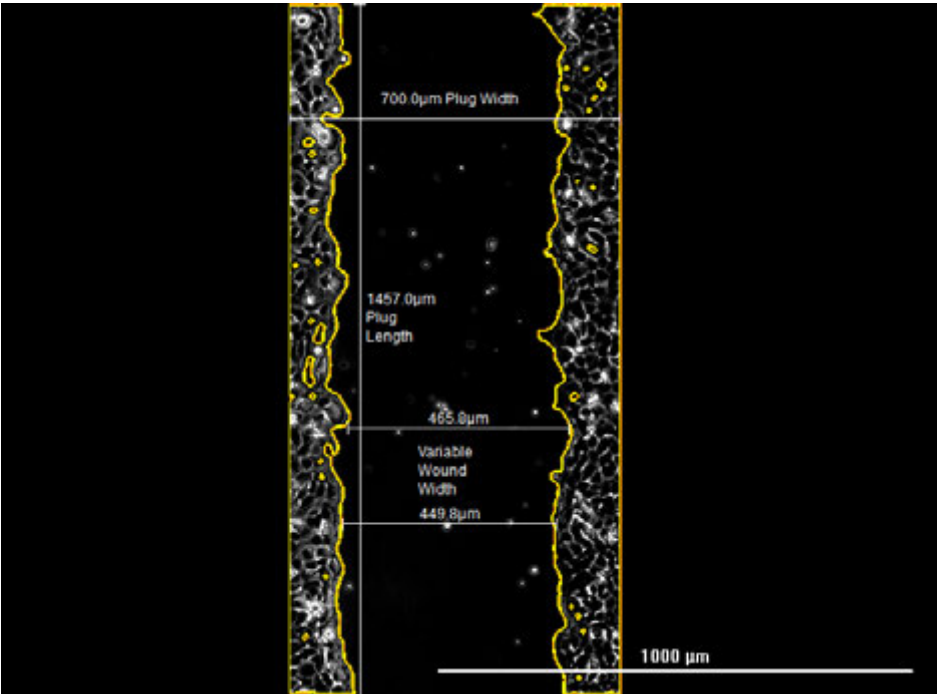
**Image-based analysis of cell movement:** Cellular analysis was performed on all 4x phase contrast images captured over the complete incubation period. Using the cellular analysis parameters in Table 1, Gen5 automatically places masks around the objects meeting the designated criteria.

**Table 1.** Agilent BioTek Gen5 cellular analysis parameters for object mask placement. Objects meeting the designated criteria will be marked with a yellow line.

Cellular Analysis Parameters	
Channel	Phase contrast
Threshold	500
Background	Dark
Split Touching Objects	Unchecked
Fill Holes in Masks	Checked
Minimum Object Size	100 µm
Maximum Object Size	10,000 µm
Include Primary Edge Objects	Checked
Analyze the Entire Image	Unchecked
Plug	Width: 700 µm Length: 1,457 µm
Advanced Options	
Evaluate Background On	1%
Image Smoothing Strength	20
Background Flattening Size	25 µm (rolling ball diameter)

In addition to the cellular analysis parameters, advanced detection options were used. Evaluation background was lowered to 1% to allow accurate identification of areas containing cells and empty areas within the image plug. Background flattening was applied to enhance the contrast between the object and background and reduce subtle changes in the signal within migrating cell areas. Finally, image smoothing was used to further smooth out variations in the phase contrast signal within cellular areas to accurately mask the collective cell monolayer. Instead of analyzing the entire field of view, an image plug with 700 µm width and 1,457 µm in length was applied to the entire 4x image (Figure 2). A width of 700 µm was chosen to include not only the complete wound area but also a small cell area on each side to use and compare phase contrast for cell analysis. The plug length of 1,457 µm equaled the 4x field of view in that dimension.

**Calculation of wound healing metrics:** The cellular analysis parameters listed in Table 1 were used to define a primary mask that quantified the area in the image plug that contained cells. This mask is depicted in Figure 2. Gen5 Calculated Metric of (Sum Area)<sub>t</sub> obtained from the mask can be used to quantify progression of cell migration as a function of time in the kinetic assay using a series of data reduction steps outlined below.



**Figure 2.** Used metrics for determination of wound healing results. Width of the plug is 700 µm, the length is 1,457 µm. Two representative lines were drawn to show the variable wound width across the entire wound length. The yellow line shows the objects masked placed by Agilent BioTek Gen5 using the parameters defined in Table 1.

**Average wound width as function of time:** While the Culture-Inserts provide a simple method to create a cell free gap, an intra and interexperiment variability in its width remains. This can easily be accounted for using for Gen5 image analysis by calculating the average wound width per time point.

**Formula 1.** Average wound width as function of time.

$$W_t = \frac{P_A - (\text{Sum Area})_t}{P_L}$$

Where  $W_t$  is the average wound width ( $\mu\text{m}$ ) as a function of time;  $P_A$  is the image plug area ( $\mu\text{m}^2$ );  $P_L$  is the image plug length ( $\mu\text{m}$ ), and the  $(\text{Sum Area})_t$  is the Calculated Metric from Gen5 ( $\mu\text{m}^2$ ).

**Percent confluency within the wound area:** The average wound width at the first kinetic time point was then also used to calculate the % confluency within the wound area according to the following formula.

**Formula 2.** % Confluency within the wound area.

$$\% \text{ Confluency} = \frac{(\text{Sum Area})_t - P_L(P_w - W_{t=0})}{P_L W_{t=0}}$$

Where  $W_{t=0}$  is the average wound width ( $\mu\text{m}$ ) at the first time point,  $P_w$  is the image plug width ( $\mu\text{m}$ ),  $P_L$  is the image plug length ( $\mu\text{m}$ ), and the  $(\text{Sum Area})_t$  is the Calculated Metric from Gen5 ( $\mu\text{m}^2$ ).

**Percent confluency in a fixed wound width gap:** In addition, % confluency in a fixed wound area ( $500 \mu\text{m}$ ) and the maximum wound-healing rate were calculated for the HT-1080 cell line according to the following formula.

**Formula 3.** % Confluency in a fixed wound gap.

$$\% \text{ Confluency}_{\text{Fixed}} = \frac{(\text{Sum Area})_t - P_L(P_w - W_L W_w)}{W_L W_w}$$

Where  $W_L$  is the fixed wound length,  $W_w$  is the fixed wound width,  $P_w$  is the image plug width ( $\mu\text{m}$ ),  $P_L$  is the image plug length ( $\mu\text{m}$ ), and the  $(\text{Sum Area})_t$  is the Calculated Metric from Gen5 ( $\mu\text{m}^2$ ).

Finally, the Kinetic Analysis tool of Gen5 3.0 calculated the highest slope of the actual wound area curve to provide the maximum wound healing rate in  $\mu\text{m}^2$  per minute.

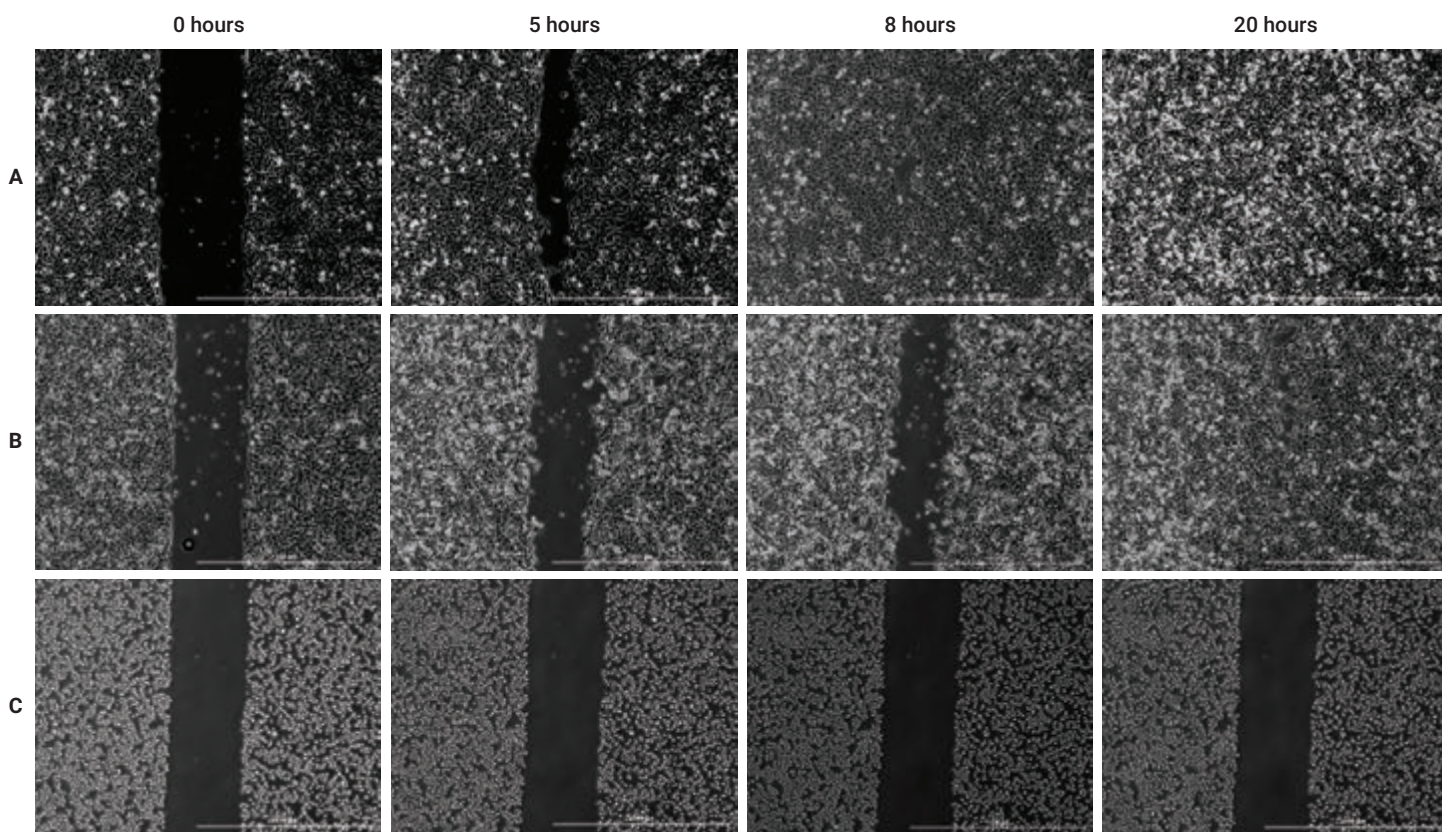


## Results and discussion

### Label-free monitoring of wound closure over time and its inhibition with Cytochalasin D

Figures 3 to 5 illustrate the ability of the Lionheart FX to accurately image the process of gap closure with three different cell types. The cells types were treated with 10,000, 100, or 0 (control) nM Cytochalasin D, and imaging was performed over a 20-hour period at 30-minute intervals.

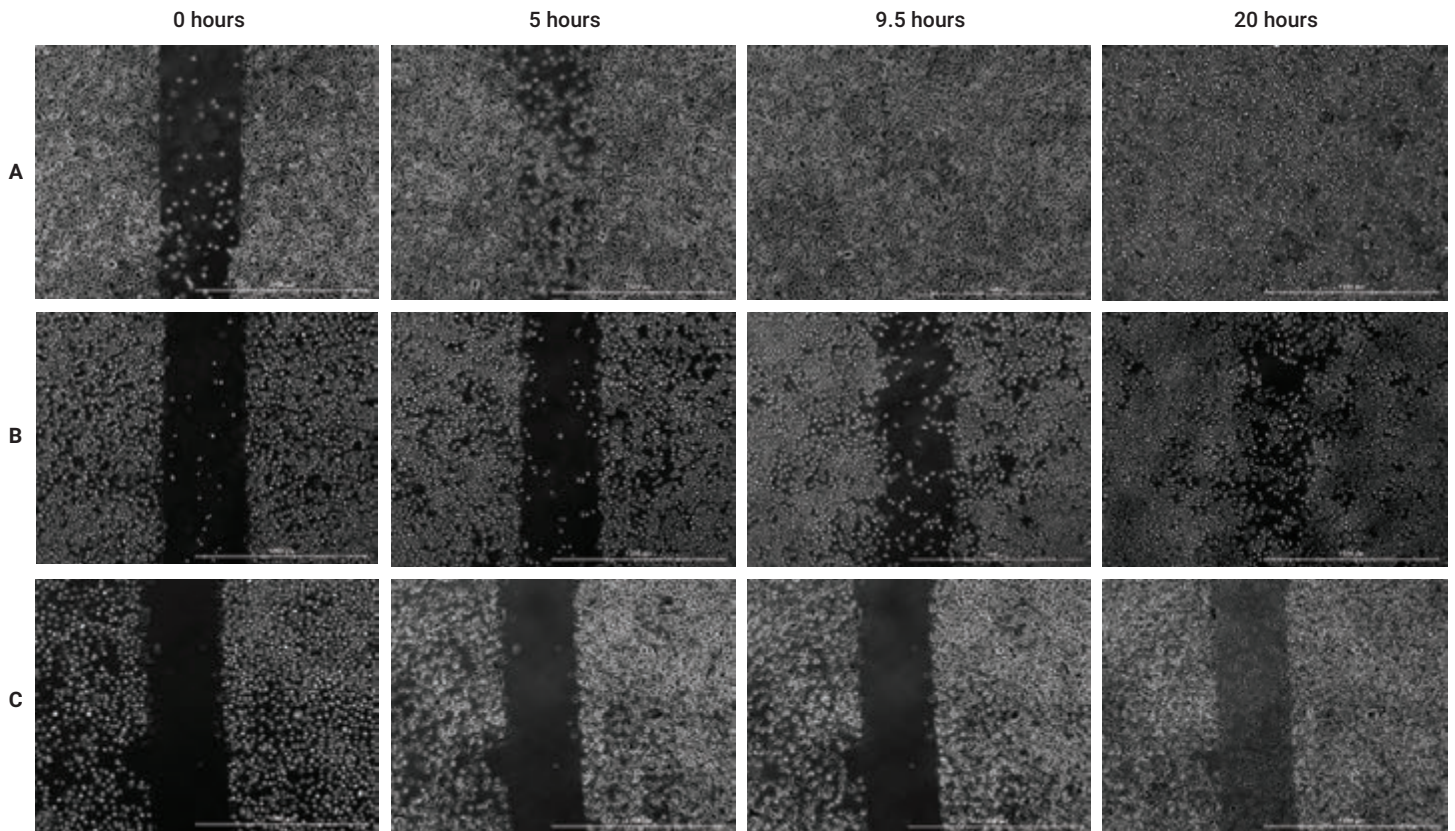
The HT-1080 cells showed a collective phenotypic movement, causing the gap to close after 8 hours (Figure 3A). While wound healing was mildly inhibited with 100 nM Cytochalasin D (Figure 3B), the higher concentration of 10,000 nM Cytochalasin D completely prevented cell migration compared to the control (Figure 3C).



**Figure 3.** Image-based monitoring of HT-1080 migration. Imaging was performed over 20 hours using phase contrast and 4x magnification. Kinetic images shown for cells treated with (A) 0, (B) 100, or (C) 10,000 nM Cytochalasin D.

Keratinocytes, in comparison, started migrating as single cells and ended up closing the wound after 9.5 hours (Figure 4A). Compared to the control, the lower concentration of Cytochalasin D had an impact on total gap closure and even

after 20 hours, the artificially created wound was not closed (Figure 4B). Keratinocytes treated with 10,000 nM compound showed no cell movement compared to the control, and therefore no migration (Figure 4C).



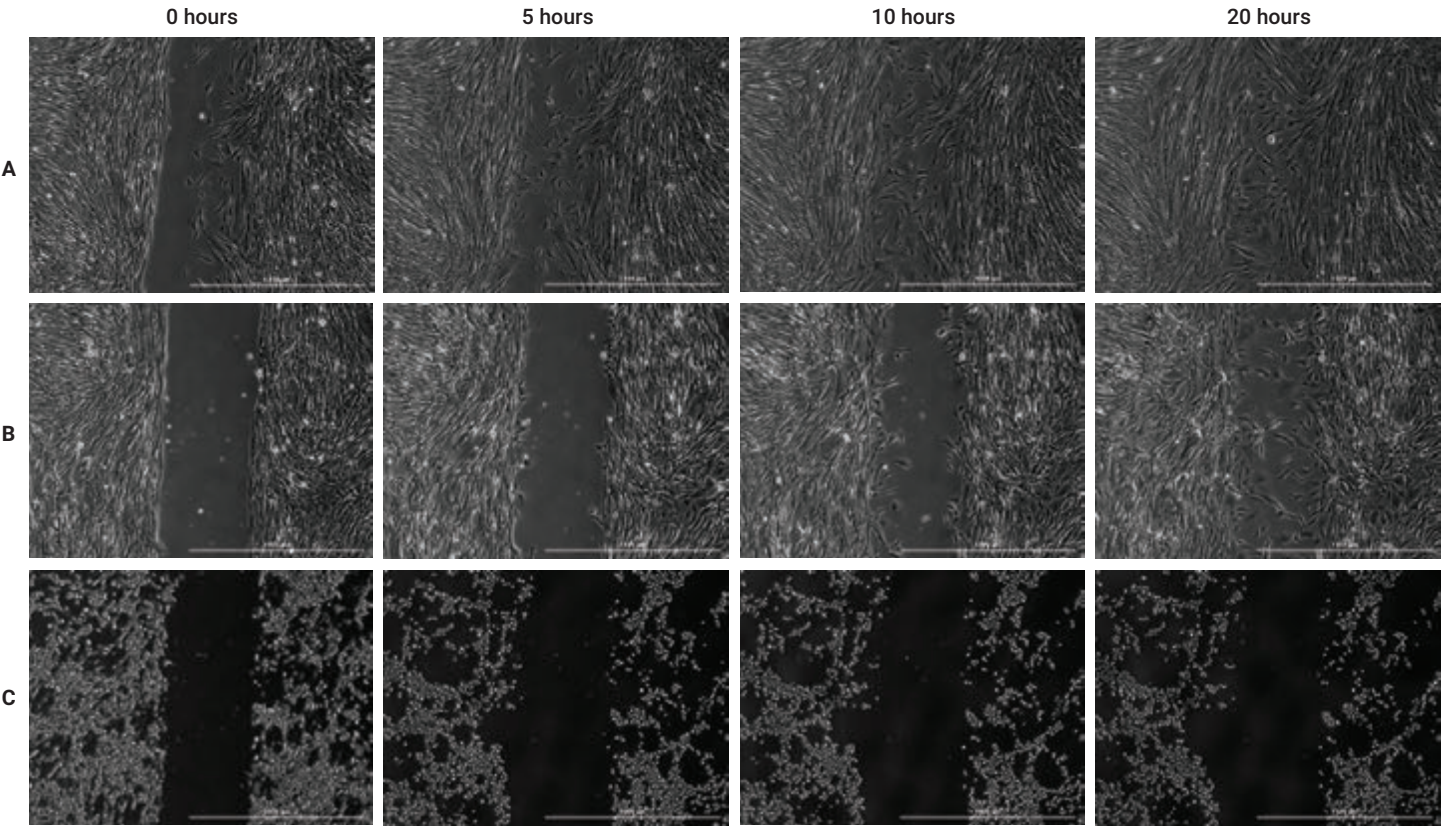
**Figure 4.** Visual confirmation of keratinocyte wound healing. Imaging was performed over 20 hours using phase contrast and 4x magnification. Kinetic images shown for cells treated with (A) 0, (B) 100, or (C) 10,000 nM Cytochalasin D.



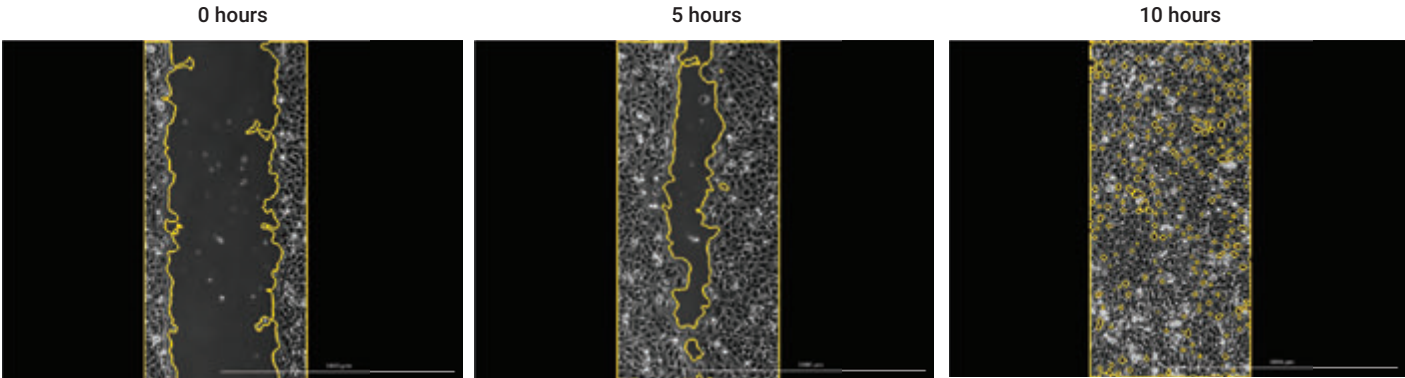
Finally, it was also possible to monitor the larger, filamentous fibroblasts moving across the gap (Figure 5A). Within the 20-hour incubation period, the fibroblasts were not able to reach confluency within the wound area. Adding 100 nM Cytochalasin D to the medium slowed down wound healing compared to the control (Figure 5B). The highest concentration of Cytochalasin D had the greatest observed impact of all tested cell models. As illustrated in Figure 5C, not only cell migration but also cell adherence is affected, indicating a potential necrotic effect on the fibroblasts.

**Quantification of wound healing and its inhibition using image-based cellular analysis**

As previously described, the cellular analysis parameters used (Table 1) were able to accurately quantify the area occupied by cells in the plug image using a primary mask (Sum Area)<sub>t</sub>. This is depicted in Figure 6 over three time points for the HT-1080 cell model. Then, using Formulas 1 and 2, the average wound width,  $W_t$  and % Confluency can be determined in the data reduction tool in Gen5.



**Figure 5.** Image-based monitoring of fibroblasts migration. Imaging was performed over 20 hours using phase contrast and 4x magnification. Kinetic images shown for cells treated with (A) 0, (B) 100, or (C) 10,000 nM Cytochalasin D.

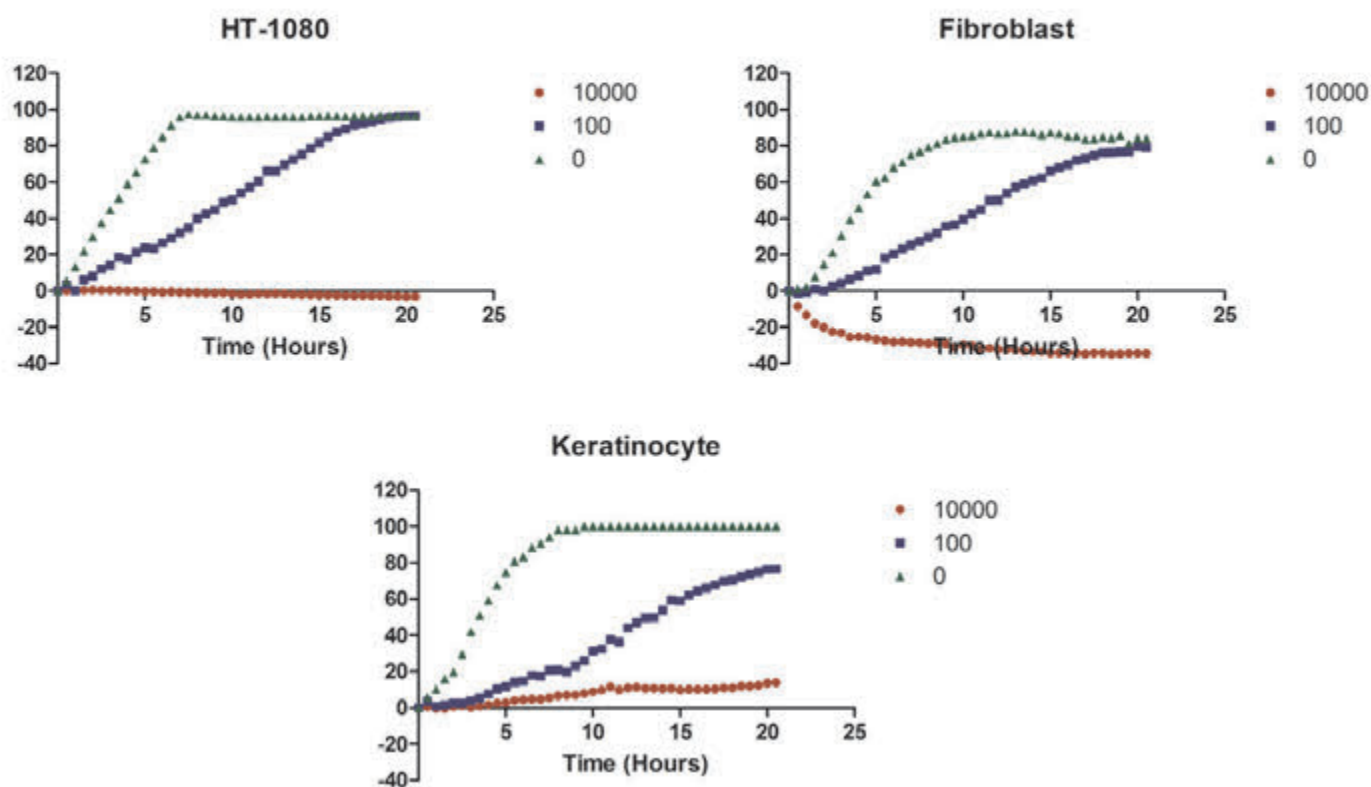


**Figure 6.** Representative images at different time points showing the applied object mask of the cellular analysis tool provided by Gen5 3.0.

Figure 7 shows kinetic plots spanning a 21-hour period for the three cell models using the three separate Cytochalasin D concentrations. Cell migration is somewhat different between the three cell lines, although in each case, Cytochalasin D demonstrates the ability to inhibit cell migration in a dose-dependent manner, albeit to different extents. The cell-free gap using the HT-1080 cells was closed after

7 to 10 hours, while for the keratinocytes it was delayed somewhat to 8 to 10 hours. The fibroblasts, by comparison, only reach 80% confluency after approximately 10 hours of incubation and demonstrated a loss of adherence at the highest inhibitor concentration, noted by negative % confluency.

## % Confluency in the Wound Area



**Figure 7.** Quantification of % confluency in the wound area to show cell migration. Formula 2 was used to calculate the values.



To better explain observed phenotypic results, Gen5 also provides user-friendly methods to monitor wound healing using alternative metrics. These are demonstrated using results from the HT-1080 cell model in Figure 8. Calculations with a fixed wound area, as shown in the upper right graph (Figure 8), can, for example, help explain variances seen amongst replicates. This is important especially when cells start to migrate before initial imaging takes place (Figures 4A and 5A). To show actual wound closure, plotting the average wound width over time (lower left graph in Figure 8) provides information about cell movement itself. In the absence of Cytochalasin D, HT-1080 cells move together, narrowing the

gap to a value approaching 0. In comparison, HT-1080 cells treated with the highest concentration of inhibitor showed a constant average wound width. Finally, the kinetic analysis tool included in Gen5 evaluates each curve to determine the maximum wound healing rate. As shown in the lower right graph of Figure 8, Cytochalasin D slows wound healing until the value reaches 0 with the highest concentration tested. Together, these metrics, which are automatically determined by Gen5, provide high information content and give a more complete understanding of wound healing and how the process can be altered by adding an inhibiting compound.

### Wound Healing Analysis Methods

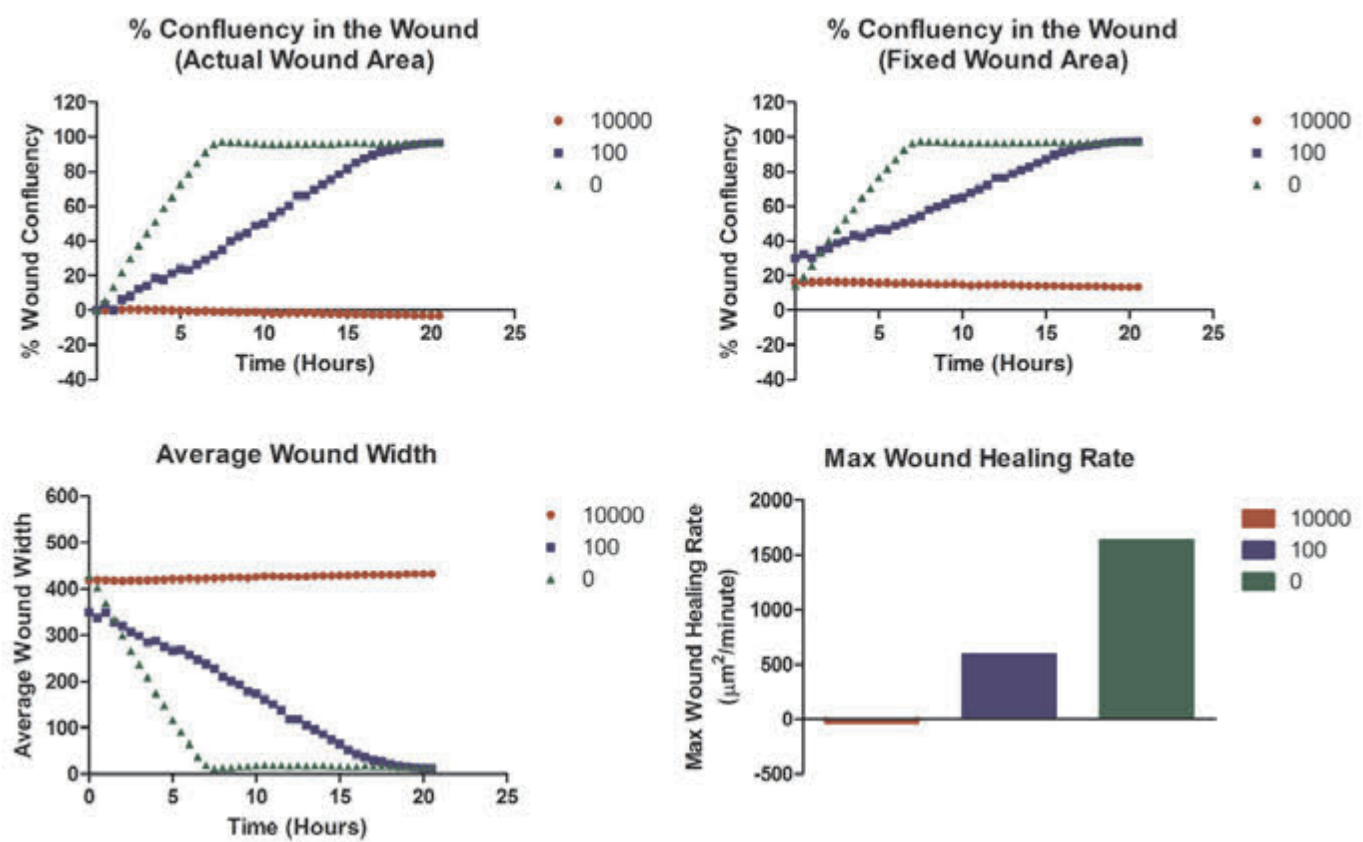


Figure 8. Variable determinations of wound healing and cell migration. Formulas 1, 2, and 3 were used to calculate the values.

To validate the analysis performed with Gen5, a second independent software provided by MetaVi Labs was also used. Similar analysis results were shown in both analysis software.

The graph showing the Percent Gap Closed All Positions in Figure 9 is comparable with the graph showing the % Confluency in the fixed Wound Area presented in Figure 8.

The Scratch Open Area All Positions can be compared to the Average Wound Width in Figure 8. While the curve shapes agree, the values represent different metrics. Instead of calculating the wound width in  $\mu\text{m}$ , the wound area in  $\mu\text{m}^2$  was calculated. Both metrics show the same progress over time. Finally, the software also calculates the Gap Closure Speed (Figure 9), which agrees with Gen5 software kinetic analysis of the Max Wound Healing Rate (Figure 8).

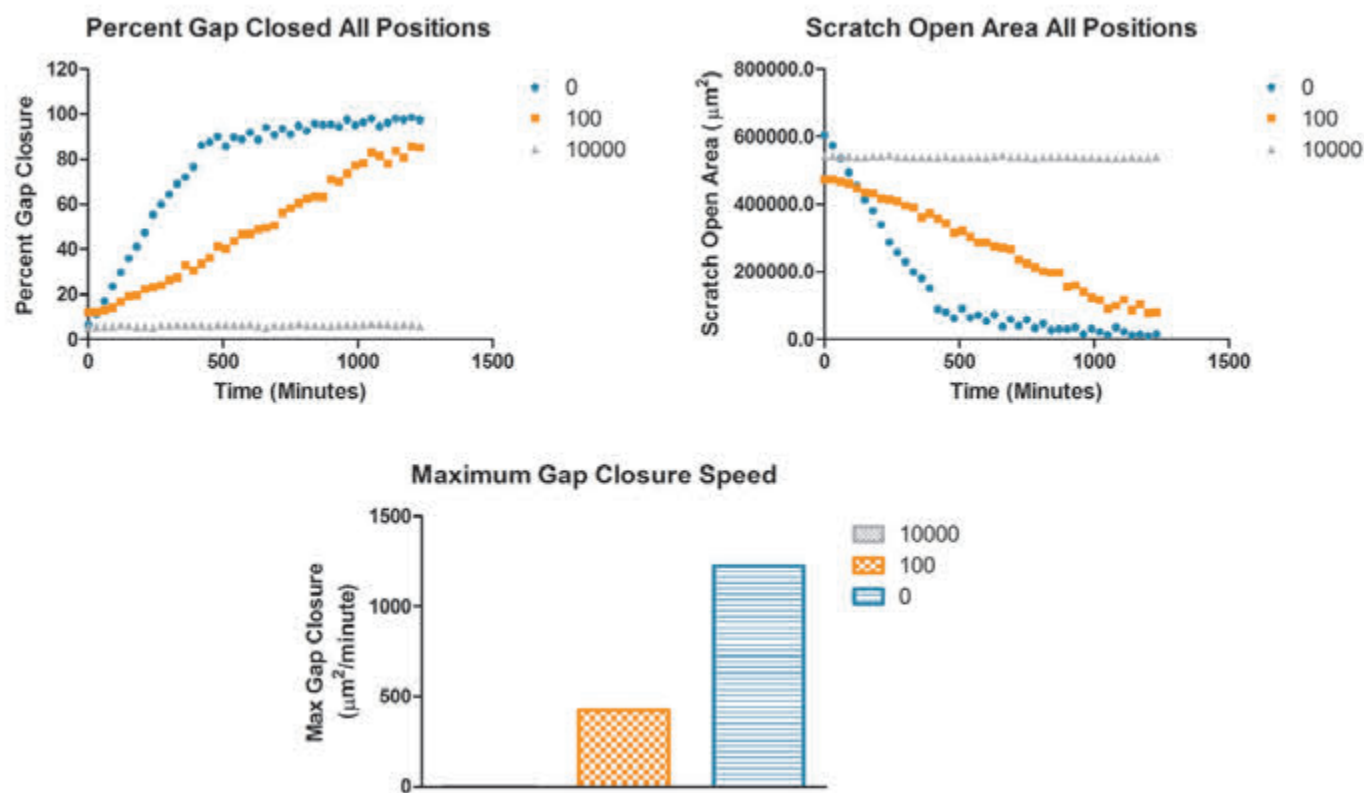


Figure 9. Determination of wound healing using an independent cell analysis software.

## Conclusion

The wound healing assay is a powerful method to investigate cell migration. The culture insert and the 35 mm  $\mu$ -Dish provided by ibidi simplifies the performance and allows easy compound addition. The results seen with the combined experimental procedure delivered results that agree with what has been shown previously in the literature. The Agilent Biotek Lionheart FX automated microscope is perfectly suitable to perform long-term kinetic imaging of the wound healing process. The Agilent Biotek Cytation 5 cell imaging multimode reader is also capable of performing the imaging needed for the experimental setup. Proper cell health can be ensured through the incorporation of appropriate environmental settings over the incubation time. Integrated Augmented Microscopy combines accurate time-lapse imaging with precise data analysis. Percent confluency was calculated for each cell type, while various determinations of gap closure can also be automatically calculated using the Agilent Biotek Gen5 microplate reader and imager software. The results were verified by using an independent second analyzing software provided by MetaVi Labs. The combination of an easy experimental setup, imaging, and precise data analysis creates a powerful tool to investigate cell migration.

## References

1. <http://www.nature.com/subjects/cell-migration>
2. Ilina, O.; Friedl, P. Mechanisms of Cell Migration at a Glance. *Journal of Cell Science* **2009**, *122*, 3203–3208.
3. Henson, J. H. *et al.* Wound Closure in the Lamellipodia of Single Cells: Mediation by Actin Polymerization in the Absence of an Actomyosin Purse String. *Molecular Biology of the Cell* **2002**, *13*, 1001–1014.

## Acknowledgements

The authors would like to thank Michael Wyler, Joel Mailliet, and Gary Prescott for their contributions to the final analysis methods included in the application note.

[www.agilent.com/lifesciences/biotek](http://www.agilent.com/lifesciences/biotek)

For Research Use Only. Not for use in diagnostic procedures.

RA44200.500787037

This information is subject to change without notice.

© Agilent Technologies, Inc. 2016, 2021  
Printed in the USA, February 1, 2021  
5994-2582EN  
AN062716\_12