

Kinetic Live-Cell Visualization and Quantification of Cell Migration

A. Ward*, J. Trigg, K. Barnes, G. Lovell, T. Dale

Sartorius UK Ltd., Royston, Hertfordshire, UK
*Corresponding author: Amber.Ward@Sartorius.com

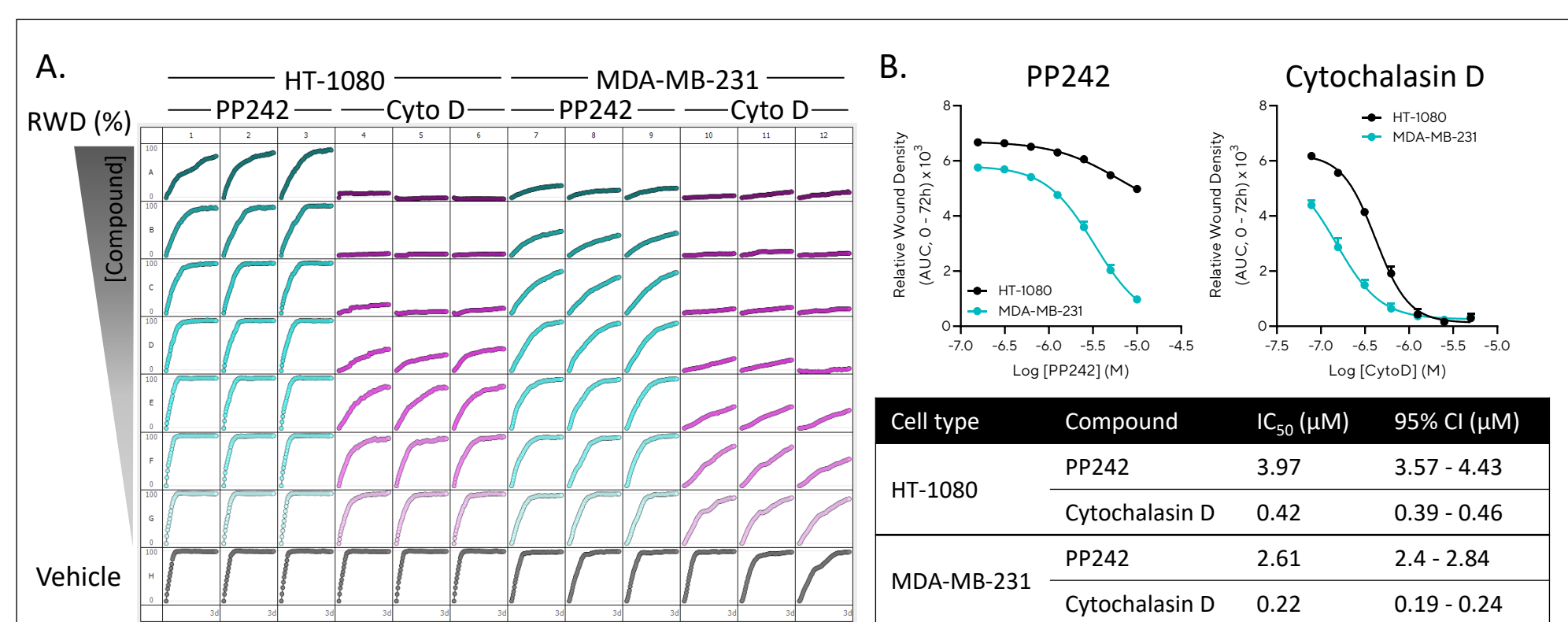
Summary & Impact

- Cell migration is a multistep process that enables cell movement in response to an environmental stimulus, this plays a role in physiological and pathological processes, including wound healing and tumor metastasis.
- The rapid rearrangement of actin filaments leads to a cycle of leading edge protrusion and lagging edge retraction, summing in the associated cellular morphological changes.
- Here we exemplify an assay for the visualization and kinetic assessment of cell migration and its modulation in a 96-well plate. Incucyte® Live-Cell Analysis enables automated image-

based measurements of cell migration *in vitro* via label-free or dual fluorescence readouts.

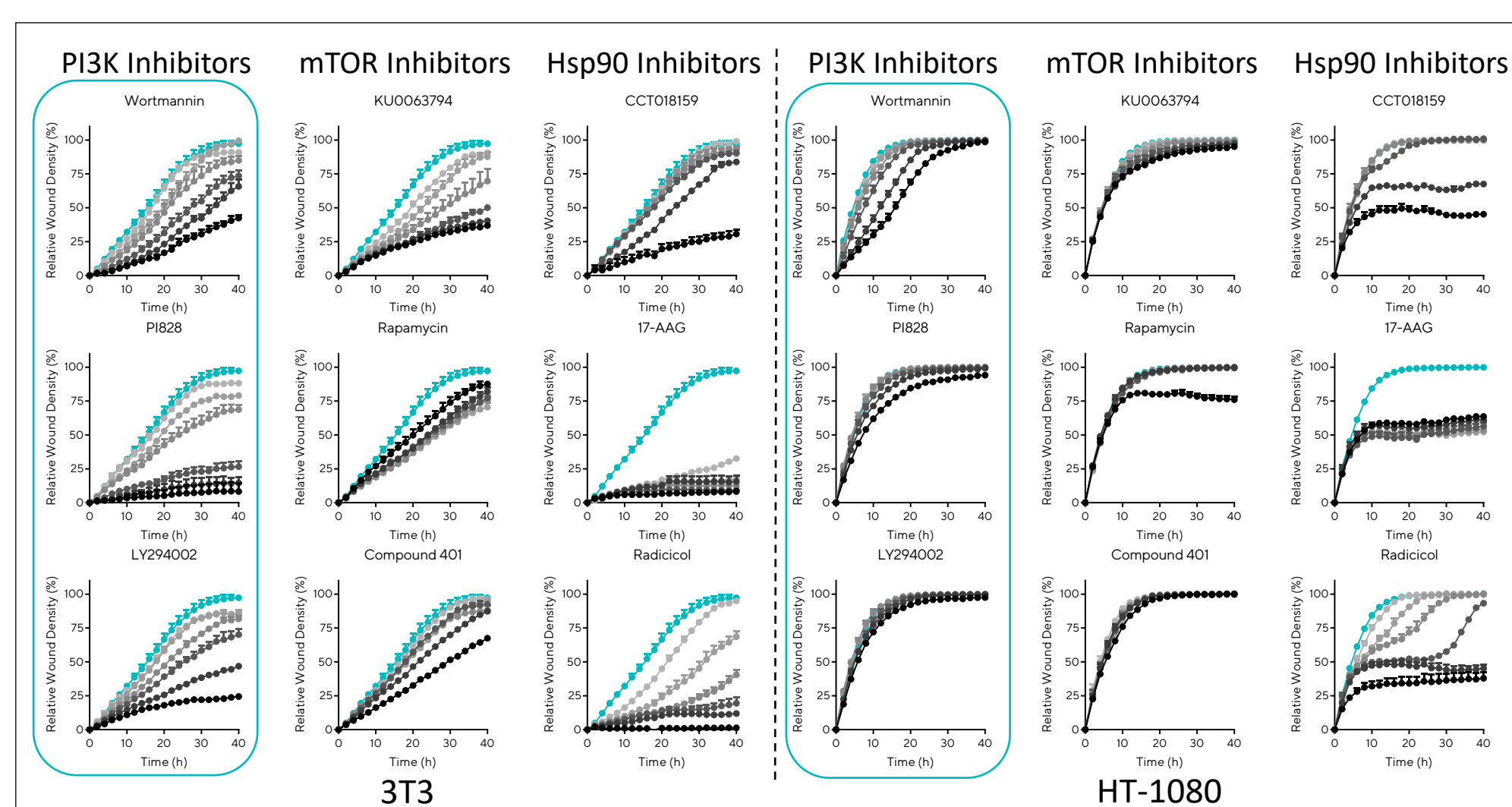
- To simultaneously monitor cell cycle dynamics, cells expressing the Incucyte® Cell Cycle Lentivirus Reagent were used. Images were acquired using the Incucyte® Live-Cell Analysis System and analyzed using integrated software.
- These data demonstrate that live-cell analysis methods provide a solution to robustly measure cell migration in real time without relying on end point analysis.

Pharmacological assessment of inhibitors of cell migration



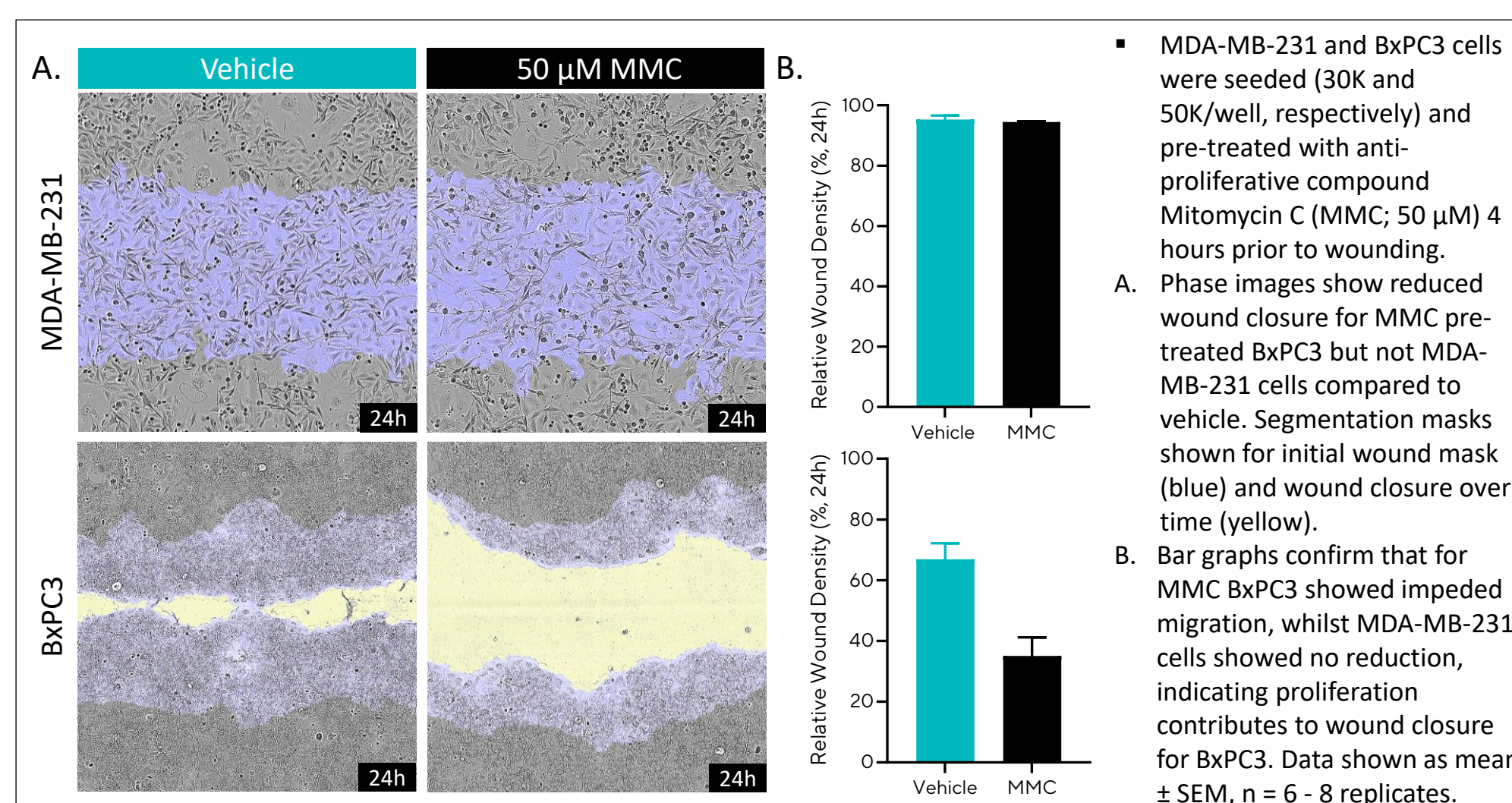
- HT-1080 and MDA-MB-231 cells were seeded at 30K/well and left for 24 hours to form confluent monolayers. Cells were wounded using the Woundmaker Tool and treated with compounds.

Cell- and molecule-dependent cell migration profiles

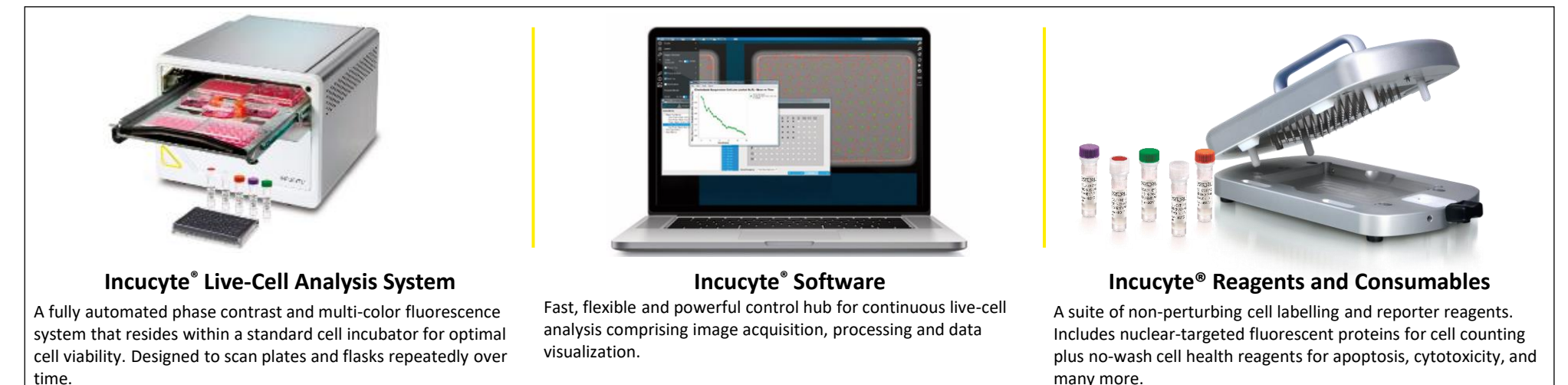


- 3T3 and HT-1080 cells (30K/well) were wounded and treated with a concentration-range (0.1 – 30 µM) of 9 inhibitors of PI3K, mTOR (both associated with Akt pathway), and Hsp90 (chaperone protein for stability).
- PI3K inhibitors showed a significant difference in efficacy with HT-1080 appearing more resistant than 3T3 cells. HT-1080 have an activated n-RAS oncogene (upstream activator of PI3K-Akt pathway) which via sustained signaling, may contribute to observed resistance. Data shown as mean ± SEM, n = 3 replicates.

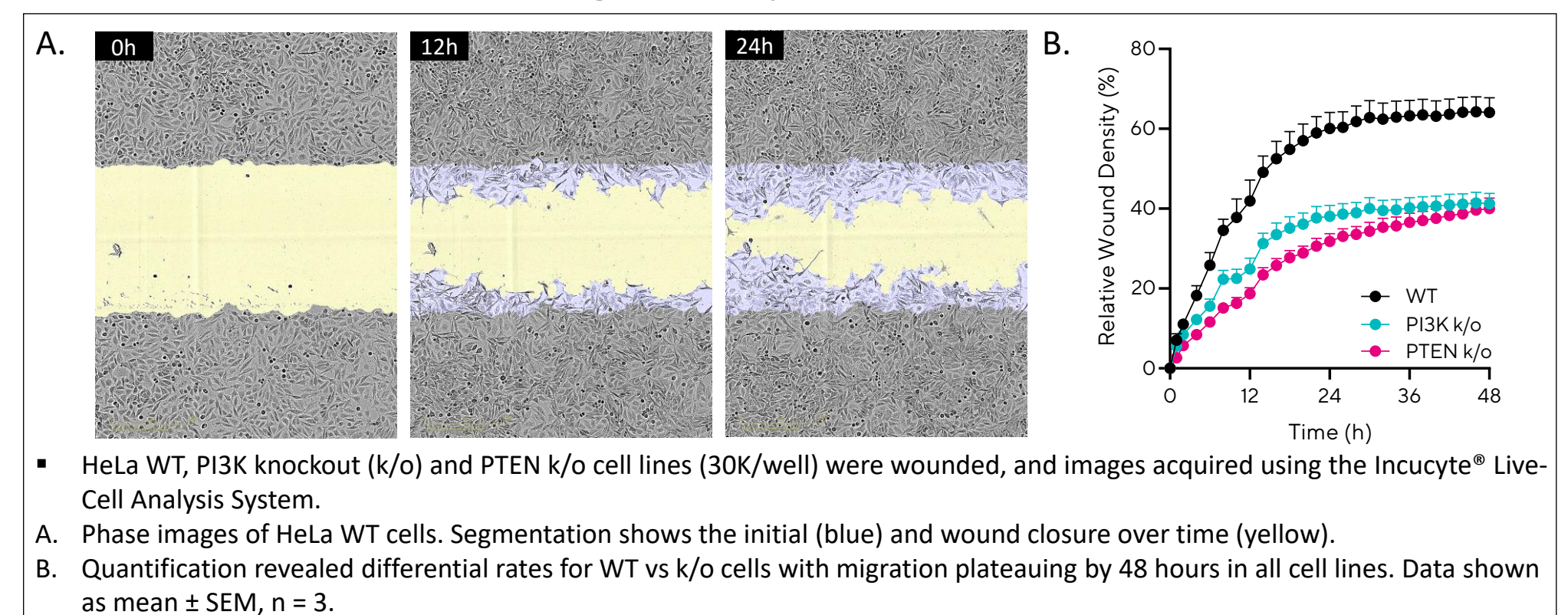
Anti-proliferative effects of mitomycin C during cell migration



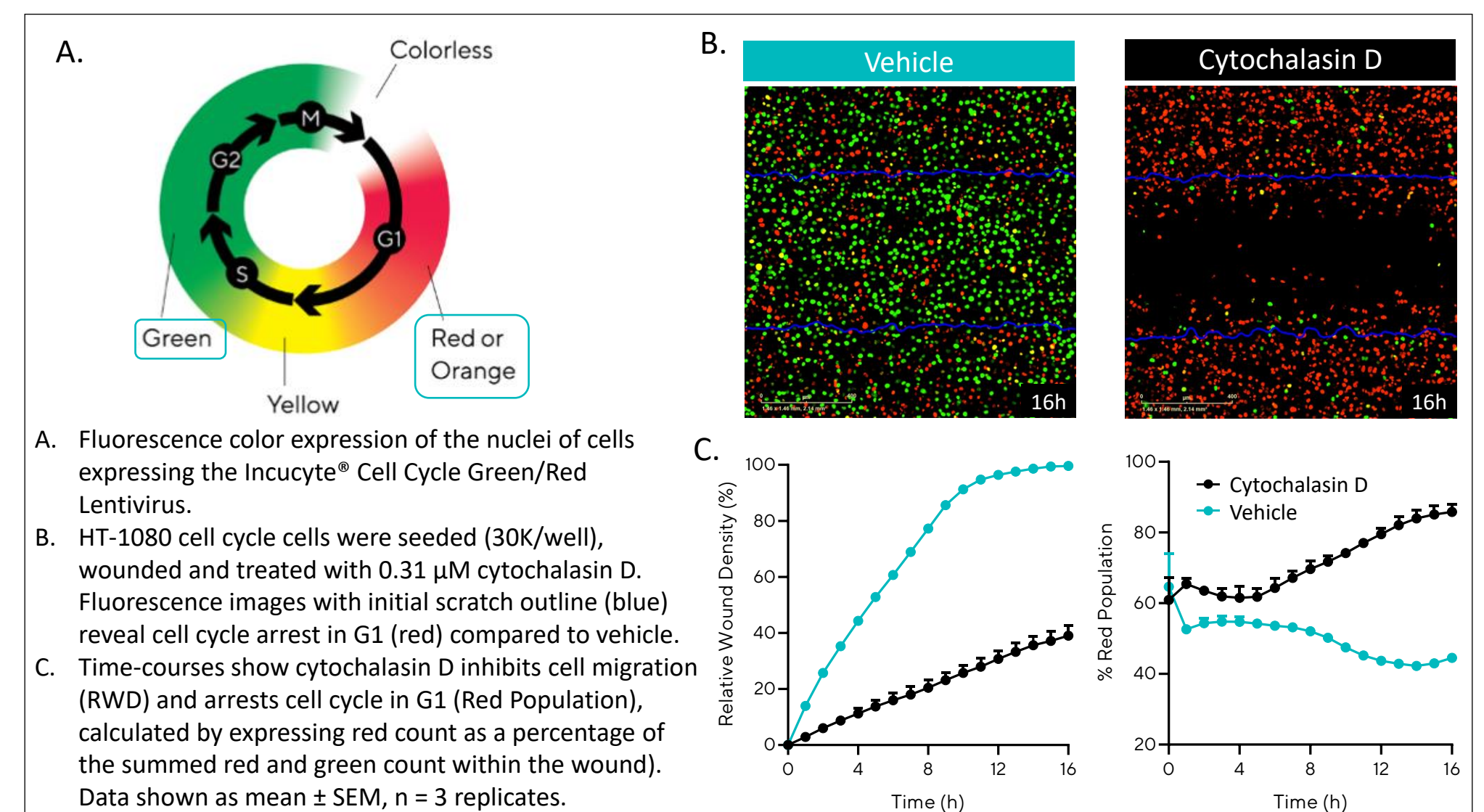
Incucyte® Live-Cell Imaging and Analysis Solutions



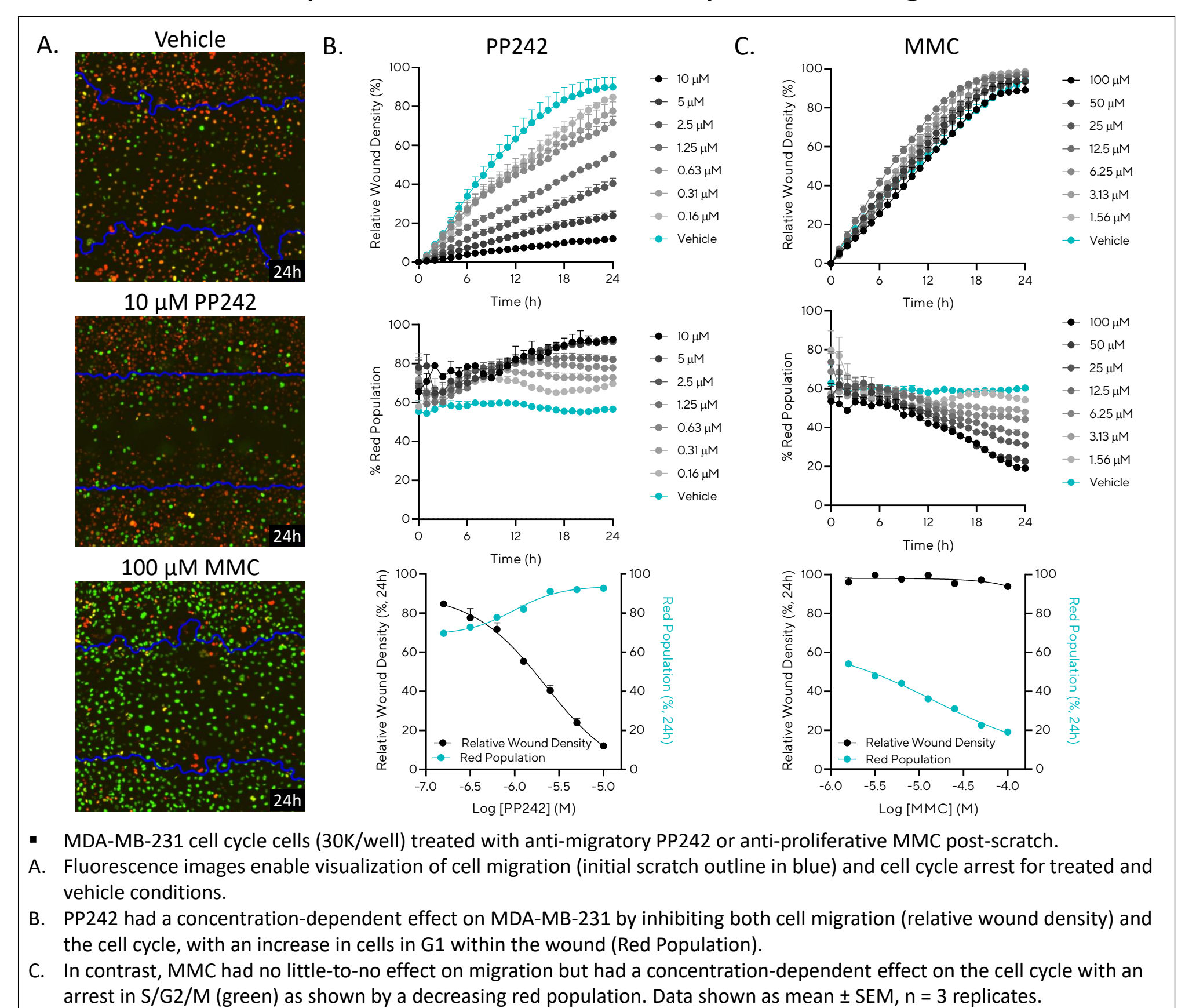
Quantification of cell migration profiles in real time



Cytochalasin D inhibits cell cycle and migration



Differential compound effects on cell cycle and migration



- MDA-MB-231 cell cycle cells (30K/well) treated with anti-migratory PP242 or anti-proliferative MMC post-scratch.
- Fluorescence images enable visualization of cell migration (initial scratch outline in blue) and cell cycle arrest for treated and vehicle conditions.
- PP242 had a concentration-dependent effect on MDA-MB-231 by inhibiting both cell migration (relative wound density) and the cell cycle, with an increase in cells in G1 within the wound (Red Population).
- In contrast, MMC had no little-to-no effect on migration but had a concentration-dependent effect on the cell cycle with an arrest in S/G2/M (green) as shown by a decreasing red population. Data shown as mean ± SEM, n = 3 replicates.