

CloneSelect Imager

Objective quantification of cell migration

- Rapid scanning to monitor cell migration
- High quality imaging and label-free detection of living cells using white light

Background

Cell migration is a key process in normal development, immune function and wound healing. Abnormal cell migration is a feature of conditions such as cardiovascular disease and cancer. The CloneSelect Imager system reveals the *in vitro* effects of therapeutic candidates and cell culture matrices on cell migration as well as the study of cell-to-cell interactions. Migration application software in CloneSelect Imager enables measurement of cell migration based on objective cell detection.

Method

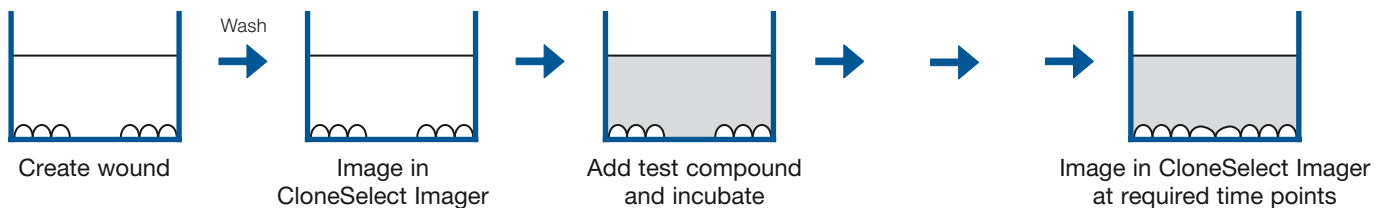
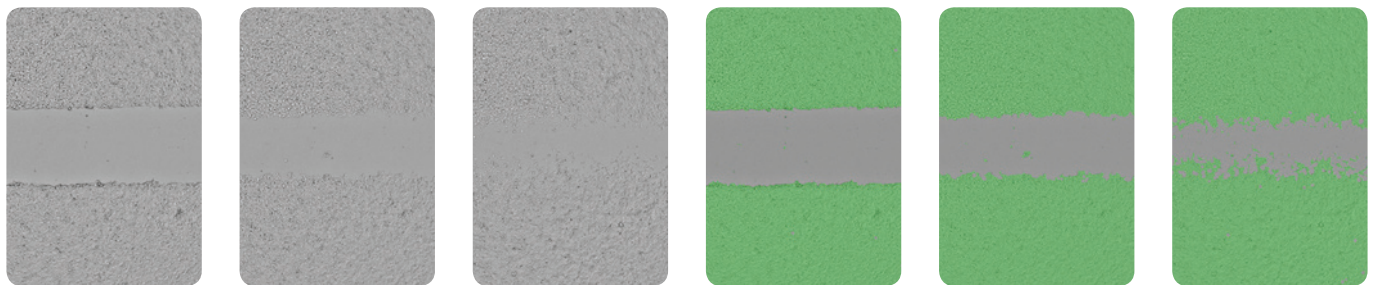


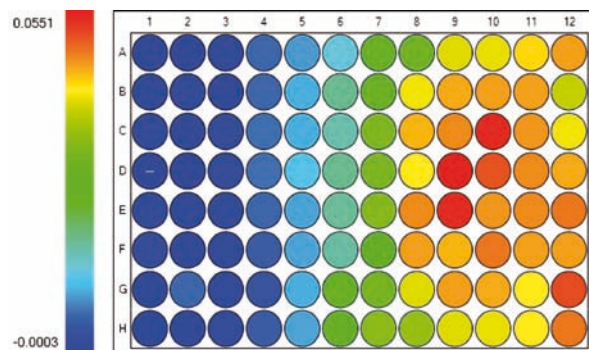
Image capture and analysis



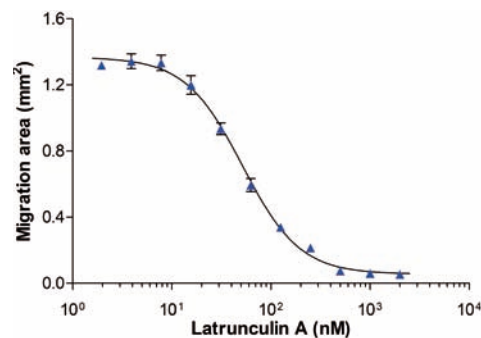
Label-free imaging of cells over time

Detection of cell area by migration application software

High quality data generation



Heat-map representation of cell migration across plate



Dose-response curve generated from exported image analysis data

CloneSelect Imager – objective, quantitative assessment of cell growth

The CloneSelect™ Imager system utilizes non-invasive, white light imaging to enable rapid, quantitative measurement of cell confluence and generation of growth curves well by well.

Fast results

- Replace time-consuming manual inspection
- Make confident, image-driven decisions
- Obtain consistent results in under three minutes per 96-well plate

For applications including:

- Rapid confluence measurement
- Growth curve generation
- Cell number estimation
- Verification of monoclonality
- Colony forming assays
- Cell migration assays

Key parameters

- Accepts multi-plate formats from single well to 384 well
- 4x objective lens
- Resolution to 1.8 μm
- Integrated barcode reader

Objective analysis

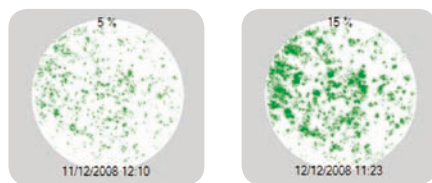


Productivity easily increased

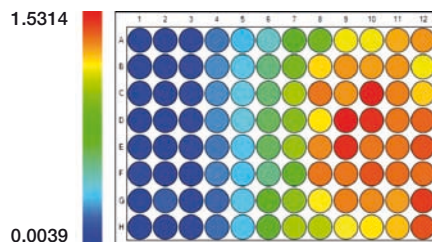


Compatible with a range of robotic systems

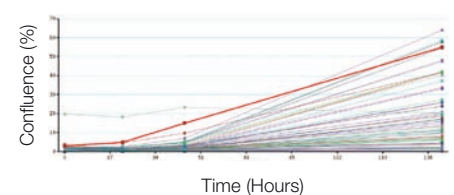
Rapid image generation and clear data analysis



Simple overlay reveals cell confluence



Heat maps facilitate analysis



Growth rates for every well viewed at every time point

ClonePix, CloneSelect, CellReporter are trademarks of Molecular Devices (New Milton) Ltd. All third party trademarks are the property of their respective owners. For a listing of trademark owners, visit www.moleculardevices.com/genetix

Application Highlight



www.moleculardevices.com/genetix