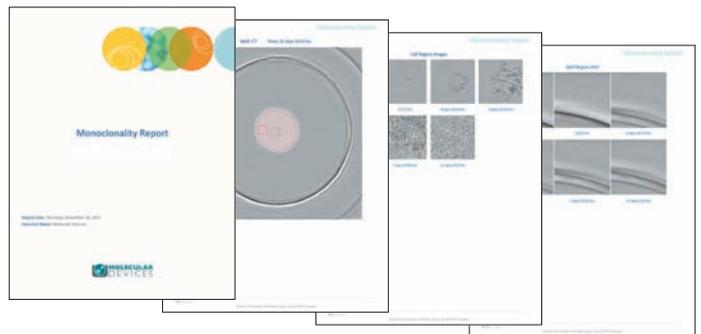




CloneSelect Imager Monoclonality Report feature

Providing image evidence of monoclonality in the cell line development process is not as simple as exporting an image of a single cell. For example, high-resolution images of the entire well should also be inspected to ensure the absence of a second cell.

With a few simple clicks, the Monoclonality Report feature on the CloneSelect® Imager (CSI) objectively organizes the supporting image evidence needed to establish clonality into an easily shareable report, saving researchers hours typically required to do the same process manually.



Example pages from the monoclonality report.

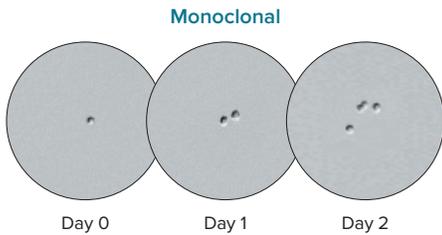
Key features

- Easily identify and select single-cell and artifact regions to include in report
- Export high-resolution images of single cells, artifacts, and entire well (optional)
- Automatically identify single cells with fluorescence version of CloneSelect Imager
- Publish monoclonality report in PDF or Word Format

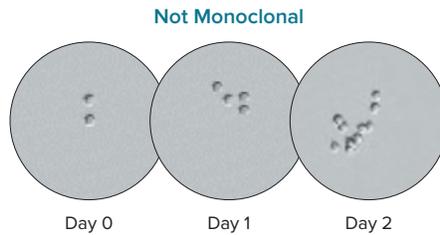
The reporting feature automatically generates the following data in a presentable format:

Monitor cell line development over time

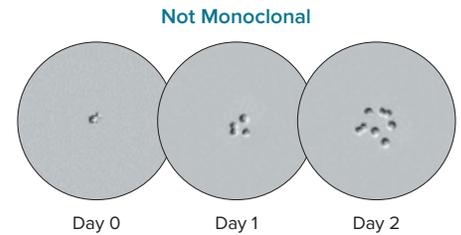
To characterize the growth from a single cell to a colony, cell regions can be designated and adjusted for each time point in a series.



A single cell on Day 0 and two cells on Day 1 confirm monoclonality.



Two cells are visible on Day 0, confirming that the cell line is not monoclonal.



Four cells have developed on Day 1, confirming that the cell line is not monoclonal.

View entire well at final time point

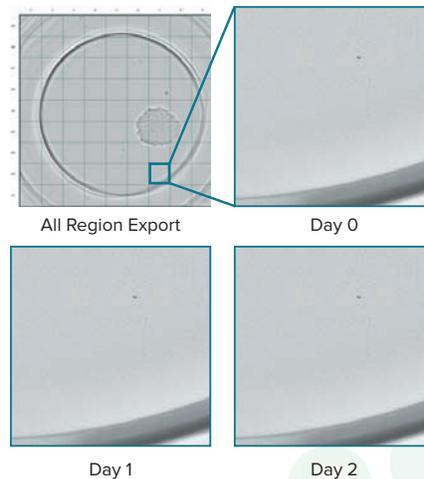
Quickly determine clonality of a cell line by visually inspecting the presence of multiple colonies in a single well.



Wells captured on Day 10. The well on the left shows one colony while the well on the right shows two colonies.

Divide an entire well into individual images

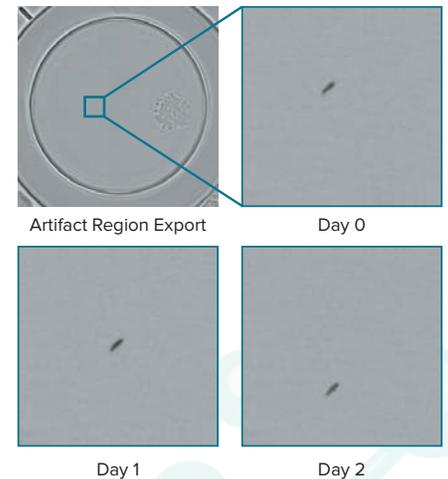
Export an entire well into 81 separate images to objectively confirm the absence of another cell.



An image of a well is divided into regions, and a selected region is displayed over time.

Highlight regions to review non-cell objects

Selectively highlight parts of a well to differentiate cells from ambiguous objects.



A selected artifact region and its corresponding images over time is shown here.

Additional data included in report

- Report Date
- Well Name
- Date Imaged
- Region Locations
- Operator Name
- % Single Colony Wells
- Barcodes
- Experimental Notes

Contact Us

Phone: +1.800.635.5577
Web: www.moleculardevices.com
Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+91.73.8661.1198

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

