



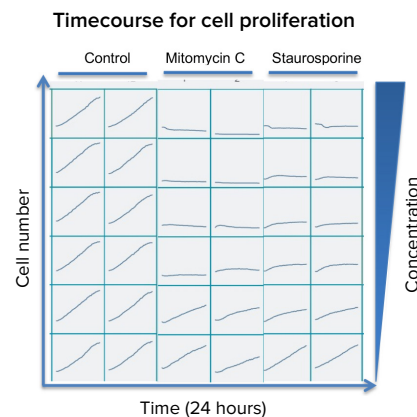
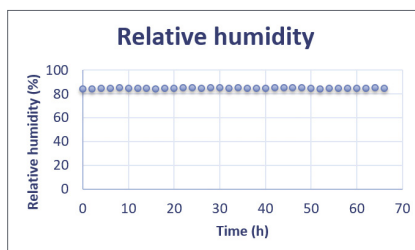
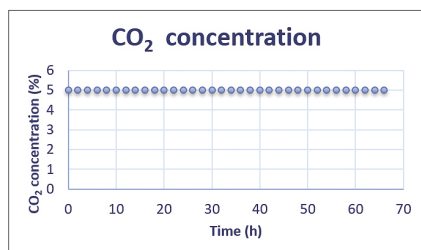
ImageXpress Pico system with environmental control

Perform extended time-lapse acquisitions on living samples while maintaining optimal conditions

Enable live cell imaging and real-time environmental control

Environmental control mimics the cell environment and enables you to run multi-day studies, time-lapse, and live cell assays. With CellReporterXpress® software, the ImageXpress® Pico Automated Cell Imaging System provides visibility for environmental control during acquisition to ensure that the system is running at peak performance during your assay.

Reproduce and maintain environmental conditions

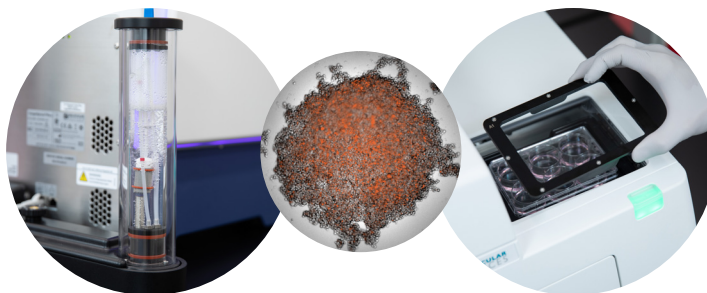


Monitor and measure changes in cell proliferation and cell phenotype over time with reliable time-lapse imaging and on-the-fly image analysis.

Key features

- Enable live cell imaging and 3D acquisition in an easy-to-use system to better identify phenotypic changes
- Generate sharp images with a greater depth of field
- Minimize evaporation with humidity controls
- Maintain healthy cell growth environment with user-defined CO₂ settings
- Store the results of control parametrics with data for review even after the experiment ends
- Study the effects of hypoxic conditions over time
- Precise temperature regulation to mitigate thermal Z-drift throughout the entirety of the time-lapse acquisition
- Ensure optimal environmental conditions at the plate level without the need to flood the entire system

The external **humidity chamber** integrates via the rear of the ImageXpress Pico system to maintain a small surface area for small lab spaces.

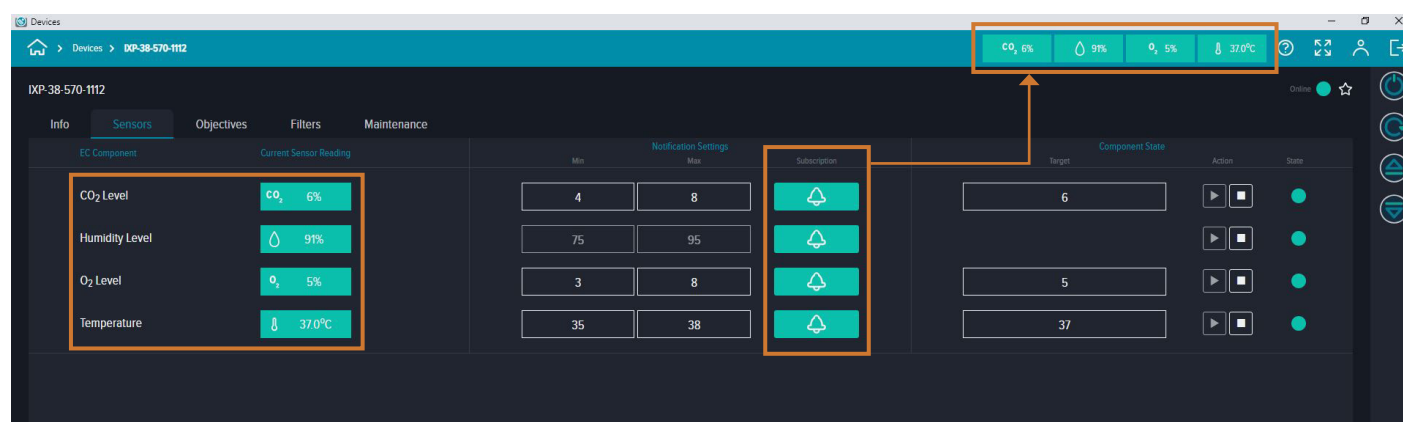


Microplates are placed in an **environmental control cassette** which introduces and maintains gas and humidity while reducing gas usage as compared to flooding the entire system. Cassettes can also be sterilized through UV.

Environmental control settings

Settings can be configured through CellReporterXpress software and monitored during acquisition

- Monitor readings in real time
- Acquired sensor data is linked to the experiment
- Subscribe to alerts to watch multiple sensors and get notifications while working within CellReporterXpress software



Environmental control specifications

Temperature control	Ambient +8°C to 40°C
Temperature control homogeneity	37°C ± 0.5°C at 23°C ambient
Gas control	O ₂ control, 1–15% and ambient, CO ₂ control, ambient to 15%
Humidity control	Active humidity control. Sample compartment controlled to 85% nominal humidity.

For more information, please visit moleculardevices.com/pico

The ImageXpress Pico system features optics by Leica Microsystems.

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
 United Kingdom +44.118.944.8000
 Europe* 00800.665.32860
 China +86.4008203586
 Taiwan/Hong Kong +886.2.2656.7585
 Japan +81.3.6362.9109
 South Korea +82.2.3471.9531
 India +91.73.8661.1198
 *Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland