

# CLIMATE CHANGE IN YOUR LAB CAN CAUSE POOR INSTRUMENT PERFORMANCE

Keeping your lab environment at the right climate is crucial to get the best out of your instruments and your results.

## Signs It's Stormy in Your Lab



### Temperature Changes

(18 °C to 25 °C) Lab temperatures outside recommended requirements affect instrument performance. Peaks can jump around with temperature swings.



### Dry Conditions

(< 20% RH) Instruments are susceptible to electrostatic discharge resulting in damage to electronic equipment.

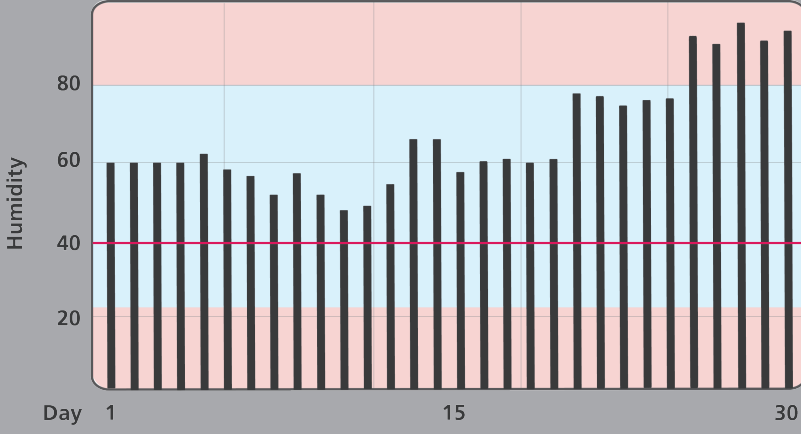


### High Humidity

(>80% RH) Moisture vapor settles on component resulting in corrosion and component failure.

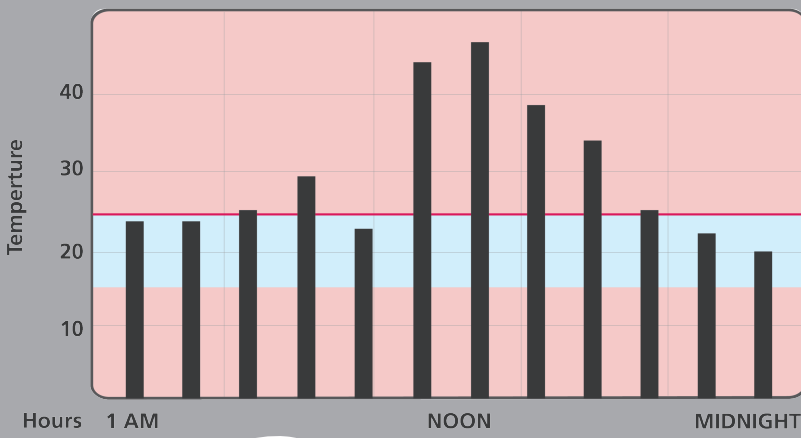
## MONITORING TEMPERATURE AND HUMIDITY

Lab climate conditions can vary dramatically, and these swings can have adverse effects on productivity and results. Here are some temperatures and humidity readings from a typical lab.



### 1 MONTH HUMIDITY VARIATIONS

In summer, high relative humidity can be common in lab environments. With high humidity, moisture becomes a problem and can impact instrument operations.



### 24 HOURS TEMPERATURE VARIATIONS

Temperatures intraday can vary widely. In this lab, a spot near the window in late afternoon causes a serious heat spike.

Out of Range! Ideal Range!  
\*PerkinElmer client: Boston, MA.

## Blue Skies, Nothing but Blue Skies for You and Your Lab!

With our lab-environment and asset-monitoring solution, you can ensure data integrity with automatic data capture and remote monitoring, giving you peace of mind that your instruments are functioning at peak performance.

### OneSource® Asset Genius™ Monitoring

#### 1 Wireless Monitoring

Monitors temperature, humidity, air pressure, and light with minimal IT support.

#### 2 Connected Data Platform

Turnkey solution for connecting, collecting, and analyzing data from your lab.

#### 3 Reporting and Notification

Reports environmental lab conditions and sends alerts if outside recommended parameters.