



Smart-View Pro monitoring case study

Challenge

Finding the right monitoring system to better maintain sample integrity in cold storage can be a challenge. In research, samples in cold storage are precious and often irreplaceable. Many labs store over \$250,000 of product in ultra-low temperature (ULT) freezers. Maintaining a cold storage environment is vital to the future of their research.

Universities and research institutes often insure the contents of their cold storage units. One American university lost over ten million dollars of scientific products due to failures in their sample storage and monitoring process. Their deductible rose from \$5,000 in 2017 to \$750,000 in 2019 due to the number of claims. A different university had to move to a self-insurance model for their research due to a lack of proactive monitoring.

Traditional laboratory refrigerators and freezers are equipped with local alarms, but they require independent monitoring with configurable notification methods to alert remotely. In many academic settings, there is not a consistent process for effectively monitoring these high-value samples. In some instances, labs depend on a security guard hearing a local alarm and contacting the appropriate personnel.

Solution

This American university wanted to standardize its remote monitoring system for cold storage equipment across campus. They chose the Thermo Scientific™ Smart-View™ Pro remote monitoring solution after testing the system and evaluating multiple monitoring solutions.

Here are some of the reasons why this institution chose the **Smart-View Pro system:**

1. Monitor critical parameters
2. Secure and compliant cloud
3. Extended wireless range
4. Scalable design
5. Easy to use and maintain
6. Customized solutions
7. Minimize hardware investment
8. End-to-end support and service

Monitors the key parameters

The Smart-View Pro device is designed to monitor critical parameters of various laboratory equipment types (ULTs, lab refrigerators and freezers, incubators, etc.). It utilizes redundant probes for measuring temperature and other parameters, with battery back-ups.

Notifies the customer remotely if alarm sounds

Customers receive advanced notice through email, text, and/or phone so they can react before sample damage occurs. Through advanced alarms and notifications, users can reduce the risk of sample loss due to environmental excursions.

Promotes lab responsibility for sample protection

The Smart-View Pro system has full traceability to support 21 CFR Part 11 compliance. Performance history, including environmental conditions and alarms, can be used for training and auditing purposes.

Turnkey solutions and support

Thermo Fisher Scientific, through their Unity Lab Services solutions, is here to help by providing presales, project planning, installation, compliance services, and technical support.

Results

The Smart-View Pro monitoring system can provide peace of mind for researchers and university leaders through expert installation services, robust user training, and ongoing remote support services. The system can also help save on costs through reduced insurance deductibles. In this case, the American university will save \$745,000 annually.



Find out more at thermofisher.com/smartvuepro

ThermoFisher
SCIENTIFIC