

Smart Notes



Why should scientists invest in a remote monitoring solution as part of their risk mitigation strategy?

A system like the Thermo Scientific™ Smart-Vue™ Pro Remote Monitoring System will support sample security, workflow efficiency, asset management and operational cost savings.

Your work represents hope – for an answer, for a breakthrough, for a therapy, or maybe a cure. Sample integrity is paramount to your success. Confidence that your one-of-a-kind patient samples are at the proper temperature is necessary for peace of mind. Regulatory bodies require validation of the conditions of your clinical products and materials. The reality is that many factors can affect instrument performance, including routine maintenance, usage behavior, power fluctuations, and mechanical stress over time. If the instruments fail to maintain settings accurately, then your scientific products could be at risk. You want to be prepared for the unexpected.



Thermo Scientific™ Smart-Vue™ Pro Remote Monitoring System

Why should scientists invest in a remote monitoring as part of their risk mitigation strategy?

Sample safety and security

- **a. Situational awareness:** Instant alerts via email/SMS/phone for out-of-range sample conditions enable urgent proactive response to manage your scientific products before they are compromised. Confidence in knowing the conditions of the scientific products anytime, anywhere allows the scientific team to focus on their next scientific breakthrough.
- **b. Compliance:** Automatic and continuous data logging creates a permanent record of readings, alarms, and acknowledgements, which assists with compliance to SOP, cGxP and supports 21 CFR part 11 among other standards. Choosing a monitoring system that is “future-ready” and scalable is important as your science evolves from research to production.

Workflow efficiency

- **a. Downtime prevention:** Proactive early notification can mitigate downtime by minimizing the number of people and hours spent managing excursions.
- **b. Automation:** Automated monitoring systems log data and create reports instantly, making the process effortless. Flexible data transmission, data storage options, and data security are all considerations in choosing the best monitoring solution for your lab.
- **c. Redundancy:** Should on-board sensor technology be compromised; a third-party remote monitoring sensor can automatically provide a redundant data log. Don't let storage conditions be a variable in your experimental data; keeping a reference log will support consistent and reproducible results.

Asset management and cost savings

A remote monitoring system can serve as part of your asset management strategy, providing an opportunity for capital and operational cost avoidance. An early warning system promotes proactive response to an alarm event that can save hundreds of thousands of dollars in sample and product replacement. Additional measurement parameters can also provide early warning to equipment changes and signal preventative maintenance, prolonging the life and optimizing performance of the equipment investment. Well-maintained equipment can also reduce the burden on building operations – reducing the demand for energy and HVAC resources, resulting in operational cost efficiencies. Finally, a remote monitoring system can contribute to minimizing cost for insurance policies that cover replacement of equipment and products. These factors combined can free up budget for the company or research institute to fund additional scientific programs and activities.

Find out more at thermofisher.com/wirelessmonitoring