

Cold storage

Enabling global leadership in vaccine manufacturing

South Korean governing authority quickly deploys cold storage equipment and solutions to support fast-track national vaccine development program

Summary

With robust biopharmaceutical manufacturing capabilities and \$5.1 billion (US) of exported biopharmaceutical goods in 2020, South Korea is set to become one of the top five global vaccine manufacturers by 2025 [1].

A local governing authority needed to run clinical trials as part of the nation's effort to provide local pharma and biopharma companies with sample data and results to aid in the development of homegrown vaccines. With clinical trials conducted in various stages for the development of four COVID-19 vaccines, the governing authority needed reliable ultra-low temperature (ULT) freezers to help ensure that vaccines were stored properly without compromising on sample integrity, data, and results.

“We are honored to be selected as the partner and provider of vaccine storage equipment and solutions to help fast-track the nationwide vaccine development initiative to become one of top global leading vaccine manufacturers.”

—Seok Soojin, Vice President and General Manager,
Thermo Fisher Scientific South Korea

The challenge: a race against time while protecting the integrity of clinical trial data and results




Cold storage equipment and solutions required at short notice to support the development of homegrown vaccines for COVID-19

Integrity of vaccine development throughout workflow must not be compromised, as data and results were vital to aid local biopharma companies in the development of vaccines

Prompted by a nationwide initiative to become one of the top global vaccine manufacturers by 2025 [1], a vaccine development governing body in South Korea needed cold storage equipment and solutions to support local vaccine-developing pharma and biopharma companies in the research and development of homegrown vaccines for COVID-19—and they needed them fast.

Acting as both the consultant for the development of homegrown vaccines and the final approver for their commercial use, the governing authority needed to run clinical trials as part of their effort to provide sample data and results to aid in the development of vaccines. If sample analysis results and data from the clinical trials were compromised, not only would it set the local companies back in terms of resources, they might also slip off the required timeline to fast-track the national vaccine development program.

The solution: timely delivery of high-performance vaccine storage equipment and solutions

	17 units of Thermo Scientific™ TSX Series Ultra-Low Temperature Freezers
	14 units of Thermo Scientific™ TSX Series High-Performance Manual Defrost Freezers
	17 units of Thermo Scientific™ TSX Series High-Performance Lab Refrigerators

Working with our channel partner, our cold storage equipment was delivered and installed with a 300% reduction in lead time, meeting the governing authority’s requirements. The equipment will be used to store vaccine candidates of various types for stability testing and research in terms of time and responsiveness.

“We are proud to be a part of a national project that contributes to the public health where our partner trusted us to provide the right expertise, equipment, and solutions.”

—Edward Park, Product Management Specialist, Controlled Temperature Technologies, Thermo Fisher Scientific South Korea

The impact: sample analysis results and data integrity contributing to seamless vaccine development

	Effective, efficient, and uninterrupted vaccine research and development
	Accuracy and integrity in data and results from clinical trials that help local pharma and biopharma companies in their vaccine development initiatives

With our cold storage solutions, the governing authority was able to establish a systematic workflow of vaccine development, providing guidelines to the local pharma and biopharma companies in the development of homegrown vaccines.

Additionally, the governing authority was able to provide accurate sample analysis data and results from their clinical trials, supporting local companies with the information required to fast-track vaccine development. This contributed to the national initiative of becoming one of the top global vaccine manufacturers.

“We were able to meet our partner’s expectations and are honored that we were able to support them with our corporate mission of enabling our customers to make the world healthier, cleaner and safer.”

—Jee Bonggyu, Account Manager, Thermo Fisher Scientific South Korea

Reference:

1. <https://www.straitstimes.com/asia/east-asia/south-korea-aims-to-be-among-top-5-global-vaccine-producers-by-2025>

Learn more at thermofisher.com/vaccinestorage

thermo scientific