



Cardinal Meat Specialists Uses NetPlus Reports to Ensure Product Quality

Cardinal Meat Specialists is one of Canada's leading meat processors and sealed-environment cooked protein ("sous-vide") manufacturers. Operating two plants in Ontario, Cardinal supplies top restaurants and retailers, emphasizing food quality, consistency, and safety.

Cardinal produces a wide range of meat products, and quality and consistency are key priorities. So the company invested in a DA 7250 SD NIR analyzer to test raw-material blends, optimize formulas, and verify finished-product quality. The system has become a workhorse for Cardinal, analyzing more than 100 samples per day.

In addition to monitoring results from the instrument's main screen, Cardinal implemented our NetPlus Reports web-based software for additional verification. "The NIR instrument is located on the busy production floor," says Jerzy Karwowski, Cardinal's QA manager. "Normally this would present a challenge: one cannot easily access the results history while the machine is in heavy use by operators. NetPlus Reports solves this by allowing remote access to the database. Our quality assurance team can remotely verify product parameters, specifically fat content, used for product release."

NetPlus Reports enables Cardinal to monitor trends from their office computers and helps during audits by facilitating retrieval of stored data. Results are quickly sorted and viewed based on sample analysis data and identification.

The reports generated by the software help Cardinal's operations and quality assurance teams analyze data and make fact-based decisions. As a result, overall process efficiency is improved, enabling Cardinal to deliver high-quality products to their customers.

Meeting the Challenge

"Since the introduction of NetPlus Reports to the DA 7250 NIR instrument, accessing critical data remotely and in real time has helped Cardinal strengthen our commitment to product quality. In a short period of time, NetPlus Reports – which is easily accessible, fast, and reliable – has become an essential tool for our quality assurance department."

Jerzy Karwowski
Quality Assurance Manager
Cardinal Meat Specialists