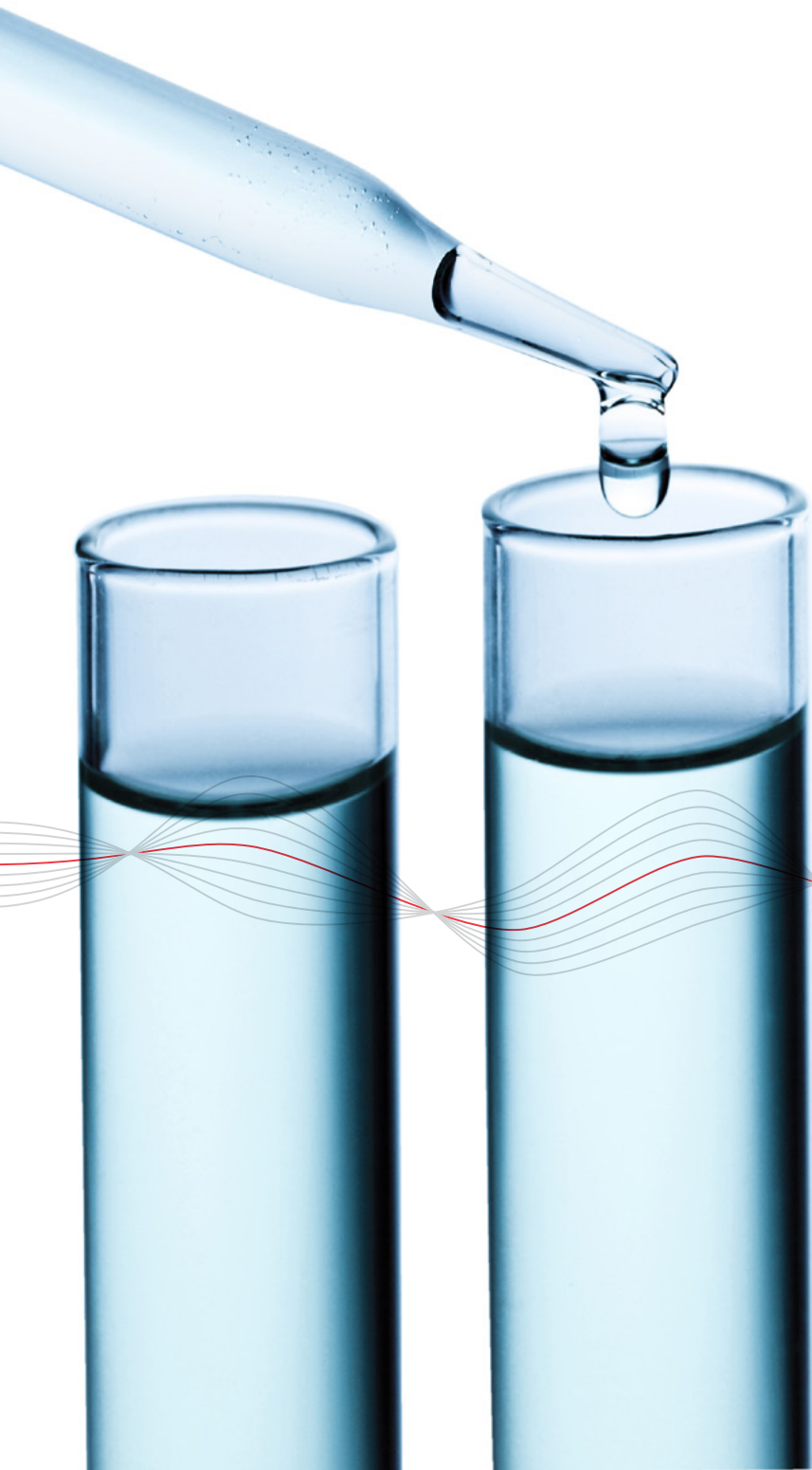


Environmental & industrial applications compendium

Gallery Discrete Analyzer



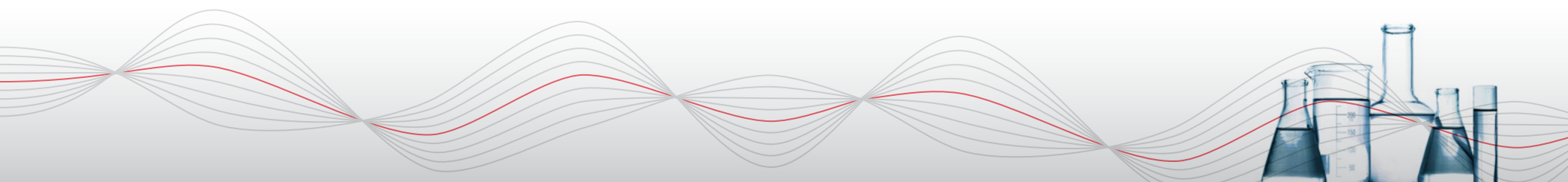


Table of contents

General technology and product overview

Smart note: What are automated discrete analyzers? How do automated discrete analyzers work?

Smart note: What is better for automating wet chemical analysis? Integrated discrete analyzer or flow analyzers?

Brochure: Easy, fast, and simultaneous pH and conductivity measurements

Industrial water analysis applications

Brochure: Solution for routine and comprehensive water analysis

C&EN webinar executive summary: Process-water characterization using discrete analyzers and ion chromatography

Application note: Determination of water pollutants using photometric analysis

Application note: Enzymatic analysis of urea in swimming pool water

Nutrient analysis of drinking water, wastewater, and soil samples

Brochure: Fast, safe, and easy nutrient analysis

Smart note: Wastewater surveillance of COVID-19 with Thermo Scientific™ Gallery™ Discrete Analyzers

Application note: Automated total oxidized nitrogen method using vanadium as reductant with correlation to cadmium and hydrazine reductant methods in sea, natural, and wastewaters

Case study: Increasing analytical efficiencies in the city of Atlanta

Case study: Faster and better analysis of complex samples in a geochemical laboratory

Case study: Easy soil testing at a South African fertilizer manufacturer

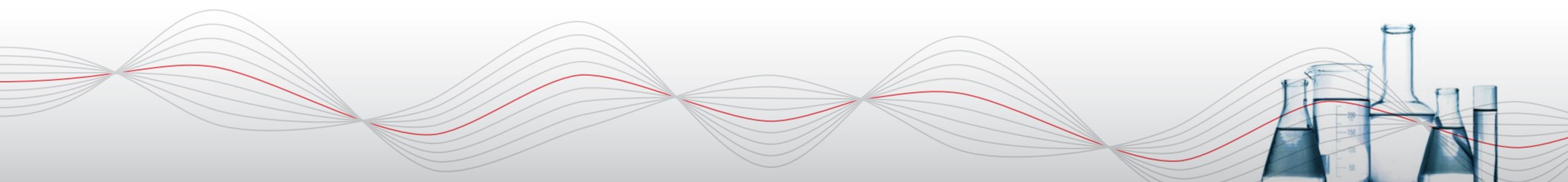
Case study: Precise analysis for inorganic fertilizers at the Finnish Food Authority

Case study: National Laboratory Service achieves high-throughput biomarker testing for SARS-CoV-2 wastewater surveillance using Gallery Plus discrete analyzers

Case study: How the Eurofins Nutrition Lab in Nantes increased capacity by 180% and reduced runtimes by 60%—Optimizing efficiency, productivity, and sensitivity through collaboration

Process water and nutrient analysis

Analytical guide: Automated nutrient analysis and water quality monitoring



What are automated discrete analyzers? How do automated discrete analyzers work?

Overview

Automated discrete analyzers utilize colorimetric and enzymatic measurements from a single sample through photometric analysis. The discrete analyzer imitates the lab chemists' operation sequence of dispensing samples, mixing reagents, incubation, and photometric measurement. The discrete analyzer provides fast and reproducible results. In discrete analysis, each individual reaction cell is isolated and the temperature is stabilized, enabling highly controlled reaction conditions.

This smart note introduces you to automated discrete analyzer technology and how it functions.

thermo scientific

DISCRETE ANALYZER SELECTION GUIDE Gallery automated discrete analyzers

Smart Note

QA

What are automated discrete analyzers? How do automated discrete analyzers work?

Automated discrete analyzers utilize colorimetric and enzymatic measurements—of several analytes simultaneously—from a single sample through photometric analysis. The discrete analyzer imitates the lab chemists' operation sequence of dispensing samples, mixing reagents, incubation, and photometric measurement; however, the discrete analyzer provides fast and reproducible results. Discrete analyzers consist of four components: a photometer with a specific number of filter positions; dispensing probes; an incubator to control the reaction temperature; and a mixer. In discrete analysis, each individual reaction cell is isolated and the temperature is stabilized, enabling highly controlled reaction conditions.

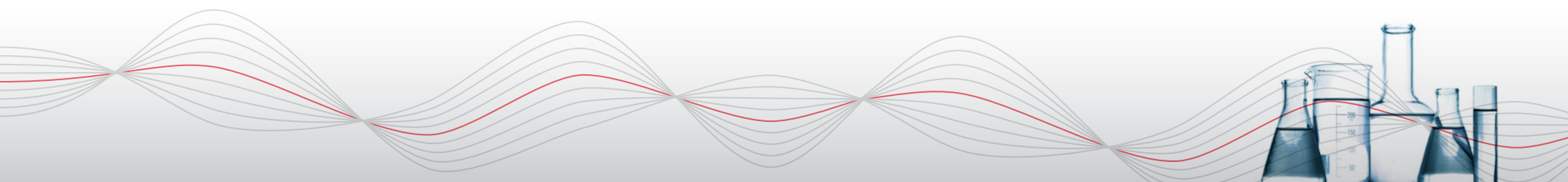
Dispensing → Mixing → Incubation → Photometric measurement → Result interpretation

Figure 1. Discrete analyzer workflow.

After the reagents and samples are prepared, they are loaded onto the instrument. Next, the individual cuvettes are loaded into the incubation chamber and the samples and reagents are dispensed to the individual cuvettes and then mixed. Finally, the combined samples and reagents undergo photometric detection, depending on the absorbance of specific wavelengths of light. Each measurement is done using single discrete cuvettes and this data is then interpreted through integrated software platforms.

ThermoFisher SCIENTIFIC

[View the full smart note](#)



What is better for automating wet chemical analysis? Integrated discrete analyzer or flow analyzers?

Overview

In general, Flow Injection Analyzers (FIA), Segmented Flow Analyzers (SFA) or Continuous Flow Analyzers (CFA) are batch analyzers, meaning they are particularly suitable for analyzing a few parameters for a large number of samples. Flow systems use specific detector modules, reagents delivery and mixing, which limits the number of parameters that they can test per sample. Typically, flow systems test 2 to 6 maximum different parameters per sample. If a laboratory is looking for an easy-to-use, high throughput, expandable, multiparameter, wet chemistry analyzer for large numbers of samples, then integrated discrete analyzers are better suited than FIA, SFA or CFA.

This smart note provides detailed comparisons between the different technologies, and walks you through the advantages of each technology.

thermo scientific

SMART NOTE Gallery discrete analyzers

What is better for automating wet chemical analysis? Integrated discrete analyzer or flow analyzers?


QA

Technology selection considerations should be based on various factors:

- Current and future sample analysis load
- Additional costs incurred by adding additional tests
- Number of parameters per sample
- Complexity of wet chemistry parameters testing for each sample
- Detection limits
- Reagent consumption
- Waste generation
- Cost per analysis
- User's skill level
- Regulatory requirements
- Maintenance and bench space requirement of equipment
- Total cost of ownership and Return on Investment (ROI)

In general, Flow Injection Analyzers (FIA), Segmented Flow Analyzers (SFA) or Continuous Flow Analyzers (CFA) are batch analyzers, meaning they are particularly suitable for analyzing few parameters for large number of samples. Flow systems use specific detector modules, reagents delivery and mixing, that limits the number of parameters that they can test per sample. Typically, flow systems test 2 to max 6 different parameters per samples. If a laboratory is looking for an easy to use high throughput, expandable, multiparameter, wet chemistry analyzer for large numbers of samples, then integrated discrete analyzers are better suited than FIA, SFA or CFA.

ThermoFisher
SCIENTIFIC

 [View the full smart note](#)

[General technology
and product overview](#)

[Industrial water
analysis applications](#)

[Nutrient analysis of drinking water,
wastewater, and soil samples](#)

[Process water and
nutrient analysis](#)

 [Back to contents](#)

Easy, fast, and simultaneous pH and conductivity measurements

Overview

pH and conductivity measurements provide crucial insights for a variety of industries, including the food and beverage, industrial process, enzyme kinetics and water analysis sectors. Fast, accurate and cost-effective pH and conductivity measurement workflows create the stream of regular, meaningful data that drives important decisions. By regularly testing all parameters of a manufacturing line, process problems are detected early, enabling intervention and improvements that protect equipment, product consistency and quality standards. In turn, the collection of accurate and timely data provides evidence for regulatory approval and audit submissions.

This brochure discusses how the integrated electrochemical measurement (ECM) module in Gallery discrete analyzers provides parallel, automated electrochemical measurement of pH and conductivity along with complete photometric testing.



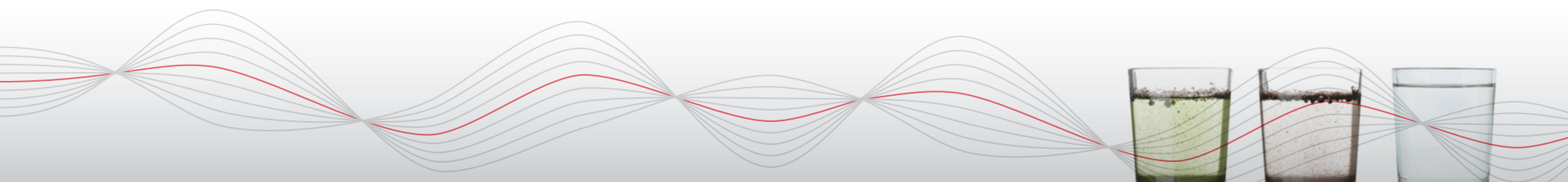
[View the full brochure](#)

[General technology
and product overview](#)

[Industrial water
analysis applications](#)

[Nutrient analysis of drinking water,
wastewater, and soil samples](#)

[Process water and
nutrient analysis](#)



Solution for routine and comprehensive water analysis

Overview

Process water quality is an integral part of the industrial manufacturing process. Multi-component analysis of corrosive ions, scaling ions, corrosion inhibitors, and corrosion indicators provides detailed information for successful and predictable unit operations. Routine testing of water samples for multiple analytes increases the overall sample testing demand on utilities and central laboratories.


This brochure summarizes how [Thermo Scientific™ discrete analyzers and ion chromatography \(IC\) systems](#) can offer rapid multiparameter testing, unattended high throughput analysis of large sample numbers, and cover a wide concentration range for sample types of varying ionic strengths for industrial process water quality monitoring. Together, discrete analysis and IC solve a majority of water analysis needs for improved productivity and reliability.

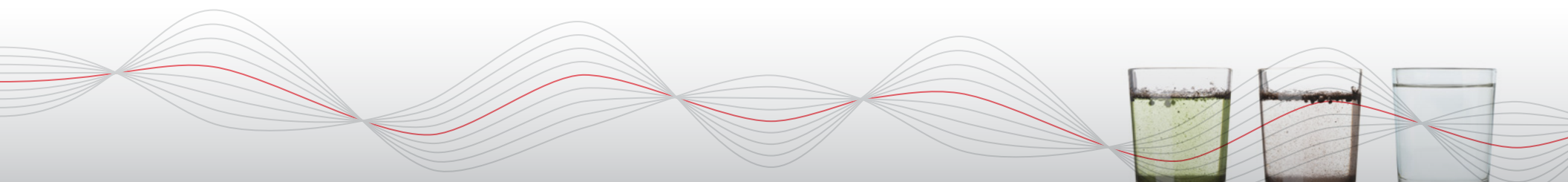
thermo scientific

Solution for routine and comprehensive water analysis

Thermo Scientific Disc-IC Systems—
a fully-automated high throughput solution

ThermoFisher
SCIENTIFIC

 [View the full brochure](#)



C&EN webinar: Process-water characterization using discrete analyzers and ion chromatography

Overview

Process water, the non-drinkable water that many companies rely on for a variety of industrial applications, undergoes extensive treatment. This water must be analyzed and processed to meet the needs of specific industries.

Process-water analysis can be challenging. This C&EN executive summary provides an overview of the appropriate instruments and methods to address those challenges. [The Gallery discrete analyzer](#) provides routine, high-throughput analysis for many parameters of interest, including pH, conductivity, and concentration, and can achieve up to ppb levels of detection and does not require skilled operators. It is a “walkaway” solution for many water analysis needs.

While the discrete analyzer platform is suitable for a wide range of user needs, those who require a comprehensive solution for water analyses should consider also incorporating a Thermo Scientific™ Dionex™ ion chromatography (IC) system into their lab, which can characterize a much broader range of anions and cations and achieve up to ppt levels of detection. IC can also help minimize signal interference from a sample matrix.

c&en | WEBINAR

PROCESS-WATER CHARACTERIZATION USING DISCRETE ANALYZERS AND ION CHROMATOGRAPHY

OVERVIEW
Process water, the non-drinkable water that many companies rely on for a variety of industrial applications, undergoes extensive treatment. This water must be analyzed and processed to meet the needs of specific industries. Several technologies can address a variety of water analysis needs and can increase the speed, quality, and consistency of process-water data analysis.

INTRODUCTION
Maintaining high-quality water is critical for many industries that use water as part of their manufacturing processes. Poor-quality water can corrode costly equipment, waste resources, and compromise the end product.

To meet the needs of a specific application, source water—which includes groundwater and raw water—must first be converted to deionized water using a series of treatment steps. Each water sample must be tested for multiple parameters, to ensure that the produced and reused process water is free of contamination. Companies must test the water samples at the point of use and multiple times throughout the industrial process, as the water quality changes as it passes through the various process stages (Figure 1).

Figure 1: Sample types and testing parameters involved in process-water analysis. “DM” stands for “demineralized” in this figure.

Image credit: Thermo Fisher Scientific

AMERICAN CHEMICAL SOCIETY | Brought to you by **ThermoFisher SCIENTIFIC**

[View the full executive summary](#)

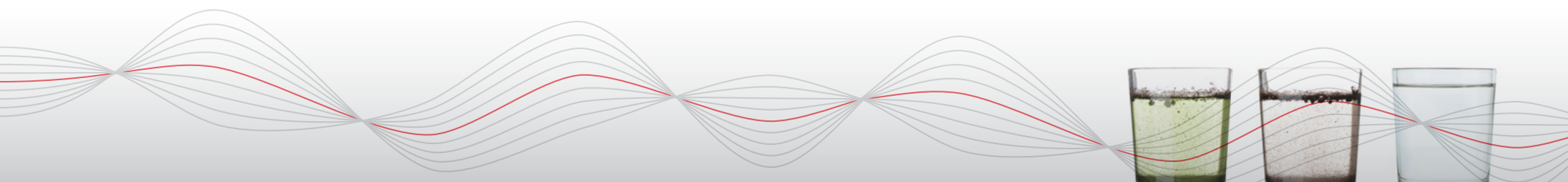
General technology and product overview

Industrial water analysis applications

Nutrient analysis of drinking water, wastewater, and soil samples

Process water and nutrient analysis

[Back to contents](#)



Determination of water pollutants using photometric analysis

Overview

The International Organization of Standardization (ISO) has published more than 550 standards to address water issues, which provide a common language for water sampling, reporting, and monitoring in order to ensure purity and other desired characteristics that apply to industrial processes as well as natural water.

In this study, the methods and performance of [Thermo Scientific™ discrete analyzers](#) are described according to the analysis guidelines of ISO 15923-1. All data is generated using the automated Gallery discrete analyzers and [Thermo Scientific™ system reagents](#) for applications via photometric and enzymatic analysis.

thermoscientific

APPLICATION NOTE 71728

Determination of water pollutants using photometric analysis

Authors
Marco Rastetter,¹ Jo Gawler,¹
Mari Kiviluoma,² and
Annu Suoniemi-Kähärä³
¹Thermo Fisher Scientific, UK
²Thermo Fisher Scientific, Germany
³Thermo Fisher Scientific, Finland


Introduction
Clean water in sufficient amounts is an ongoing global challenge. Water resources are experiencing increasing pressure in many parts of the world, requiring countries to improve the management and protection of water ecosystems. In 2000, the European Union's (EU) Water Framework Directive (WFD) announced an integrated approach to protecting water quality, quantity, and the role of habitat. As a result, governmental bodies of different EU member states are required to work with one another to ensure that the requirements of the directive are met and water quality is maintained. Quality criteria include nutritional composition as well as chemical composition, i.e. the level of pollutants.

Keywords
DA, Automated analysis,
ISO 15923-1, water analysis,
discrete analysis

Goal
To demonstrate compliance with the ISO 15923-1 standard for the determination of water pollutants using an automated photometric procedure.

The International Organization of Standardization (ISO) develops standards to benefit state authorities, regulatory bodies, and industry for the purpose of equitably and durably managing shared water resources. As a network composed of 163 countries, ISO published more than 19,500 international standards covering almost every industry from technology and food safety to agriculture and healthcare. More than 550 standards currently exist to address water issues and these quality standards provide a common language for water sampling, reporting, and monitoring in order to ensure purity and other desired characteristics that apply to industrial processes as well as natural water.

ThermoFisher
SCIENTIFIC


 [View the full application note](#)

General technology
and product overview

**Industrial water
analysis applications**

Nutrient analysis of drinking water,
wastewater, and soil samples

Process water and
nutrient analysis

 [Back to contents](#)

Enzymatic analysis of urea in swimming pool water

Overview

Nitrogen-containing impurities such as urea, ammonia, amino-acids, creatinine, and uric acid introduced to swimming pool water by bathers react with free chlorine to form chlorine-containing compounds. It is important to control the level of urea in swimming pool water, as urea is a potential source of hazardous ammonia chloramines and a possible nutrient for bacteria and algae, all of which pose a hygienic risk.

Urea is typically measured by the Koroleff method, which is based on persulphate digestion. Urea can also be measured using an enzymatic method using the [Thermo Scientific discrete analyzers](#). In the SYKE (Finnish Environment Institute) swimming pool water report, six of the proficiency test participants used the Koroleff method and three laboratories used the enzymatic method.

This application note demonstrates that the Koroleff method results were generally lower than results obtained from the enzymatic method. Results of the enzymatic method were closer to calculated results of the proficiency test samples. By switching from the Koroleff method to the enzymatic method, the result levels are expected to be more accurate and therefore higher than previously reported results.

thermoscientific

APPLICATION NOTE 71461

Enzymatic analysis of urea in swimming pool water

Authors
Kirsti Nikkila,¹ Marita Jokinen,²
Tuula Jumppanen,¹ Mari Klemm,²
Juhanii Airo,¹ and Annu
Suonemi-Kähärä²
¹Metropolilab Oy, Helsinki, Finland
²Thermo Fisher Scientific,
Vantaa, Finland

Introduction
Nitrogen-containing impurities such as urea, ammonia, amino-acids, creatinine, and uric acid introduced to swimming pool water by bathers react with free chlorine to form chlorine-containing compounds. It is important to control the level of urea in swimming pool water because urea is a potential source of hazardous ammonia chloramines and a possible nutrient for bacteria and algae, all of which pose a hygienic risk.

In Finland, urea concentration in swimming pools is regulated by Valvira (National Supervisory Authority for Welfare and Health). Based on current guidelines, urea concentration must be less than 0.8 mg/L.

Urea is typically measured by the Koroleff method,² which is based on persulphate digestion. Urea can also be measured using an enzymatic method. In the SYKE (Finnish Environment Institute) swimming pool water report,³ six of the proficiency test participants used the Koroleff method and three laboratories used the enzymatic method. Based on the study, the Koroleff method results were generally lower than results obtained from the enzymatic method. Results of the enzymatic method were closer to

Keywords
Discrete Analyzer, Photometric Analyzer, Aquakem

Goal
To use an enzymatic method for determination of urea in the water of swimming pools

ThermoFisher
SCIENTIFIC

[View the full application note](#)





Fast, safe, and easy nutrient analysis

Overview

Determining nutrient levels in drinking water, wastewater, and soil samples is critical for protecting aquatic habitats and maintaining clean and safe drinking water supplies. Utility companies and environmental laboratories must regularly measure elemental phosphorus and nitrogen in sewage water, along with a range of other pollution indicators, to ensure discharge streams are compliant with regulatory standards. Wastewater nutrient analysis can also be used for the assessment of population-level infection, including SARS-CoV-2 surveillance, by providing important biomarker indications for population size.


Traditional wet chemistry techniques, including titrations, flow injection analysis, and other colorimetric techniques, have been used for many years to undertake nutrient analysis of drinking water, wastewater, and soil samples. However, these approaches are slow, labor-intensive and often unreliable, involving hazardous reagents that add substantial costs for waste disposal.

This brochure summarizes how [Thermo Scientific discrete analyzers](#) can provide a faster, safer, and more reliable replacement to traditional wet chemistry methods for nutrient analysis.



Fast, safe, and easy nutrient analysis

Thermo Scientific Gallery and Gallery Plus discrete analyzers – for complete and simultaneous nutrient analysis of drinking water, wastewater, and soil samples

 [View the full brochure](#)

Wastewater surveillance of COVID-19 with Gallery discrete analyzers

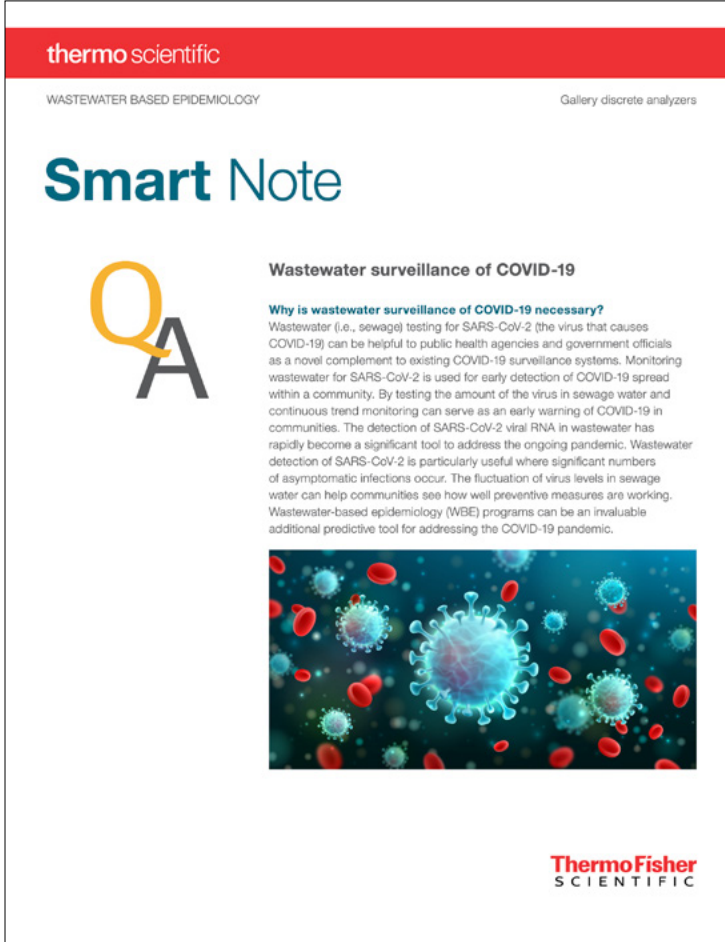
Overview

The detection of SARS-CoV-2 viral RNA in wastewater has rapidly become a significant tool to address the ongoing pandemic. Wastewater detection of SARS-CoV-2 is particularly useful where significant numbers of asymptomatic infections occur.

Determining the various nutrient parameters such as nitrogen, phosphorus, ammoniacal nitrogen, urea and creatinine, pH, and conductivity for a large number of samples on a daily basis is highly challenging. A discrete analyzer consolidates these parameters and offers a high throughput solution.

In this smart note, learn about strategies to consolidate chemical reference parameters and population biomarker analysis to support the wastewater-based epidemiology.

For more information, [watch this interactive webinar with Reg Godwin, from the Environment Agency UK](#), as he discusses sewage water monitoring by discrete analyzer for population biomarkers and shares their early study results.



thermo scientific

WASTEWATER BASED EPIDEMIOLOGY Gallery discrete analyzers

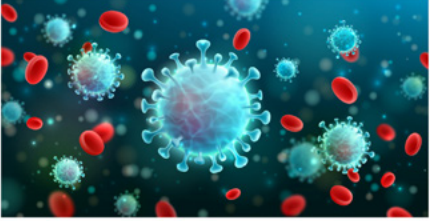
Smart Note

QA


Wastewater surveillance of COVID-19

Why is wastewater surveillance of COVID-19 necessary?

Wastewater (i.e., sewage) testing for SARS-CoV-2 (the virus that causes COVID-19) can be helpful to public health agencies and government officials as a novel complement to existing COVID-19 surveillance systems. Monitoring wastewater for SARS-CoV-2 is used for early detection of COVID-19 spread within a community. By testing the amount of the virus in sewage water and continuous trend monitoring can serve as an early warning of COVID-19 in communities. The detection of SARS-CoV-2 viral RNA in wastewater has rapidly become a significant tool to address the ongoing pandemic. Wastewater detection of SARS-CoV-2 is particularly useful where significant numbers of asymptomatic infections occur. The fluctuation of virus levels in sewage water can help communities see how well preventive measures are working. Wastewater-based epidemiology (WBE) programs can be an invaluable additional predictive tool for addressing the COVID-19 pandemic.



ThermoFisher
SCIENTIFIC

 View the full case study

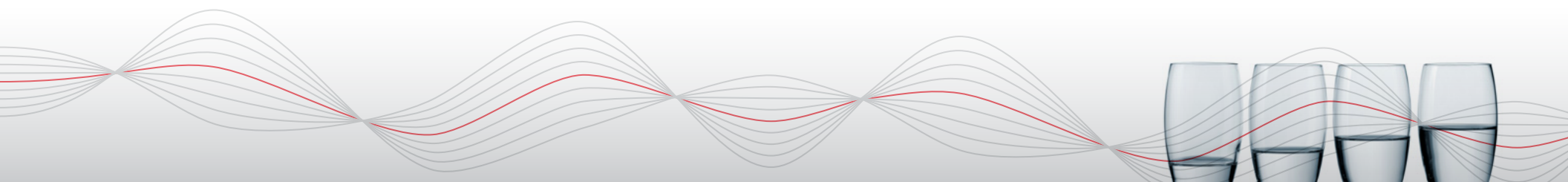
General technology
and product overview

Industrial water
analysis applications

Nutrient analysis of drinking water,
wastewater, and soil samples

Process water and
nutrient analysis

 Back to contents



Automated total oxidized nitrogen method using vanadium as reductant with correlation to cadmium and hydrazine reductant methods in sea, natural, and wastewaters

Overview

Measurement of total oxidizable nitrogen (TON) is required for accurate determinations of nitrate concentrations in drinking water and wastewater. Traditional TON measurement methods based on cadmium reduction coils have several limitations, such as costly waste disposal and carcinogenic health risks. [Thermo Scientific discrete analyzers](#) can be used as a safer alternative for TON measurement, as they support and automate new EPA-approved enzymatic reduction methods.

This application note details the automated TON method using the Gallery discrete analyzer.

thermoscientific

APPLICATION NOTE 71395

Automated total oxidized nitrogen method using vanadium as reductant with correlation to cadmium and hydrazine reductant methods in sea, natural, and waste waters

Authors
Tuula Juntippanen,¹ Marita Jokinen,¹ Juhani Aho,² Mari Klemm,³ Daniel Höger,⁴ and Annu Suoniemi-Kihäri⁵
¹Thermo Fisher Scientific, Vantaa, Finland; ²Metropolilab Oy, Helsinki, Finland; ³Thermo Fisher Scientific, Sydney, Australia


Introduction
Total oxidized nitrogen (TON) or the sum of nitrate and nitrite in water samples can be determined using a variety of techniques. Hydrazine is often used in automated analysis as the reducing agent, converting nitrate to nitrite.¹ Unfortunately the hydrazine method cannot be used to analyze a seawater sample limiting its use as a general purpose method.² The Cadmium (Cd) Reduction method, therefore, is still the reference method used in many countries. The Cd Reduction method has several drawbacks which include the personal risks to users when handling the Cd-column and the column itself may be easily damaged by using an oily sample matrix.³ Enzymatic reduction is a very good alternate method^{4,5} however it is not widely known to most laboratories, despite recently having been described as an alternate method by the United States Environmental Protection Agency (US EPA).⁶

Keywords
Discrete Analyzer, Aquasem 250, Gallery Plus

Goal
Correlation of a Nitrate method (TON Vanadium) to TON Cadmium and TON Hydrazine methods using automated discrete analysis

In this application note an automated, two reagent method using vanadium chloride as a reducing agent was studied and compared to both the hydrazine and Cd reduction methods. Vanadium chloride as reductant has routinely been used in nitrate analysis. In one application, the reduction step in the method is performed at a temperature greater than 80 °C and nitrous oxide (NO) is measured by chemiluminescent detection.^{6,7} When the

ThermoFisher
SCIENTIFIC

 [View the full application note](#)

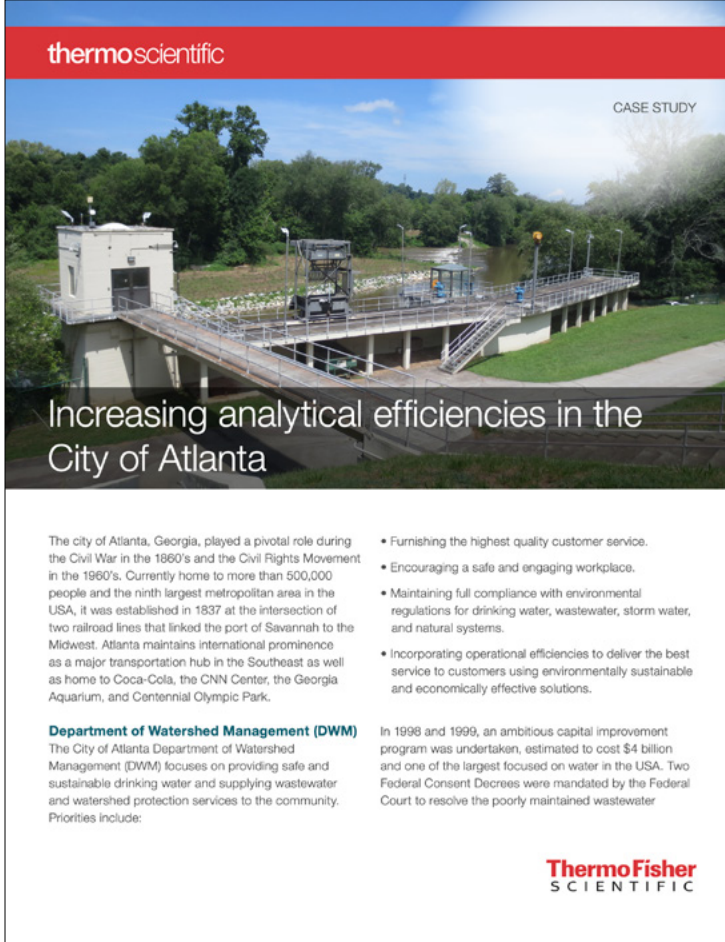
Increasing analytical efficiencies in the City of Atlanta

Overview

The City of Atlanta Department of Watershed Management (DWM) focuses on providing safe and sustainable drinking water and supplying wastewater and watershed protection services to the community. Affectionately dubbed “Gertrude,” their [Thermo Scientific™ Gallery™ Plus discrete analyzer](#) has an attached electrochemical (ECM) unit that allows them to test pH and conductivity in addition to the desired alkalinity and hardness tests.

“Using the Gallery Plus analyzer freed up a technician. We are actually saving manpower and everyone can go home on time”, says Carolyn Duncan, Senior Chemist in the City of Atlanta Watershed Protection Department.

Read more about the department’s experience with their discrete analyzer and automated test workflow in this case study.



The case study cover features a photograph of a wastewater treatment facility with a red header containing the 'thermoscientific' logo and the text 'CASE STUDY'. The main title is 'Increasing analytical efficiencies in the City of Atlanta'. The text on the cover includes:

The city of Atlanta, Georgia, played a pivotal role during the Civil War in the 1860's and the Civil Rights Movement in the 1960's. Currently home to more than 500,000 people and the ninth largest metropolitan area in the USA, it was established in 1837 at the intersection of two railroad lines that linked the port of Savannah to the Midwest. Atlanta maintains international prominence as a major transportation hub in the Southeast as well as home to Coca-Cola, the CNN Center, the Georgia Aquarium, and Centennial Olympic Park.

Department of Watershed Management (DWM)
The City of Atlanta Department of Watershed Management (DWM) focuses on providing safe and sustainable drinking water and supplying wastewater and watershed protection services to the community. Priorities include:

- Furnishing the highest quality customer service.
- Encouraging a safe and engaging workplace.
- Maintaining full compliance with environmental regulations for drinking water, wastewater, storm water, and natural systems.
- Incorporating operational efficiencies to deliver the best service to customers using environmentally sustainable and economically effective solutions.

In 1998 and 1999, an ambitious capital improvement program was undertaken, estimated to cost \$4 billion and one of the largest focused on water in the USA. Two Federal Consent Decrees were mandated by the Federal Court to resolve the poorly maintained wastewater

ThermoFisher
SCIENTIFIC

[View the full case study](#)

Faster and better analysis of complex samples in a geochemical laboratory

Overview

Murphy and Son Limited, The Council for Geoscience (CGS), located in Pretoria, South Africa, offers a wide range of geoscientific and analytical services. Previously, the CGS laboratory used a relatively uncomplicated traditional method for determination of the following anions: fluoride (F), chloride (Cl), nitrite (NO₂), bromide (Br), nitrate (NO₃), phosphate (PO₄), and sulphate (SO₄). They were particularly interested in examining anion concentrations in water as well as leachable anion concentrations in rock and soil samples.

CGS's investment in the [Gallery Plus discrete analyzer](#) enabled them to achieve a huge improvement in anion analysis capability. The lab is now able to examine additional parameters such as nitrate and phosphate at low concentrations, hexavalent chromium (Cr (VI)), ammonia, and cyanide with high throughput testing capacity and minimal maintenance.

Read more about the CGS story and their testing capability improvement for environmental analyses, drinking water quality monitoring, mine water evaluations and geochemical and hydrological mapping in this case study.



thermoscientific

CASE STUDY

Faster and better analysis of complex samples in a geochemical laboratory

In 1887 Murphy and Son Limited, The Council for Geoscience (CGS), located in Pretoria, South Africa, offers a wide range of geoscientific and analytical services. For the past fifteen years the CGS laboratory used a relatively uncomplicated traditional method for determination of the following anions: fluoride (F), chloride (Cl), nitrite (NO₂), bromide (Br), nitrate (NO₃), phosphate (PO₄), and sulphate (SO₄). They are particularly interested in examining anion concentrations in water as well as leachable anion concentrations in rock and soil samples.

Their fairly rigid existing method could not handle high concentrations of Cl and SO₄, in addition to very low concentrations of NO₂ and PO₄, within the same analytical run, a typical requirement of environmental or mine water analysis. Often, samples had to be repeatedly diluted to achieve the desired analytical range. Only 45–50 samples (without dilutions) could be handled in a standard overnight run and the instrument was continually susceptible to damage from high acid or metal matrix concentrations.

Cover image – typical samples at the CGS laboratory

ThermoFisher
SCIENTIFIC

[View the full case study](#)

General technology
and product overview

Industrial water
analysis applications

Nutrient analysis of drinking water,
wastewater, and soil samples

Process water and
nutrient analysis

[Back to contents](#)

Easy soil testing at a South African fertilizer manufacturer

Overview

Omnia Fertilizer is part of the Agriculture Division of Omnia and a market leader in its field in South Africa. In the past, the R&D lab used a Segmented Flow Analyzer (SFA) to measure soil samples. With 4,000 samples to test each day during the growing season, operating a SFA was labor intensive, with high maintenance required for troubleshooting and high operation expertise required.

To solve these issues, the managers of the lab purchased a [Gallery discrete analyzer](#). The company discovered it is operator friendly, simple to use, and capable of saving significant amounts of time. The test results were demonstrated with good repeatability and the LIMS interface was an easy integration.

“Maintenance is substantially less than with our older test methods”, says Edna Laubscher, Chemtech Analytical Services Manager

Read more about the fertilizer manufacturer’s experience with discrete analyzers in this case study.



thermoscientific


CASE STUDY

Easy soil testing at a South African fertilizer manufacturer

Omnia Fertilizer is part of the Agriculture Division of Omnia and a market leader in its field in South Africa. This strong agronomic unit plays an important role in advising and providing tools so farmers can improve their practices. They produce and sell ammoniumnitrate based, chemically granulated compounds, liquids, and specialty fertilizers that optimize the quality and yield of crops while also reducing risk to farmers and the environment. Customers include farmers' cooperatives and wholesalers in East Africa, Australia, New Zealand, and Brazil. They plan to expand into the Europe, South America, and Asia in the near future. With their world class laboratories and production facilities, clients are provided with products that are in compliance with strict standards.

ISO 17025 Standards
ISO 17025 standards specify general requirements for competence in performing tests or calibrations. This directive encompasses sampling, testing, and calibrations performed using standard methods, nonstandard methods, and laboratory developed methods and is recommended for use by laboratories following good practices in quality and technical management.

ThermoFisher
SCIENTIFIC

 [View the full case study](#)

[General technology and product overview](#)

[Industrial water analysis applications](#)

[Nutrient analysis of drinking water, wastewater, and soil samples](#)

[Process water and nutrient analysis](#)

 [Back to contents](#)

Precise analysis for inorganic fertilizers at the Finnish Food Authority

Overview

Evira, the Finnish Food Authority located in Helsinki, Finland is charged with providing effective, efficient, consistent, and risk-based control of the entire food chain. The lab at Evira routinely measures potassium (K) and phosphate (PO_4) for approximately 1,000 inorganic fertilizer samples per year. In just two hours, they are able to complete all the required analyses using the automated [Gallery discrete analyzer](#). The same analysis used to consume two days with old titration and gravimetric methods.

“This is the best method (with discrete analyzers) for measuring PO_4 in inorganic fertilizers. It is much more precise since there is no interference from other analytes”, says Aija Pelkonen, Research Chemist at Evira.

Read more about Evira's experience with discrete analyzers for inorganic fertilizer analysis in this case study.



The banner features a photograph of the Evira building in Helsinki, Finland. The text is overlaid on the image, with a red header containing the Thermo Scientific logo and the words 'CASE STUDY'. The main title of the case study is 'Precise analysis for inorganic fertilizers at the Finnish Food Authority'.

thermoscientific CASE STUDY

Precise analysis for inorganic fertilizers at the Finnish Food Authority

Evira, the Finnish Food Authority located in Helsinki, Finland is charged with providing effective, efficient, consistent, and risk-based control of the entire food chain. This includes foodstuffs and any products used in primary production or agriculture. Ensuring a high level of hygiene throughout the entire production chain, their control begins with seeds and plants, continues through field inspections, and monitors finished products in the market. Control is specifically targeted at fertilizers, pesticides, animal feeds, genetically modified, and organic products. The quality of operations is assured by the Finnish Accreditation Service (FINAS) and seed analytics are determined by the International Seed Testing Association (ISTA).

The lab at Evira routinely measures potassium (K) and phosphate (PO_4) for approximately 1000 inorganic fertilizer samples per year. In two hours they are able to complete all the required analyses using the Thermo Scientific™ Gallery™ automated discrete analyzer they purchased in 2011. The same analysis used to consume two days with old titration and gravimetric methods.


The lab at Evira routinely measures potassium (K) and phosphate (PO_4) for approximately 1000 inorganic fertilizer samples per year. In two hours they are able to complete all the required analyses using the Thermo Scientific™ Gallery™ automated discrete analyzer they purchased in 2011. The same analysis used to consume two days with old titration and gravimetric methods.

The lab at Evira routinely measures potassium (K) and phosphate (PO_4) for approximately 1000 inorganic fertilizer samples per year. In two hours they are able to complete all the required analyses using the Thermo Scientific™ Gallery™ automated discrete analyzer they purchased in 2011. The same analysis used to consume two days with old titration and gravimetric methods.

For validation, Mrs. Pelkonen compared the PO_4 test to a titration method. She found it was good for measuring a high concentration of PO_4 and resulted in an acceptable uncertainty under 8%.

“Phosphate samples are not an issue with the Gallery analyzer and there is no glassware to clean. I hate cleaning glassware,” Mrs. Pelkonen admitted with a smile.

ThermoFisher
SCIENTIFIC


 [View the full case study](#)

General technology
and product overview

Industrial water
analysis applications

Nutrient analysis of drinking water,
wastewater, and soil samples

Process water and
nutrient analysis

 [Back to contents](#)

National Laboratory Service achieves high-throughput biomarker testing for SARS-CoV-2 wastewater surveillance using Gallery Plus discrete analyzers

Overview

Wastewater testing for SARS-CoV-2 can complement public health agencies' existing COVID-19 surveillance systems with early detection and prediction of COVID-19 spread in a community. The National Laboratory Service (NLS) of Environment Agency UK is one of the laboratories leading the way in the emerging field of wastewater epidemiology.

Determining the chemical reference parameters such as pH, conductivity, and population biomarkers for a large number of samples on a daily basis is extremely challenging, and until now not practical. The automated Gallery discrete analyzer overcomes these challenges. Automated features and parallel determination of multiple analytes from a single sample increase sample throughput and analytical certainty to support public health agencies. In addition, the analyzer's automation features provide true walkaway time for operators to perform other tasks or to monitor multiple systems.

Learn about how NLS consolidates chemical reference parameters and population biomarker analysis to support the wastewater-based epidemiology in this case study.



thermoscientific

CASE STUDY 000263

National Laboratory Service achieves high-throughput biomarker testing for SARS-CoV-2 wastewater surveillance using Gallery Plus discrete analyzers

"If you can warn local or national governments of a COVID outbreak about to occur, that's hugely powerful. Now we can see three to five days ahead if there's going to be a lot of cases presenting at hospitals. Having that early warning via quick, 24-hour turnaround on our biomarker testing, together with the molecular biology results, is something that we're really proud to have implemented. Without the Gallery Plus discrete analyzer, testing wastewater biomarkers at the scale and throughput needed wouldn't be feasible."

—Reg Godwin, Technical Specialist, National Laboratory Service, Environment Agency UK

Several Thermo Scientific™ Gallery™ Plus Discrete Analyzers are in operation providing high-throughput wastewater biomarker analysis at the NLS's Exeter laboratory. Photo courtesy of the NLS.

Gallery Plus discrete analyzer enables wastewater epidemiology program

Wastewater testing for SARS-CoV-2 can complement public health agencies' existing COVID-19 surveillance systems with early detection and prediction of COVID-19 spread in a community. Monitoring SARS-CoV-2 in wastewater is particularly useful where significant numbers of asymptomatic infections occur. In addition, the fluctuation of virus levels in sewage water can help agencies determine the effectiveness of preventive measures. The National Laboratory Service (NLS) of Environment Agency UK is one of the laboratories leading the way in the emerging field of wastewater epidemiology.

ThermoFisher SCIENTIFIC

[View the full case study](#)

General technology and product overview

Industrial water analysis applications

Nutrient analysis of drinking water, wastewater, and soil samples

Process water and nutrient analysis

[Back to contents](#)

How the Eurofins Nutrition Lab in Nantes increased capacity by 180% and reduced runtimes by 60%—Optimizing efficiency, productivity, and sensitivity through collaboration

Overview

The Eurofins Nutrition Lab in Nantes sought to increase the capacity of their equipment, improve turnaround times, and maximize the sensitivity of their testing—without placing an additional burden on their technicians. The result is the lab has an upgraded workflow that has boosted capacity by 180%, reduced runtimes by 60%, and improved and streamlined the global workflow for sugar profile analysis.

“Thermo Fisher Scientific is already highly recommended as a supplier for equipment for [chromatography](#), inorganic analysis, [colorimetric and enzymatic analysis](#), sample preparation, and more—and we would recommend them to any Eurofins laboratories looking to improve their performance.”, says Romain Rivoallan, Production Manager in the Eurofins Nutrition Lab.

In this case study, we expand upon the lessons learned and benefits experienced by the Eurofins team in Nantes from the collaboration and demonstrate how advanced analytical equipment can benefit not only nutritional analysis but Good Manufacturing Practice workflows for any product.

thermoscientific

CASE STUDY 73674

How the Eurofins Nutrition Lab in Nantes increased capacity by 180% and reduced runtimes by 60%

Optimizing efficiency, productivity and sensitivity through collaboration

The result? An upgraded workflow that has:

- Boosted capacity by 180%
- Reduced runtimes by 60%
- Improved and streamlined the global workflow for sugar profile analysis

Why upgrade?

The Eurofins Nutrition Lab in Nantes was facing a number of challenges in its carbohydrate analysis, a prominent one being productivity. “Sugar profile analysis

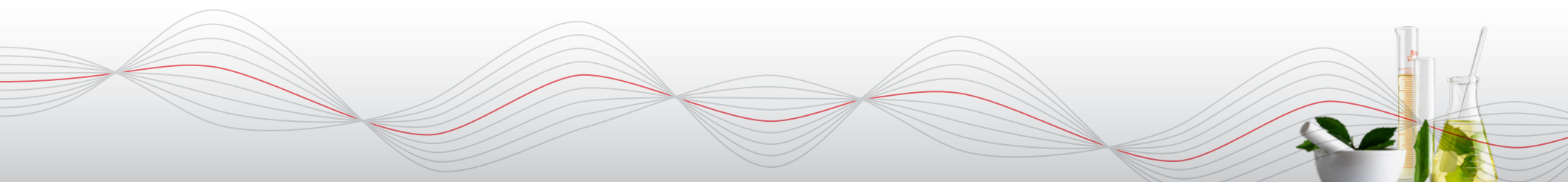
The Eurofins Nantes analytical campus hosts one of the largest nutritional analysis laboratories in Europe. To maintain their high testing standards, they sought a way to improve the efficiency, sensitivity, safety and capacity of their workflows, which rely upon advanced ion chromatography systems (IC/ICS). They opted to collaborate with Thermo Fisher Scientific and, in December 2017, deployed an optimized method of carbohydrate analysis, implemented eluent generation on their existing Thermo Scientific™ ICS-3000 and Thermo Scientific™ ICS-5000+ systems (which run on Thermo Scientific™ Chromleon™ Chromatography

Data System (CDS) software, and installed a new IC column: the Thermo Scientific™ Dionex™ CarboPac™ SA10. The laboratory has continued this collaboration since, installing additional eluent generation in 2019 and a Thermo Scientific™ ICS-6000 in 2020.

In this case study, we expand upon the lessons learned and benefits experienced by the Eurofins team in Nantes from the collaboration and demonstrate how advanced analytical equipment can benefit not only nutritional analysis, but Good Manufacturing Practice workflows for any product.

ThermoFisher SCIENTIFIC

[View the full case study](#)




Automated nutrient analysis and water quality monitoring

Overview

This analytical guide summarizes the use of Gallery discrete analyzers to detect the various nutrient and water analytes of interest, their respective testing chemistries, applicable regulatory method references, reagents required, sample matrixes, calibration curves, method detection limits, precision summaries, and method performance linearities.

The image shows the cover of an analytical guide. On the left is a grey sidebar with a white 'Table of Contents' section listing various analytes: Water analysis, Alkalinity, Ammonia, Calcium, Chloride, Chromium (VI), Fluoride, Iron (Ferrous), Magnesium, Nitrite, Phosphate, Silica, Sulfate, Total Hardness, Total Oxidized Nitrogen (TON-Hydraxine), Total Oxidized Nitrogen (TON-Vinadum), Total Oxidized Nitrogen (TON-Enzymatic), and Gallery analyzers. The main cover area features a red header with the 'thermoscientific' logo. Below the header is a photograph of a hand in a blue nitrile glove pouring liquid from a glass beaker into a larger glass container. The text 'Automated nutrient analysis and water quality monitoring' is overlaid on the bottom of the photo. Below the photo, it says 'Fast and accurate Thermo Scientific Gallery discrete industrial analyzers' and the 'ThermoFisher SCIENTIFIC' logo is in the bottom right corner.

 [View the analytical guide](#)

Discrete analyzer products

Thermo Scientific Gallery discrete analyzers with ready-to-use system reagents are optimized for speed, flexibility, and precision for wine, beer, malt, beverages, enzymes, soil, process water, ground water, wastewater, and drinking water analysis that enables improved quality control through consolidated testing.



Find out more at thermofisher.com/discreteanalysis

thermo scientific