

Customer-focused solutions

We've got your application covered!

Applications include:

- Fossil-Fuel, Steam-Generation Units
- Waste Incinerators
- Portland Cement Plants
- Petroleum Refineries
- Gas Turbines
- Basic Oxygen Furnaces
- Sewage-Treatment Plants
- Fertilizer Plants
- Ethanol Plants
- Pulp and Paper Mills
- Glass Manufacturing
- Grain Elevators
- Pharmaceutical Manufacturing
- Metal Smelters
- Chemical Plants
- Sulfur-Recovery Units
- Landfill Gas Combustions
- Air-Quality Networks
- Area/Fenceline Monitoring
- Munitions Disposal

Compliance

As your partner in compliance, our systems are designed with the necessary protocols to ensure your monitoring reports are accurate and accessible. Our system solutions can be designed to help you comply with U.S. EPA Part 60 or Part 75 performance standards, as well as other U.S and International ambient and source air monitoring standards.

Proven designs

We are dedicated to providing our customers with a cost-effective, high-performance solution that exceeds expectations and sets the standard for system instrumentation. Traditional, field-proven platforms, such as our iSeries instruments platform, are known in the industry for advanced electronics and ease of maintenance. Advanced platforms for ambient and particulate monitoring, like our 1405 Series of monitors and the new Partisol iSeries samplers, prove that traditional methodologies, combined with today's technology, can make monitoring easier than ever.

Technologies

Our global team of scientists investigates the latest technologies to drive product advancement and partner with all other internal teams to ensure that only quality products, with the best price and performance, are available to our customers. Utilizing our own proprietary, field-proven technologies, in addition to industry preferred methods, we offer innovative and customer-focused solutions.

Technologies include:

- Chemiluminescence
- Pulsed UV Fluorescence
- NDIR
- UV Absorption
- FID
- Cold-Vapor Atomic Fluorescence
- Beta Attenuation
- Light Scattering (Nephelometry)
- TEOM

Methods

We recognize that preferred methods for monitoring may vary by application or by geography. While a wet-basis dilution method may be particularly suitable for coal-fired utility monitoring, it is more common to use a dry-basis, full-extractive method for a gas-turbine application. We are proud to offer numerous monitoring methods, such as hot-wet or cold-dry full extractive, dry and wet-basis dilution, as well as continuous or manual particulate sampling, ensuring there is a Thermo Scientific monitoring method that's right for your application.

Methods include:

- Cold-dry Full Extractive (needs chiller)
- Wet-basis Dilution
- Dry-basis Dilution
- Continuous Sampling
- Manual Sampling

Configurations

In addition to offering traditional monitoring solutions, we also provide flexible system designs to meet your specific facility needs. Through our comprehensive portfolio of probe and Data Acquisition Handling Systems (DAHS) offerings, enhanced by our exclusive third party relationships, and a wide range of systems components, we are able to offer a superior choice of system options that are customizable to your application needs.

Our instruments and systems

Are designed for multiple applications, leveraging platform commonalities and providing field-proven performance.

For example, our Thermo Scientific Mercury Freedom System was specifically designed for the utility industry to comply with CAMR and led the way with more than 500 shipments of systems for installation in coal-fired utilities, using a simple design that combines performance with ease of use. Since then, we've successfully adapted the Mercury Freedom System for monitoring mercury in other applications, such as cement kilns, waste incinerators, munitions disposal plants and process control.

USA

27 Forge Parkway
Franklin, MA 02038
Ph: (508) 520-0430
Fax: (508) 520-2800
orders.aqi@thermofisher.com

India

C/327, TTC Industrial Area
MIDC Pawane
New Mumbai 400 705, India
Ph: +91 22 4157 8800
india@thermofisher.com

China

+Units 702-715, 7th Floor
Tower West, Yonghe
Beijing, China 100007
Ph: +86 10 84193588
info.eid.china@thermofisher.com

Europe

Ion Path, Road Three,
Winsford, Cheshire CW73GA
UK Ph: +44 1606 548700
Fax: +44 1606 548711
sales.epm.uk@thermofisher.com

Find out more at thermofisher.com/cems