

Mass spectrometry

Purpose-built performance for complex contaminant analysis

Orbitrap Exploris EFOX Mass Detector



thermo scientific

Built-for-purpose mass detection



The Orbitrap Exploris EFOX Mass Detector coupled to the Thermo Scientific™ Vanquish™ Flex Duo dual channel system delivers built-for-purpose mass monitoring.

This high-resolution accurate mass (HRAM) system is designed to satisfy market industry demands with the ability to perform targeted analysis and precise quantitation, suspect screening, as well as sample profiling.

Welcome to the Thermo Scientific™ Orbitrap Exploris™ EFOX (Environmental Food Organic Xenobiotics) Mass Detector, the perfect fit for environmental and food safety testing. Beyond the renowned high-resolution accurate mass detection provided by Thermo Scientific™ Orbitrap™ mass analyzers, this built-for-purpose system is simple to operate and compliance ready with Thermo Scientific™ Chromeleon™ software. HRAM data provides utmost confidence for enabling retrospective data analysis and the ability to easily add new compounds of interest without redeveloping methods and the need to reinject your sample, unlike conventional systems that require time-consuming MS/MS development and optimization, in addition to retention time adjustments.

Ease of use | Operational simplicity | Reliability

- **Simplicity for environmental and food safety laboratories**

The Chromeleon software offers built in methods, control, data processing and management that simplify tasks and reduce errors for more 'right first time' analyses. A suite of smart tools make work faster and easier, while ensuring reproducible, high-quality mass monitoring.

- **Productivity to accelerate commercialization**

High uptime boosts productivity. Rapid "set and forget" calibration procedures provide consistent mass stability for at least four weeks at prescribed conditions. Instrument status monitoring and optimized pressure control alert users when maintenance is needed, avoiding unnecessary downtime and repeat analyses.

- **Compliance for competitive advantage**

Designed for data integrity, data security, and compliance, fully scalable Chromeleon software gives you a competitive edge in meeting global regulatory requirements.

- **Consistent results deliver confidence**

Robust and reliable performance from system to system and from site to site ensure high-confidence results used to make critical decisions. Industry-proven Orbitrap technology, now purposely designed for mass monitoring in future-proof lab operations.

- **Complexities in environmental testing**

The Orbitrap Exploris EFOX Mass Detector simplifies environmental testing workflows, enabling faster and more reliable detection of a wide range of pollutants, including ubiquitous contaminants like per- and polyfluoroalkyl substances (PFAS). Its high-resolution capabilities and streamlined data processing reduce turnaround time and operational burden, helping labs achieve and maintain regulatory compliance adherence more efficiently.

- **Challenges in food safety testing**

With out-of-the-box workflows for key environmental and food safety applications, including PFAS and pesticide testing, the high-throughput Orbitrap Exploris EFOX Mass Detector is ready for compliance, supporting labs in meeting stringent regulatory requirements with confidence and efficiency.





Technical and online support: peak performance for your instruments

Helping you keep your instruments running at peak performance is our goal. Whether you're looking for an instrument manual or spare parts, want to submit a repair request, or check on the status of your warranty or service contract, we have every support option you're looking for.

thermofisher.com/technicalresources



Services and support

Unity™ Lab Services provides comprehensive service solutions for your instruments—everything from installation and maintenance to repair and compliance support. Keep focused on what matters. We've got your back.

thermofisher.com/unitylabservices



One online service platform that does it all

Managing service for your instruments and equipment can be time-consuming and frustrating. Now it doesn't have to be. You and your team can easily manage service for your Thermo Fisher Scientific assets on one convenient online platform.

Use Services Central to request service, view service request status and service history, check service plan or warranty coverage, access manuals, and more.*

thermofisher.com/servicescentral



Targeted MS quantitation—controlled and accessed from anywhere

Streamline targeted MS quantitation by combining Chromeleon software with Thermo Scientific™ MS systems for a superior solution. Chromeleon software offers data processing tailored for targeted MSⁿ quantitation and screening workflows.

thermofisher.com/chromeleonms

*Supported instruments and equipment vary by region. Visit thermofisher.com/servicescentral to view what is supported in your location.

 Learn more at thermofisher.com/efox

For Research Use Only. Not for use in diagnostic procedures. © 2025 Thermo Fisher Scientific Inc. All rights reserved.
 All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. BR-004182-EM-EN 1025

thermo scientific