# Complete Sample Preparation Workflow Solutions

#### Sample Preparation, Gas Chromatography-Mass Spectrometry, and Data Analysis

When performing analyses, the combination of Thermo Scientific Dionex sample preparation systems with Thermo Scientific GC-MS/MS products enable unstoppable productivity and higher efficiency; helping chemists save time, labor and money by simplifying routine workflows and providing rapid pathways from sample to results.



SCIENTIFIC

# Complete Sample Preparation Workflow Solutions

### **Sample Preparation**

The automated **Dionex ASE 150 and 350 Accelerated Solvent Extractor** systems enables extraction of solid and semisolid samples using organic solvents and water at elevated temperatures and pressures.

- Unattended, automated extraction of up to 24 samples (up to 100 mL); requiring less time and solvent, and reducing costs
- Performs automated extractions within 12 to 20 minutes

The automated **Dionex AutoTrace 280 SPE** instrument extracts large-volume samples (20 mL – 20 L) for the isolation of trace organics in water or aqueous matrices more quickly with better analyte recovery than manual liquid-liquid extraction (LLE) methods.

- Reduces solvent usage and time with the SPE technology
- Increase productivity or sample throughput with unattended operation

The **Rocket Evaporator** is a revolutionary solvent evaporator that concentrates or dries up to 18 ASE<sup>™</sup> tubes or 6 large-volume flasks unattended.

- Effective bumping and cross-contamination protection with precise sample temperature regulation
- Proven to yield fast, unattended operation that significantly improves laboratory productivity

For more information on these Sample Preparation products, visit thermoscientific.com/samplepreparation

#### **Separation and Detection**

The high-performance **TRACE 1300 Series GC** features a userexchangeable miniaturized, instant connect injectors and detectors that reduce maintenance downtime and enable the user to quickly tailor instrument capability to specific applications and daily workload.

- Instant connect, user-installable injectors and detectors
- Easy implementation of existing methods with ultimate productivity in routine analysis

For more information on the TRACE 1300 Series GC, visit thermoscientific.com/trace1300

The **Thermo Scientific<sup>™</sup> ISQ<sup>™</sup> Series** single quadrupole GC-MS system is an affordable and robust solution for dependable, routine GC-MS analyses that offers operational simplicity and unstoppable productivity.

- GC-MS venting not required to exchange the ion source, for continuous operations
- Inert sample path for highest chromatographic integrity

For more information on the ISQ Series GC-MS, visit thermoscientific.com/ISQSeries

The **TSQ 8000 Triple Quadrupole GC-MS/MS** is a reliable, easyto-use system that enables faster, more precise, error-free analyses, saving time and reducing laboratory costs.

- Maximum source robustness for high throughput analysis
- MS/MS simplicity for effortless method development and operation

For more information on the TSQ 8000 GC-MS/MS, visit thermoscientific.com/tsq8000

**TraceGOLD GC Columns** offer you a leap forward in column performance, delivering ultra low bleed, superior inertness, and a high level of reproducibility, guaranteed.

- High levels of reproducibility both run-to-run and column-tocolumn. You can expect consistent high-level performance from every column
- Superior inertness ensuring excellent peak shape and sensitivity, especially for highly active or difficult compounds

## **Data Analysis**

**Chromeleon 7.2 CDS** software is the first CDS that combines separation (GC/IC/LC) and MS in an enterprise (client/server) environment.

- One set of methods, one set of reports and one central location for chromatography and MS data
- Comprehensive suite of Chromeleon CDS productivity tools from dynamic data linking, instant processing, to SmartPeaks<sup>™</sup> detection - now available for MS

For more information on Chromeleon CDS, visit thermoscientific.com/chromeleon

**TraceFinder 3.0 Software** is an easy-to-use software for routine GC, GC-MS, LC, and LC-MS quantitation, and targeted screening, that increases productivity with powerful method development, simplified acquisition and comprehensive data review.

- Develops complex, processing method in minutes
- Runs the chromatography, processes the data, and produces final results

