

THE SMART BEHIND SEPARATIONS



GC 2400™ Platform with Detachable Touchscreen
Gas Chromatography

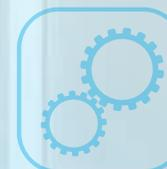
The New Way to GC

Every day, analytical laboratories are challenged with balancing high productivity with efficient operations. And with the growing trend of hybrid work models, gas chromatography operators seek remote access to their GC – whether they're working in a different part of the lab or at a completely different location.

Discover the **GC 2400™ Platform**, offering innovative technology that enables access to real-time information on the go. Efficiently monitor the status of your sample runs with a detachable, intuitively designed touchscreen, helping you make faster decisions from anywhere – in or out of the lab.

Maximize GC productivity and increase automation of lab operations with easy-to-use **SimplicityChrom™ Chromatography Data System (CDS) Software**. It enables the integration of every step in the GC workflow, from instrument control to data processing.

Experience innovative GC workflows with the **smart, simplified, and sustainable** GC 2400 Platform.



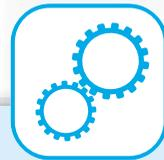
Smart. Simplified. Sustainable.

Bring value to your lab. Discover how the GC 2400 Platform helps your lab driving productivity with innovative user-experience.



SMART CONNECTIVITY

- Comprehensive and automated workflows maximize productivity
- Offers remote connectivity functionality supported by live system status
- SimplicityChrom CDS Software is designed for all experience levels to expedite results



SIMPLIFIED OPERATIONS

- Optimizes operations with smart sampling capabilities, status notifications, and easy-to-maintain hardware
- Software-driven tasks and automated, comprehensive workflows so labs can focus on results.



SUSTAINABLE SOLUTIONS

- Smart gas management to reduce gas consumption
- Alternative use of hydrogen as a carrier gas with built-in sensor enables lab and operator safety
- Supports sustainability throughout the growth of the lab with flexible and smart sampling capabilities – built to grow with your lab's needs

What Innovation Looks Like

Smart connectivity, simplified operations, and sustainable design come together to create complete, efficient workflow solutions for gas chromatography. The GC 2400 Platform is backed by a portfolio of separation chemistries and solutions that offer labs an efficient workflow experience – bringing the same smart technology we use in everyday life into the lab environment.

AS 2400™ Liquid Sampler

Plug-and-play functionality, driving uptime and growing with labs' productivity.

MS 2400™ SQ Detector

Dependability, efficiency, and intuitive user experience (UX) drives productivity.

GC 2400 System

Small footprint, advanced gas management, and information on-the-go.

HS 2400™ Headspace Sampler

Unparalleled precision with unique pressure balanced sampling technology.

Simplicity Vision™ App

Real-time information, direct operations, system updates, and status of analysis via the detachable touchscreen or any PC – no matter where you are.

SimplicityChrom Chromatography Data System (CDS) Software

Novel user experience with intuitive workflow and iconographic style.



Welcome to a New GC Experience

Our innovative **GC 2400 System** brings a new experience to GC operators. The touchscreen, complete with customizable LEDs and sounds, provides status updates and notifications that deliver crucial information at a glance – inside and outside the lab – helping you make better decisions, faster.

The automation capabilities are powered by easy-to-use SimplicityChrom CDS Software. Plus, with simplified maintenance features, the GC 2400 System supports lab productivity and increases instrument uptime.

With the Asset Genius™ solution for laboratory monitoring, you can return time to your scientists by automating the collection, visualization, and reporting of key performance data for insights into your most critical lab assets, including the GC 2400 Platform. And with 24/7 monitoring and alert notifications, you can have total confidence in your assets – and your lab environment.



GC 2400 System at a Glance

- Detachable touchscreen for faster decisions on the go
- Customizable LEDs and sounds for real-time status
- SimplicityChrom CDS Software control
- Small footprint
- Fast, efficient oven
- Easy maintenance
- Asset Genius

Sampling Productivity Maximized

To maximize productivity and profit, labs need autosamplers that enable high throughput while also delivering simplified operations.

With single- and dual-tower capabilities, the **AS 2400 Liquid Sampler** features plug-and-play towers that can be easily moved from one analytical channel to another, or from one GC to another, without the need for manual reconfigurations. Downtime is significantly minimized, enhancing your lab's overall output and ROI. It's the perfect balance between investment and profit, making your lab economically sustainable.

AS 2400 Liquid Sampler at a Glance

- Single and dual tower configuration
- Plug-and-play towers for easy reconfigurations
- Multiple tray capacities



Engineered for Perfection

Boasting unparalleled sensitivity, pinpoint precision, and ultimate ease-of-use, the robust **HS 2400 Headspace Sampler** features unique pressure-balanced sampling technology that introduces samples into the column without the use of a gas syringe or valve and loop.

By eliminating the valve and loop, many sources of variability and contamination are decreased, allowing the HS 2400 Headspace Sampler to avoid dead volumes, carryover, and peak distortion. That leads to sharper peaks, exceptional reproducibility, and ensures more accurate results.

HS 2400 System at a Glance

- Detachable touchscreen for faster decisions on the go
- Pressure-balanced technology for fast GC method development and exceptional reproducibility
- Supports constant pressure and constant flow modes
- Intuitive design and ease of use
- Injection volume adjustment for rapid method development
- Integrated into the GC workflow
- Optional sensor for safe use of H₂ as a carrier gas



Increased Productivity and Confidence in Your Results

Fully integrated into the GC 2400 Platform, the **MS 2400 SQ Detector** is a powerhouse. Offering accurate detections in full-scan mass spectra, selected ion full ion (SIFI), and selected ion monitoring (SIM) modes, our single-quad system gives your lab the flexibility it needs to maximize sample throughput and operate smoothly and efficiently. The exclusive Marathon™ Filament technology boasts exceptional durability and extended lifetime.

SMARTsource™ Technology provides tool-free source access, cleaning, and maintenance – giving your lab greater instrument uptime and sample throughput.

MS 2400 SQ Detector at a Glance

- Maximize productivity with full-scan mass spectra, SIFI, and SIM modes
- SMARTsource Technology (Simplified Maintenance and Removal Technology)
- Marathon Filament – for ultimate robustness



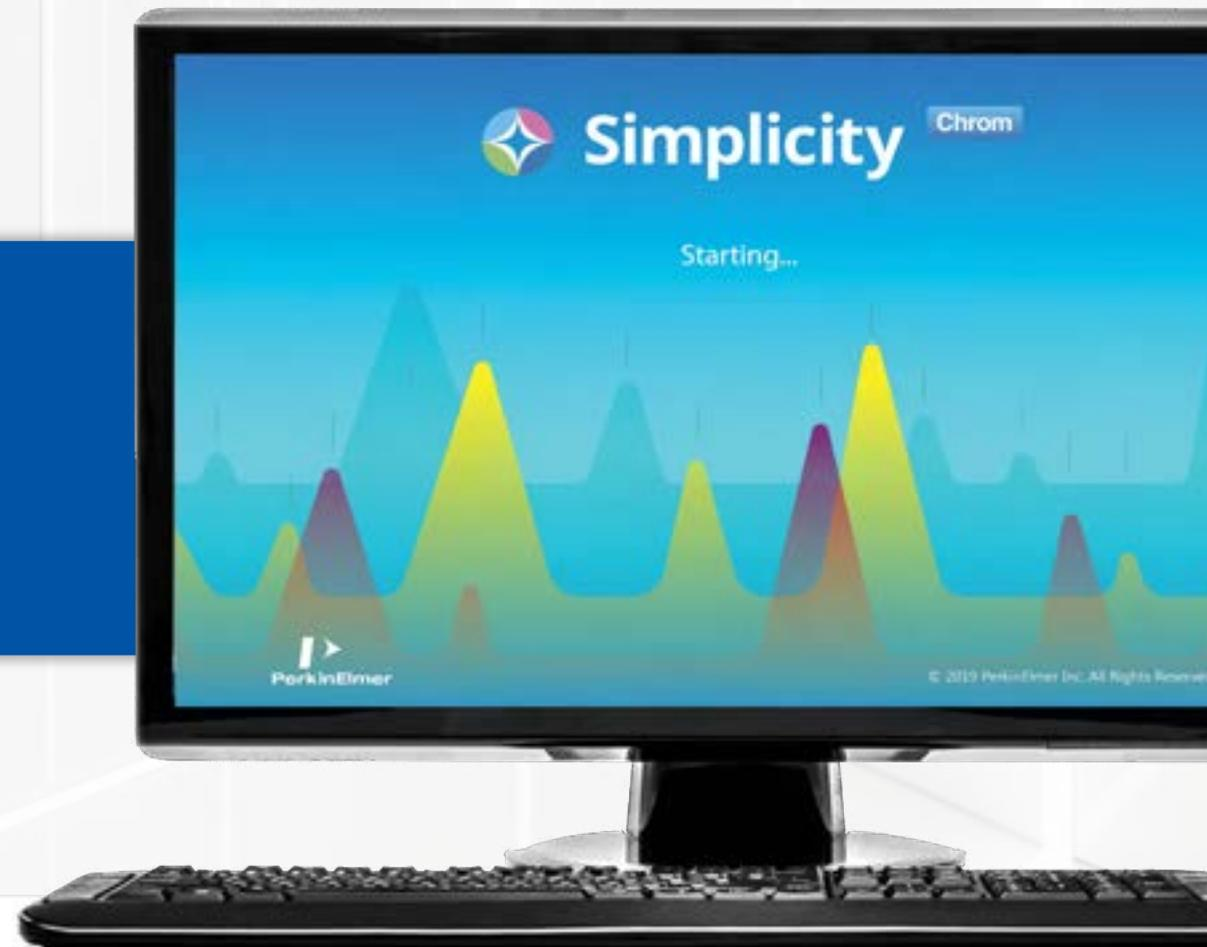
Designed with You in Mind

The GC 2400 Platform's smart features, automated processes, and flexibility throughout the day is driven by **SimplicityChrom CDS Software**.

SimplicityChrom CDS Software truly connects and integrates all steps of the GC workflow in a way that simplifies the processes. It's intuitive and flexible, enabling you to achieve your analytical goals and get results quickly. For labs that test highly regulated samples such as pharmaceuticals and food, SimplicityChrom CDS Software supports compliance with 21 CFR Part 11.

SimplicityChrom CDS Software at a Glance

- Modern design, user-friendly iconography
- Easy-to-learn workflows for any skill level
- User account and profile management
- Powerful search and filter features
- Supports and facilitates 21 CFR Part 11 compliance



Visualize Better Efficiency

Simplicity Vision is a browser interface that provides flexible access to the GC 2400 Platform. The technology allows for real-time status updates and monitoring of every step of the GC workflow as well as instrument setup and prerun manual operations. Used as a single or multi-instrument interface, operators can benefit from working remotely and making decisions on the go.

Simplicity Vision App at a Glance

- Assess instrument readiness status
- Check GC methods and sequences status
- Direct instrument operations
- Browse consumables and accessories



Complete, Integrated Workflows

The GC 2400 Platform can be used in many different analytical sectors and applications, from industrial and environmental analysis to food and pharmaceutical quality and safety testing. The platform offers the power of connectivity and workflow integration throughout the entire process.



OneSource
Laboratory Services Support, from beginning to end.

Complete Workflow Solutions



For additional information on workflows, visit perkinelmer.com/GC2400workflows

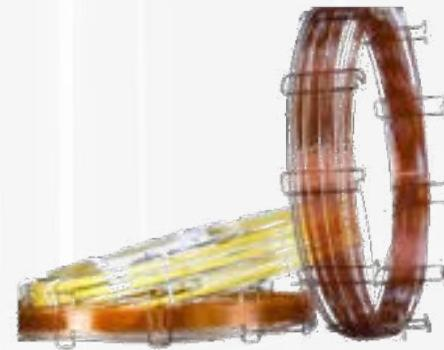
The Right Consumables for Your Analysis

The innovative design and smart features of the GC 2400 Platform make it easier than ever for cutting-edge labs to access, operate, and monitor their gas chromatographic separations – separations made from the reliable, reproducible, and rugged columns that are the heart of any GC instrument.

Whether your analytes are volatile or semivolatile, polar or nonpolar, at high concentrations or at ultratrace levels, we'll help you select the best column to separate and accurately quantitate your sample – injection to injection, column to column. Plus, we offer a variety of stationary phases and column dimensions to fit your unique analytical needs.

Starting with our most popular GC column part numbers, we've invested in original in-house column manufacturing capabilities to bring you a new level of reliability and performance. Each new GC 2400 is shipped with a PerkinElmer-made column to make sure your instrument – and your analysis – start off right at the very beginning.

But we don't stop at columns. We offer a complete line of vital consumables to support your analytical workflow. Each consumable is carefully chosen and engineered to give you the best results.



Consumables at a Glance

- A complete line of GC columns for a wide range of applications
- Instrument consumables specifically engineered for optimal performance
- Consumables and accessories for preventative instrument maintenance
- Consumables to support the entire workflow, from sample preparation to detection



At Your Service

A true analytical workflow cannot exist without dedicated support – every step of the way. We're here to help maximize your GC workflow, from beginning to end.



Trust in Our Expertise

Instrument Services

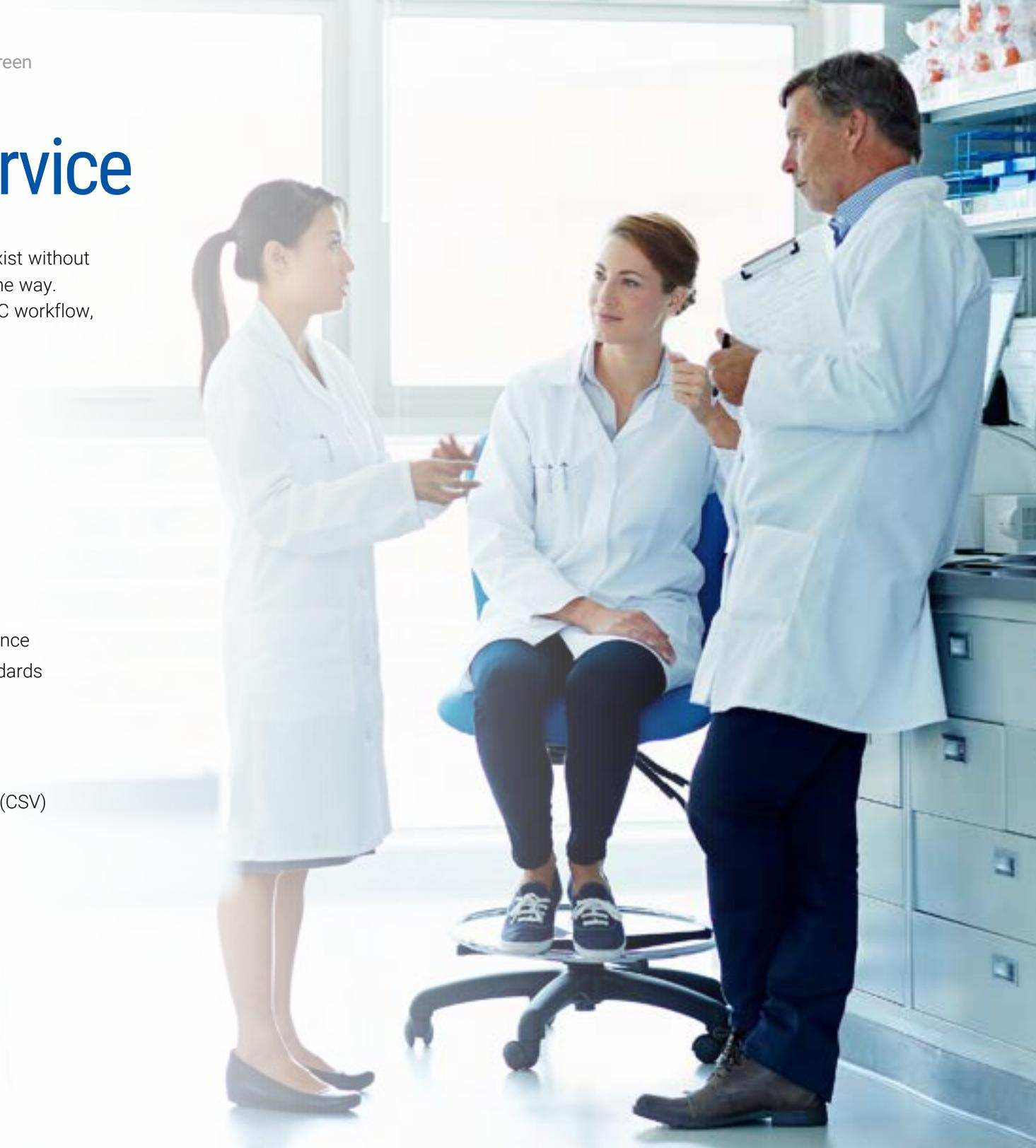
- Optimize for uptime
- Enhance productivity
- Routine and scheduled maintenance
- Compliance with regulatory standards

Compliance Services

- Instrument qualification
- Data integrity
- Computerized system validation (CSV)

Education Services

- Classroom and online training
- Basic on-site training
- Advanced on-site training



[For more information, visit perkinelmer.com/GC2400](https://www.perkinelmer.com/GC2400)

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

581533

PKI