

HUMAN HEALTH

ENVIRONMENTAL HEALTH

OneSource
Laboratory Services

IMPROVE YOUR
WORKFLOW WITH
CDS SERVICES

OneSource
Chromatography Data System Services


PerkinElmer
For the Better



PerkinElmer’s OneSource Chromatography Data System (CDS) services span the entire software lifecycle from requirements analysis to production support. Our CDS service offerings are focused on the strategic deployment of software applications, developing solutions to enable optimal use, and achieving and maintaining a solid, validated state while minimizing operational downtime.

ENTERPRISE CDS DEPLOYMENT

As a recognized leader in chromatography solutions, PerkinElmer® can leverage its practical experience in the deployment of a CDS as well as its experience and knowledge of other informatics solutions such as Laboratory Information Management Systems (LIMS), Electronic Laboratory Notebooks (ELNs), Laboratory Execution Systems (LES), and Chemistry and Biology Informatics. This deep expertise enables us to develop a CDS strategy that is appropriate for your current landscape and aligns with your long-term informatics vision. We know that the deployment of an enterprise CDS requires an effective planning phase that yields a roadmap that can be executed on-time, in scope, and on budget with minimal disruption to your laboratory operations. PerkinElmer possesses strong knowledge of laboratory workflows and operation as well as the demands of many different laboratory types and associated regulations. Through our long history with chromatographic solutions, we also possess the expertise for the successful integration of laboratory instrumentation with a CDS. PerkinElmer will inject learned best practices to yield an effective strategy and plan so as to ensure successful implementation as required by the client’s policy. Our planning phase deliverables include:

Project Plan	✓
User Requirements Specification	✓
Business Process Diagrams	✓
Technology Selection Evaluation Matrix	✓
Functional Requirements Specification	✓
Design Specification	✓
Quality Assurance Plan	✓
Data Strategy	✓
Systems Integration Strategy	✓

CDS CONFIGURATION

Chromatography represents over 90% of the data generated by analytical laboratories, including highly regulated quality control laboratories. As a result, the chromatography workflow, including the use of the CDS, warrants critical analysis for effectiveness, efficiency, and compliance gaps. In addition, PerkinElmer’s experience is that the CDS requires focused configuration to yield optimal laboratory workflow. As an example, many laboratories rely on standalone business tools such as spreadsheets to generate intermediate or final chromatographic results and word processing software to generate reports. Additionally, laboratories often do not leverage electronic signatures to realize a paperless solution. This omission creates the need for a hybrid process marred by inefficiencies, errors and expensive paper retention costs. PerkinElmer leverages our strong knowledge of laboratory workflow analysis and CDS technical expertise to develop configuration solutions that minimize this risk. These configurations include:

Custom Fields to Capture Metadata	✓
Custom Calculations	✓
Custom Reports	✓
Electronic Signature Configuration	✓
Methods Standardization	✓
Integration Configuration	✓

CDS VALIDATION AND CHANGE MANAGEMENT

Chromatography, more than any other technique, yields data that establishes and supports product integrity. The CDS requires a solid validation and change management program that can withstand regulatory scrutiny. PerkinElmer leverages expertise in chromatographic solutions, instrument qualification, computer system validation, and the CDS to contribute in all phases of validation from the development of the validation strategy to the development of the summary report to withstand this scrutiny. PerkinElmer has experience with the risk-based approach to validation and will leverage already established vendor qualification as part of the strategy. PerkinElmer will execute validation as required by the client's policy while injecting learned best practices.

Change management is a critical component of maintaining the validated state of the CDS. Changes include CDS configuration changes as well as application patches and changes to supporting systems such as the operating system. PerkinElmer possesses the validation and technical knowledge to develop a CDS specific change management program that is consistent with the client change control policy. The change management program may include:

Change Management Plan	✓
Change Control Procedures	✓
Updates to Configuration Specification	✓
Asset Inventory	✓

CDS APPLICATION MANAGEMENT

The support and management of the CDS requires a unique blend of skills – domain specific scientific, compliance expertise and laboratory knowledge coupled with the inherent technical capability in the CDS. PerkinElmer offers CDS management solutions to perform production support and implement application enhancements while maintaining the validated state.

Production support is provided using a structured support process driven by accountability against a defined Service Level Agreement (SLA) using scalable onsite, onshore, near-shore and offshore delivery options. Support responsibilities may include:

Application Administration	✓
End-User Problem Resolution	✓
Data Management	✓
Instrument Connection / Removal	✓
Upgrade / Patch Management	✓
Training	✓

BENEFITS

PerkinElmer's Chromatography Data System (CDS) services are an integral part of our OneSource® Laboratory Services program and extend our ability to enable laboratories to concentrate on science. Benefits of PerkinElmer's CDS service offerings include:

- Reliable delivery through program and project management capabilities
- Complete system lifecycle capabilities
- CDS expertise
- Chromatography and laboratory domain knowledge
- Regulatory and compliance knowledge
- Global reach
- Onsite, onshore, near-shore and offshore delivery options

Application enhancements deliver functionality or process improvements that can increase the effectiveness of the scientist and/or increase compliance. Enhancements can be directed at the laboratory workflow or supporting processes. This may include:

Custom Calculations to Eliminate Use of Off-Line Tools	✓
Custom Reports to Facilitate Data Review and Decision Making	✓
Streamlined Validation Processes for Common Changes	✓
Configuration of Part 11 Application Settings	✓
Synchronization of Laboratory CDS Workflows	✓

OneSource is the Only One You Need

From preventative maintenance and repair to qualification, calibration and laboratory relocation, and Lab IT, we're the only source – the ONE source – you need to help optimize your operations and cost-effectively manage laboratory assets throughout their entire lifecycle.

We pride ourselves on offering a complete portfolio of laboratory services. So, no matter what the name on the front of an instrument and no matter what the technology inside, we have the knowledge and expertise to take care of it.

OneSource Services

- Instrument Service and Repair
- Qualification and Validation
- Analytical Method Services
- Asset Procurement & Disposition
- Laboratory Relocation
- Lab IT
- Business Intelligence Solutions
- Chromatography Data Systems Services

To learn more about Lab IT and additional services available from OneSource Laboratory Services, please visit www.perkinelmer.com/labIT or contact your PerkinElmer OneSource Representative.

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

OneSource

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 ©2012-2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.