

# The iLAB™ Laboratory Execution System (LES)

## Efficient, accurate and compliant laboratory execution

The PerkinElmer iLAB™ Laboratory Execution System (LES) module provides QA/QC laboratories with a structured platform that reduces dependence on paper based processes while automating and controlling testing procedures.

The iLAB™ LES module enables QA/QC laboratories to:

- Reduce risk by creating a right first time environment for procedural execution.
- Improve efficiency and productivity by reducing workload for lab personnel and management.
- Use industry leading technology to reduce the time for deployment and achieve benefits sooner.
- Ease the burden of regulatory compliance.

### Get it Right the First Time

The iLAB™ LES module creates a right-first-time environment - at the bench level, as tests are being executed – that eliminates many of the opportunities for error that are inherent in manual systems.

- Testing procedures are controlled on a step-by-step basis to ensure compliance with SOP conditions and requirements.
- Data capture, calculations and status checks on reagents, instruments and training records can be automated to reduce errors.
- Product and result specifications are enforced at the time of analysis so analysts can immediately take appropriate action when results are out of specification.

Creating a right-first-time environment helps QA/QC laboratories avoid situations that can lead to costly and time consuming rework or in worst cases release of products or raw materials that don't meet corporate or brand standards.

### Work More Efficiently

The iLAB™ LES module frees analysts from many of the manual tasks that need to be carried out before, during and after test execution. Instrument logs and inventory records can be automatically accessed and updated so analysts spend less time looking for information and manually updating systems. Automatic data collection, calculations and reporting of results to LIMS and other systems help to further reduce analyst's workload. Analysts are able to work more efficiently without compromising the quality of their work.

**3 Preparation of HPLC Reference Standard**

**3.1 HPLC Standard 1 Overview:**  
This is a standard stock solution containing 100 mg/L of XYZ. The standard stock solution will be diluted with ultrapure water and stored in the dark between runs, at +4°C.

**3.2 Weighing Steps:**  
Weigh out 10 mg of XYZ.

Standard ID: 1000036

Standard Status: Active

Standard Expiry Date: 2012-11-23 00:00:00

Balance ID: Balance 1234

Balance Status: Calibrated

Standard Weight: mg

Standard Amt Remaining: 100.00 mg

Figure 1: iLAB LES worksheets provide step by step automation and control over testing procedures and generation of testing documents.

Laboratory management can have increased confidence in the quality of test execution allowing them to spend less time monitoring and reviewing day to day activities. This enables them to focus more time and attention on reviewing and analyzing laboratory operations in order to achieve a higher level of efficiency and productivity.

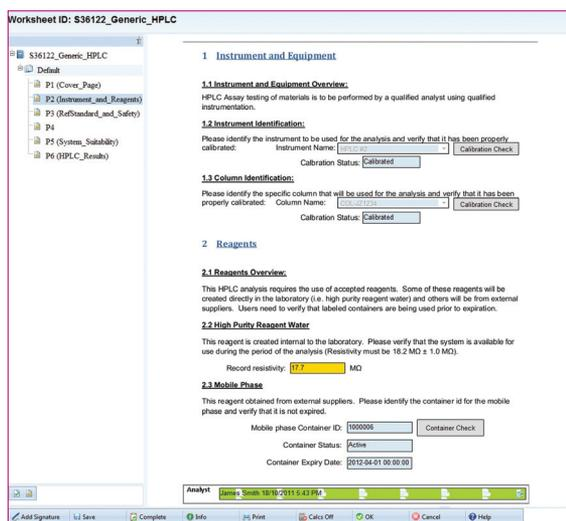


Figure 2: Quickly convert paper forms, lab notebooks, log books, sign out sheets etc. into electronic worksheets.

## Leading Technology for Rapid Deployment

One of the roadblocks to deployment of an LES can be the process of translating laboratory workflows and processes from paper into the electronic environment.

The iLAB™ LES module utilizes industry leading ReDI™ technology to overcome this challenge by quickly converting paper forms, lab notebooks, log books, sign out sheets, etc. into electronic worksheets.

Right away, analysts and technicians are able to work with electronic worksheets that provide them with a familiar and comfortable interface.

- Processes don't need to change in order to conform to the software.
- Training requirements are reduced, system acceptance is increased.
- Advanced automation can be quickly added to worksheets using simple 'drag and drop' ReDI™ technology.

For more information, please visit [www.perkinelmer.com/informatics](http://www.perkinelmer.com/informatics) or call 1-800-315-7300

PerkinElmer  
100 CambridgePark Drive  
Cambridge, MA 02140  
P: (800) 315-7300 or  
(+1) 617-588-9300  
[www.perkinelmer.com/informatics](http://www.perkinelmer.com/informatics)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2012 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. The iLAB™ platform application software and the iLAB™ Laboratory Execution System software module are provided by CambridgeSoft Corporation, a wholly owned subsidiary of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

## Ease the Burden of Regulatory and GxP Compliance

A Laboratory Execution System can help laboratories reduce the cost and effort of meeting regulatory and GxP compliance requirements.

Electronic worksheets, data files, methods and processes are automatically archived to a secure server reducing or eliminating the risk of lost or damaged paper forms or notebooks and reducing the costs of long term paper storage and retrieval. Electronic documents can be quickly retrieved when needed for an audit or problem investigation.

In addition to controlling procedural execution the security features, audit trail and electronic signature challenges in the iLAB™ LES module also control access to the application and document analyst activity during test execution.

## A Standardized Informatics Platform for QA/QC

The iLAB™ LES module is part of a suite of integrated modules for the iLAB™ platform, a highly configurable solution for QA/QC laboratories that can be deployed and supported with minimal impact on IT operations. The iLAB™ platform enables organizations to:

- Quickly meet end user needs with configurable forms and templates.
- Utilize industry leading technology to simplify integration with instruments and with other informatics systems.
- Leverage PerkinElmer Professional Services for implementation and support assistance.