

HUMAN HEALTH

ENVIRONMENTAL HEALTH

# RUNNING A LAB CAN BE A WALK IN THE PARK

## AxION eDoor Open-access software for streamlined operations.

From running an individual sample to managing a complete network of instruments and users, AxION® eDoor™ has the power to simplify every aspect of your laboratory operation. A robust yet intuitive open-access software, eDoor works with the Flexar™ SQ 300 MS to deliver reliable, high quality MS results no matter how complex your samples or workflows.

### Greater Simplicity. Greater Control.

With real-time, web-based access and live updates, AxION eDoor allows lab managers to control every aspect of their operation—from sample submission to final results—more efficiently than ever. Data fields and final reports can be customized and tailored to suit a particular application or user group. The software also simplifies the lives of scientists, making it easy to log in samples and receive results anywhere, anytime.

#### AxION eDoor at a Glance.

- Open-access system and intuitive interface for effortless operation.
- Rapid, first-available instrument assignment.
- Simple web-based sample submission process and easy-to-use, walk-up, touch screen sample drop off.
- Fast, convenient access to results via web or email on any PDA.
- Predefined methods and reporting enable any user to access mass spec data.
- Compatibility with workflows in all types of laboratories.



A modern, tile-based interface quickly and easily guides users through each step of an analysis and gives managers instant access to all the tools and information they need.

## Take Control of Your Instruments. Take Control of Your Lab.

Managing analytical instrumentation in a multi-user environment is no longer a complicated, time-consuming task. AxION eDoor's robust features and remote access capabilities enable lab managers to perform a variety of functions for unprecedented control and simplicity:

- Define instrument acquisition parameters based on specific methods.
- Assign an analysis to a particular instrument, group or person to optimize productivity and results.
- View sample queues on instruments to manage lab activity.
- Monitor group/individual activity and track instrument usage for effective preventive maintenance scheduling.
- Archive results and reports with assigned access privileges for specific users or groups.

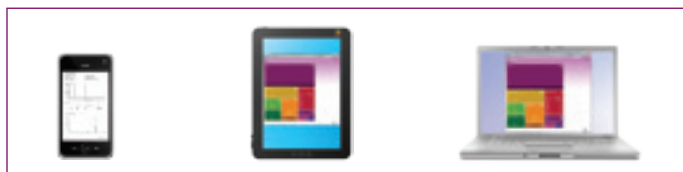
## Let Your Scientists Focus on Science.

With its simple sample-submission process and remote access capabilities, AxION eDoor minimizes time away from the bench for scientists, optimizing laboratory efficiency and productivity.

Step 1: Log in sample from any web-enabled device.

Step 2: Place sample on assigned instrument.

Step 3: Retrieve results via web or email.



## A Better Solution for Everyone on Every Level.

Discover how a real-time software can offer real-world benefits throughout the lab. Simplify every analysis for your chemists. Optimize the efficiency of your workflows. Streamline operations for your business.

AxION eDoor open-access software. Let yourself in on a whole new level of laboratory insight, control and productivity.

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

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. All other trademarks are the property of their respective owners.

010119\_01

Printed in USA

## Let Yourself in on the Power of MS.

No matter what your application or sample type, AxION eDoor makes it easy to obtain complete sample insight through mass spectrometry. Couple the software with the advanced Flexar SQ 300 MS and you can take advantage of:

- Grounded ESI and Field-Free APCI sources to analyze any compound.
- Capillary exit CID for detailed structural information.
- Multi-stage ion path ensures superior transmission for exceptional sensitivity.
- Fast MS detection capabilities (ideal for HPLC and UHPLC applications).
- Easy cleaning and maintenance—no tools required.
- Compatible with any PerkinElmer LC or Agilent LC running ChemStation® under Windows XP®.
- Reliable long-term performance.

The Flexar SQ 300 MS with AxION eDoor software—identifying, quantifying and confirming compounds has never been done with such ease, speed and confidence.



PerkinElmer's unique AxION Separation Probe can be attached to the interchangeable, grounded ionization probes, allowing chromatography to be performed directly on the source for minimized dead volumes.

The innovative design of the snap-in ionization probes allows them to be dedicated to specific users or applications in an open-access environment, minimizing cross contamination while delivering greater control, speed and productivity.

  
**PerkinElmer®**