

BRILLIANT
PERFORMANCE
FOR ROUTINE
QUANTIFICATION



QSight® 210 MD Screening System

The complete solution for high throughput routine quantification

QSight® 210 MD Screening may not be available in all countries. Please check out availability with your local PerkinElmer representative


PerkinElmer
For the Better

WHEN
MAXIMUM
PERFORMANCE
AND MINIMUM
DOWNTIME
MATTER



RUN THOUSANDS
OF SAMPLES PER DAY



The QM1000® 210 MD Screening system is designed for routine high throughput quantitation. With easy to use instrument control software, the ability to automatically download sample and control information, a mass spectrometer designed to reduce maintenance and repair and data analysis software designed to make it quick and easy to review results, the QM1000® 210 MD Screening system makes it easy to run thousand samples per day.

WHAT MAKES THE QM1000® 210 MD SCREENING SYSTEM SO EXCEPTIONAL?

- ✓ A simple three step process to set up and run a method
- ✓ Ability to download worklists, internal standard & calibration information directly into the software to eliminate transcription errors
- ✓ Remote instrument diagnostics software to quickly identify problems before they occur
- ✓ Unique StayClean™ source design to eliminate daily maintenance
- ✓ Unique Laminar Flow Ion Guide™ to reduce the need for re-tuning
- ✓ Easy-to-use, visual based database software to simplify data review and analysis
- ✓ Unique Unifield Detector™ design to detect more analytes within a run
- ✓ A modular mass spectrometer design to simplify maintenance and repair

DESIGNED TO MAXIMIZE PERFORMANCE AND THROUGHPUT

Higher Sensitivity and Productivity

Unique StayClean™ technology employs hot surface induced solvation (HSID™), and an orthogonal sampling interface to increase uptime. A continuous flow of hot gas is used to de-solvate and guide ions of interest into the instrument to increase signal and reduce chemical noise while forming a barrier that prevents unwanted analytes from dirtying the source.

Higher Performance

Ions are transferred from the HSID™ Interface into the system's Laminar Flow Ion Guide™ and through to the analyzer by a flow of background gas, without the need for an axial electrical field or the associated field fluctuations, delivering more consistently higher levels of performance with reduced drift or the need for frequent re-optimization.

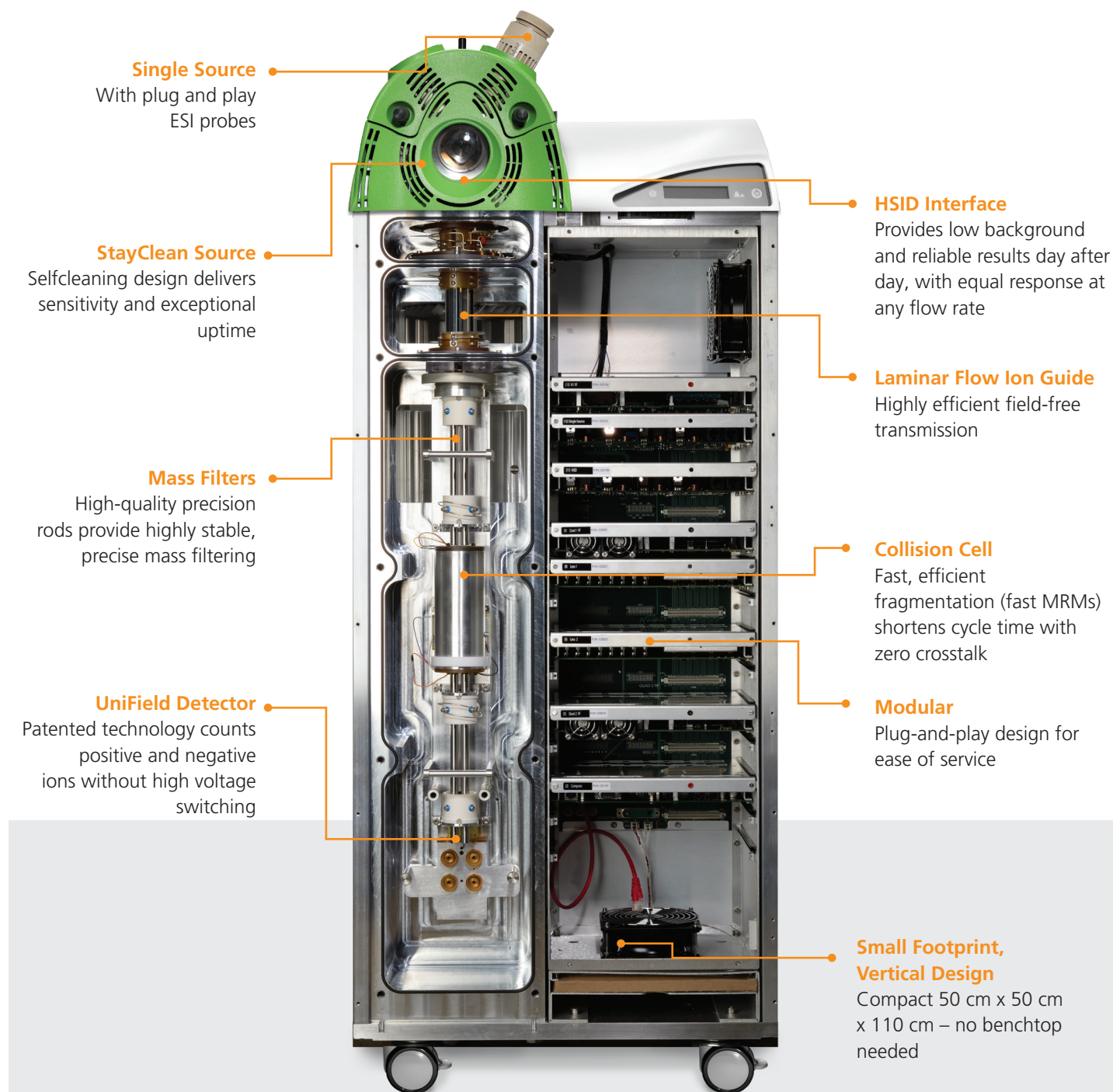
More analytes per run

With its high energy dynode to attract positive ions, the QSight system's Unifield Detector™ causes positive ions to collide with the dynode to form electrons that cascade into the detector while negative ions are detected as usual. This provides nearly simultaneous detection of positive and negative ions without the need for high voltage switching. Polarity switching happens in microseconds, limited only by the ion source and path, making it easier to detect a wider range of analytes in a single run.



IT'S SMALL, IT'S SIMPLE, YET SENSITIVE AND HIGHLY SPECIFIC

HIGH THROUGHPUT LC TECHNOLOGY DESIGNED TO DELIVER CONSISTENT RESULTS



Single Source

With plug and play
ESI probes

StayClean Source

Selfcleaning design delivers
sensitivity and exceptional
uptime

Mass Filters

High-quality precision
rods provide highly stable,
precise mass filtering

UniField Detector

Patented technology counts
positive and negative
ions without high voltage
switching

HSID Interface

Provides low background
and reliable results day after
day, with equal response at
any flow rate

Laminar Flow Ion Guide

Highly efficient field-free
transmission

Collision Cell

Fast, efficient
fragmentation (fast MRMs)
shortens cycle time with
zero crosstalk

Modular

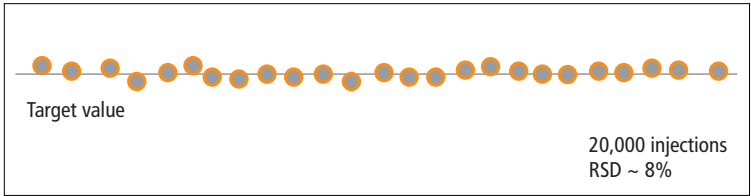
Plug-and-play design for
ease of service

Small Footprint, Vertical Design

Compact 50 cm x 50 cm
x 110 cm – no benchtop
needed

PROVEN INSTRUMENT RELIABILITY

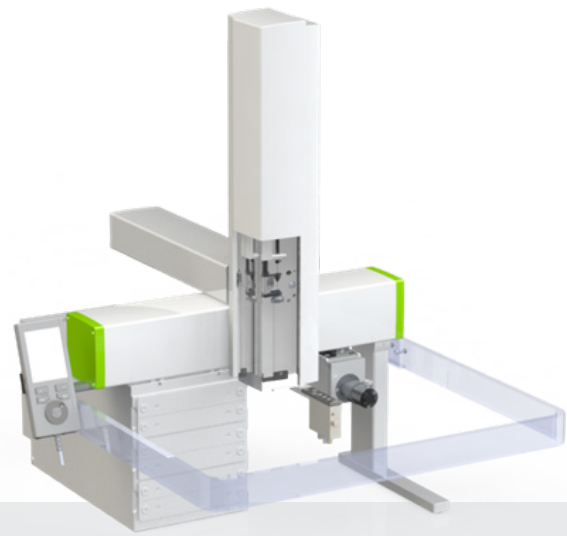
Recognizing that consistent results are essential to proper data interpretation, the QSight™ system, with unique HSID™ interface, Laminar Flow Ion Guide™ and Unified Detector™, have been designed to make it easier to achieve consistent, reproducible results across one or multiple instruments, week-in and week-out, without the need for frequent cleaning or re-optimization.



FOR EXCEPTIONAL SENSITIVITY, THROUGHPUT AND PRODUCTIVITY

MORE RESULTS WITH BETTER ACCURACY IN LESS TIME

The QSight system includes a high capacity autosampler, precision binary pump, best in class MS detector, and easy to use instrument control and data analysis software that have been optimized to run hundreds of samples per day, for weeks at a time, to generate consistent results with minimum maintenance or downtime.



QSIGHT® HC AUTOSAMPLER MD

- Sample Injector Module with loop injector design
- Injection volume up to 100 µl
- Holds 12 x 96 or 384 position well plates
- Foil cutter to prevent clogging of the needle
- Sample compartment settable from 4 to 40 °C
- Hand-held control module for manual operation



QSIGHT® BINARY PUMP MD

- Run up to two solvents in any proportion
- Flow range from 10 to 5000 µl/min
- Operation from 0 – 6,000 psi operating pressure
- Re-settable log counter to automatically record pump seal usage and operation time



SIMPLICITY™ 3Q MD SOFTWARE

- Control of all instrument components
- Method set-up
- Data acquisition, calculations and flagging
- Sample based result processing
- Instrument calibration and tuning

QSIGHT® 210 MD

- 5 – 1,500 m/z Mass Range
- 27,000 Da/Sec Scan Speed
- Mass drift < .05 amu / 24 hrs
- Polarity Switching < 10 ms
- Minimum dwell time 1 ms



OPTIONAL: MSMS WORKSTATION SOFTWARE*

- Stand alone IVD software for result handling
- Capture worklist directly via interfaced laboratory equipment
- Import ISTD info from barcoded source document
- Consolidate data generated from one or more instruments
- Program thresholds, cut-offs by sample type
- Colored coded results view
- Flag based upon results vs user defined cut-offs
- View sample results & spectra
- View results grouped by disorder or other formats

* MSMS workstation software may not be licenced in accordance with the laws in all countries, such as the United States and Canada. Please check with your local representative for availability



SERVICE AND SUPPORT

- Standard one year warranty including:
 - Twice yearly preventative maintenance
 - Phone support
 - Remote diagnostics
 - 48 hours on-site response
 - Firmware updates
 - Parts, travel and labor

(Service options may vary by location)

DELIVERING SOFTWARE THAT IS SIMPLICITY ITSELF

SIMPLICITY™ 3Q MD SOFTWARE was designed with ease of use and productivity in mind. With a graphical user interface and guided three step workflow, the software makes it faster and easier to get up and running generating results. The software can also automatically capture internal standard and calibration target values used during a run, eliminating manual transcription errors. Capable of real-time acquisition of thousands of MRMs per run, with results shown in real time, the software provides the ability to monitor data acquisition as it occurs as well as review results after each run, making it easier to check instrument performance prior to leaving the instrument to run overnight or access priority results.

OPTIONAL: MSMS WORKSTATION SOFTWARE* collects data from one or more instruments into a database that can format flagged concentration or enzyme activity based results into a well plate map, grid, disorder based groupings or other user definable formats to quickly and easily process hundreds of samples. The software can also be used to perform quality control trend analysis to monitor instrument performance.

* MSMS workstation software may not be licenced in accordance with the laws in all countries, such as the United States and Canada. Please check with your local representative for availability

3 STEP WORKFLOW

STEP 1: DOWNLOAD SAMPLE WORKLIST

Download sample worklist from a dried blood spot puncher, plate map or other software

STEP 2: CHECK INSTRUMENT STATUS

Check the color coded instrument status bar to confirm all components are ready

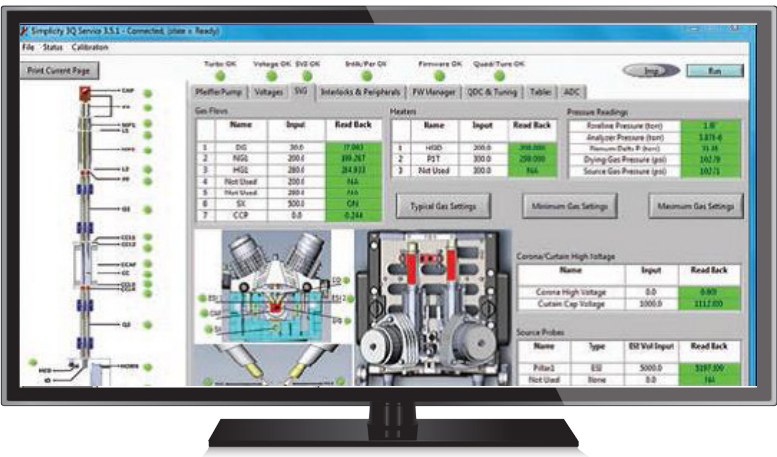
STEP 3: START A RUN

Click to start acquiring data



A COMMUNITY OF SUPPORT

PerkinElmer's global team of clinical Diagnostic service and support engineers are dedicated to provide the training and support required to get customers up and running as quickly as possible as well as providing the maintenance and repair services required to ensure maximum productivity. The Simplicity™ 3Q Service Module supports this effort by allowing service engineers to login to a system remotely to review instrument status real-time in order to quickly identify potential issues and initiate repairs to reduce service costs and maximize instrument uptime.



QSite® 210 MD Screening may not be available in all countries. Please check out availability with your local PerkinElmer representative

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 ©2017, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

1599-9816-2

BRO QSite 210 MD Screening system , August 2017