

The NBS solution
that's as unique as
they are.



revvity

MSMS Newborn Screening Solutions

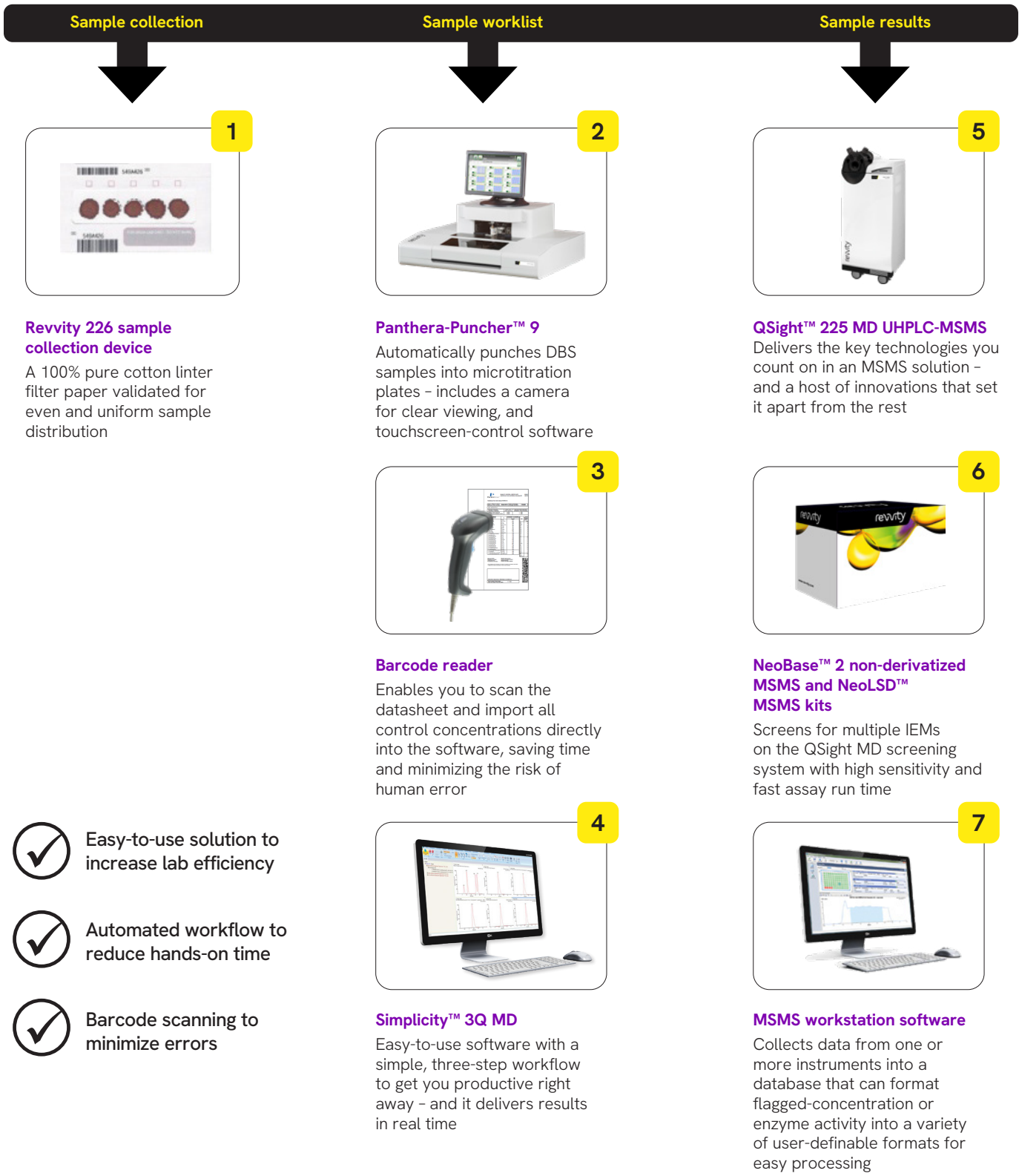
Everything comes together to deliver more screening

Early detection of metabolic and other inherited and rare disorders can improve the quality of life for at-risk babies and their families. With the **QSight™ 225 MD UHPLC-MSMS**, we're enabling expanded high-throughput screening for a variety of inborn errors of metabolism (IEMs) with a wide analyte panel, and offering access to tier 2 chromatography.

Our solution delivers a unique combination of self-cleaning MSMS instrumentation that provides exceptional system uptime, a simple three-step method setup, and easy-to-use database software that simplifies review and analysis.



A fluent workflow, from sample to results



Designed for your workflow

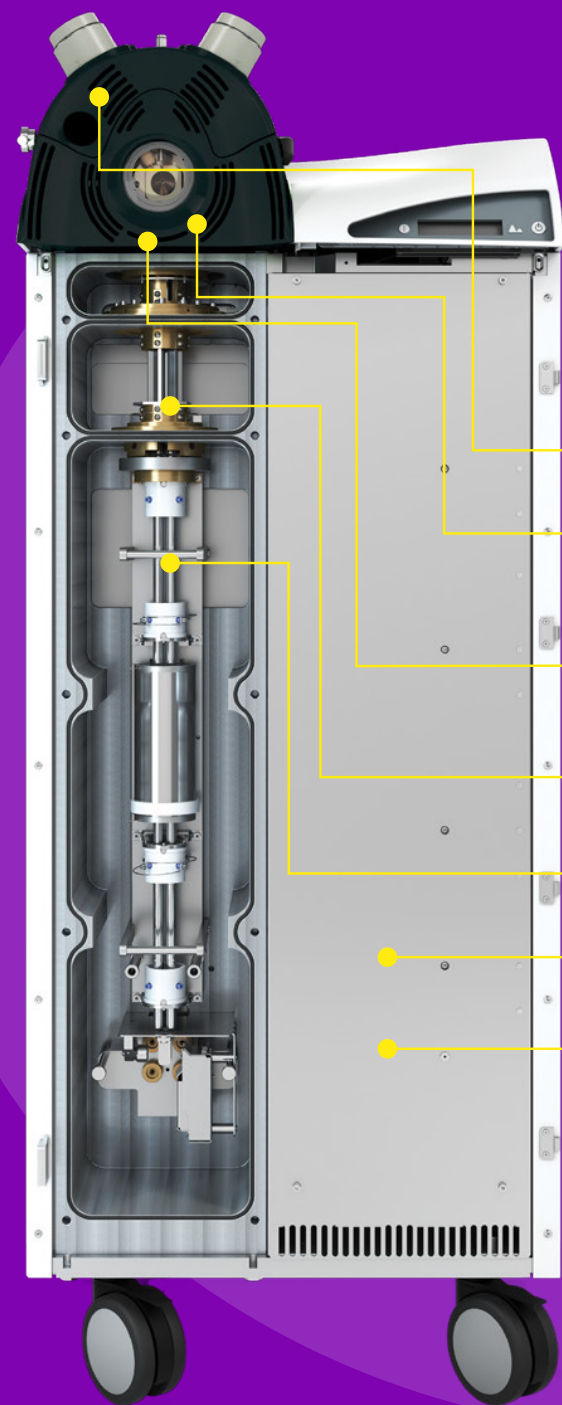
QSight™ 225 MD UHPLC-MSMS

Take your MSMS workflow to the next level with tier 2 chromatography

- Dual source to run multiple applications without reconfiguring the system
- Designed with class-leading sample capacity for high-throughput newborn screening as well as tier 2 chromatography
- Optimized UHPLC workflows: Integrated column switcher to easily switch between tier 1 and tier 2 runs without replumbing
- Intended for screening and confirmation workflows requiring chromatography



Inside and out, it's MSMS done right



- **Dual source (QSight 225)** - Dual source with automatic switching - no need to stop to change plumbing
- **StayClean™ source** - Self-cleaning design delivers exceptional sensitivity and uptime
- **HSID interface** - Provides low background and reliable results day after day, with equal response at any flow rate
- **Laminar Flow Ion Guide™** - Delivers highly efficient field-free high transmission
- **Mass filters** - Feature high-quality precision rods that deliver highly stable, precise mass filtering
- **Small footprint, vertical design** - Compact 50 cm x 50 cm x 110 cm form factor - no benchtop needed
- **Modular** - Plug-and-play design for ease of service

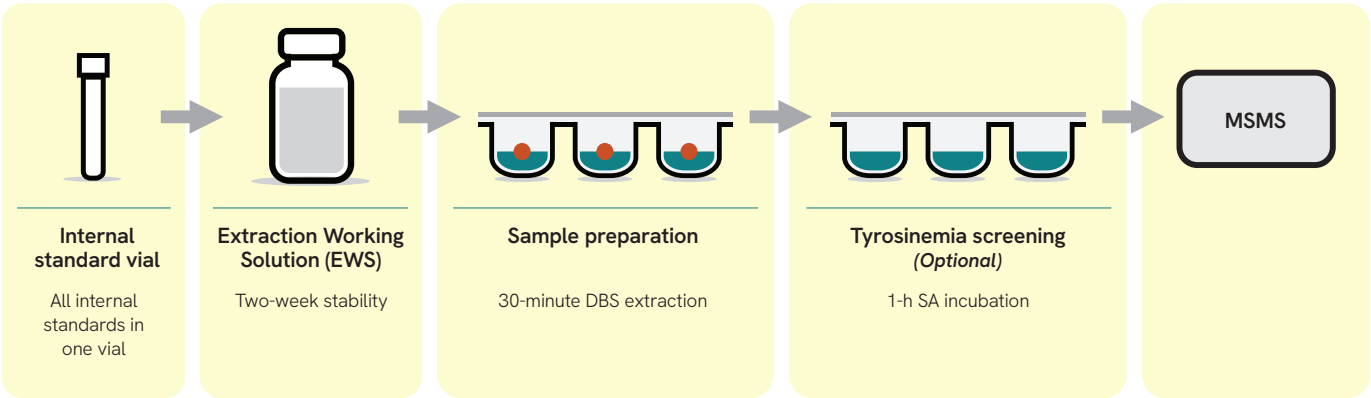
- ✓ High sensitivity and productivity
- ✓ High throughput, low maintenance
- ✓ Compatibility with Revvity assays

Next generation expanded newborn screening

NeoBase™ 2 Non-Derivatized MSMS Kit

- Test for up to 57 analytes including markers for X-ALD and ADA-SCID
- Short sample preparation time
- Automated import of QC targets and IS concentrations by a barcode reader
- Incorporated data analysis software for results analysis and management
- Automatic switch between NeoBase™ 2 non-derivatized MSMS kit and NeoLSD™ MSMS kit
- Pioneer in AAAC screening with the highest number of reference sites globally using NeoBase™ 2 non-derivatized MSMS Kit and NeoLSD™ MSMS Kit

NeoBase™ 2 Non-Derivatized MSMS Kit Workflow



Disorders

- Amino Acidopathies
- Organic Acidemia
- Fatty Acid Oxidation Disorders
- Purine Metabolism Disorder
- Peroxisomal Disorder

Kit contents:

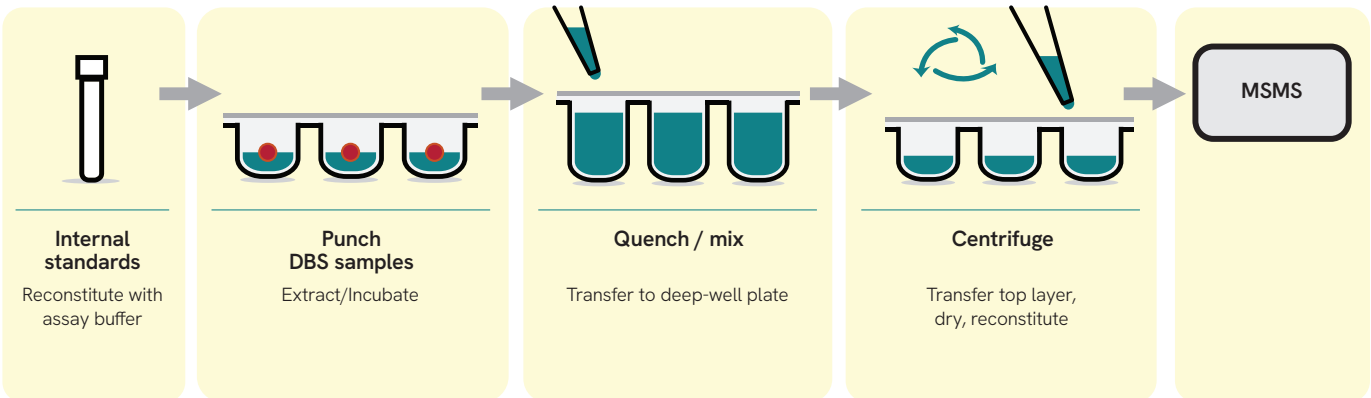
- NeoBase 2 non-derivatized assay solution
- NeoBase 2 succinylacetone assay solution
- NeoBase 2 internal standards
- NeoBase 2 controls
- U-bottomed microplates
- Adhesive microplate covers
- Barcode labels for plates
- Lot-specific quality control certificate

Easy and efficient LSD screening

NeoLSD™ MSMS Kit

- Screen for up to six LSDs from a single DBS punch and single incubation
- Less manual pipetting steps by incorporating automated liquid handler
- High analytical precision near screen cutoff
- Wide dynamic range and low retest rate
- Up to 1000 samples per day on a single QSight™ MD Screening System
- Flexibility for the expansion of NBS programs by MSMS to include additional LSDs

NeoLSD™ MSMS Kit Workflow



Disorders

- | | |
|--------------------|-----------|
| ▪ Pompe | ▪ Krabbe |
| ▪ MPS-1 | ▪ Gaucher |
| ▪ Niemann-Pick A/B | ▪ Fabry |

Kit contents:

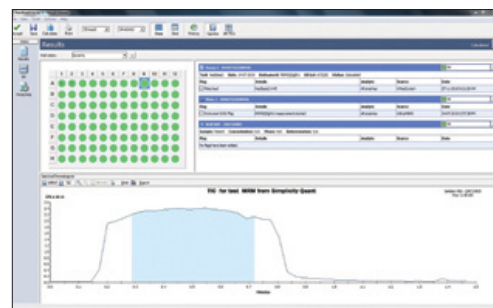
- NeoLSD extraction solution
- Neo MSMS flow solvent
- NeoLSD assay buffer
- NeoLSD kit controls
- NeoLSD substrates and internal standards
- Deep-well microplates
- U-bottom microplates
- Adhesive aluminum foil microplate covers
- Adhesive microplate covers
- Barcode labels for plates
- Lot-specific quality control certificate

Process more samples quickly and accurately

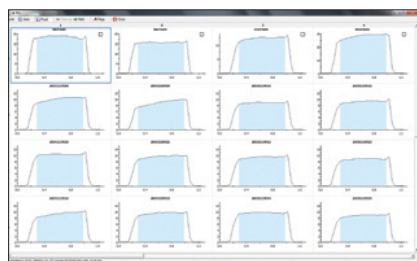
MSMS Workstation software (IVD) provides a user-friendly interface to store data, set limits, review results, and perform QC trend analysis for monitoring laboratory performance. The software enables you to process thousands of results per day, with maximum efficiency, while minimizing transcription or operator errors.

Key functionalities

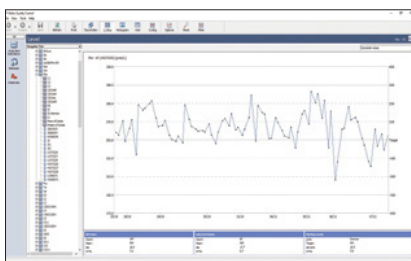
- Download a sample worklist, internal standard, and control target values from kit inserts, directly into the system, to avoid manual transcription errors
- Define flagging rules and cutoffs, with continuous adjustments to cutoff values, to reduce false positive results
- Analyze plate data, review flagged results, and group results by disorder to simplify data review and analysis
- Archive results from one or more instruments into a central database, and/or export results to CLIR or other LIMS, for easy review and retrieval



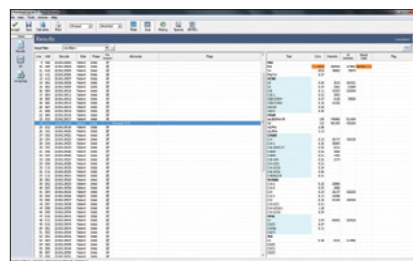
Set sample, internal standard, control limits and ratios to automatically flag results



Quickly view and interpret TIC data for the entire plate



Perform QC trend analysis across one or more instruments to monitor performance



Interpretation and grouping of results is simple and effective



revvity

www.revvity.com