

One powerful kit,
seven screens.
Now includes MPS II.



revvity

NeoLSD™ 7plex LC-MSMS kit

Expanding your screening capabilities

NeoLSD™ 7plex LC-MSMS kit enables efficient, accurate screening for **seven lysosomal storage disorders (LSDs)**, all in a single injection from two DBS punches.

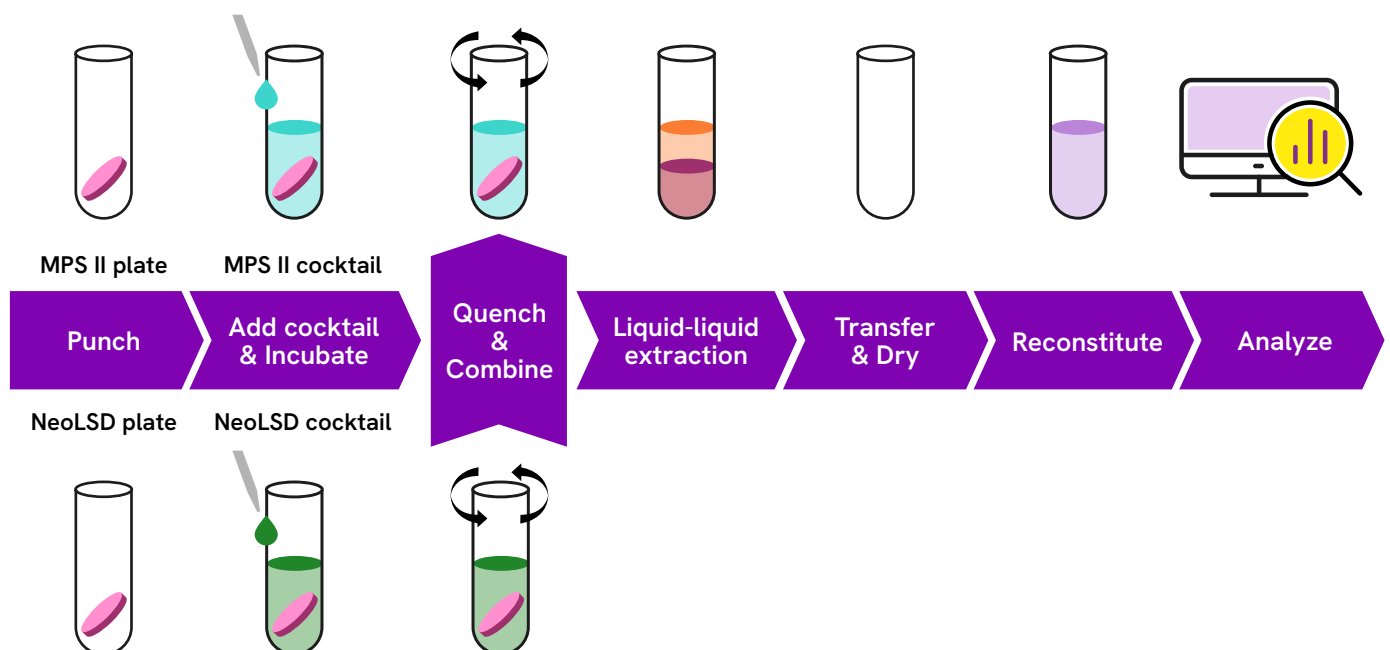
Screen for:

- Gaucher
- Niemann-Pick A/B
- Pompe
- Krabbe
- Fabry
- MPS I
- MPS II (New)

Key benefits:

- Multiplex 7 LSDs for increased efficiency
- Specifically designed software for streamlined, easy processing of LC-MSMS produced data
- LC-MSMS-based assay ensures accuracy and enhanced separation
- 2 punches consolidated to 1 injection
– driving efficient use of the MSMS instrumentation
- Trusted, proven technology – now expanded

NeoLSD 7plex LC-MSMS kit workflow



7 lysosomal storage disorders from **2** punches,
combined to **1** dataset

Smarter screening with MSMS Workstation software

Our **IVDR compliant MSMS Workstation** software is tailored for high-throughput newborn screening labs:

Seamless workflow integration that can...

- Track samples through multi-punch workflows
- Combine data from separate preparation pathways
- No manual data entry - Enables the import of IS concentrations, QC targets and SD from QC certificate with a barcode reader
- Easy optimization of cut-offs
- Intuitive and familiar Revvity user interface
- Easy result acceptance and customizable exporting and reporting tools

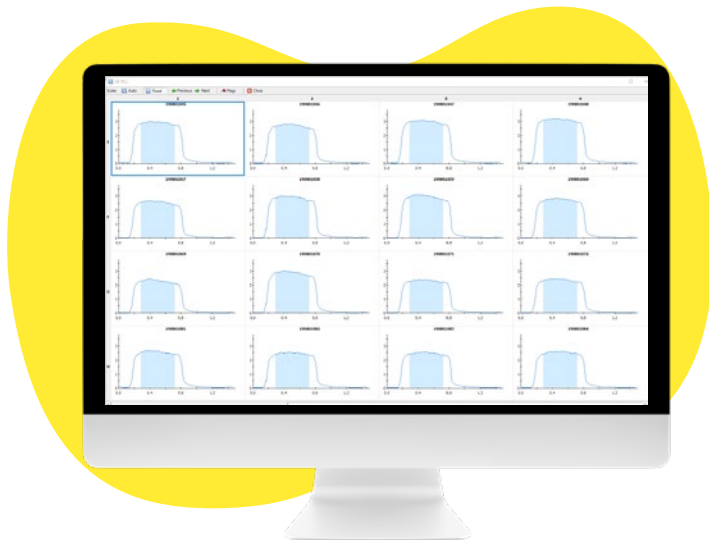
MSMS Workstation modules:

- Result Viewer
- Plate Generator
- Kitlot Editor
- Cutoff Analyzer
- Quality Control



QSight® 225MD UHPLC screening system

MSMS Workstation software at the center of your workflow:



- ✓ Intuitive
- ✓ Adaptable
- ✓ Accurate

Certified confidence: CE-IVDR-compliant

The NeoLSD 7plex LC-MSMS kit is CE-IVDR-certified, ensuring the highest standards of diagnostic accuracy and safety.

Why it matters:

- Provides peace of mind, knowing you can move forward with complete confidence
- Enhances lab compliance
- Supports the best outcomes in newborn care



Complete workflow integration

The NeoLSD 7plex LC-MSMS kit is optimized for use with the QSight® 225MD UHPLC Screening System, delivering a streamlined workflow with the new Simplicity 2.3 software version optimized for high-throughput LC assays:

- Automated peak detection and integration
- Peak quality flagging for improved reliability
- Injection counter for individual LC-columns

Built on success with a globally trusted solution.

- NeoLSD MSMS Kit: Over 12 million babies screened
- Lower false positive rate vs. other methods*
- 20+ years of Revvity leadership in MSMS innovation

*Viall S, Held P. Five years of newborn screening for Pompe, Mucopolysaccharidosis type I, Gaucher, and Fabry diseases in Oregon. Mol Genet Metab Rep. 2025 Apr 15;43:101221



revvity

www.revvity.com

NeoLSD™ 7plex LC-MSMS kit is a CE marked *in vitro* diagnostic medical device, compliant with IVD Regulation (EU) 2017/746, certified by Notified Body number 0537, Eurofins Electric & Electronics Finland Oy.

QSight® 225MD UHPLC Screening System is a CE marked *in vitro* diagnostic medical device, compliant with IVD Regulation (EU) 2017/746.

MSMS Workstation is a CE marked *in vitro* diagnostic medical device, compliant with IVD Regulation (EU) 2017/746.

Copyright ©2025, Revvity, Inc. All rights reserved.
Revvity is a trademark of Revvity, Inc.
All other trademarks are the property of their respective owners.

For *in vitro* diagnostic use. This product is only available where licensed in accordance with law. Please contact your local representative for availability.