



# Newborn health: Screening solutions, services, and research.

## How you can work with Revvity

Revvity is pleased to offer a complete portfolio of services and solutions to meet your laboratory, hospital, or health center's unique needs. See how you can partner with the global leaders in newborn screening today.



### Screening solutions

#### Supporting the entire NBS workflow.

- Instruments for labs: Ideal for startups or high-volume automation across MSMS, molecular, Immunoassay, and IEF
- Kits designed for over 50 different disorders, all integrating smoothly with Revvity instruments
- Software that is cloud based, secure, and built to deliver results analysis fast and efficiently



### Testing services

#### Ensuring a healthier start for every newborn.

- Newborn screening Panels- covering metabolic disorders
- StepOne Test: Comprehensive biochemical profiling for over 70 disorders in newborns
- Newborn genetic panel: Next Gen Sequencing for Newborns



### Research solutions

#### A complete and customizable NGS workflow.

- An end-to-end NGS workflow for newborn research
- Compatible with existing lab instrumentation (modular option)
- Customizable to your laboratory
- 390+ gene panel
- 87K+ variant database
- Genomic software suite