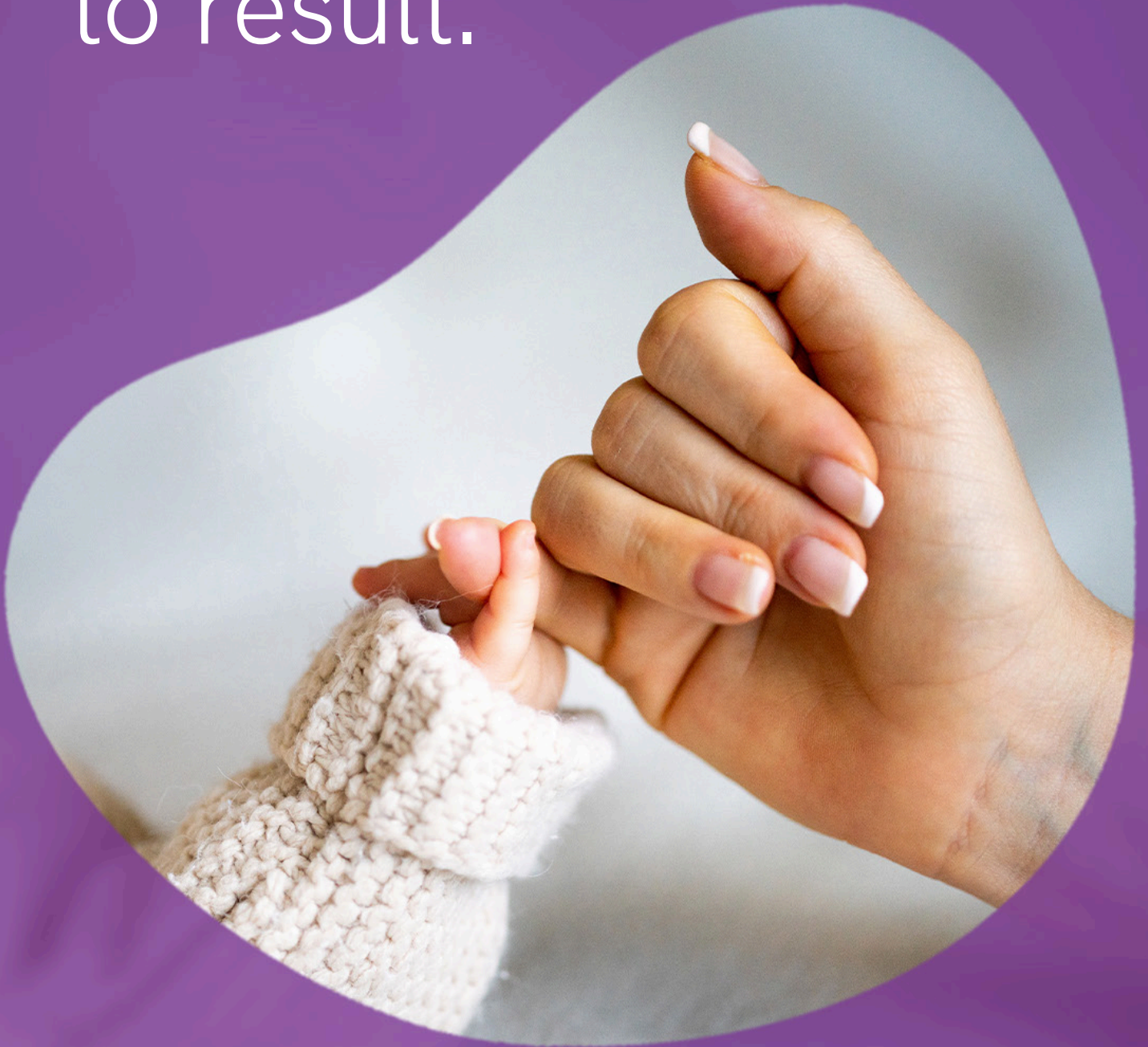


Complete solutions  
from sample  
to result.



revvity

Complete solutions for newborn screening



# Helping babies survive and thrive.

Newborn screening can prevent hundreds of thousands of instances of physical and mental disabilities and even deaths. As the global leader in newborn screening, over 800 million babies have been screened with our products over the last 25 years.

## **140 million**

babies are born worldwide annually.

## **1 in 3 babies**

receive screening of any kind, and many are only screened for one or two conditions.

## **50+ conditions**

are known to have good outcomes when detected and diagnosed early.

## **85 newborns**

get a healthier start to life everyday thanks to screening using our products.

## **Over 110 countries**

have newborn screening programs supported by our products.

# Earlier detection with a complete range of screening solutions for NBS disorders

Revvity's solutions support healthcare professionals in identifying affected newborns early, enabling timely medical intervention and care decisions.

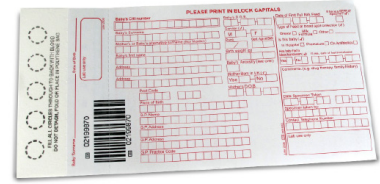
Disorders	Analytes	Instruments				
		GSP™	VICTOR2™ D	QSight™ MD	Eonis™ System	Migele™ Gel Electrophoresis Unit
Congenital hypothyroidism	hTSH	✓	✓			
	T4	✓	✓			
Congenital adrenal hyperplasia	17-OHP	✓	✓			
Cystic fibrosis	IRT	✓	✓			
Phenylketonuria	Phenylalanine	✓	✓			
Galactosemia	GALT	✓	✓			
	TGAL	✓	✓			
G6PD deficiency	G6PD	✓	✓			
Biotinidase deficiency	Biotinidase	✓	✓			
Duchenne Muscular Dystrophy	CK-MM	✓	✓			
Hemoglobinopathies	Hb Variants					✓
In-born errors of metabolism	Multiple			✓		
ADA-SCID	Ado, dAdo			✓		
X-ALD	Multiple LPCs			✓		
7 types of Lysosomal disorders	Multiple			✓		
Severe combined immunodeficiency	TREC				✓	
Spinal Muscular Atrophy	SMN1				✓	
X-linked agammaglobulinemia	KREC				✓	
Toxoplasmosis	IgM	✓				

Products may not be licensed in accordance with the laws in all countries, such as the United States and Canada. Please check with your local representative for availability.

# From sample collection to result, we cover the entire screening process

## 1 - Sample collection

**Revvity 226 Sample Collection Device** is a dried blood spot collection card comprised of 100% pure cotton linter filter paper validated for even and uniform sample distribution. Revvity can custom-print and manufacture a device format that meets your specialized newborn screening requirements.



## 2 - Punching

**DBS Puncher™ Instrument** is a semi-automatic device for punching dried blood spot samples into microtitration plates. The 2-plate capacity puncher is an easy-to-use, robust option for lower capacity screening.



**Panthera-Puncher™9 Instrument** is the next-generation semi-automatic dried sample punching device with an automatically adjusting punch pattern. Panthera-Puncher instrument allows you to process sample materials with optimum efficiency and performance.



## 3 - Sample preparation

**DELFIA™ Trio** is a 3-in-1 instrument developed for sample processing of DELFIA manual kits. It combines disk removing, washing, and dispensing into one instrument offering a streamlined workflow for manual DELFIA users.



**DELFIA Plateshake** for sample processing with two preset shaking speeds optimized for DELFIA assays.



**TriNEST™ incubator shaker** is a solution intended for use in the Victor2D and MSMS workflows.



## 4 - Analysis and measurement platforms

**VICTOR2™D instrument** is the ideal measurement platform for low volume laboratories. Easy to use and reliable, Victor2D provides start-up screening programs with the right balance of cost-effectiveness and control.

**Disorders:** Congenital Adrenal Hyperplasia, Congenital Hypothyroidism, Galactosemia, Phenylketonuria, G6PD Deficiency, Phenylalanine, Cystic Fibrosis, Biotinidase Deficiency



**GSP™ instrument** for neonatal dried blood spot assays is the system of choice for medium to large laboratories. Performing all assay stages automatically with superior accuracy and traceability, GSP instrument supports our complete range of analytes and employs several measurement techniques for screening of core NBS disorders. From plate loading to results output, GSP instrument offers laboratories outstanding efficiency and confidence throughout the screening process.

**Disorders:** Congenital Adrenal Hyperplasia, Congenital Hypothyroidism, Galactosemia, Phenylketonuria, Cystic Fibrosis, Biotinidase Deficiency, G6pd Deficiency, Duchenne Muscular Dystrophy, Toxoplasmosis



## 4 - Analysis and measurement platforms

**QSight™ MD systems** enable expanded high-throughput mass spectrometry screening and, with the QSight 225 MD UHPLC Screening System, the option for tier 2 testing – using just a single heel-prick dried blood spot (DBS) card. The QSight system delivers a unique combination of self-cleaning MSMS instrumentation that provides exceptional system uptime, a simple setup process, and easy-to-use database software that simplifies review and analysis.

**Disorders:** Metabolic Disorders, ADA-SCID, X-ALD, and Lysosomal Storage Disorders



**Eonis™ System** - Multiplexing of SMA, SCID and XLA screening with one assay utilizing Real-Time PCR technology, without increasing daily hands-on workload and complexity. The design of the EONIS assay and system enable automation without compromising the sample traceability from punch to result.

**Disorders:** Severe Combined Immunodeficiency, X-linked Agammaglobulinemia, Spinal Muscular Atrophy



**Migele™ Gel Electrophoresis instrument** is a robust, scalable and easy to use instrument for detection of crucial lifesaving disorders. It works with the RESOLVE™ hemoglobin kit to detect hemoglobinopathies in newborns and adults, including sickle cell disease and other hemoglobin variants and thalassemias.

**Disorders:** Hemoglobinopathies



# EVOYA software platform

An interface layer that connects the entire workflow, the EVOYA™ software platform has been designed uniquely for newborn screening workflows while managing your data securely.

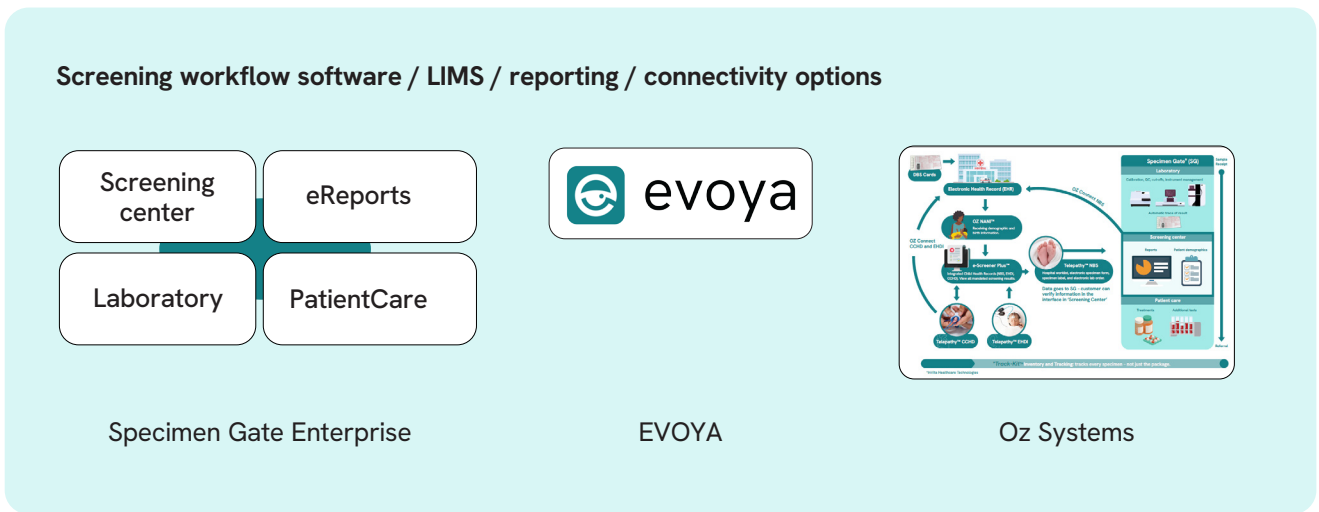
For those involved in newborn screening, keeping that data secure is paramount. Information about newborns needs to be provided quickly and securely without compromising your service.

In a system dealing with such sensitive data, security must be a key component and compliant with the highest standards.

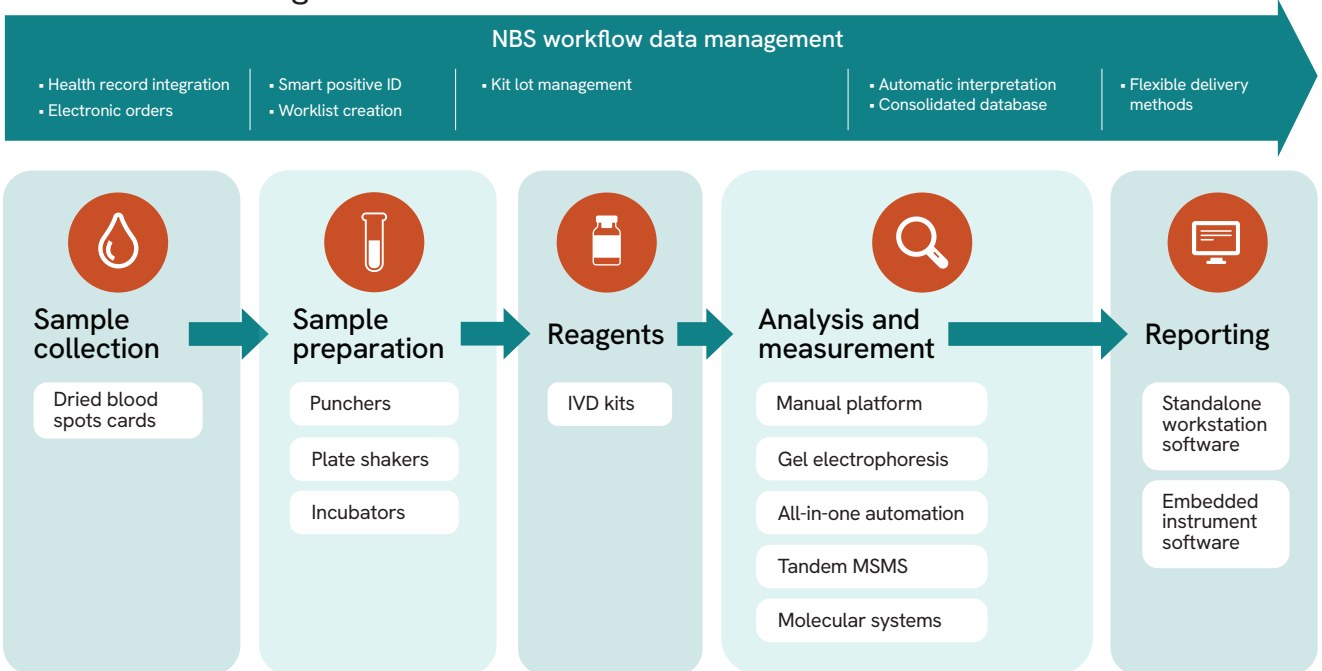


# Data management systems: Enhancing efficiency, accuracy, and patient care

Control and monitor the entire newborn screening process - including patient reporting and follow-up -- in one comprehensive information management system. Our informatics software for neonatal screening is the result of years of continuous development based on our experience as the recognized leader in newborn health and laboratory information management technology. The range of products are widely used in all parts of the world, and represents the state-of-the-art in data organization, storage, analysis and security.



## Newborn screening workflow



# Elevated support, delivering efficient transformation

Through our experts in newborn screening, we can help you to create service programs that meet the standards of even the most demanding screening laboratories. Our global presence enables Revvity experts to share both innovation and deployment expertise worldwide.



**Consultation** - Revvity develops, manufactures and supports a market leading range of instruments, reagents, and consumables to the screening community - providing unparalleled knowledge and depth of expertise with over 75 years of experience.

**Training** - Our training team offers diagnostics product training courses for customers. Training is conducted by product managers and specialists from customer and technical support, software services and R & D. We aspire to meet your training expectations and to give you the best possible training experience.

**Service excellence** - Revvity's dedicated diagnostics service team comprises trained and certified engineers who have an average of 15 years of experience maintaining Revvity and partner equipment.

**Customer support** - Our product specialist and technical support teams provide country and regional support, ensuring you have the necessary touchpoints when you need them and where you need them.

# Scalability is key



## Startup or small laboratories:

### Cost-effective solutions:

Affordable, manual solutions ideal for small or startup labs.

### Scalability:

Expand easily by adding parameters or increasing sample volume.

### Flexible setup:

Start small, grow as your lab's needs increase.



## Growing mid-size laboratories:

### Full automation:

Automated solutions streamline workflows for neonatal screening.

### Efficiency focus:

Automation frees up time for other lab management tasks.

### Enhanced performance:

Transition to mid-level automation for greater convenience.



## Large laboratories with high-capacity:

### Expanded screening:

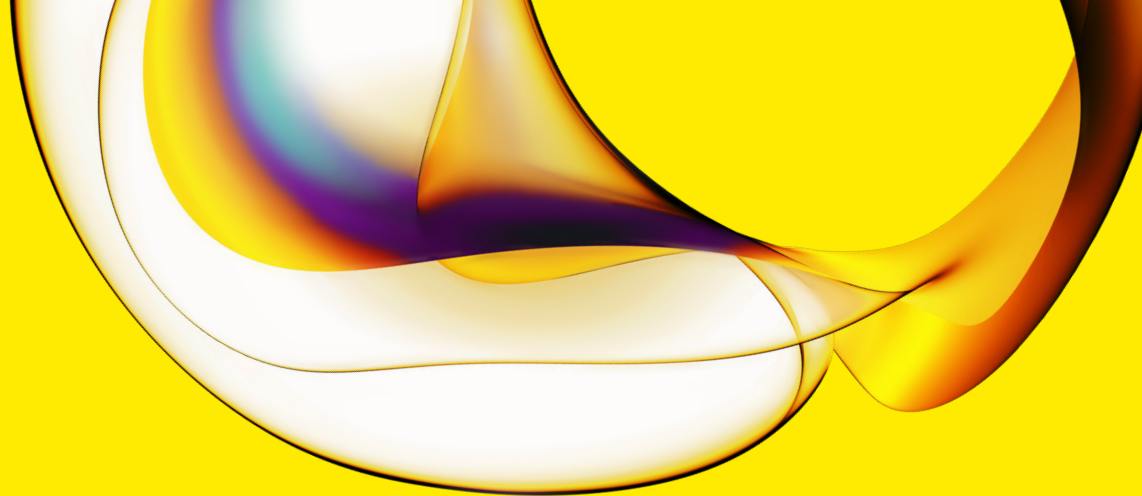
Add more disorders to your screening menu with ease.

### High throughput:

Handle large volumes with automated, high-throughput quantitation.

### Seamless integration:

Fully integrate your systems for maximum efficiency.



[www.revivity.com](http://www.revivity.com)

revvity

[www.revivity.com](http://www.revivity.com)

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.

Revvity does not endorse or make recommendations with respect to research, medications or treatments. All information presented is for informational purposes only and is not intended as medical advice