

Patient safety.
Quality of care.
Intervention
timeliness.



revvity

Oz newborn screening suite
End-to-end newborn blood spot
screening management



- Introduction
- Blood spot screening
- Oz NANI
- Oz telepathy NBS
- Oz connect NBS
- Integrated bundle

Automate timeliness

An end-to-end solution with real time, verifiable birth information, efficient automated specimen collection, smart tracking for effective alerts and timely results reporting.

With all newborn screening, there is a window of opportunity to change a baby's future and prevent life-lasting disability or death. We are baby-centric because we know you hold that baby's future in your hands. We are specimen-centric, offering Track-Kit™ because in your world, accurate, timely specimens are the first step in changing the future.



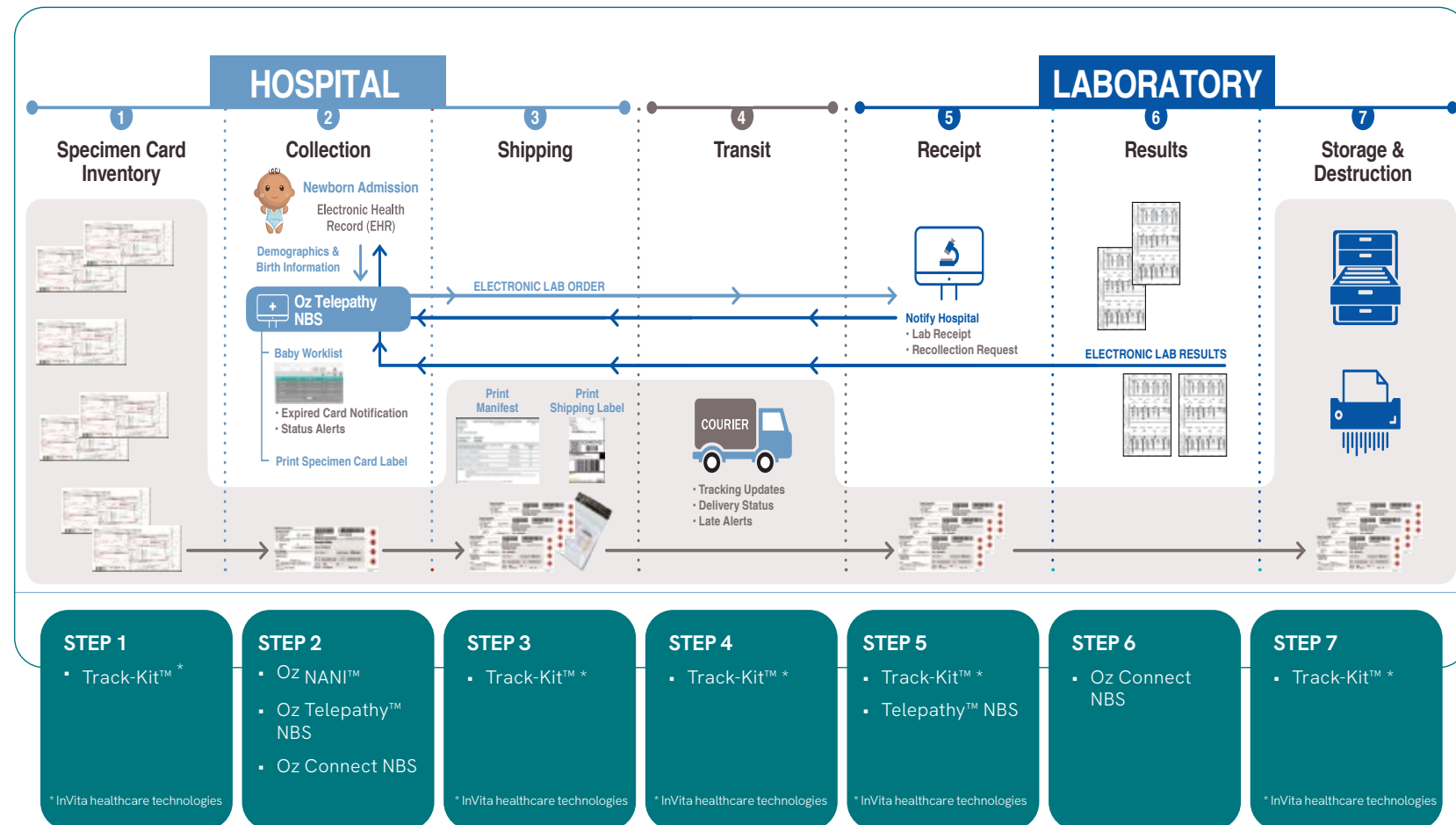
**Blood Spot Screening
(NBS)**





End-to-end newborn blood spot screening management

- Introduction
- Blood spot screening
- Oz NANI
- Oz telepathy NBS
- Oz connect NBS
- Integrated bundle





- Introduction
- Blood spot screening
- Oz NANI
- Oz telepathy NBS
- Oz connect NBS
- Integrated bundle

End-to-end newborn blood spot screening management

Oz NANI™

Integrates patient demographics from the hospital's EHR (Electronic Health Record) to all Oz software solutions, giving you an accurate denominator of hospital births in real time. You know who needs to be screened.*

Oz Telepathy™ NBS

Oz Telepathy Newborn Blood Spot (NBS) prepopulates an electronic form with the state laboratory's required data elements, which are printed on a label and affixed to the blood spot specimen card. Sends an electronic lab order with complete information to the state lab.

- Manages the care of each baby by providing a worklist of babies needing screening and status for each specimen collected.
- Eliminates handwritten specimen cards to improve information content, legibility, and accuracy
- Improves accuracy by scanning the specimen barcode
- Simplifies hospital workflow by streamlining the process
- Displays age in hours in order to collect specimen at appropriate time
- Captures multiple specimens for an individual patient

- Improves timeliness for reporting back to the primary care physician and parent
- Provides specimen shipping and delivery notifications
- Captures the submitted information and stores it for future audit

*The hospital EHR and interface engine will dictate which fields can be electronically sent.



Global view of quality assurance metrics

- Specimen card inventory
- Accurate demographics
- Age at collection
- Percentage of newborns screened
- Percentage of newborns needing rescreen
- Incoming specimen cards and shipments
- Transit time

Our integrated collection-to-reporting solution simplifies workflow in the hospital and state lab, eliminates handoffs to deliver timely results so your Newborn Blood Spot (NBS) screening program is more cost-effective, improves patient safety, and reduces human error using information technology standards.



- Introduction
- Blood spot screening
- Oz NANI
- Oz telepathy NBS
- Oz connect NBS
- **Integrated bundle**

End-to-end newborn blood spot screening management

Oz Telepathy NBS

Prior to the lab receiving the specimen, Oz Telepathy NBS generates an electronic lab order (ELO) to the state laboratory information management system (LIMS) in an HL7 standards-based message with required newborn blood spot data elements. ELO provides the highest quality demographic data when coupled with Oz NANI.

- Informs lab staff that the specimen has been collected
- Saves time by eliminating the re-keying of data elements at the lab
- Reduces missing or incorrect data received at the lab
- Allows for a cross check to confirm a specimen is received

Oz Connect™ NBS

Lab results are sent electronically to the EHR using HL7 standards-based messaging. Oz Connect NBS improves accuracy of the data reported to the newborn's medical home. Includes message monitoring to ensure messages are consumed.

- Saves time by eliminating printing and faxing at the lab as well as scanning results at the hospital
- Improves timeliness of data returning to the hospital and ultimately to the newborn's medical home



Track-Kit™ *

Track-Kit, from InVita Healthcare Technologies, reduces late specimens. Specimen tracking from collection to storage gives you an automatic end-to-end audit trail.

- Inventory management ensures never running out of viable specimen cards and identifies card wastage
- Notifies collectors of specimen card expiration (via Oz Telepathy NBS) which increases specimen quality
- Increases shipping efficiency of specimen cards by generating the shipping manifest and label
- Tracks packages AND specimens during shipping process
- Provides delivery notification to hospitals (via Oz Telepathy NBS) to meet CAP/CLIA regulations
- Alerts the lab and hospital when specimen delivery is missing or late
- Tracks specimen storage and destruction
- Provides global real-time view of state-wide activity



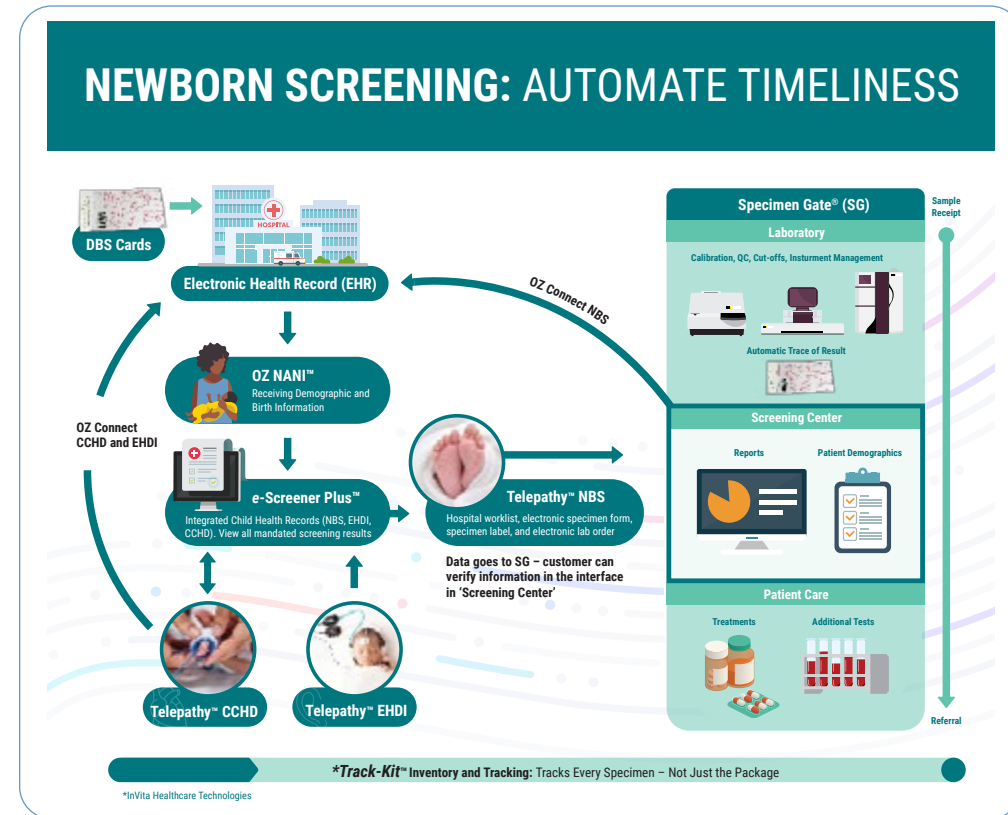
An integrated bundle

- Introduction
- Blood spot screening
- Oz NANI
- Oz telepathy NBS
- Oz connect NBS
- Integrated bundle

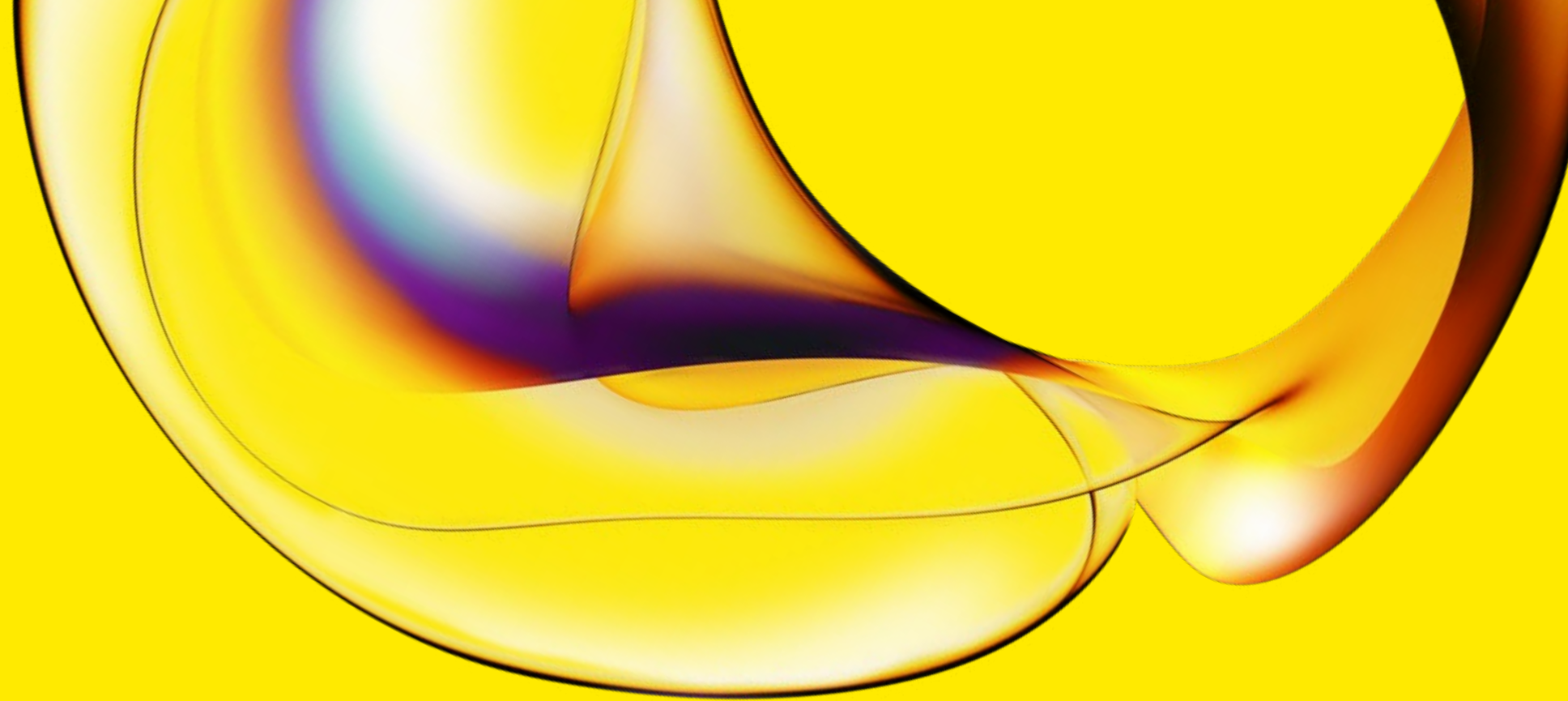
Revvity gives caregivers the tools – and the time – to deliver the personal care and responsiveness patients experienced decades ago. Our end-to-end newborn screening solutions reduce human error and improve patient safety, quality of care and intervention timeliness.

Revvity newborn screening

We are a global market leader in neonatal screening, currently serving customers in more than 100 countries. We are a total solution provider offering complete systems based on a broad range of high quality, validated products, including newborn screening kits, consumables, instruments and software. Since 1985, over 700 million babies have been screened with our products. Every day more than 70 babies around the world get a healthier start to life thanks to the early detection of a serious disorder.



* Telepathy™ EHD and Telepathy™ CCHD are required for results reporting back into the EHR.



www.revivity.com

revivity

Revvity, Inc.
940 Winter Street, Waltham, MA 02451 USA
www.revivity.com

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2023, Revvity, Inc. All rights reserved.

1200800