

We seek out
and introduce
new innovative
data sharing
solutions.



revvity

Oz e-Screener Plus™
Newborn screening
information management



- Introduction
- For healthcare providers
- Hospital interfaces
- Data acquisition
- Case management
- Continuum of care
- Contact us

Oz e-Screener plus

We stay close to the pulse of newborn screening to learn and understand the changing needs of newborns and all who care for them.

Oz eSP™ provides a baby-centric solution that allows you to manage all newborn screening data in one integrated child health record that stores birth admission, newborn screening information, and risk factors. It is the foundational infrastructure for all implementations. It has been widely used in state screening programs for two decades to monitor quality and assure care. Oz eSP meets or exceeds all HIPAA and HITECH requirements and adheres to NIST 800 guidelines.



Blood spot screening
(NBS)



Hearing screening
(EHI)



Heart screening
(CCHD)



Birth conditions and
Birth defects monitoring

Products may not be licensed in accordance with the laws in all countries. Please check with your local representative for availability.



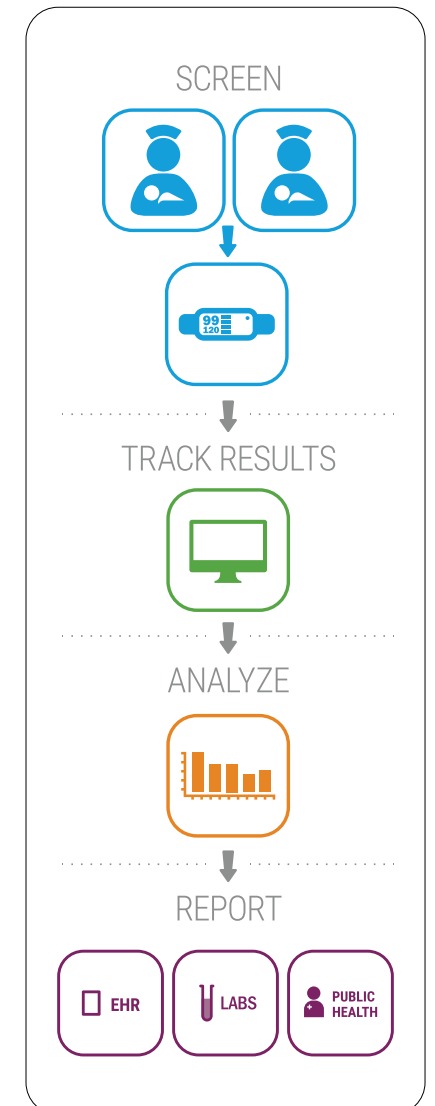
For healthcare providers

- Introduction
- For healthcare providers
- Hospital interfaces
- Data acquisition
- Case management
- Continuum of care
- Contact us

- Integrated continuum of care for all stages of the newborn screening program: inpatient and outpatient screening, diagnostic follow-up and intervention
- Fully secure web-based data management accessible anywhere, anytime, with no hardware or software required; complete audit trail of user and system changes protects programs from negligence and improves risk management and quality assurance functions
- Rule-based care paths that support next steps and best practices with clinical decision support
- Real Time Results—less time finding results and determining next steps means more time caring for babies
- Error reduction in reporting and data collection
- Verifiable data—not only aggregate data—for reporting and monitoring program success
- Analytics across individual hospitals and entire state programs
- Automated audit trails of user and system changes
- Provides data for risk management and quality improvement.
- Privacy and security of patient-sensitive information through rule-based and role-based access rights
- Standardized reporting mechanisms from state public health to federal agencies—HL7 electronic messaging tested and IHE proven
- National Quality Forum (NQF) measures and new Stage 2 Meaningful Use support for reporting from hospitals to state programs

Additional modules

- **CDC Dashboard** makes reporting data to the CDC easier and complete. Reports quality data based on verifiable data as well as aggregate regional reporting.
- **Vital Records Match** makes sure every baby in your state is in eSP regardless of where they are born - home or hospital. Compares patient records in eSP to Vital Records.
- **Early Intervention Data Linkage** provides data to your early intervention agencies for tracking babies who need further care.





- Introduction
- For healthcare providers
- Hospital interfaces
- Data acquisition
- Case management
- Continuum of care
- Contact us

Hospital interfaces benefiting public health

Revvity provides interfaces to hospitals that also benefit the goals of public health.

Oz NANI™

Interface between a hospital's EHR (Electronic Health Record) and all software solutions including our Oz eSP™ database and Oz Telepathy tools. Oz NANI gives you an accurate denominator of hospital births in real time. Revvity assists the hospital in setting up the ADT (Admission, Discharge and Transfer) message using IHE and HL7 standards-based protocols.*

Oz Telepathy™ NBS (electronic lab order)

An electronic lab order (ELO) is generated and sent to the state laboratory information management system (LIMS) in HL7 standards-based messaging with required newborn blood spot data elements (Oz NANI required).

- Saves time by eliminating the re-keying of data elements at the lab
- Reduces missing or incorrect data received at the lab
- Allows for cross check on whether a specimen is received
- Assists the laboratory in understanding the workload and allowing for better allocation of staffing
- Prints a label that is affixed to the specimen card providing legible, complete demographics.

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

Oz Telepathy™ CCHD & EHDI results into eSP™

- Eliminates duplicate data entry for hospital staff
- Reduces errors by eliminating manual data entry
- Improves timeliness for hospital reporting results to the state

Oz Connect NBS (electronic lab results)

Interface between the laboratory and hospital's EHR (Electronic Health Record) making electronic consumption of lab results possible. Includes message monitoring to ensure messages are consumed.

- Eliminates printing and faxing of results to the hospitals
- Increases timeliness of data returning to the hospital and ultimately to the newborn's medical home
- Improves accuracy of the data reported to the newborn's medical home





- Introduction
- For healthcare providers
- Hospital interfaces
- Data acquisition
- Case management
- Continuum of care
- Contact us

Direct acquisition of data from screening hardware: critical to quality and accuracy.

Revvity eliminates manual entry and human interpretation of the results. When a human has to transpose results from one piece of equipment to the next, errors do occur. A passed screening result that is accidentally marked as a refer causes the family undue stress. But worse, a referred result that is mistakenly marked as a pass can delay intervention for the baby. In the case of CCHD wrong screening results could result in death. Eliminating manual entry of results and electronically capturing them greatly increases accuracy and ultimately improves patient care. Revvity provides the best solution by interfacing to many manufacturers' equipment.

Oz Telepathy EHDl solution

- Collects results directly from screening devices and into Oz eSP™

Oz Telepathy CCHD solution

- Captures multiple screening values directly from the pulse oximeter
- Provides clinical decision support increasing accuracy
- Monitors outcomes of the clinical decision support algorithm chosen by your state
- Transfers screening results into Oz eSP database





Case management and care coordination

- Introduction
- For healthcare providers
- Hospital interfaces
- Data acquisition
- Case management
- Continuum of care
- Contact us

Optimal care and delivery of services are paramount for every state agency or department working with the public. Case management provides a view of systems and services that impact targeted populations — merging information between health settings while informing the entire care team and public health in a meaningful way. Care coordination assures that infants who do not pass a screening receive necessary follow-up including rescreening, diagnostic evaluation, treatment, and referral to early intervention service agencies. Care coordination supplies information to families and providers so they can more effectively advocate and receive access to appropriate care... to detect and connect, one baby at a time.

Case management

- Secure online, web-based tool for managing caseloads of children needing follow-up care after newborn screening
- Customizes workflow and care needed based on user's tasks/access
- Case level tracking of follow-up care, including detailed reports and communications with family and providers
- Allows users to document care, create reminders and enter case notes at each step in the case management process
- Applicable for public health reporting and diseases surveillance





- Introduction
- For healthcare providers
- Hospital interfaces
- Data acquisition
- Case management
- Continuum of care
- Contact us

Integrated continuum of care

The whole reason for newborn screening is identifying the babies who need follow-up. Accurate and timely results are essential to prompt newborn care. With all newborn screenings, there is a critical window of opportunity to change a baby's future and prevent life-lasting disability or death. Revvity creates solutions to ensure caregivers know what care is needed within that window of opportunity.

Diagnostic audiological assessment

Revvity is proud of their partnership with several states that have a high report rate for diagnostic assessment results. Oz eSP provides access to pediatric audiologists to enter diagnostic audiological results. A diagnostic summary as well as full diagnostic assessment data for each ear including amplification and intervention information can be entered into the child's health record.

Early intervention

Oz eSP provides data to your early intervention agencies for tracking babies who need further care. Based on the care path, once a referral to early intervention is needed Oz eSP will trigger a referral to early intervention and mark the child's record as requiring early intervention.





www.revivity.com

revivity

Products may not be licensed in accordance with the laws in all countries.
Please check with your local representative for availability.

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.

1200821