



Screening solutions improving new lives.

Oz NANI™ (Newborn admission notification information)

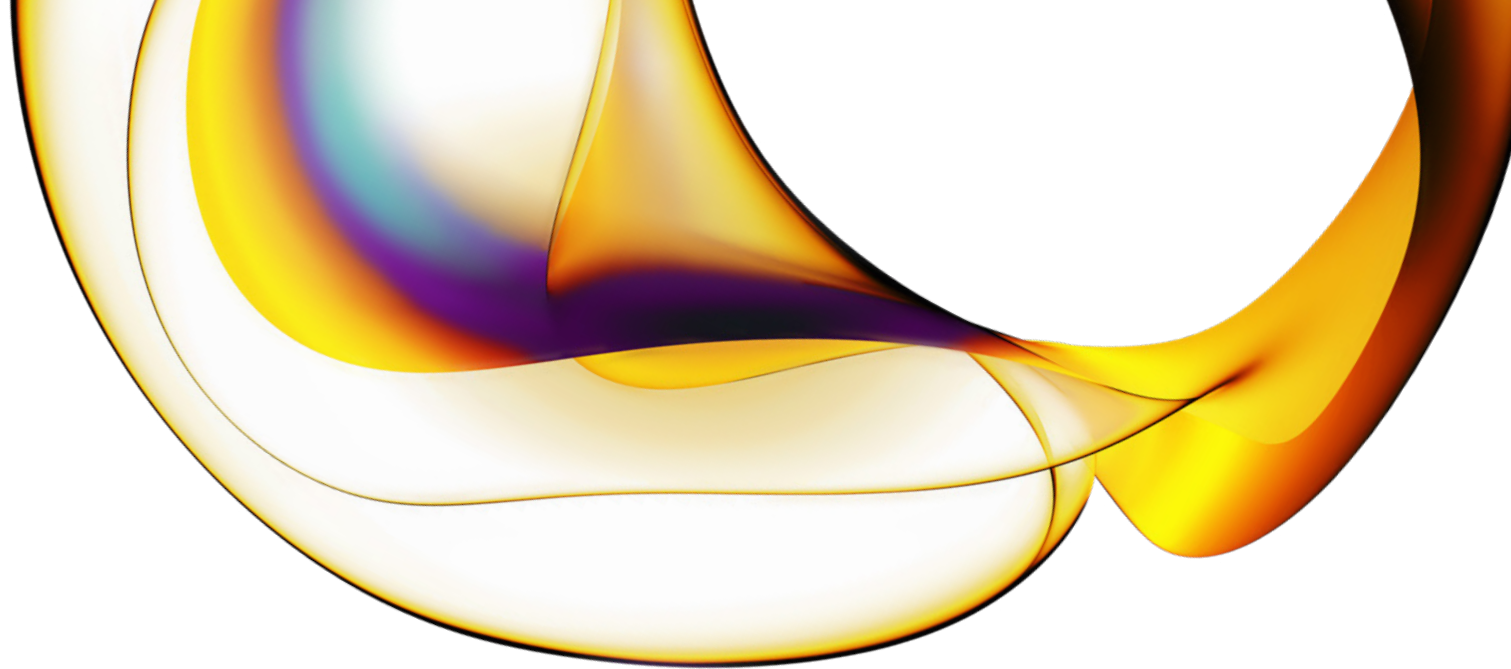
Eliminate the time, repetition and opportunity for error created with manual data entry. Let your staff focus on patient care and let Oz NANI ensure complete, accurate, up-to-date patient demographics and birth information. Oz NANI is the interface between a hospital's EHR (Electronic Health Record) and all Revvity software solutions.*

Oz NANI improves quality and reduces risk by providing:

- Automated newborn admission demographics
- Automated updates including discharge and internal transfer information in real time
- Bolsters patient safety by providing secure patient demographics
- Accurate, automated denominator of births for screening workflow optimization
- Eliminates duplicate data entry and other manual entry errors
- Updates the patient name in real time so follow-up is more efficient

Revvity assists your hospital in setting up the ADT (Admission, Discharge and Transfer) message using Integrating the Healthcare Enterprise (IHE) and HL7 standards-based protocols.

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



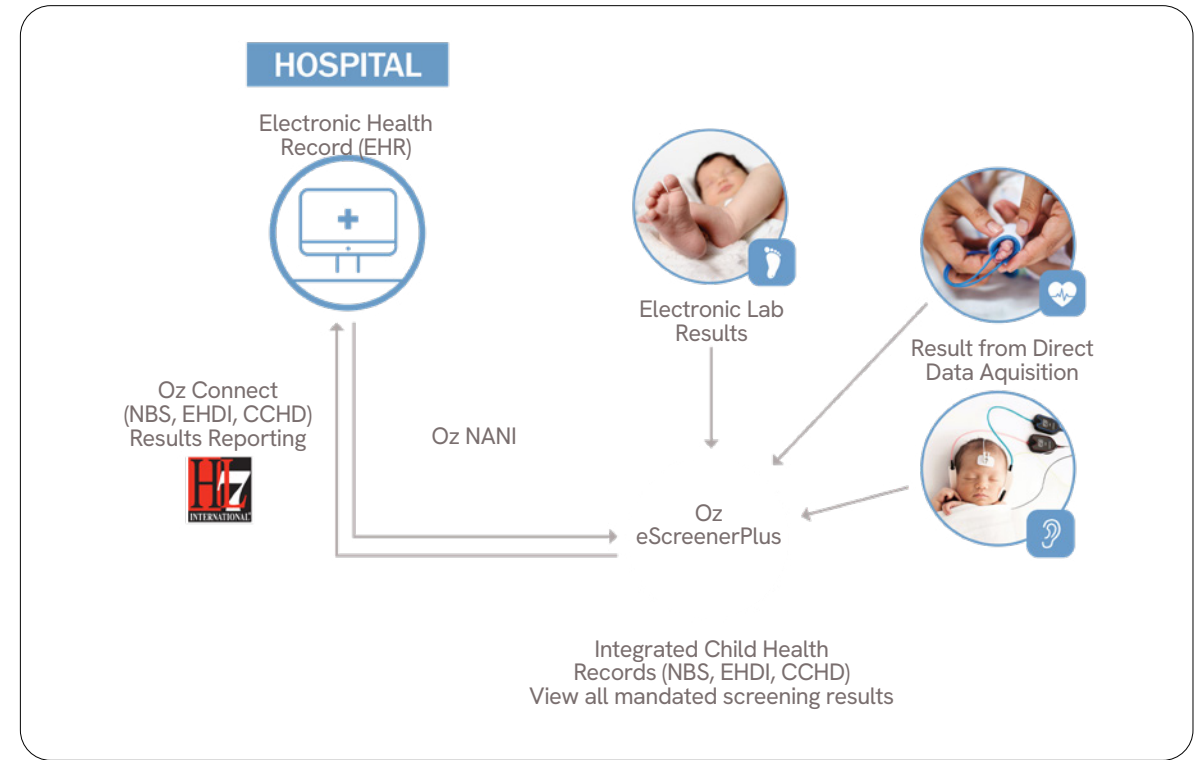
Tracey Kunkel, Secretary, UnityPoint Health-St. Luke's, Sioux City, IA

"Oz NANI has made my work more efficient, plus it saves time. Oz NANI operates in real time and it automatically updates information. The baby's given name gets automatically updated even though the baby is not named at birth. Plus, other data that was not available at birth is added later. There is less room for human error; the records are completed much sooner; the doctors, the state, and I are all much happier with the results."



Revvity – an 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.



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

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.

revvity