

revvity

# The EVOYA™ Platform.

## 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.

## EVOYA™ Platform Newborn Screening Software

EVOYA newborn screening management software has been designed to protect cloud-based applications, data, and infrastructure from unauthorized access, data breaches, and other security threats.

EVOYA security encompasses a range of technologies, protocols, and best practices designed to ensure your data's confidentiality, integrity, and availability.



## Key security features

With EVOYA, you can rest assured that your data is held to the highest security standards available, and data security will continue to evolve as any potential threat arises.



### Centralized security management

A dedicated team to check and keep EVOYA safe and secure



### Certified to the latest standards

Certified under the ISO/IEC 27001 information security management principles enabling HIPAA and GDPR information security standards



### Authentication and Access Control

Controlling who can access the software based on their roles and responsibilities, enforcing multi-factor authentication



### Data Encryption

Encrypting data in transit (outside and inside the cloud), between each of the application components (in the cloud), and in the database and storage



### Data Loss Prevention (DLP)

Preventing accidental or intentional exposure of sensitive information, detecting and blocking the transmission of sensitive data outside the organization, or encrypting data before it leaves the cloud environment



### Regular testing

External companies routinely test EVOYA's security to keep it robust

Speak to your local representative or visit [Revvity.com](https://www.revivity.com) for more information.

