



Data management, simply done.

The EVOYA Platform version 2.5 is now available

Making data easier to manage

The EVOYA™ platform v2.5 enables you to manage your laboratory's newborn screening data through Cloud-based solutions. The intuitive interface is designed to help you manage your data easily. Automatic import of test result data to the cloud is now available along with a redesign of the user interface. The latest version also allows for follow-up cases to be managed within the application

Interpret your results

Result files are automatically imported into the EVOYA system where interpretation reviews take place. Demographic information is available to be used in the interpretation logic. You can approve, adjust, or reject interpretation outcomes as well as provide additional text to be included on final reports — all with the touch of a button. Manual result entry is also available to manage test results which do not come from result file imports.

Generate reports

Once the workflow is complete, you can generate reports from easy-to-use templates (normal, abnormal, unsatisfactory). Additional flexibility with roll-up and roll-out logic for your panel of tests and results is now available. Your reports may be printed, downloaded, or retrieved remotely through the EVOYA platform, allowing you to manage your data flexibly. Email notifications may be set up to alert your remote users of reports available for download.

Review your data

Searches, patient summary and patient journey views help facilitate data review related to specimens, test orders, results, and reports. In addition, the ability to link specimens from the same patient together is now included. External users may also access EVOYA to enter demographics into the system remotely.



Technical information

- The EVOYA platform is designed as a microservice-based solution, supported by Cloud via Amazon Web Services (AWS). The EVOYA platform is compatible with Chrome, Firefox, and Edge chromium-based browsers.
- In addition, version 2.5 is available in English, Brazilian Portuguese, French, Italian, Canadian French, Arabic, German and Mexican Spanish languages.
- Design and development of the EVOYA platform is certified under ISO/IEC 27001 information security management principles.