

One informatics suite,
purpose-built for
genomic analysis.



revvity

Revvity Genomics Software

The gold-standard in NGS software.

To date, genomic software has helped labs to unravel the mysteries behind the DNA helix. However, the high computational demand inherent in DNA sequencing has led laboratories to make tradeoffs in their software selection, often choosing between ease of use, ease of implementation, or genomic specialization.

Revvity offers a suite of configurable, NGS-specific modules that can address all needs. With LIMS, DNA analysis, and interpretation capabilities, our software can amplify your current NGS research efforts or can jumpstart your lab's sequencing ambitions. Grounded in genomics and packed with a host of lab-agnostic features to synchronize with your current laboratory setup, we're here to simplify your science.

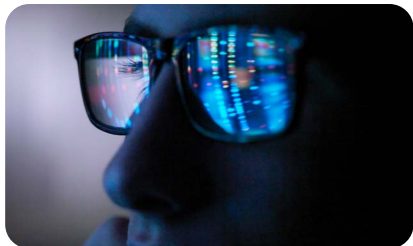
Discover how you can elevate your research with Revvity Genomics Software today.



Why select our software?



Engineered by laboratory specialists
Accessible for labs of all backgrounds, designed for secure genomic analysis.



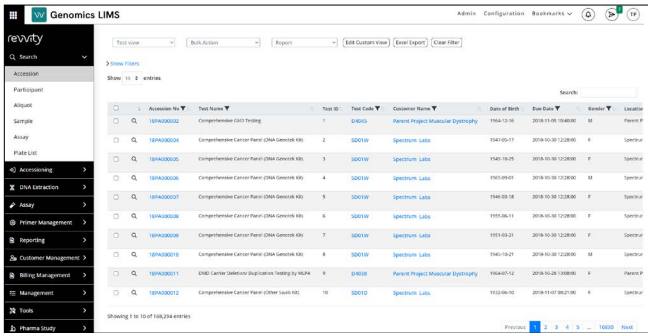
Primed for deep genomic insight
Carefully curated genetic panel of 390+ genes and 100K+ variants.



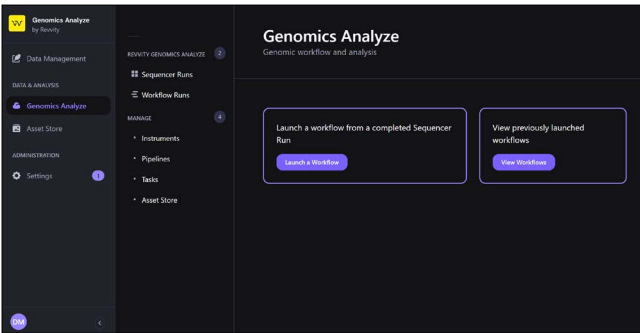
Configurable and adaptable for laboratories
Compatible with lab's existing software, multiple assays and application areas.*

* As 3rd party software will vary in orientation, please speak with your local Revvity representative to confirm its compatibility.

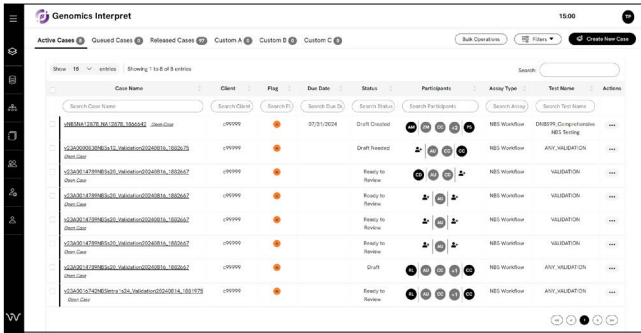
Key components



Revvity Genomics LIMS



Revvity Genomics Analyze



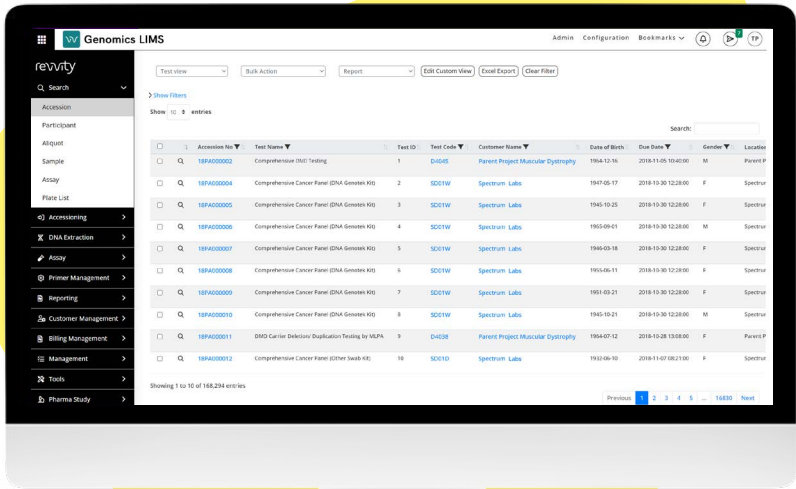
Revvity Genomics Interpret

Revvity Genomics LIMS

Create efficiency inside your lab.

Key features

- Workflow and assay management
- Sample tracking
- Aliquot management
- Tag and ticket management
- Index management
- Global accessibility
- Billing functionality

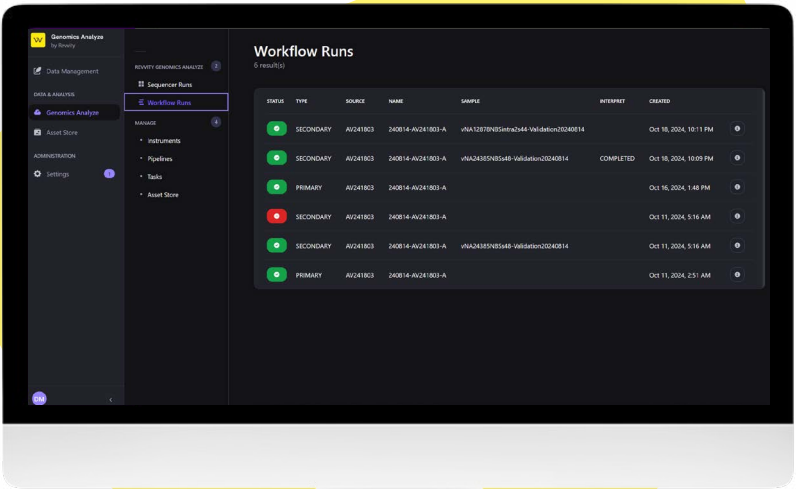


Revvity Genomics Analyze

Centralized hub for primary and secondary analysis.

Key features

- Primed for high data volumes
- Robust QC steps for assurance
- Simplified user interface
- Accommodates multiple pipeline templates
- Global accessibility



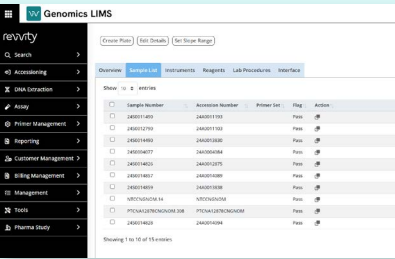
Customize all parameters for primary and secondary analysis, facilitating easier management and categorization across your research cases. Backed with a precise quality control system, bringing confidence to your calculations.

A robust LIMS system designed specifically for genetic analysis. Refined over years of use across the diverse Revvity Omics laboratory network, reimaged for laboratories across the world.

End-to-end connectivity for assurance

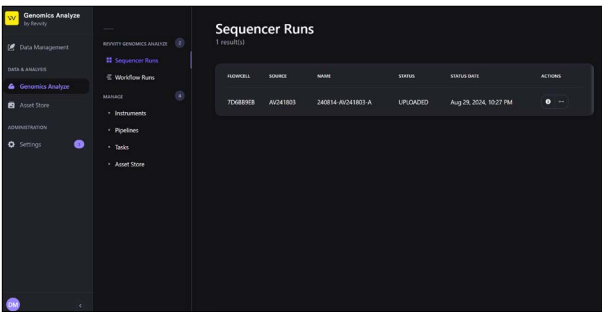
- ✓ Accessioning
- ✓ DNA extraction
- ✓ Sequencing & assay
- ✓ Data capture
- ✓ Final documentation
- Delivery of findings

Centralized hub for lab activities

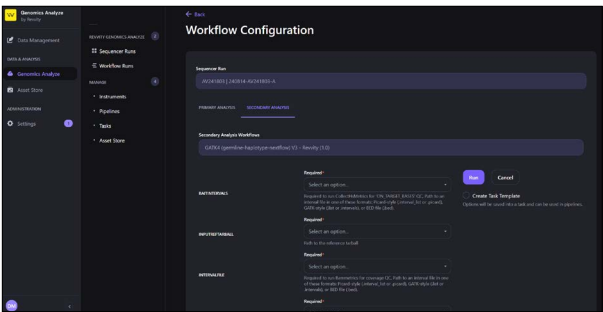


Built to maximize compatibility

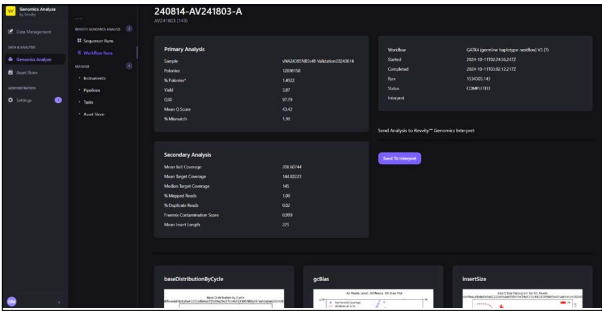
- Support multi-assay analysis
- Analysis workflow management
- Batch processing
- Extensive plate management
- Sample journey across test, assay, batch, plate and aliquot levels



Consolidated overview of sequencing and workflow runs



Create templates for running workflows



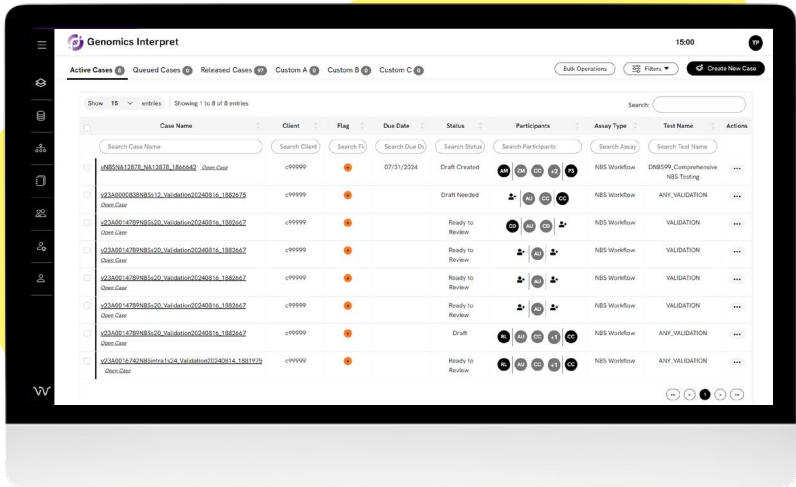
Robust QC steps for assurance

Revvity Genomics Interpret

Deeper insight, actionable findings.

Key features

- Automatic variant annotation
- Case ownership management
- Documentation templates
- Direct data input from Analyze
- Integration with LIMS
- Global accessibility



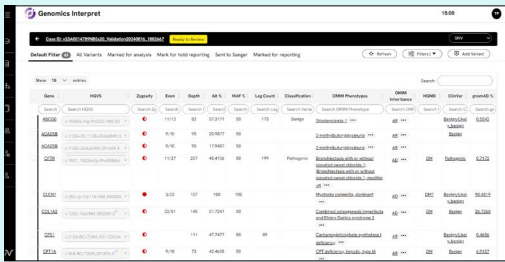
Introducing Revvity’s end-to-end NGS workflow

If your laboratory is also seeking an NGS workflow, we’d be thrilled to share our configurable and complete NGS solution - compatible with both targeted sequencing and WGS methods. Built on the same underlying principles as our software suite, see how you can amplify your sequencing aspirations alongside the global leaders in newborn screening today.

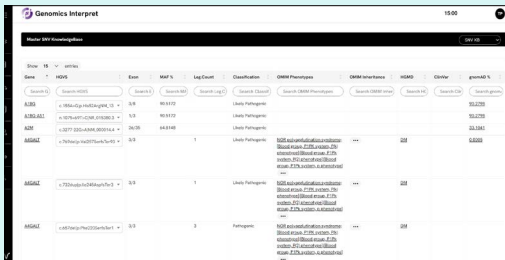
Optional connection

Required connection

The crown jewel of Revvity’s NGS software suite. Features an expert-curated panel of 390+ genes for labs of all backgrounds, spanning 17 research areas (e.g. endocrinology). Backed by an ever-growing database of 100K+ variants, updated with new findings from 3rd party and Revvity laboratories. Become a pioneer in your field of science.



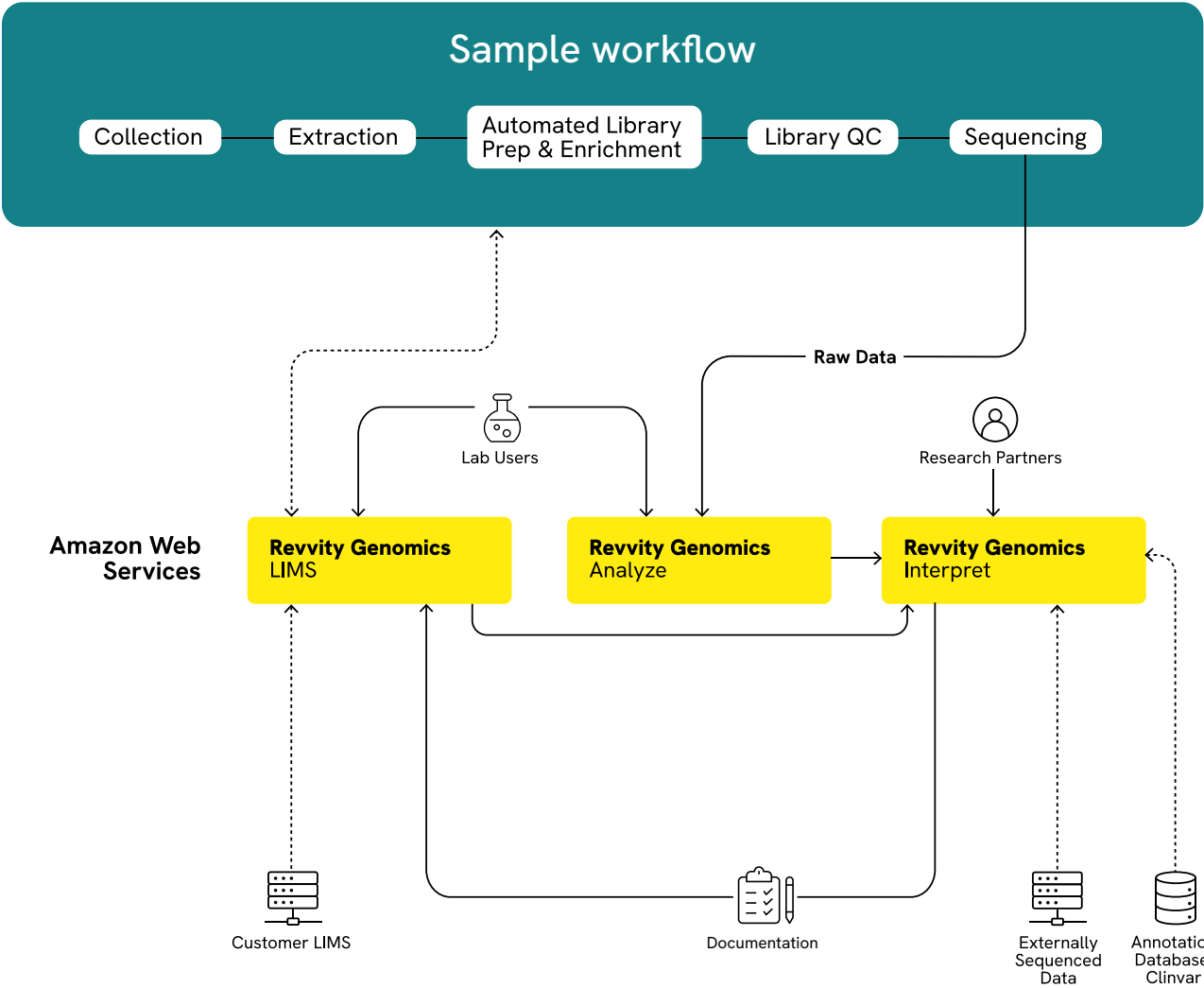
Simplified case management



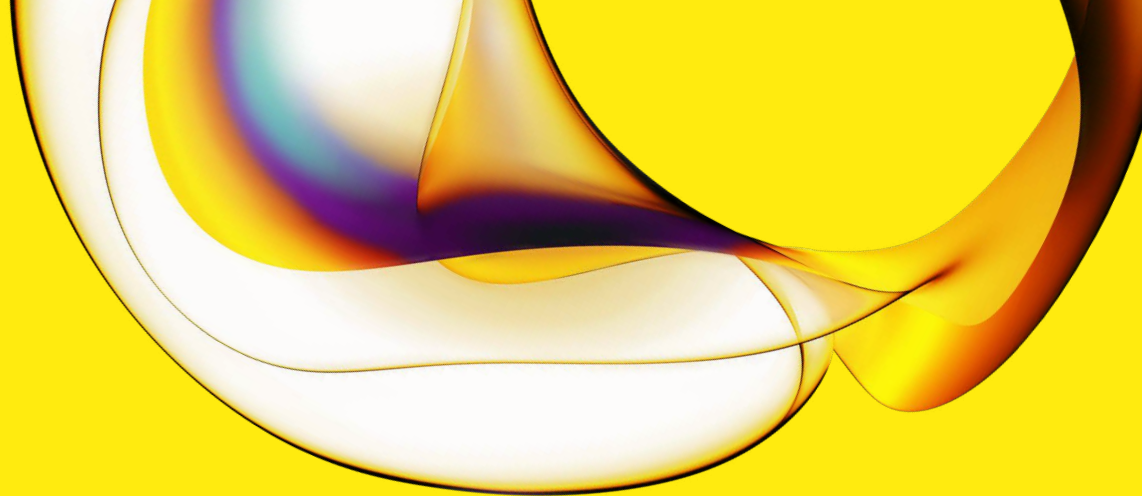
Variant database enables faster assessment



Customizable set of templates for exploring



The above diagram reflects a full implementation of Revvity Genomics Software. Please reach out to your local representative to discuss alternative arrangements.



[Learn more](#)

Scan or click
to learn more.



[Contact us](#)

Scan or click
to contact us.

www.revvity.com

revvity