

The future of NGS software



revvity

**A modular & customizable software
suite, designed for genomic analysis.**

For research use only. Not for use in diagnostic procedures.

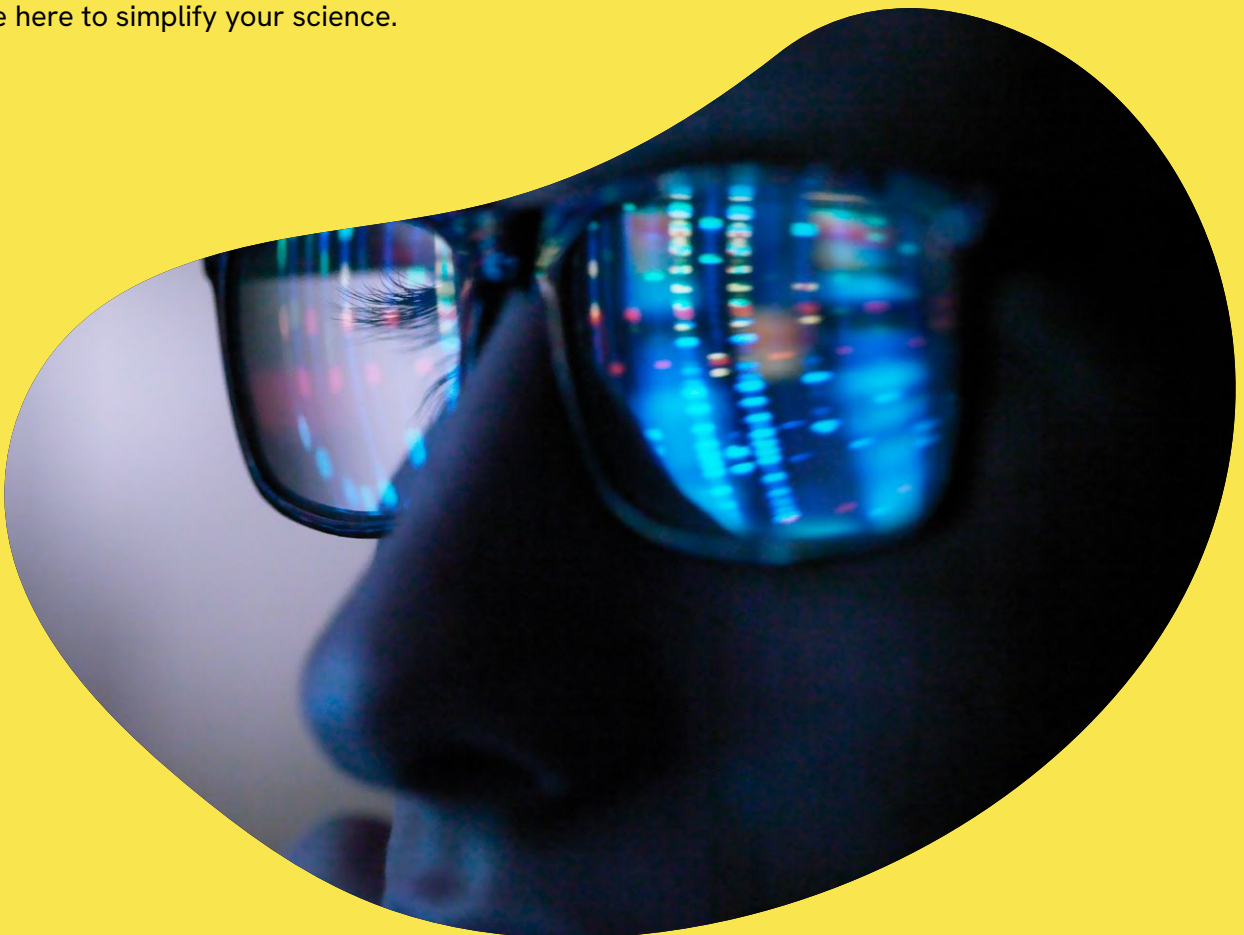
The gold-standard in NGS software.

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.

Introducing a suite of configurable, NGS-specific software that covers all steps. 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's
NGS software suite today.**

[Contact Us](#)



Why select our software?



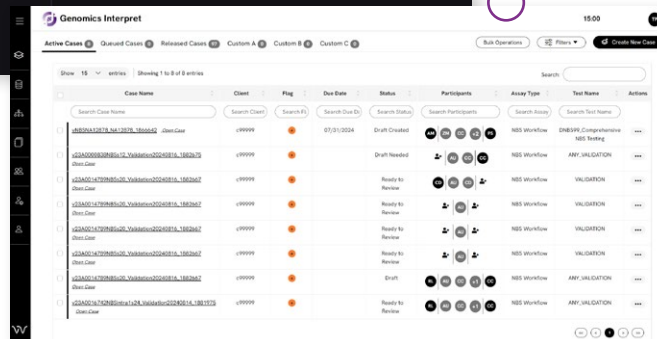
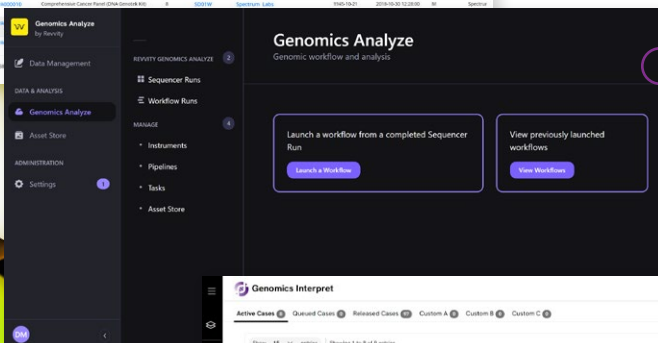
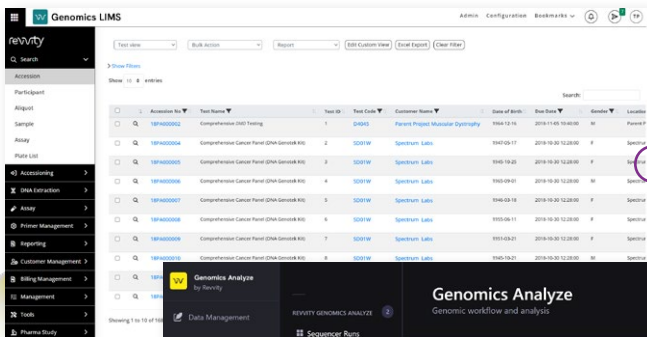
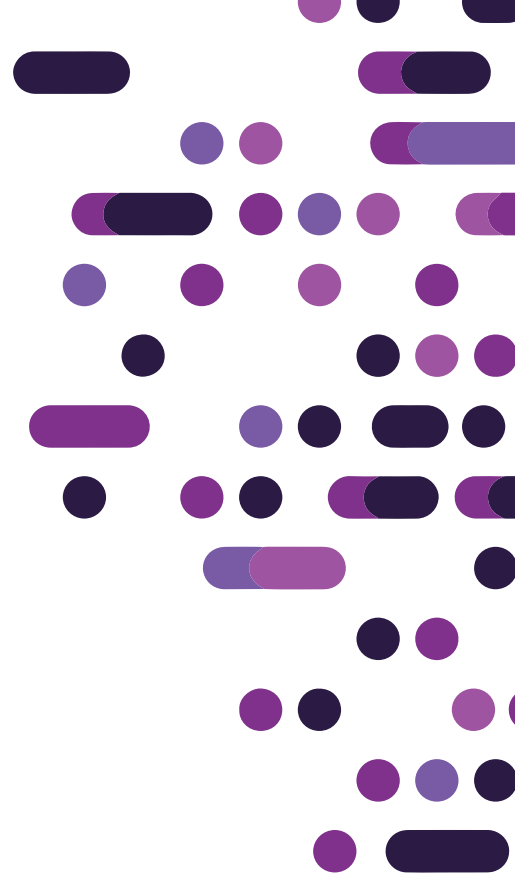
Engineered by laboratory specialists
Accessible for labs of all backgrounds,
designed for secure genomic analysis.



Primed for deep genomic insight
Carefully curated genetic panel of 390+ genes
and 87K+ variants.



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



Key Components

Revvity Genomics LIMS

Revvity Genomics Analyze

Revvity Genomics Interpret

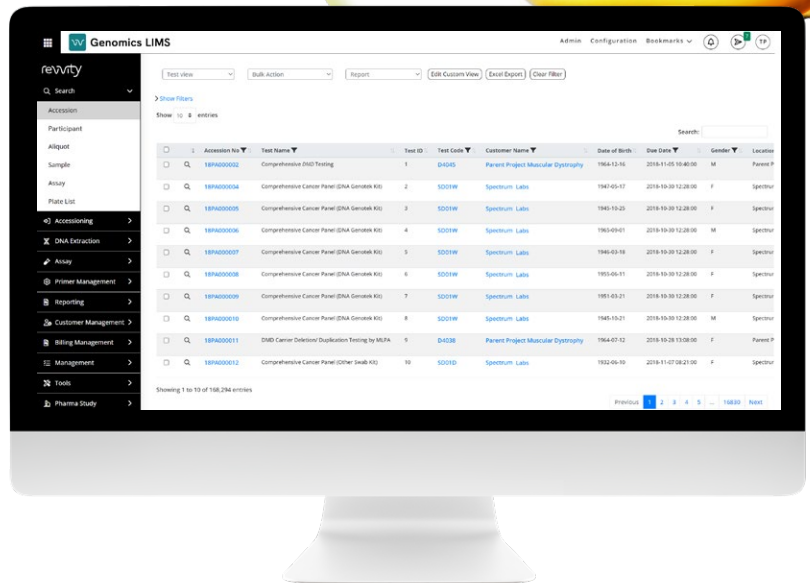
* As 3rd party software will vary in orientation, please speak with your local Revvity representative to confirm its compatibility.

Revvity™ Genomics LIMS Software

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

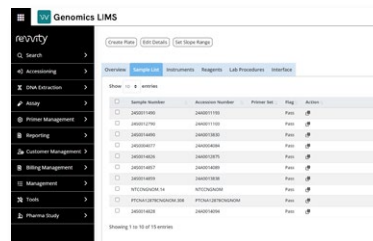


A robust LIMS system designed specifically for genetic analysis. Refined over years of use across the diverse Revvity Omics laboratory network, reimaged for laboratories across the world.

END-TO-END CONNECTIVITY FOR ASSURANCE

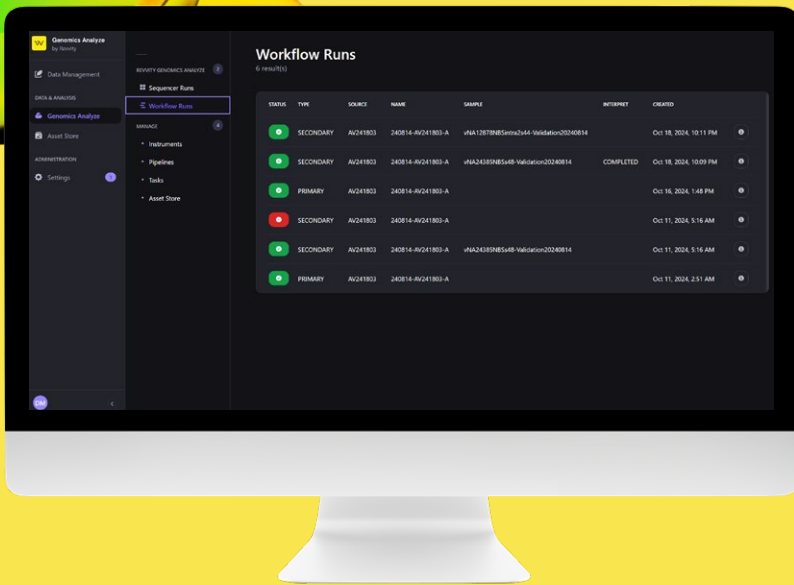
- ✓ Accessioning
- ✓ DNA Extraction
- ✓ Sequencing & Assay
- ✓ Results Capture
- ✓ Report Generation
- Report Delivery

CENTRALIZED HUB FOR LAB ACTIVITIES



BUILT TO MAXIMIZE COMPATIBILITY

- Support Multi-Assay testing
- Flexible workflow management
- Batch processing
- Extensive plate management
- Sample journey across test, assay, batch, plate and aliquot levels



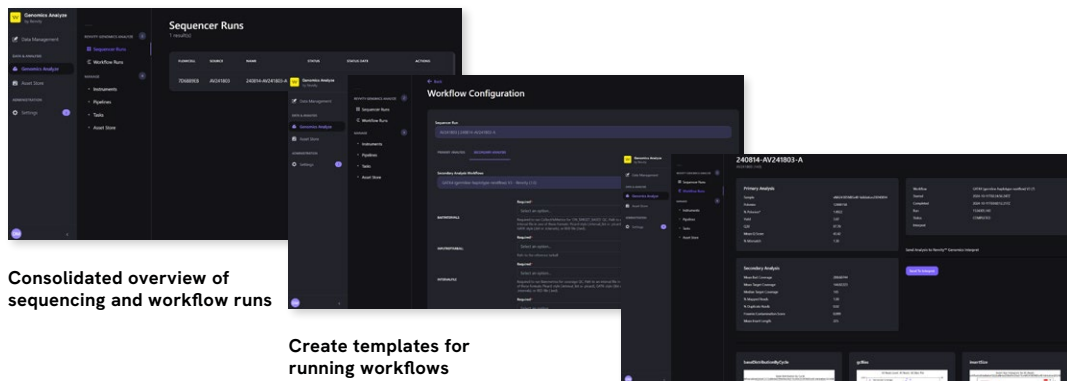
Revvity™ Genomics Analyze Software

Centralized hub for primary & 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.



Consolidated overview of sequencing and workflow runs

Create templates for running workflows

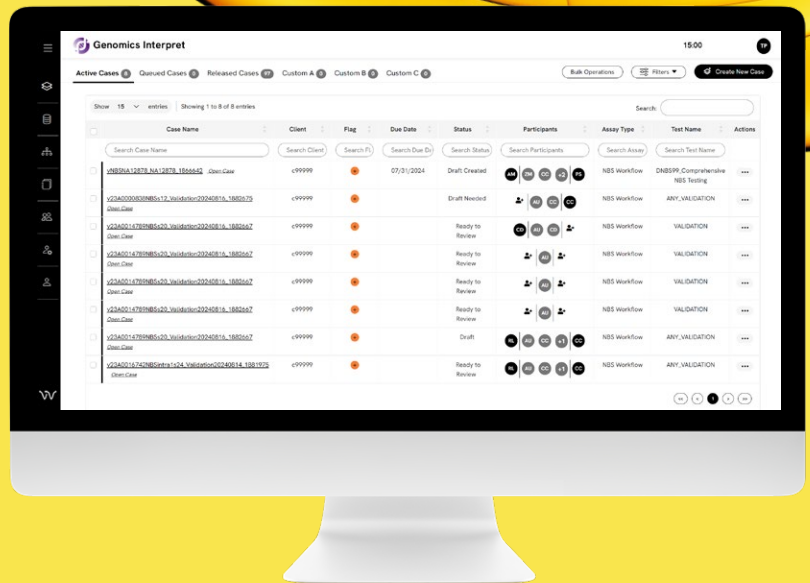
Robust QC steps for assurance

Revvity™ Genomics Interpret Software

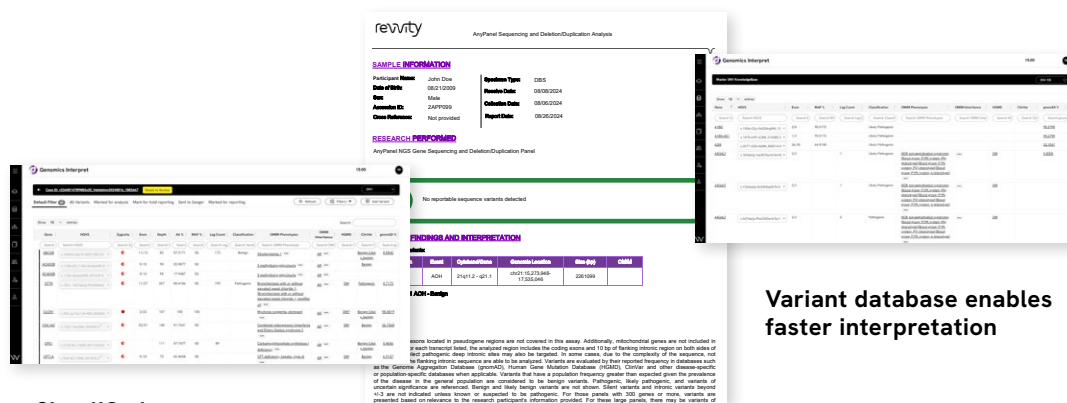
Deeper insight, actionable reporting.

Key Features

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



The crown jewel of Revvity's NGS software suite. Features a panel of 390+ genes for labs of all backgrounds, spanning 17 research areas (e.g. endocrinology). Backed by an ever-growing database of 87K+ variants, updated with new findings from 3rd party and Revvity laboratories. Become a pioneer in your field of science.



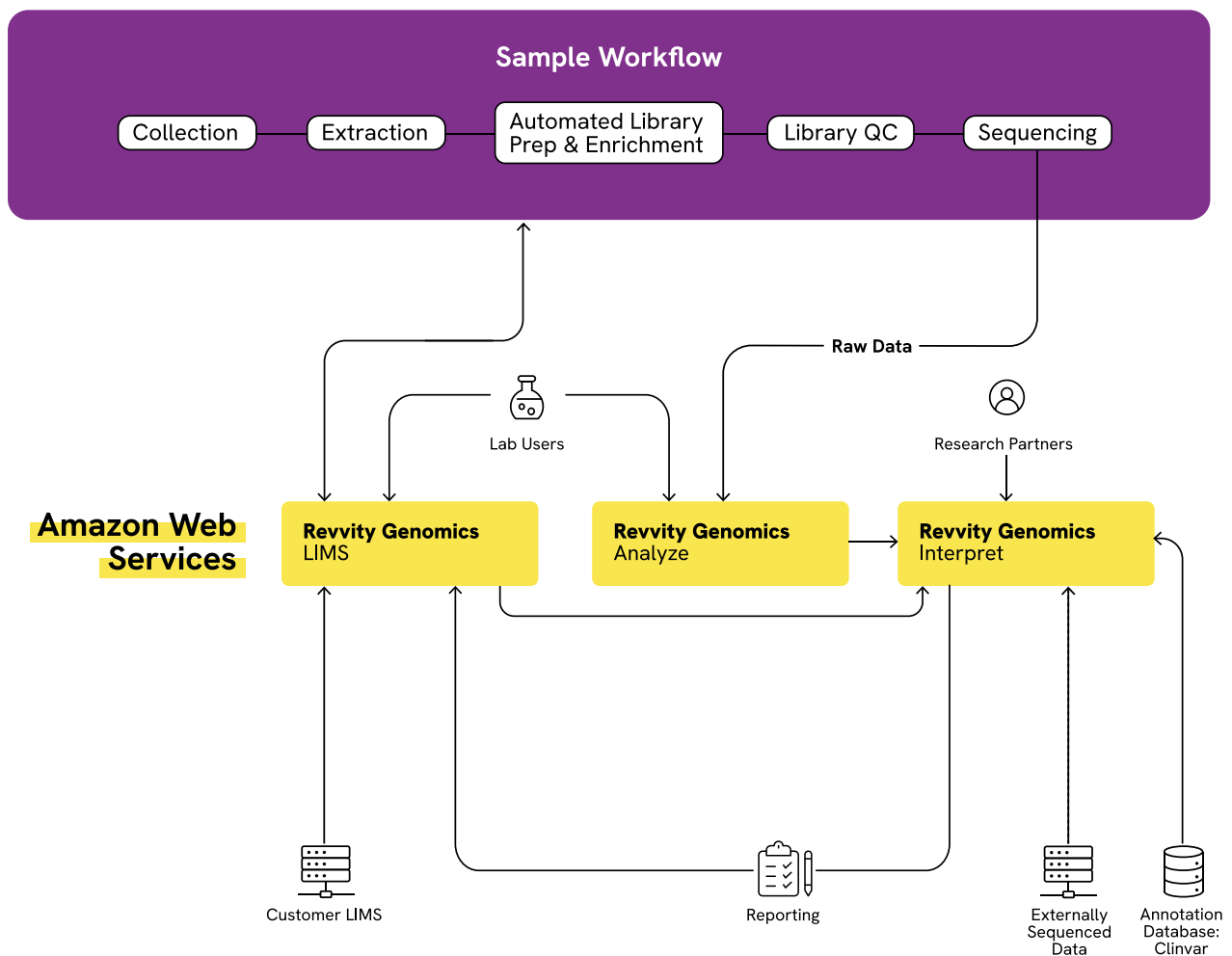
Simplified case management

Customizable set of templates for exploring

Variant database enables faster interpretation

Introducing Revvity's end-to-end NGS workflow

If your laboratory is also seeking an NGS workflow, we'd be thrilled to share our customizable and complete NGS solution. 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.



The above diagram reflects a full implementation of Revvity's software suite. Please reach out to your local representative to discuss alternative arrangements.



Want to learn more?
Request a software
demonstration from
your local Revvity
representative.

Contact Us

revvity

Revvity, Inc.
940 Winter Street, Waltham, MA 02451 USA
(800) 762-4000 | www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2024, Revvity, Inc. All rights reserved.

For research use only. Not for use in diagnostic procedures.

www.revvity.com