

The NGS software designed by a lab, for your lab.

A complete and customizable software suite, built for genetic research.



Ignite your lab research.

Introducing the modular NGS software that covers all bases. Complete with LIMS, DNA analysis and interpretation capabilities, our software can elevate your existing NGS workflows or kickstart your sequencing aspirations. No matter where you begin, Revvity is here to support.

Discover how you can achieve your laboratory sequencing goals with Revvity today.

[Learn More](#)

Why select our software?



Engineered by lab specialists

- Designed specifically for genomic analysis
- Caters to labs of all backgrounds
- Backed by the global leaders in newborn screening



Primed for deep genomic insight

- 390+ carefully curated genes
- Backed by a database of 87K+ variants
- Scalable for targeted & whole genome research



Customizable & adaptable NGS solution

- User-defined, lab-friendly configuration
- Compatible with existing software arrangements
- Supports multiple assays and workflows

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

revvity