



A complete and configurable informatics suite, built for accessible genetic research.

Revvity Genomics Software.

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

Why select our software?

Engineered by lab specialists

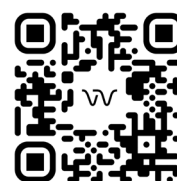
- Designed specifically for genomic analysis
- Caters to labs of all backgrounds
- Backed by the global leaders in newborn screening

Primed for deep genomic insight

- 390+ carefully curated genes
- Backed by a database of 100K+ variants
- Scalable for targeted & whole genome research

Adaptable NGS solution

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



[Learn more](#)

Scan or click
for more information



[Contact us](#)

Scan or click
for more information