



A streamlined approach to genotyping with low-pass WGS.

NEXTFLEX HT Agrigenomics Low-Pass WGS kit

Low-pass whole genome sequencing (LP-WGS) offers a cost-effective approach to sequencing entire genomes at coverages ranging from 0.1x to 10x. This technique, followed by imputation, is becoming the preferred method for genotyping among plant and animal breeders.

Revvity and CURIO™ are now offering a comprehensive solution including reagents, automation, and analysis for your WGS data, addressing some of the main challenges found by breeders.

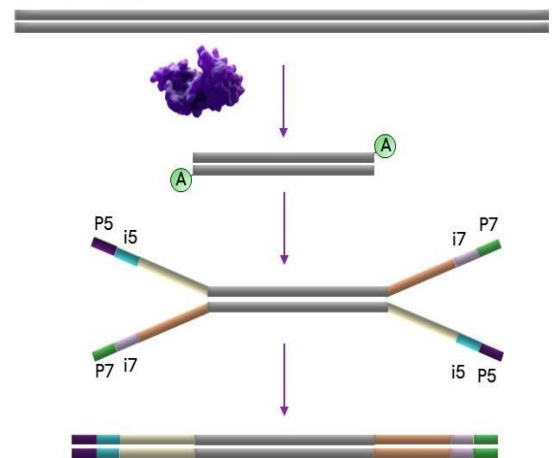
The NEXTFLEX™ HT Agrigenomics Low-Pass WGS Kit is a new, streamlined workflow compatible with Illumina® and Element® Biosciences platforms. With a wide input range, it is suitable for all kind of plant samples and can be run either manually or in automated workstations. The libraries were homogenized and extracted with Revvity solutions.

Key features

- **Robust** - fast and easy protocol
- **Wide input range** - from 1 ng to 500 ng
- **PCR-free** - compatible workflow
- **High throughput** - up to 1,536 UDI barcodes



From early research to large-scale daily operations of breeders, CURIO provides best-in-class imputation analysis for plant and animal breeders with speed and scalability found in no other agrigenomics platform.





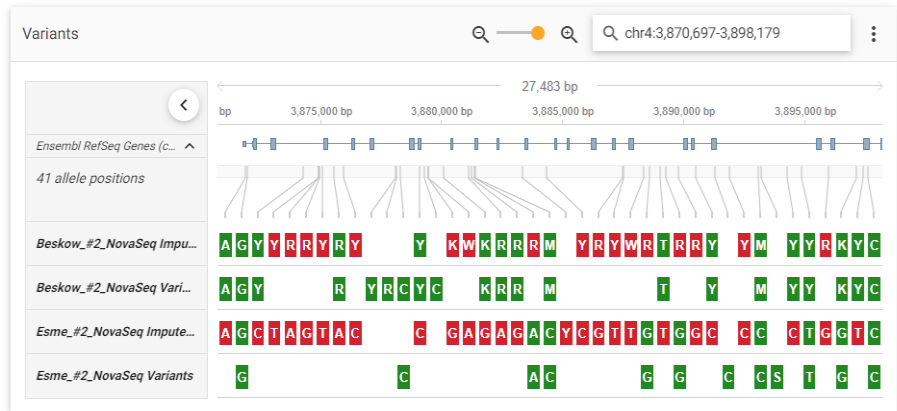
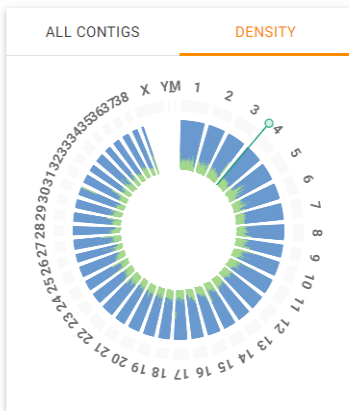
- **Speed** - Better results in minutes, not hours or days
- **Versatility** - Works with any species with a reference
- **Flexibility** - Massively scalable and cost-effective cloud-based solution
- **Transparency** - Fully transparent and validated best-in-class algorithms
- **Convenience** - Build your own reference panel using a simple interface
- **Efficiency** - Simple and easy integration for fully automated workflows
- **Approach** - Partner-based approach ensures successful outcomes

Next-generation bioinformatics: LP-WGS + imputation

Curio Genomics is a bioinformatics software company focused on accelerating agriculture genomics research.

Using proprietary ultra-parallelized cloud architecture, the CURIO platform enables unprecedented data processing speed of large and complex genomic data sets.

Through the platform's intuitive user interface, breeders can process, organize, visualize, and interpret their data independently, resulting in faster, more informed breeding decisions.



From sample to insights, Revvity + CURIO enables simpler, more efficient agrigenomics with better results for more informed data-driven breeding decisions.

Scan or click to learn more about this solution for Low-Pass WGS for agrigenomics today.



Scan or click to learn more about Revvity + CURIO today.

