



End-to-end whole genome sequencing metagenomics solutions

Power your research

Whole genome sequencing (WGS) provides a complete picture of the genetic material in a microbial community, allowing for the identification of all microorganisms present. By analyzing the entire genome, researchers can gain insights into the functional capabilities of the microbial community, such as metabolic pathways, resistance genes, and potential virulence factors. This information is difficult to obtain with other approaches such as 16S rRNA or targeted sequencing.

Revvity and CosmosID® are now offering a comprehensive solution including reagents, automation, and analysis for your WGS data.

The NEXTFLEX® Rapid XP V2 DNA-seq Kit for metagenomics is a streamlined workflow for WGS on Illumina® and Element® Biosciences platforms. With a wide input range, it is suitable for all kind of samples. Libraries produced using this kit are normalized during library construction using proprietary NEXTFLEX® normalization beads, delivering consistent mass and sequencing cluster density for all samples in a library pool. Up to 1,536 barcodes for multiplexing available for purchase separately.

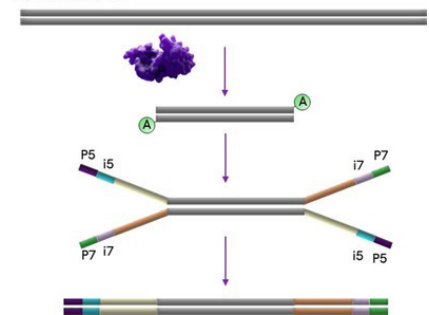
The combination of the robust, fast and easy library preparation workflow from Revvity with the speed and superior performance of the AI-driven CosmosID® platform will provide a head start to your next metagenomic project.



Key features

- Access to KEPLER™ & CHAMP™ profilers
- Quantifies species abundance
- Superior strain level resolution
- Compare microbiome data in minutes
- Easy to interpret charts and visualizations
- No coding experience required

Genomic DNA



Two cutting-edge pipelines for best results

CHAMP - Human Microbiome Profiler



- Employs a database of > 30,000 microbiome samples obtained from humans
- Identifies low-abundance species within samples.
- Quantify species abundance
- Phage and virome profiling
- High resolution annotation at the species level

KEPLER - Host Agnostic Profiler



- Designed for analysis of environmental, animal, soil, food sample and many others
- Best-in-class low abundance taxa detection
- Quantify species abundance
- Annotates microorganisms at the species level
- >600 peer-reviewed publications to date

Explore the next generation of library prep solutions

NEXTFLEX® Rapid XP v2

- Robust, fast and easy protocol
- Wide input range - from 100 pg to 1 µg
- Possibility of PCR-free workflow
- Proprietary bead-based normalization
- Up to 1,536 UDI barcodes

Kit information

[NEXTFLEX Rapid XP v2 Kit for Metagenomics](#)



<https://cosmosidhub.com/>

revvity