



End-to-end metagenomics solutions.

Power your research

By utilizing Next-Generation Sequencing (NGS), entire communities within a sample can be efficiently analyzed from minimal quantities of DNA. Revvity and CosmosID® are now offering a comprehensive solution including reagents, automation, and analysis for your 16S data.

Revvity offers a robust NEXTFLEX™ portfolio for 16S rRNA applications. We offer three distinct 16S rRNA amplicon panels that have been designed to target the 16S rRNA V1-V3, V3-V4 and V4 hypervariable regions, respectively. Maximize the economy and speed of targeted sequencing for the metagenomic needs of your laboratory.

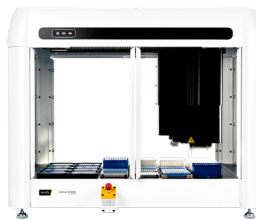
Analysis powered by CosmosID-HUB®

These panels are now bundled with access to CosmosID-HUB®, an online software solution that enables fast and easy analysis of complex microbiome data. CosmosID-HUB® gives scientists user-friendly access to version-controlled and validated 16S pipelines. The software also enables rapid data interpretation through a comparative analysis software with features such as tables, heatmaps, bar charts, multiple Alpha and Beta Diversity indexes, abundance distribution plots, differential abundance testing, as well as comprehensive statistics between groups.



Key features

- Characterize any sample type.
- Protocols automated on Revvity's suite of liquid handlers to increase your throughput.
- DADA2 & PICRUST2 bioinformatics pipelines for taxonomic and functional analysis
- Advanced machine learning to differentiate signal from noise
- Create your own reports and publication-ready visualization.



Automated on the Revvity Sciclone™ G3 NGSx workstation

Let us help automate your metagenomics

NEXTFLEX 16S rRNA Amplicon-Seq kits

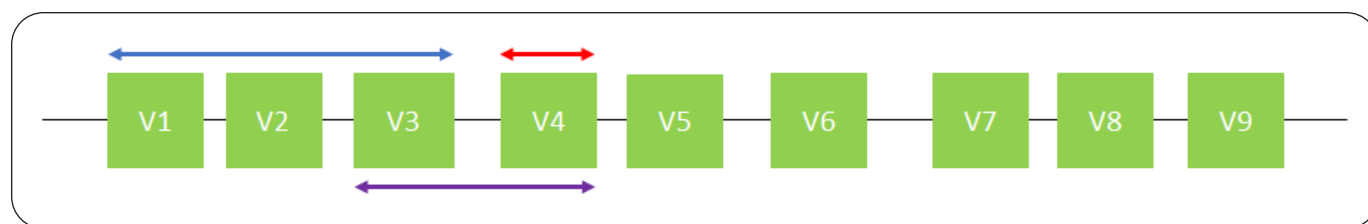


Figure 1: Schematic representation of targeted regions covered by the NEXTFLEX 16S series of amplicon-seq kits.

CosmosID-HUB®: Automated taxonomic and functional profiling with interactive visualizations and no bioinformatics experience needed.



Kit information

Kit name
NEXTFLEX 16S V1-V3 amplicon-seq kit
NEXTFLEX 16S V4 amplicon-seq kit 2.0
NEXTFLEX 16S V3-V4 amplicon-seq kit



cosmosidhub.com

revvity