



revvity



Quick Guide to CosmosID-HUB WGSData Upload

Redeeming your CosmosID-HUB Credits

1. Go to www.cosmosidhub.com/revvity
2. Complete the form, including your unique Kit ID (you may enter multiple).
3. A member of the CosmosID-HUB team will reach out to complete your onboarding.
4. Once onboarding is completed, HUB credits will be issued to your account.
5. Upload your shotgun data and run your microbiome study for up to 60 days.

WGS Data Upload Requirements

- WGS profiling workflow supports *paired-end* sequencing data only.
- Sequencing data files should not exceed a size of 50 GB per sample, in fastq.gz format.
- For each sample ID, two paired-end fastq files must be uploaded with "R1_001 " or "R2_001 " followed by the sequencing suffix (e.g., *Sample0123_R1_001.fastq.gz* + *Sample0123_R2_001.fastq.gz*)
- File names should not have any spaces and special characters in them.
- Maintaining an average base quality score >PHRED 20 is recommended across all individual reads inside a fastq file.

If you have any questions, please reach out at ngs@revvity.com