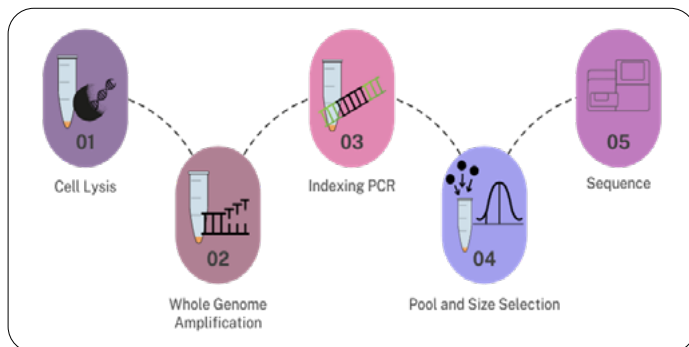




End to end solution for PGT-A.

PG-Seq™ Rapid v2 Kit was developed to detect whole chromosome aneuploidies and structural arrangements (such as unbalanced translocations and segmental errors) down to 7 Mb in ~3 hours using a robust and streamlined workflow.



| Figure 1: PG-Seq™ Rapid v2 workflow.

- Ready-mixed reagents. Minimal tube openings and pipetting step.
- With up to 384 barcodes available, suitable for labs from any size, from low to high-throughput.
- Detection of sample swapping and external DNA contamination with no additional effort.

- Compatible with all Illumina sequencers and also with Element Biosciences. Full control of your throughput by selecting the flow cell that fits best your need.
- Only 250,000 reads per sample required to achieve 10Mb of resolution. Run more samples in the same batch.
- Automated in the Sciclone G3 NGSx workstation.
- PG-Find software included in the kit:
 - **Chromosome copy number detection:** it identifies chromosome copy number changes in the human genome, which is crucial for detecting aneuploidies.
 - **Compatibility:** the software is compatible with Illumina® and Element Biosciences® sequencing platforms.
 - **Automatic analysis:** it provides automatic calling of aneuploidy and copy number variants, streamlining the analysis process.
 - **User-friendly interface:** the software includes features for easy data loading, analysis, and visualization.
 - **High accuracy:** it has been extensively tested for accuracy in detecting whole chromosome aneuploidies and sub-chromosomal abnormalities.

- mtDetect software included in the kit:
 - **Mitochondrial DNA analysis:** it extracts and analyzes mtDNA information to monitor external DNA contamination and detect sample swapping or mislabeling.
 - **Integration with PG-Seq™ Rapid v2:** the software works seamlessly with the PG-Seq™ Rapid v2 workflow.
 - **Sample grouping:** the software creates identity matrices and sample groupings to ensure accurate embryo classification.

Part number	Description
4340-0248	PG-Seq Rapid kit v2 (48 rxn)
4341-IL48A	PG-Seq Indexing Primers, Illumina (BC1-48)
4341-IL48B	PG-Seq Indexing Primers, Illumina (BC 49-96)

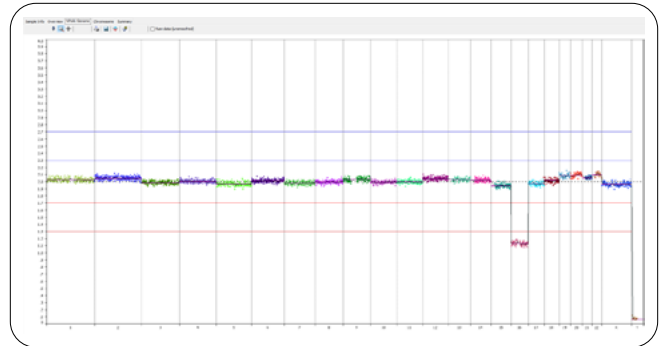


Figure 2a: D5 embryo biopsy showing chr16 monosomy.



Figure 2b: Zoom into the affected chromosome.



revvity