

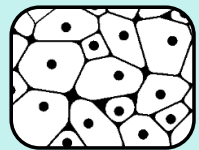


# Reveal the secrets hidden in complex samples.

## Get the answers you need with single-cell sequencing

There are 40 trillion cells in the human body. Each cell type provides a different function, and each cell has its own evolving agenda. Single-cell analysis unravels these cell communities and identifies critical information often overlooked by bulk approaches. Ultra revealing single-cell sequencing data is becoming more accessible due to technology advances, which are simplifying experiments and reducing costs.

**Single-cell multiomics enables transcriptome-wide gene expression analysis to be combined with protein detection at single-cell resolution. This allows researchers to:**



Investigate sample heterogeneity



Examine different development states



Identify rare cells



Investigate existing treatments



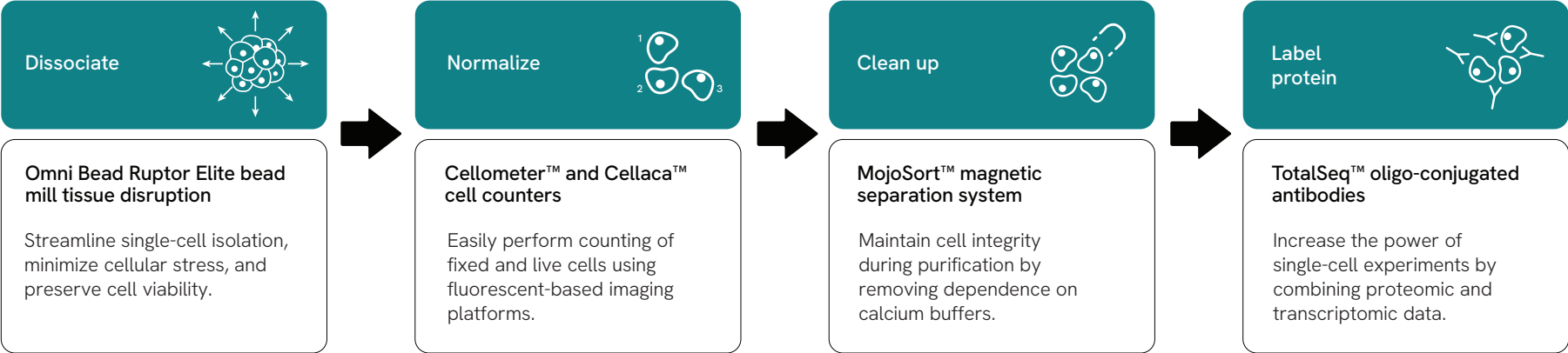
Discover new targets and biomarkers

Keep reading for solutions at every step of your experimental process ➡

# Revvity solutions for key steps of your single-cell sequencing journey.

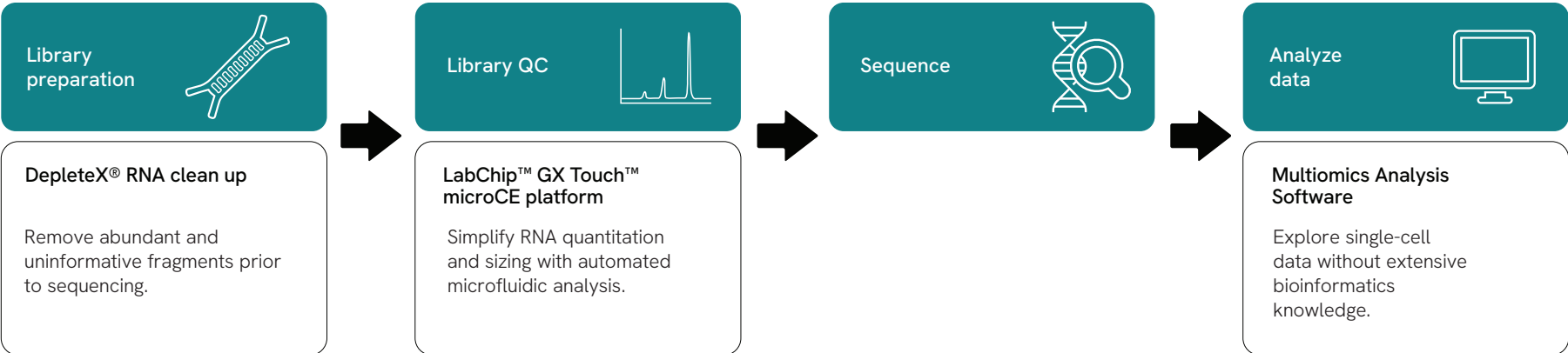
Specifically designed to work alongside your existing processing steps, making implementation simple.

## Upstream sample preparation



## Partition and barcode

## Downstream fragment analysis



Scan to learn more.

