

# Elevate your single-cell sequencing: Beat the stats.

Despite single-cell sequencing demonstrating potential as a valuable technique to identify new biomarkers and therapeutic targets, wide-scale adoption may be limited by high costs and noisy data.

Discover some simple strategies to help overcome these obstacles.



## Expand tissue preparation throughput by 12x

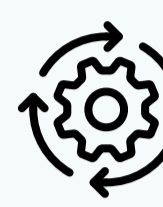
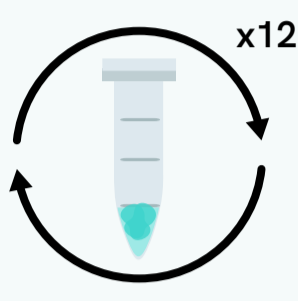
Dissociate multiple samples simultaneously to scale studies and expedite time to results.



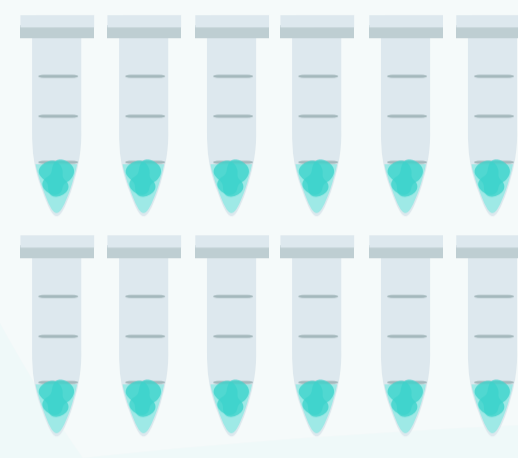
**Did you know?**  
It typically takes ~2 hrs to dissociate each tissue sample manually.



Manual repeated processing

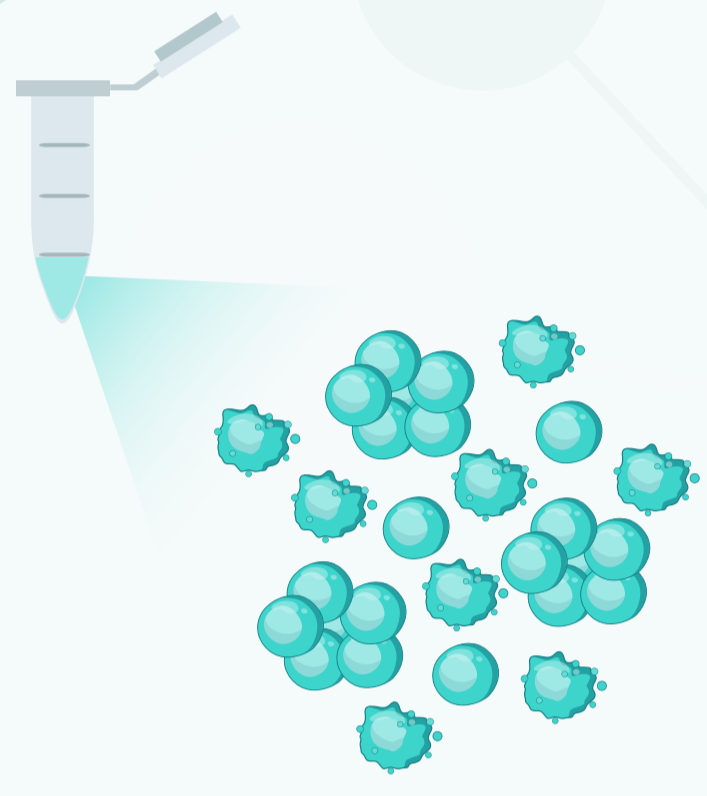


Parallel processing

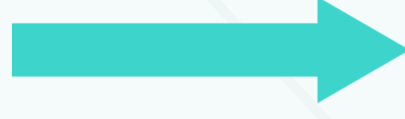


## Confirm >90% cell viability

Enrich for your desired cell populations to generate high-quality data.



Remove unwanted cells from suspension



Identify and filter cell clumps and debris



Count viable cells to confirm sample quality



**Did you know?**  
Cell viability can be as low as 20% post dissociation.

## Lower single-cell sequencing costs by ~40%

Increase value in sequencing runs by multiplexing samples into fewer lanes.

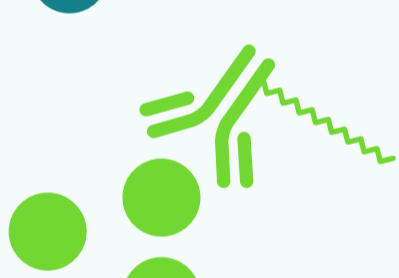


**Did you know?**  
Depending on technology used, the cost of a single-cell sequencing kit can range from \$1.5k - \$17k per run.

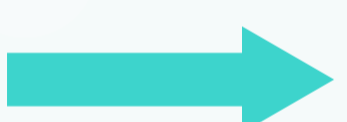
Sample 1



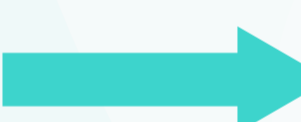
Sample 2



Hashtag



Pool



Sequence

## Increase usable mapped reads by 100%

Reduce noise in genomic libraries prior to sequencing by removing many uninformative fragments.



**Did you know?**  
Genomic library contains just **25%** usable reads.



2-fold increase in usable reads



Remove uninformative fragments:

- Ribosomal
- Mitochondrial
- Non-variable
- Non-transcriptomic

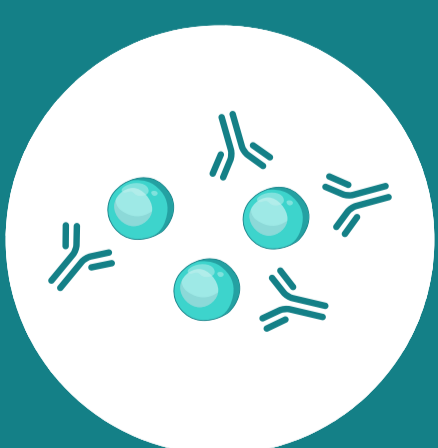


## Products to turn your bottlenecks into breakthroughs



### Sample preparation

- Omni Bead Ruptor Elite™ sample homogenizer
- Cellometer™ cell counters
- MojoSort™ dead cell removal



### Multiplex analysis

- TotalSeq™ oligo labeled antibodies for multiplex and multiomic analysis



### Data resolution

- DepleteX® RNA boost technology

For more information, visit: [www.revvy.com/single-cell-sequencing](http://www.revvy.com/single-cell-sequencing)

For research use only. Not for use in diagnostic procedures.

Copyright ©2024, Revvity. All rights reserved. Revvity is a registered trademark of Revvity.

1485044

revvity