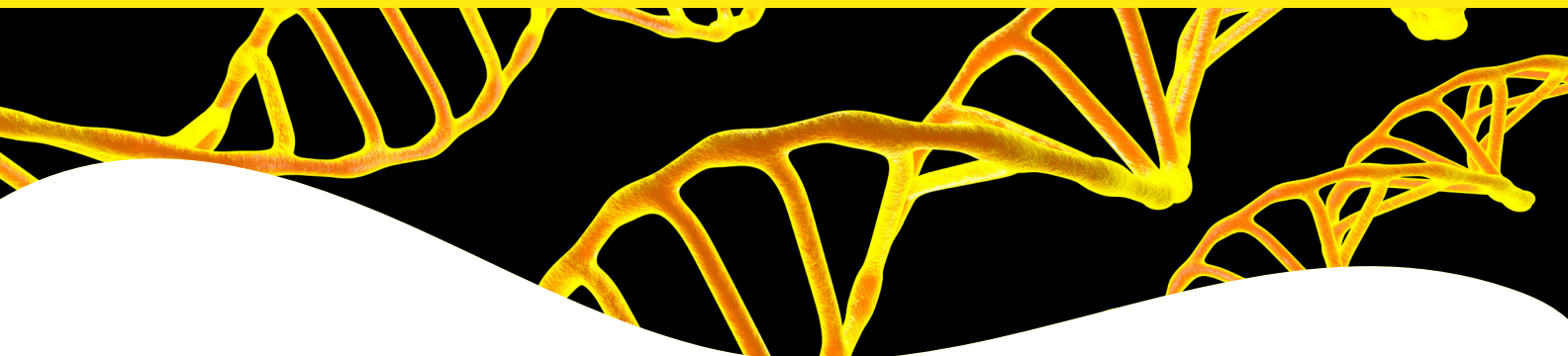


Sample homogenization solutions



More processing. Less programming.

LH96

- Fully automated homogenization with integrated liquid handling
- Automate liquid handling, pipetting, liquid transfer, weighing, diluting, and sample transfer
- Process up to 96 samples in a batch
- Rapid, repeatable, reliable results



Bead Ruptor 96+

- Multipurpose homogenization, offering a blend of traditional milling and high-throughput capabilities.
- Sealed jar milling, well plate or tube homogenizing in one unit
- Variable speeds up to 30 Hz
- Digital controls ensure reproducibility
- Homogenize in tubes ranging from 2 - 50 mL
- Large processing swath ensures high impact forces



Bead Ruptor Elite

- Fastest and broadest range of multi-sample processing up to 8 m/s
- No cool-down between runs
- Touch screen interface
- Homogenize in 0.5, 2, 7, 15, 30 or 50 mL tubes
- Minimize thermal degradation with Cryo cooling



Bead Ruptor 24

- Fixed 2 mL tube carriage processing
- Touch screen interface
- Minimize thermal degradation with Cryo cooling
- Speeds up to 6.5 m/s
- Built on the tried and tested Elite platform



Bead Ruptor 12

- Process up to 16 tubes at a time
- Variable speeds up to 6 m/s
- Homogenize in 0.5, 2 or 7 mL tubes
- No cool down between runs
- Near-linear motion for the highest processing efficiency



OM500

- 500 Watt Ultrasonic Homogenizer
- Include pulse mode, timer, amplitude control, and a digital display of both wattage and joules
- Process samples up to 1 L in volume
- Six different probes available
- Sound abatement chamber available to mitigate processing noise



Bead Ruptor 4

- Homogenize in 0.5, 1.5, 2 or 7 mL tubes
- Unique inclined carriage
- Small footprint
- Variable speeds up to 5 m/s



OM100

- 100 Watt Ultrasonic Homogenizer
- Include pulse mode, timer, amplitude control, and a digital display of both wattage and joules
- Programmable operation - Set time and amplitude for hands free operation
- Ideal for small samples up to 50 mL
- Three different probe sizes available



Cryo Unit

- Compatible with the Bead Ruptor Elite and Bead Ruptor 24
- Minimizes thermal degradation by cooling the processing chamber
- Protect heat sensitive analytes such as RNA and proteins
- Operates with liquid nitrogen or dry ice
- Compatible with 0.5, 1.5, 2 x 24 and 7 mL Tube Carriages Only



OM50

- 50 Watt Ultrasonic Homogenizer
- Process sample volumes from 200 µL up to 50 mL
- Automatic tuning for easy interchange of probe tips
- Three different probe sizes available
- Minimized footprint



THQ

- Integrated digital speed display and control
- Whisper Drive Technology ensures ultra-quiet operation
- Process sample volumes from 200 µL up to 100 mL with Omni Tips™, and up to 120 mL with Generator Probes
- Variable speeds from 8,000 - 32,000 rpm in increments of 500 rpm



TH

- Omni Tip™ Plastic Probes support the elimination of cross-contamination
- Designed for homogenizing hard or soft tissue, PCR, RNA/DNA, sample prep & many other applications
- Variable speed controller for speeds of 5,000 - 30,000 rpm



Macro-ES

- 1,800 watt digitally controlled motor for consistent, powerful sample processing
- Speeds variable from 1,000 - 20,000 rpm
- Suitable for processing sample volumes 250 μ L to 30 L
- 12 Programs with 12 time and speed steps, including ramping function
- Wide variety of accessories including generator probes, cutting blades, and sealed chambers



GLH850

- Powerful & versatile for a wide range of sample types & volumes
- Compatible with stainless steel probes 5 to 30 mm diameter
- Compatible with 7 and 12 mm Omni Tip™ Plastic Probes and Hybrid Probes
- Process sample volume ranging from 200 μ L - 5 L
- Variable speed from 5,000 to 25,000 rpm



CF-350

- Universal laboratory centrifuge
- Maximum rotational speed of 15,000 rpm (21,630 rcf)
- 13 different rotors available to support tubes from 2 - 200 mL
- Rotors for deep well plates and PCR tubes available
- Quiet, maintenance-free operation



CF-260

- High-speed laboratory centrifuge
- Maximum rotational speed of 18,000 rpm (31,150 rcf)
- 10 different rotors available to support tubes from 2 - 50 mL
- Rotors for deep well plates and PCR tubes available
- Store up to 100 programs



CF-220

- Ergonomic and user-friendly clinical centrifuge
- Maximum rotational speed of 15,500 rpm (23,100 rcf)
- Seven different rotors available to support tubes from 2 - 50 mL
- Large, illuminated LCD interface
- Rotor for PCR tubes is available



Omni Tips

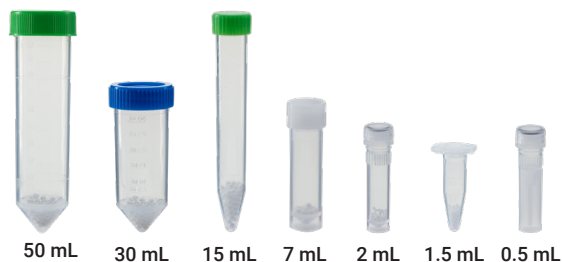
- Omni Tips™ are disposable plastic homogenizing probes. They are available in 5, 7 and 12 mm diameters and compatible with nearly all Revvity homogenizing products
- The Omni Tip™ design eliminates the hassle of disassembly and maintenance of traditional stainless-steel generator probes, eliminating the chances of cross contamination
- The 12 mm Omni Tip™ is designed for larger or harder samples where processing with a 7 mm tip is no longer efficient



Shop online

Bead Tubes

- Revvity offers a selection of Beads and Tubes for ultimate versatility
- A variety of bead media are available for different sample types. Omni's various tube sizes fit a range of sample volumes.
- Various tube sizes accommodate a range of sample volumes. Our selection includes 0.5 mL, 1.5 mL, 2 mL, 7 mL, 15 mL, 30 mL and 50 mL tubes.



Visit website



Learn more

Contact us



Learn more



revvity