



Omni Bead Ruptor 12

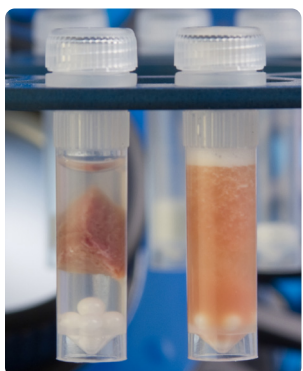
Bead Mill Homogenizer

Part number: 19-050A

Ideally suited for nucleic acid, drug, protein, enzyme, and virus extractions.



- Includes snap-in tube carriage for 2 mL and 7 mL tubes:
 - Holds 12 x 2 mL or 0.5 mL tubes and/or 4 x 7 mL tubes
 - Includes semi-permanent ring



Before (left) and after (right)

Specifications

Power rating: 115-230V, 50/60 Hz
Performance range: 0.8 - 6.0 m/s in increments of 0.15 m/s
Run time: 1 second to 9:59 minutes
Cycles: 1-10
Programs: 99
Processing volume: 25 µL to 6 mL
Dimensions: W: 15" (38.1 cm), D: 17" (43.2 cm), H: 13" (33.0 cm)
Weight: 65 lbs. (29 kg)
Warranty: One year
Certification: CE, UKCA

Features

- Optimized tube motion with speeds up to 6 m/s
- Process up to 16 samples per cycle
- Convenient front loading design
- Locking safety lid
- Reproducible extraction of proteins, nucleic acids, and small molecules

