



Part number: 25-010

Omni Bead Ruptor 4

Bead Mill Homogenizer

Built upon the innovative design and unsurpassed quality of the Omni Bead Ruptor™ family. The Omni Bead Ruptor 4 bead mill homogenizer is ideally suited for:

- Homogenizing up to 4 samples in seconds
- Reproducible extraction of proteins and nucleic acids
- Simultaneously processing of up to 4 x 2 mL, 0.5 mL tubes, 4 x 1.5 mL or 1 x 7 mL tubes.

Efficiently lyse:

- Human tissues
- Animal tissues
- Plant and food
- Micro-organisms

Specifications

Power rating: 100-230V, 50/60 Hz

Performance range: 1 - 5.0 m/s in increments of 1 m/s

Run time: 1 to 5 minutes

Cycles: 1

Programs: 1

Processing volume: 25 µL to 5 mL

Dimensions: W: 8.5" (21.5 cm), D: 10" (25.4 cm), H: 11.5" (29.2 cm)

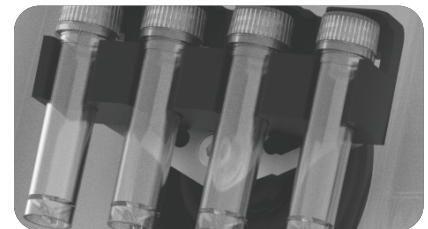
Weight: 16.5 lbs. (7.48 kg)

Warranty: One year

Certification: CE, UKCA

Features

- Broad performance range — gentle mixing to high force homogenization
- Small footprint — requires minimal bench space
- Holds 4 x 0.5 mL, 1.5 mL, 2 mL tubes or 1 x 7 mL



| Snap in tube carriage



For best results,
use with our
pre-filled bead kits

