

# Bead Ruptor™ VesselVue™

## Homogenizer



Part number: 25-030

- **Vial Size or type:** For use with VesselVue P-formulation microbubble contrast reagent, 3 mL vial, cat. # P-4007003.  
H: 1.37" (34.9 mm), Ø 0.58" (14.8 mm)
- **Quantity of vials:** 1
- **Speed range:** 1 - 5 m/s in increments of 1 m/s
- **Cycle count:** 1
- **Run time range:** 1 sec to 5 min
- **Control:** Keypad with membrane buttons
- **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)
- **Voltage requirements:** 100-230 VAC, 50/60 Hz
- **Power requirements:** 4A Max
- **Warranty:** One year
- **Standards approval:** CE and UKCA approval

## Key features

- Compact benchtop unit
- Simple intuitive user interface
- Lid safety interlock

