



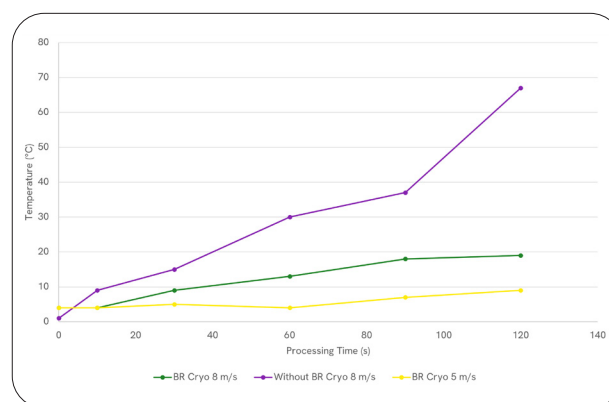
Part number: 19-8005

Omni BR Cryo Cooling Unit

The BR Cryo Cooling Unit is compatible with the Omni Bead Ruptor Elite (19-042E) and the Omni Bead Ruptor 24 (19-070) and is designed to prevent the increase of sample temperature during the homogenization process. Friction generated through the bead milling process generates heat that is imparted to the sample. In order to maintain a constant processing temperature during homogenization, the BR Cryo passes chilled air (-100 °C to -50 °C) into the sample processing chamber and across the sample tubes to dissipate heat.

Specifications and Operating Requirements

- **Dimensions:** W: 9.0" (22.9 cm), D: 13.5" (34.3 cm), H: 11.5" (29.2 cm)
- **Weight:** 16 lbs. (7.2 kg)
- **Cryo Chamber Volume:** 1 L
- **Air Source Requirements:** Compressed Dry ISO8573 Class 2 or better (water < 5 ppm, total oil < 0.1 mg/m³, vapor pressure dew point < -50 °C)
- **Minimum Air Source Pressure:** 55 psi
- **Maximum Air Source Pressure:** 120 psi
- **Recommended Air Source Pressure:** 100 psi
- **Cooling Source:** Liquid Nitrogen or Dry Ice in Ethanol
- **Liquid Nitrogen Consumption:** 1 L per 5 min of homogenization time



- **Dry Ice Consumption:** 1 L (0.5 L of ethanol + 0.5 L of dry ice) per 90 minutes of homogenization time
- **Tube Compatibility:** 0.5, 1.5, 2 x 24 and 7 mL Tube Carriages Only. Not compatible with tube sizes >7 mL or 48 Position 2 mL Carriage
- **Instrument Compatibility:** BR Cryo is compatible with Bead Ruptor Elite (19-042E) and Bead Ruptor 24 (19-070)
- **Warranty:** One Year
- **Standards Approval:** CE, UKCA Approved

