



Part number: 06-021

# Omni Prep Multi-sample Homogenizer

## Omni Prep

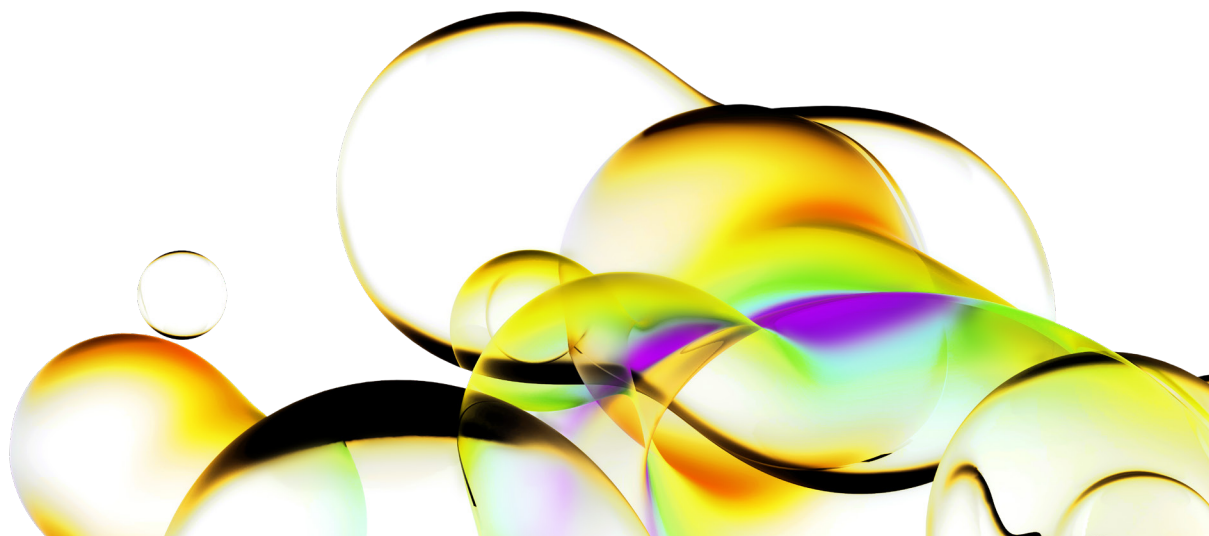
The Omni Prep Multi-Sample Homogenizer is designed to streamline the labor and time-intensive homogenization process and to make it more comfortable for the user. It features Whisper Drive Technology - a brushless motor that is virtually maintenance free and ultra quiet during operation. Unlike brush motors, Whisper Drive Technology guarantees accuracy and repeatability, since the motor holds its processing speed to +/- 1% accuracy throughout the homogenization cycle, regardless of fluctuations in sample size or sample viscosity.

An oscillating sample rack holds up to 6 tubes from 1.5 mL to 50 mL in size. Microprocessor control allows the user to control processing time and speed (500 - 30,000 rpm).

The Omni Prep increases productivity by removing the tedious manual homogenization process and dramatically improves workflow. It is small enough to fit in most laminar flow cabinets and an optional ice tray is available to process heat sensitive samples.

## Specifications

- **Volume range:** 250  $\mu$ L to 30 mL
- **Probe Compatibility:** 5, 7, 10 mm stainless steel probes and 7 or 12 mm Omni Tip™ plastic or hybrid probes
- **Power rating:** 320 watts
- **Dimensions:** 40 cm x 20 cm x 30 cm (15.7 in x 7.8 in x 11.8 in)
- **Weight:** 9.53 kg (21 lbs)
- **Timer:** 0 - 10 minutes
- **Speed control:** 500 - 30,000 rpm
- **Sound level:** < 75 db
- **Certification:** CE and UKCA
- **Warranty:** Two Years





**Tube holders** 1.5 - 2, 5, 15, 50 mL  
Each tube holder holds two tubes.

**Note:** Tube holders are required to operate the Omni Prep.

Holder for 1.5 - 2 mL Tubes **Part number: 06-209-02**

Holder for 5 mL Tubes **Part number: 06-209-05**

Holder for 15 mL Tubes **Part number: 06-209-15**

Holder for 30 - 50 mL Tubes **Part number: 06-205-50**



**Processing rack**

Movable rack for efficient processing of samples ranging in volume from 250  $\mu$ L to 30 mL.

**Note:** At least one rack is required to operate the Omni Prep.

**Part number: 06-220**



**Cooling ice trays**

Ice trays can be added to the Omni Prep rack to help keep your samples cold during processing.

Ice Tray for 1.5 - 5 mL Tubes **Part number: 06-207**

Ice Tray for 15 - 50 mL Tubes **Part number: 06-210**

[Contact us](#) for more information or to get a quote.

