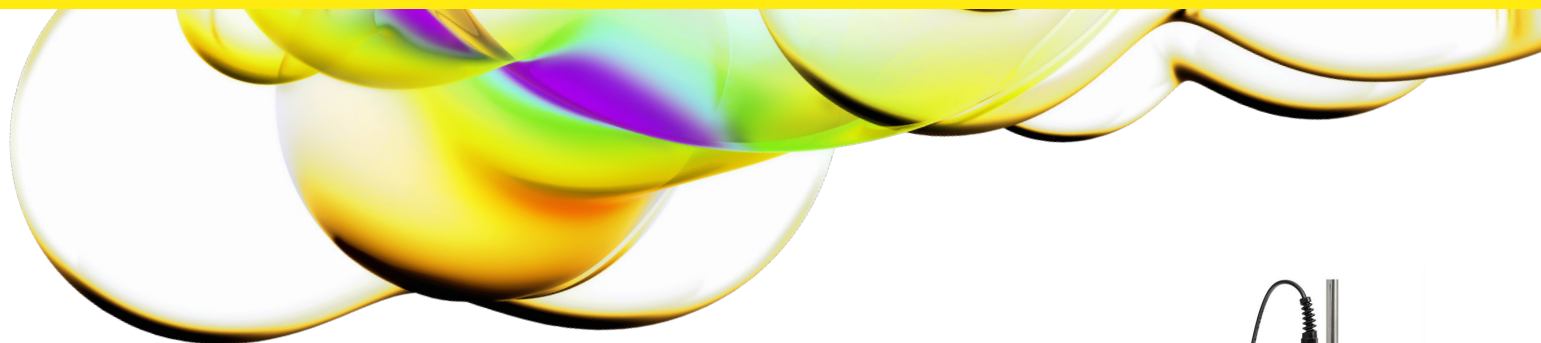
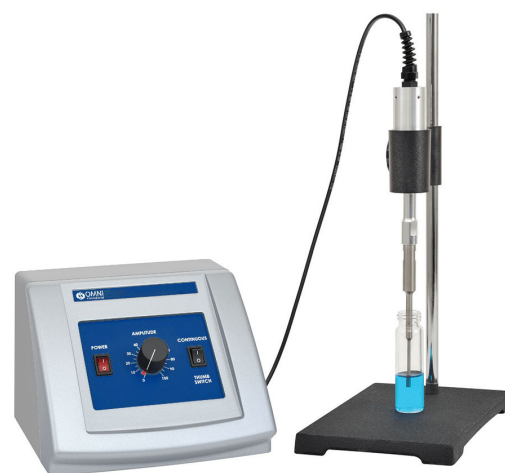


OM50 Ultrasonic Homogenizer (50W)



The OM50 is a compact and cost effective 50 watt ultrasonic homogenizer that will occupy less bench space than any unit on the market.

This model is effective for standard cell disruption and many other small volume applications. Probes are available in three different sizes.



Stand sold separately.

Technical specifications

Specify desired voltage when ordering

Power rating	50 Watts
Frequency	20 kHz
Dimensions	W: 8" (20.0 cm), D: 7.5" (19.0 cm), H: 5.75" (14.6 cm)
Voltage	110 or 220V 50/60 Hz

OM50 Includes power supply, convertor, cables and wrench set

Features

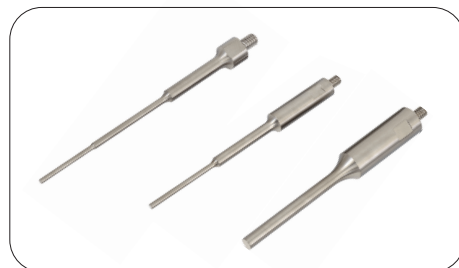
- Smallest unit available
- Thumb-switch or continuous operation
- Simple and effective operation
 - Power on/off switch
 - Amplitude control
 - Continuous or thumb switch-activated processing



OM50 accessories

Probes

Probe part #	Processing volume	Tip diameter	Amplitude
060-078	200 µL to 5 mL	5/64" (1.6 mm)	200 µm
060-125	500 µL to 15 mL	1/8" (3.2 mm)	180 µm
060-250	10 mL to 50 mL	1/4" (6.4 mm)	120 µm



Sound enclosure with converter holder

W: 12" (30.5 cm), D: 12" (30.5 cm),
H: 20" (50.8 cm)
(Part number: 060-12)



Replacement converter

(Part number: 060-03)



Support stand with converter holder

(Part number: 060-10)

