



Mixer Homogenizer

Macro Homogenizer



Macro ES Homogenizer

Homogenizer Mixers

Mixer Homogenizer

Part no: [HMG17105](#)
[HMG17106](#)

Processing range: 0.25 mL to 10 L

Generator probes: 10, 20, 30 mm

Blade assemblies:

1" and 2" in sealed chambers

Power rating: 600 watt

Speed: Variable from 500 to 18,000 rpm

Dimensions (with stand): H: 24" (61 cm),
W: 10" (25 cm)

Weight: 21 lbs. (9.5 kg)

Sound level: <68 db

Warranty: 3 year motor warranty

Macro Homogenizer

Part no: [17505](#)
[17506](#)

Processing range: 0.25 mL to 30 L

Generator probes: 10, 20, 30, 45, 55 mm

Blade assemblies:

1" and 2" in sealed chambers

Power rating: 1,800 watt

Speed: Variable from 1,000 to 20,000 rpm

Dimensions (with stand): H: 24" (61 cm),
W: 10" (25 cm)

Weight: 21 lbs. (9.5 kg)

Sound level: <68 db

Warranty: 3 year motor warranty

Macro ES Homogenizer

Part no: [MES-115 \(115V\)](#)
[MES-220 \(220V\)](#)

Processing range: 0.25 mL to 30 L

Generator probes: 10, 20, 30, 45, 55 mm

Blade assemblies:

1" and 2" in sealed chambers

Power rating: 1,800 watt

Speed: Variable from 1,000 to 20,000 rpm,
soft start, $\pm 5\%$ accuracy

Programs: 12 programs with 12 time and
speed steps, including ramping function

Interface: RS-232 and Windows software

Dimensions: W: 14.2" (36 cm),
D: 16.9" (43 cm), H: 30.7" (78 cm)

Weight: 32 lbs. (14.5 kg)

Sound level: <68 db

Warranty: 3 year motor warranty

Compliance: CE Approved