

Omni GLH 850

General laboratory homogenizer

The Omni GLH 850 homogenizer is an affordable, high performance instrument capable of processing virtually any laboratory sample in seconds. A powerful 850 watt motor provides maximum versatility for homogenizing, dispersing or emulsifying the largest variety of sample types and sizes in its class. The Omni GLH 850 features an advanced, integrated, push button digital control system that ensures accuracy and repeatability. The unit's closed loop variable speed controller (5,000 - 25,000 rpm) maintains selected speeds within +/- 100 rpm of the original set point. The digital timer can be set from seconds to minutes, and includes a 30 second countdown timer, which is ideal for remote applications in fume hoods or biosafety cabinets. A soft start feature ensures smooth processing and minimizes splashing.

The Omni GLH 850 is compatible with a wide range of interchangeable stainless steel rotor stator generator probes (sold separately), from 5 to 30 mm diameter, and processes samples ranging from 200 µL to 5 liters. Stainless steel probes are constructed of quality materials for durability, and are easy to disassemble for cleaning. Omni Tips™ plastic disposable generator probes are available to completely eliminate the risk of carryover and cross contamination between samples (adapter required).

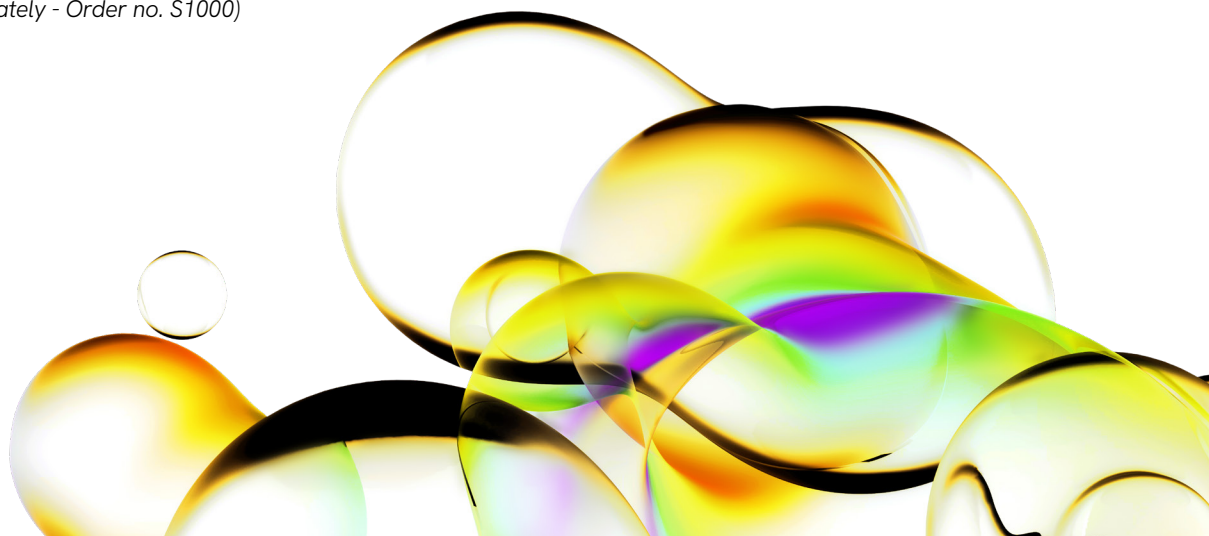
Includes


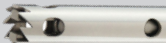



- 850 Watt Motor
(115V) - Order no. 20-010
(220V) - Order no. 20-020
- Cross rod and stand clamp
(lab stand sold separately - Order no. S1000)



Specifications

- **Processing Range:**
- **With Omni Tips:** <200 µL to 100 mL
- **With Generator Probes:** <200 µL to 5 L
- **Power Rating:** 850 Watt
- **Speed Control:** Variable from 5,000 - 25,000 rpm (increments of 500 rpm)
- **Timer:** 1 sec - 10 min with 30 sec start delay
- **Controls:** Digital
- **Dimensions:** H: 8.5" (21.6 cm), Diameter: 2.75" (7 cm)
- **Weight:** 3.5 lbs. (1.6 kg)
- **Sound Level:** < 72 db.
- **Warranty:** Two Year Motor Warranty
- **Standards Approval:** CE & UKCA Approved



| Stainless steel probes | Size | Processing range | Type | Particle size | Order no. |
|---|-------------|------------------|---------------------------|---------------|------------|
|  | 5 x 75 mm | 200 µL - 5 mL | Flat bottom | Fine | G5-75 |
| | | | Flat bottom - wide window | Coarse | G5-75W |
|  | 7 x 115 mm | 250 µL - 30 mL | Saw tooth | Fine | G7-115ST |
| | | | Saw tooth - wide window | Coarse | G7-115STW |
|  | 10 x 115 mm | 1.5 mL - 100 mL | Saw tooth | Fine | G10-115 |
| | | | Saw tooth - wide window | Coarse | G10-115W |
|  | 20 x 195 mm | 100 mL - 2 L | Saw tooth | Fine | G20-195ST |
| | | | Saw tooth - wide window | Coarse | G20-195STW |
|  | 30 x 195 mm | 200 mL - 5 L | Flat bottom | Fine | G-30NA-195 |
| | | | Flat bottom - wide window | Coarse | G-30WA-195 |

7 and 12 mm Omni Tips™ disposable homogenizing probes



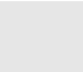


The Omni Tips' patented design eliminates the hassle of disassembly and maintenance of traditional stainless steel generators.

Proprietary Omni Tip™ Probes are engineered from durable plastics and are so economical they can be disposed of after each use to eliminate cross-contamination during homogenization.

The outer tube of Omni Tips™ is made of polycarbonate. The inner shafts are made of Ultem 1000, a polyetherimide. The Omni Tips' patented design eliminates the hassle of disassembly and maintenance of traditional stainless steel generators.

12 mm Omni Tips™ are suitable for homogenization where sample size or sample hardness dictates that processing with the 7 mm tip is prohibitive.

| A selection of our best-selling plastic probes: | | | |
|---|---|----------|-----------|
| | 7 mm Soft Tissue Omni Tip™ | | Order no. |
|  | Omni Tips™ Plastic Homogenizing Probes | 25 pack | 30750 |
| | Omni Tips™ Plastic Homogenizing Probes | 50 pack | 32750 |
| | Omni Tips™ Plastic Homogenizing Probes | 100 pack | 34750 |
| | 7 mm Hard Tissue Omni Tip™ | | Order no. |
|  | Omni Tips™ Plastic Homogenizing Probes | 25 pack | 30750H |
| | Omni Tips™ Plastic Homogenizing Probes | 50 pack | 32750H |
| | Omni Tips™ Plastic Homogenizing Probes | 100 pack | 34750H |
| | Sterile Omni Tips™ Plastic Homogenizing Probes | 25 pack | 30750H-S |
| | 12 mm Hard Tissue Omni Tip™ | | Order no. |
|  | Omni Tips™ Plastic Homogenizing Probes | 25 pack | 301250 |
| | Omni Tips™ Plastic Homogenizing Probes | 50 pack | 321250 |
| | Omni Tips™ Plastic Homogenizing Probes | 100 pack | 341250 |

GLH 850 package

The Omni GLH 850 package includes: 850 watt motor, 10 x 115 mm saw tooth stainless steel generator probe, 20 x 195 mm saw tooth stainless steel generator probe, 30 x 195 mm flat bottom stainless steel generator probe, tool kit, instruction manual, cross rod, and motor stand assembly.

