## Automated high-throughput homogenization.

revvit



## Omni Prep 96 Automated Homogenizer Workstation

For research use only. Not for use in diagnostic procedures.

# Advanced, high-throughput sample preparation

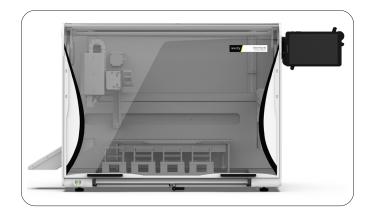
## Omni Prep 96 automated homogenizer workstation

Sample preparation is an integral step of any analytical procedure. The complexity and diversity of samples can make testing a challenge, especially at the very beginning of your process where homogeneity matters. Whatever your sample we've got you covered.

The Omni Prep 96 is a fully automated homogenization workstation, enabling true walk-away processing in your lab. As market consolidation continues and pressure mounts to keep costs low, it is necessary to rely on efficiency and predictability in your workflows. You can count on us and our technology to deliver. Our automation for sample prep enables lower costs per sample, more traceability, and better analytical reproducibility.

The Omni Prep 96 workstation was designed to provide automated homogenization using the same generator probe movement as a manual handheld homogenizer operated by a laboratory technician. During processing the motor heads are capable of both vertical and horizontal intra-tube movement to maximize sample contact with the generator probes in order to ensure complete homogenization.





#### Features

- Process volumes from 250  $\mu L$  to 40 mL in 2 mL to 50 mL tubes
- Create and save customized methods in the intuitive software
- Unique oscillating motion to ensure complete homogenization
- Variable speed from 500 28,000 rpm
- Compatible with 7 & 12 mm plastic Omni Tips<sup>™</sup>
- Robust, reliable industrial motors and controls

The small footprint fits into most bio-safety cabinets and fume hoods. User-friendly programming makes saving, recalling, or changing existing methods easy. Our patented Whisper Drive Technology<sup>™</sup> with brush-less motors delivers maintenance-free, ultra-quiet, maximum processing power.

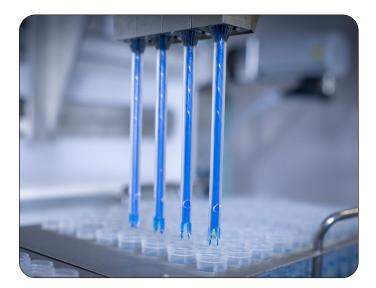
The cleaning station provides wash bath functionality to the Omni Prep 96 to support using 7 and 12-mm Hybrid Probes and 5, 7, and 10-mm Stainless Steel Generator Probes. It features flow-through, static, and ultrasonic cleaning tanks and overflow control. The flow-through tank is connected to a laboratory water supply, and both the static and ultrasonic tanks can be filled with reagents that will assist in mitigating analyte carryover. The three separate areas of the tank can be used in any order within a homogenization method.

Part Number	Description	Qty
51-02A-1	Prep 96 Automated Homogenizer (No Accessories)	1
51-BRA-1006-J02	Deck Adapter for Clean Probes/Tips Rack	1
23-HWA-22	Rack for 7 mm Clean Probes/Tips	1
23-HWA-23	Rack for 12 mm Clean Probes/Tips	1
23-HWR-11-JA	Deck Adapter for Standard Sample Racks	1
23-HWR-15	2 mL x 96 Tubes Standard Sample Rack	1
23-HWR-18	5 mL x 96 Round Bottom Tubes Standard Sample Rack	1
23-HWR-14	14/15 mL x 96 Round Bottom Tubes Standard Sample Rack	1
23-HWR-13	30 mL x 24 Tubes Standard Sample Rack	1
23-HWR-12	50 mL x 24 Tubes Standard Sample Rack	1
51-HWA-23S1	Deck Adapter for SBS Sample Racks	4
23-HWR-18SBS	SBS Sample Rack 5 mL x 24 Round Bottom Tubes	4
23-HWR-11SBS	SBS Sample Rack 14/15 mL x 12 Round Bottom Tubes	4
23-HWR-15SBS	SBS Sample Rack 14/15 mL x 24 Round Bottom Tubes	4
23-HWR-16SBS	SBS Sample Rack 30 mL x 8 Tubes	4
23-HWR-17SBS	SBS Sample Rack 50 mL x 8 Tubes	4
23-HWR-25	Passive Cooling 2 mL x 96 Tubes Sample Rack	1
23-HWR-28	Passive Cooling 5 mL x 96 Round Bottom Tubes Sample Rack	1
23-HWR-24	Passive Cooling 14/15 mL x 96 Round Bottom Tubes Sample Rack	1
23-HWR-23	Passive Cooling 30 mL x 24 Tubes Sample Rack	1
23-HWR-22	Passive Cooling 50 mL x 24 Tubes Sample Rack	1
23-HWA-13B-JA	Deck Adapter for Dirty Probes/Tips Disposal Rack	1
23-HWA-13B	Disposal Rack for Dirty Probes/Tips	1
19-6619	Replacement Disposal Tubes for Dirty Probe Tip Rack	100
23-HWA-16JA	Disposal Chute for Dirty Probes/Tips	1
51-BRA-1095	Cleaning Station - 120V	1
51-BRA-1096	Cleaning Station - 220V	1
30750H	7 mm Omni Tip™ Plastic Homogenizing Probes (7 x 110 mm)	25
32750H	7 mm Omni Tip™ Plastic Homogenizing Probes (7 x 110 mm)	50
34750H	7 mm Omni Tip™ Plastic Homogenizing Probes (7 x 110 mm)	100
35750H	7 mm Omni Tip™ Plastic Homogenizing Probes (7 x 110 mm)	500
3M750H	7 mm Omni Tip™ Plastic Homogenizing Probes (7 x 110 mm)	1000
301250	12 mm Omni Tip™ Plastic Homogenizing Probes (12 x 110 mm)	25
321250	12 mm Omni Tip™ Plastic Homogenizing Probes (12 x 110 mm)	50
341250	12 mm Omni Tip™ Plastic Homogenizing Probes (12 x 110 mm)	100
351250	12 mm Omni Tip™ Plastic Homogenizing Probes (12 x 110 mm)	500
3M1250	12 mm Omni Tip™ Plastic Homogenizing Probes (12 x 110 mm)	1000
22-007	7 mm Omni Tip™ Hybrid Probe (Stainless Steel & Plastic)	1
22-007-6	7 mm Omni Tip™ Hybrid Probe (Stainless Steel & Plastic)	6
22-012	12 mm Omni Tip™ Hybrid Probe (Stainless Steel & Plastic)	1
22-012-6	12 mm Omni Tip™ Hybrid Probe (Stainless Steel & Plastic)	6
B5-075	5 x 75 mm Stainless Steel Generator Probe	1
B7-110ST	7 x 110 mm Stainless Steel Generator Probe	1
B10-110ST	10 x 110 mm Stainless Steel Generator Probe	1

Part Number	Description	Qty
19-660-1000	2 mL Low Binding Tubes with Screw Caps	1000
19-6605-500	5 mL Round Bottom Tubes with Snap Caps	500
19-6614-500	14/15 mL Round Bottom Tubes with Screw Caps	500
19-6635	30 mL Tubes with Screw Caps	100
19-6635-500	30 mL Tubes with Screw Caps	500
19-6650	50 mL Tubes with Screw Caps	100

## Specifications

Processing Range	250 µL to 40 mL
Sample Racks	2, 5, 14/15, 30 and 50 mL
Sample Capacity	24 to 96 Samples
Programming	Local control via touch screen HMI
Motor Power Rating	100 Watts
Speed Control	500 to 28,000 rpm
Dimensions	37.5 in. (w) x 25.5 in. (d) x 28.7 in. (h), 95.2 cm (w) x 64.8 cm (d) x 72.9 cm (h)
Weight	125 lbs. (57 kg)
Sound Level	< 75 db
Certification	TUV, NRTL, CE, UKCA, RCM, FCC Part 15 Approved
Power Supply	100-230 VAC, 50/60 Hz





revvity

Revvity, Inc. 940 Winter Street Waltham, MA 02451 USA www.revvity.com

For a complete listing of our global offices, visit www.revvity.com Copyright ©2024, Revvity, Inc. All rights reserved.

1377423

### www.revvity.com