

revvity

Sample disruption and homogenization

Product catalog



Contents

Bead mill homogenizers3 - 7

Instruments and accessories

Consumables

Rotor-stator handhelds and industrial mixers8 - 15

Handheld homogenizers, bundles, accessories, and generator probes

Omni GLH850 and Omni TH

Omni THQ and Omni Micro

Omni TM

Omni Tip™ generator probes

Single-use plastic disposables

Hybrid multi-use probes

Industrial mixers/macros homogenizers and accessories

Omni Mixer

Omni Digital Programmable Macro-ES

Omni Macro

Generator probes for Mixers/Macros

Sealed chambers with stainless steel probe/blade assembly

Other portfolio solutions16

Omni automated homogenizers and workstations

Single-probe sonicators and accessories

Bead mill homogenizers

[Omni Bead Ruptor™ bead mills](#), built on over 60 years of sample prep expertise, are indispensable powerhouse units for labs looking to transition from single, manual handheld preps to multi-sample processing, whatever your lab throughput.

[Omni Bead Ruptor Elite™ \(19-042E\)](#)

- Powerful: Processing speed from 0.8 - 8 m/s. Unique motion range
- Versatile: Multiple sample tube volume compatibility
- Usability: No cool-downs between runs. Intuitive touchscreen interface
- Dimensions: 33 cm (13") H x 38.1 cm (15") W x 43.2 cm (17") D



[Accessories](#)

Part No.	Description	Capacity (max)	Volume range
19-372	1.5 mL microcentrifuge tube carriage	1.5 mL tubes x 24	200 µL -1 mL
19-373	2 mL tube carriage	2 mL screw cap tubes x 24	200 µL -1.5 mL
19-378	48 place 2 mL tube carriage w/ finger plate	2 mL screw cap tubes x 48	200 µL -1.5 mL
19-374	7 mL tube carriage	7 mL screw cap tubes x 12	0.5 mL -5.0 mL
19-375	15 mL tube carriage w/ finger plate	15 mL screw cap conical x 12	0.5 mL -10 mL
19-376-HT	30 mL tube carriage w/ finger plate	30 mL screw cap conical x 6	5.0 mL -30 mL
19-377	50 mL tube carriage	50 mL screw cap conical x 3	5.0 mL -50 mL
19-8005	Bead Ruptor™ cryo cooling unit	Tubes carriages < 7mL. Not compatible with 48 place unit	

[Omni Bead Ruptor™ 96 \(27-0001 \(115V\) and 27-0002 \(220V\) \)](#)

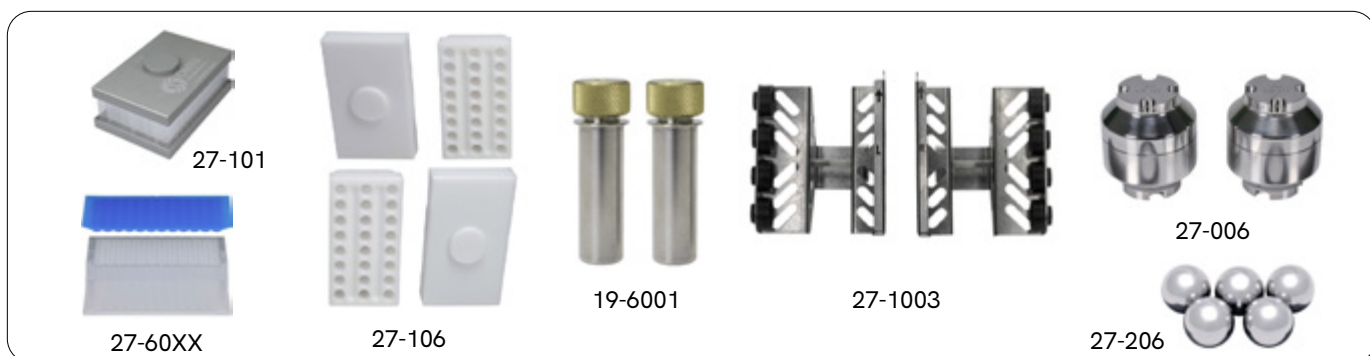
- Powerful: 30 Hz max speeds. Large, high-impact motions
- High-throughput: Up to 384 samples / cycle
- Versatile: Well-plate, tubes, milling jar (cryo-milling) compatibility
- Dimensions: 24 cm (9.5") closed / 46.5 cm (18") open H x 38.5 cm (15") W x 42 cm (16.5") D



[Tube and milling jar accessories](#)

Part No.	Description	Capacity (max)	Quantity included
27-004	25 mL milling jars and O-rings	4 -10 mL sample volumes	2 jars
27-203	10 mm milling balls	Samples used with 25 mL milling jar	10 balls
27-006	50 mL milling jars and O-rings	8 -20 mL sample volumes	2 jars
27-206	25 mm milling balls	Samples used with 50 mL milling jar	5 balls
27-1003	50 mL tube holders	< 50 mL volumes per tube	2 holders
27-106	Microtube or screw cap tube holders	1.5 mL tubes x 48 or 2 mL screw cap tubes x 48 per holder	2 holders (4 parts)
19-6001	2 mL stainless steel lysing/grinding cryo tube	< 1.5 mL volumes	2 tubes





Well-plate accessories and consumables

Part No.	Description	Capacity (max)	Quantity included
27-101	Well-plate adapters	1 x 2 mL deep well plates, 2 x 1 mL or 3 x SBS standard plates per each adapter	2 adapter set (4 parts)
27-107	96-well 2D barcoded storage tube adapter for Micronic and FluidXtubes		2 plates
27-108	96-well 2D barcoded storage tube adapter for Matrix tubes		2 plates
27-6005	2 mL 96-well deep well plates pre-filled with 5 mm stainless steel beads		5 plates
27-6006	2 mL 96-well deep well plates pre-filled with 0.1 mm ceramic beads		2 plates
27-6007	2 mL 96-well deep well plates pre-filled with 0.1 mm ceramic beads		10 plates
27-6009	2 mL 96-well deep well plates pre-filled with 1.4 mm ceramic beads		10 plates
27-6011	2 mL 96-well deep well plates pre-filled with 1.4 mm ceramic beads		2 plates
27-520	2 mL 96-well deep well plates		10 plates
27-530	Sealing mat for 2 mL 96-well deep well plates		1 sealing mat
27-7005	5 mm stainless steel beads bulk	Well-plates use recommended.	325 g (1 pack)

For labs with lower throughput, the Omni Bead Ruptor™12 or Omni Bead Ruptor™ 4 enables 12 or 4 sample simultaneous processing per run. Units require no cool-down between runs, boast a user-centric front-loading design, and come equipped with integrated lid-lock for safety.

Omni Bead Ruptor 12 (19-050A)

- Powerful: Processing speed from 0.8 - 8 m/s, unique motion range
- Throughput: Up to 12 x 2mL tubes or up to 4 x 7 mL tubes
- Balanced economics: Invest in throughput, at an attractive price
- Dimensions: 33 cm (13") H x 38.1 cm (15") W x 43.2 cm (17") D



Omni Bead Ruptor 4 (25-010)

- Powerful: Processing speed up to 5 m/s, unique motion range
- Throughput: Up to 4 x 0.5 - 2.0 mL tubes or 1 x 7 mL tube
- Compact efficiency: Budget-friendly, with small footprint
- Dimensions: 29.2 cm (11.5") H x 21.5 cm (8.5") W x 25.4 cm (10") D



Bead mill consumables

Source your homogenizing bead mixes, as well as bulk beads, tubes, and deep-well plate consumables from us. Though optimized for use together with our Omni Bead Ruptor products, they are compatible with other bead mills units.

Bead selection by sample type

	Glass		Garnet		Ceramic				Metal	
Sample type	0.1 mm	0.5 mm	0.15 mm	0.7 mm	0.1 mm	0.5 mm	1.4 mm	2.8 mm	6.5 mm	2.4 mm
Bacterial/yeast										
Spores										
Algae/plankton										
Soft soil/sediment/waste water										
Clay/feces/sludge										
Soft tissues, brain, liver, kidney, cultured mammalian cells										
Hard tissues, heart, muscle, skin										
Plant material										
Dry grinding										
Dense biological material, hair, bone, seeds										

Bulk beads and active grinding media

Part no.	Description	Part no.	Description
19-683	0.1 mm ceramic beads, 500 g	19-641	0.1 mm glass beads, 400 g
19-684	0.5 mm ceramic beads, 500 g	19-642	0.5 mm glass beads, 400 g
19-645	1.4 mm ceramic beads, 325 g	19-643	0.15 mm garnet beads, 500 g
19-646	2.8 mm ceramic beads, 325 g	19-644	0.7 mm garnet beads, 500 g
19-682	6.5 mm ceramic beads, 325 g	19-640	2.4 mm stainless steel beads, 500 g
19-900S	Active grinding media, stainless steel, 50 units	27-7005	5 mm stainless steel beads, 325 g

Tube selection by volume range

Tube size	20 µL	100 µL	1 mL	5 mL	15 mL	30 mL	50 mL
0.5 mL	20 µL - 300 µL						
1.2 mL stripwell	20 µL - 800 µL						
1.5 mL		200 µL - 1 mL					
2.0 mL		100 µL - 1.8 mL					
7.0 mL			1 mL - 5 mL				
15 mL				5 mL - 15 mL			
30 mL				2 mL - 20 mL			
50 mL				2 mL - 35 mL			

The nuances of tube and bead media selection for robust sample homogenization should not be overlooked when exploring tough tissues or complex matrix inputs upstream the extraction method for your valuable analyte.



Ready-to-use, pre-filled bead mix tubes

All pre-filled tube mixes are certified RNase and DNase free. Part numbers ending in D are also certified microbial DNA-free.

0.5 mL tubes

Part no.	Description	Popular applications
19-623	0.15 mm garnet beads, qty 50	Micro-organism lysing, soil, clay, sediment, wastewater/feces
19-625	0.25 mm carbide beads, qty 50	RNA extraction from micro-organisms, soil, and sediments
19-626 19-626D	1.4 mm ceramic beads, qty 50 Available as microbial DNA-free	Soft tissue, biopsies, tissue punches, arteries, nerves, and veins for research

1.5 mL tubes

Part no.	Description	Popular applications
19-617 19-617D	1.4 mm ceramic beads, qty 50 Available as microbial DNA-free	Soft tissues, brain, liver, and kidney
19-618 19-618D	2.8 mm ceramic beads, qty 50 Available as microbial DNA-free	Hard tissues, heart, muscle, skin, tendon, tail, whole organs, and hard plant material
19-610	2.4 mm metal beads, qty 50	Hard tissues, skin, muscle, bone, hair, roots, and seeds. Dry grind

2.0 mL tubes

Part no.	Description	Popular applications
19-621 19-621-1000	0.1 mm glass beads, qty 50 0.1 mm glass beads, qty 1000	Micro-organism lysis, cells, bacteria, gram -, gram +, and spores
19-624 19-624-1000	0.7 mm garnet beads, qty 50 0.7 mm garnet beads, qty 1000	Hard and fibrous tissues. Soil, sediment, wastewater/feces
19-620 19-620D 19-620-1000	2.4 mm metal beads, qty 50 Available as microbial DNA-free 2.4 mm metal beads, qty 1000	Hard tissues, muscle, skin, bone, hair, roots, and seeds. Dry grind
19-622 19-622D 19-622-1000	0.5 mm glass beads, qty 50 Available as microbial DNA-free 0.5 mm glass beads, qty 1000	Tough micro-organism lysis, gram +, yeast, fungi, algae, spores and plankton
19-627 19-627D 19-627-1000	1.4 mm ceramic beads, qty 50 Available as microbial DNA-free 1.4 mm ceramic beads, qty 1000	Soft tissue grinding, brain, liver, kidney, spleen, and soft plant
19-635-1000	5 mm metal beads, qty 1000	Various applications
19-631 19-631D	0.5 mm ceramic beads, qty 50 Available as microbial DNA-free	Universal microbial grinding. Larger micro-organisms, protozoans, nematodes, and parasites
19-636D	2.8 mm and 0.1 mm ceramic beads, qty 50 Microbial DNA-free	Lysis of viruses, bacteria, and fungi in gut, stool, and soil
19-628 19-628D 19-628-1000	2.8 mm ceramic beads, qty 50 Available as microbial DNA-free 2.8 mm ceramic beads, qty 1000	Hard tissues, heart, muscle, skin, tendon, tail, whole organs, and hard plant material

2.0 mL tubes (continued)

Part no.	Description	Popular applications
19-629	1.4 mm and 2.8 mm ceramic beads, qty 50	Various applications
19-632 19-632D 19-632-1000 19-632D-1000	0.1 mm ceramic beads, qty 50 Available as microbial DNA-free 0.1 mm ceramic beads, qty 1000 Available as microbial DNA-free	All micro-organisms. Gram+ and gram-bacteria, yeast, mold, algae, spores, and plankton

15 mL tubes

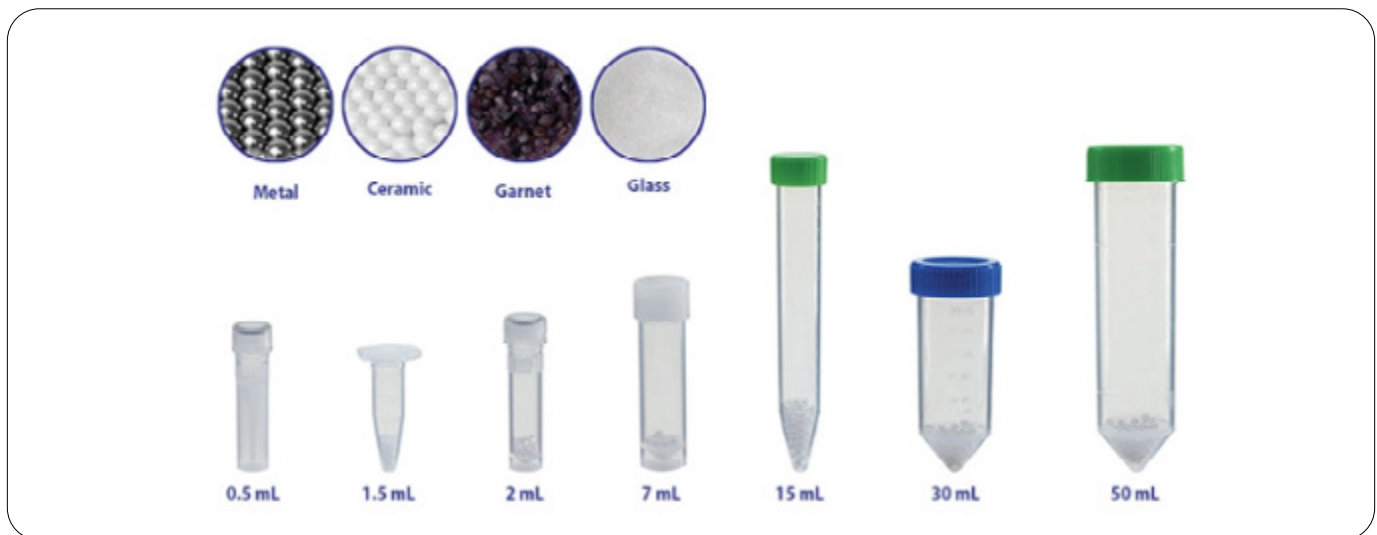
Part No.	Description	Popular applications
19-1658	2.8 mm ceramic beads, qty 50	Tough and fibrous plant material, hard tissues, seeds, and feed

30 mL tubes

Part No.	Description	Popular applications
19-6350 19-6350Z	2.4 mm metal beads, qty 50 2.4 mm metal beads with 1/4" ceramic rod, qty 50	Hard plant or animal tissues, dry grinding, roots, and seeds. Larger pieces of hard plant or animal tissues, dry grinding, roots, and seeds
19-6358 19-6358D 19-638-500 19-638Z	2.8 mm ceramic beads, qty 50 Available as microbial DNA-free 2.8 mm ceramic beads, qty 500 2.8 mm ceramic beads with 1/4" ceramic rod, qty 50	Tough and fibrous plant material, hard tissues, seeds, and feed Larger pieces of tough and fibrous plant material, hard tissues, seeds, and feed
19-6357 19-6357Z	1.4 mm ceramic beads, qty 50 1.4 mm ceramic beads with 1/4" ceramic rod, qty 50	Soft plant or animal tissues, fruit, and cereals Larger pieces of soft plant or animal tissues, fruit, and cereals

50 mL tubes

Part No.	Description	Popular applications
19-6508	2.8 mm ceramic beads, qty 50	Tough and fibrous plant material, hard tissues, seeds, and feed



[We offer a range of homogenizing beads of different composition and sizes. Pre-filled or empty bulk tubes are also available.](#)



Rotor-stator homogenizers

[Our high-shear lab homogenizers](#) are designed to process tough samples quickly. Maximum speeds vary by unit for our handheld offerings, but ranges from 25,000 rpm up to 35,000 rpm. Meanwhile, our industrial mixers and macros reach up to 20,000 rpm and offer sealed chamber processing. Handheld units are also available as stand-mounted configurations, allowing labs versatility in length of processing time. All systems are supported by a wide breadth of compatible generator probes to support a multitude of applications.

Handheld units

Omni General Lab Homogenizer - GLH 850

- Powerful: 850-watt homogenizer
- Versatile: Wide range of sample volumes from 200 μ L to 5 L. Compatible with 5 mm to 30 mm diameter stainless steel probes and 7 mm or 12 mm Omni Tips



Unit and bundles

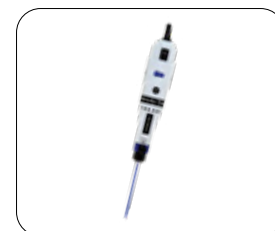
Part no.	Description
20-010 (115V)	Includes stand clamp (1), cross rod (1), and tool kit (1). Requires separate purchase of a generator probe to operate. Requires purchase of stand assembly to be stand-mounted
20-020 (220V)	
20-010-P (115V)	Includes 3 stainless steel generator probes - 10 x 115 mm saw tooth (1), 20 x 195 mm saw tooth (1), 30 x 195 mm saw tooth (1)
20-020-P (220V)	

Accessories

Part no.	Description
A1000SB	Omni Tip adapter for compatibility with 7mm or 12 mm Omni Tips (qty 1)
S1000	Stand assembly (qty 1)
S-T1001	Generator probe tool kit to remove probes knives, bearing, and a shaft holder (qty 1)

Omni Tissue Homogenizer -TH

- Powerful: 125-watt homogenizer
- Versatile: Wide range of sample volumes from 200 μ L to 100 mL. Compatible with 5 mm to 10 mm diameter stainless steel probes and 7 mm or 12 mm Omni Tips



Unit and bundles

Part no.	Description
TH115 (115V)	Includes Omni Tip adapter (1), Omni tip sample pack (1), stand clamp (1), and tool kit (1). Requires purchase of stand assembly to be stand-mounted
TH220 (220V)	
THP115 (115V)	Includes 3 stainless steel generator probes -5 x 75 mm flat bottom (1), 7 x 115 saw tooth (1), 10 x 115 mm saw tooth (1). Also, 1 motor stand assembly, 1 cross rod
THP220 (220V)	



TH bundles (continued)

Part no.	Description
TH115-K (115V) TH220-K (220V)	Includes motor stand assembly (1). Requires separate purchase of a generator probe to operate
TH115-PCR (115V) TH220-PCR (220V)	Includes Omni Tip adapter (1), 7mm Hard Tissue Omni Tips plastic probes (6), and 7 mm Soft Tissue Omni Tips plastic probes (6)
TH115-PCRD (115V) TH220-PCRD (220V)	Includes Omni Tip adapter (1) and 7 mm Soft Tissue Omni Tips plastic probes (12)
TH115-PCRH (115V) TH220-PCRH (220V)	Includes Omni Tip adapter (1) and 7 mm Hard Tissue Omni Tips plastic probes (12)
TH115-PCR5D (115V) TH220-PCR5D (220V)	Includes Omni Tip adapter (1), 5 x 75 mm standard flat bottom stainless steel probe (1), and 7 mm Soft Tissue Omni Tips plastic probes (12)
TH115-PCR5H (115V) TH220-PCR5H (220V)	Includes Omni Tip adapter (1), 5 x 75 mm standard flat bottom stainless steel probe (1), and 7 mm Hard Tissue Omni Tips plastic probes (12)
TH115-PCR7D (115V) TH220-PCR7D (220V)	Includes Omni Tip adapter (1), 7 x 115 mm standard saw tooth stainless steel probe (1), and 7 mm Soft Tissue Omni Tips plastic probes (12)
TH115-PCR7H (115V) TH220-PCR7H (220V)	Includes Omni Tip adapter (1), 7 x 115 mm standard saw tooth stainless steel probe (1), and 7 mm Hard Tissue Omni Tips plastic probes (12)
TH115-PCR10D (115V) TH220-PCR10D (220V)	Includes Omni Tip adapter (1), 10 x 115 mm standard saw tooth stainless steel probe (1), and 7 mm Hard Tissue Omni Tips plastic probes (12)
TH115-PCR10H (115V) TH220-PCR10H (220V)	Includes Omni Tip adapter (1), 10 x 115 mm standard saw tooth stainless steel probe (1), and 7 mm Hard Tissue Omni Tips plastic probes (12)

Accessories

Part no.	Description
A1000SB	Omni Tip adapter for Omni TH (black nose ring / pre-2016). For 7 mm or 12 mm Omni Tips
12-121	Omni Tip adapter for Omni TH (blue nose ring). For 7mm or 12 mm Omni Tips
S1000	Stand assembly. For units/bundles except TH115-K/TH220-K for stand mounting
S-T1001	Generator probe tool kit to remove probes knives, bearing, and a shaft holder

Stainless steel generator probes for Omni TH and Omni GLH 850 (G-style drive connect)

Part no.	Description	Volume range
G5-75	5 x 75 mm, flat bottom (standard window / fine to medium samples) (1)	200 µL - 5 mL
G5-75W	5 x 75 mm, flat bottom (wide window / medium to coarse samples) (1)	
G7-115ST	7 x 115 mm, saw tooth (standard window / fine to medium samples) (1)	250 µL - 30 mL
G7-115STW	7 x 115 mm, saw tooth (wide window / medium to coarse samples) (1)	
G10-115	10 x 115 mm, saw tooth (standard window / fine to medium samples) (1)	1.5 mL - 100 mL
G10-115W	10 x 115 mm, saw tooth (wide window / medium to coarse samples) (1)	



Large volume stainless steel probes for Omni GLH 850 (G-style drive connect)

Part No.	Description	Volume range
G20-195ST	20 x 195 mm, saw tooth (standard window / fine to medium samples) (1)	50 mL - 2.0 L
G20-195STW	20 x 195 mm, saw tooth (wide window / medium to coarse samples) (1)	
G-30NA-195	30 x 195 mm, flat bottom (standard window / fine to medium samples) (1)	75 mL - 5.0 L
G-30WA-195	30 x 195 mm, flat bottom (wide window / medium to coarse samples) (1)	



Omni Digital Tissue Homogenizer -THQ

- Energy efficient: 40-watt homogenizer
- Versatile: Wide range of sample volumes from <200 μ L to 120 mL. Compatible with 5 mm to 10 mm diameter stainless steel probes and 7 mm or 12 mm Omni Tips



Unit

Part No.	Description
12-500	Includes 7mm Hard Tissue Omni Tips plastic probes (6), 7 mm Soft Tissue Omni Tips plastic probes (6), stand clamp (1), tool kit (1), and storage box. Requires purchase of stand assembly to be stand-mounted

Accessories

Part No.	Description
S1000	Stand assembly. For units/bundles except TH115-K/TH220-K for stand mounting
S-T1001	Generator probe tool kit to remove probes knives, bearing, and a shaft holder

Omni micro homogenizer - μ Hb

- Energy efficient: 12-volt lithium ion battery powered, 35 watt motor with operating speeds up to 30,000 rpm
- Versatile: Wide range of sample volumes from <200 μ L to 100 mL, suitable for field-based studies. Compatible with 5 mm to 10 mm diameter stainless steel probes and 7 mm or 12 mm Omni Tips



Unit

Part No.	Description
UHB-115 UHB-220	Includes tool kit (1). Requires purchase of probes to operate, and stand assembly to be stand-mounted

Accessories

Part No.	Description
S1000	Stand assembly
S-T1001	Generator probe tool kit to remove probes knives, bearing, and a shaft holder



Stainless steel generator probes for Omni THQ and Omni micro (uHB) (B-style drive connect)

Part no.	Description	Volume range
B5-075 B5-075-6	5 mm x 75 mm, saw tooth (fine to medium samples) (qty 1) 5 mm x 75 mm, saw tooth (fine to medium samples) (qty 6)	200 µL - 5 mL
B7-110ST B7-110ST-6	7 mm x 110 mm, saw tooth (fine to medium samples) (qty 1) 7 mm x 110 mm, saw tooth (fine to medium samples) (qty 6)	250 µL - 30 mL
B10-110ST B10-110ST-6	10 mm x 110 mm, saw tooth (fine to medium samples) (qty 1) 10 mm x 110 mm, saw tooth (fine to medium samples) (qty 6)	1.5 mL - 100 mL



Omni tissue master homogenizer -TM

- Efficient: 125 watt motor with operating speeds up to 35,000 rpm
- Reliable: Supports sample volumes from <200 µL to 100 mL. Compatible with own set of 5 mm to 10 mm diameter stainless steel TM generator probes. Not compatible with other Omni stainless steel, hybrid or plastic probes. Preferred solution for labs with few variations in sample size



Unit and bundles

Part no.	Description
TM125-115 (115V) TM125-220 (220V)	Includes 7 mm x 114 mm stainless steel flat bottom TM probe (1)
TMP125-115 (115V) TMP125-220 (220V)	Includes 3 stainless-steel, flat-bottom TM probes -5 mm x 95 mm (1), 7 mmx 114 mm (1), and 10 mm x 114 mm (1)
TM125-115-05 (115V) TM125-220-05 (220V)	Includes 5 mm x 95 mm stainless steel flat bottom TM probe (1)
TM125-115-10 (115V) TM125-220-10 (220V)	Includes 10 mm x 114 mm stainless steel flat bottom TM probe (1)
TMP125-115-0507 (115V) TMP125-220-0507 (220V)	Includes 2 stainless-steel, flat-bottom TM probes -5 mm x 95 mm (1) and 7 mmx 114 mm (1)

Accessories

Part no.	Description
F22000	TM Stand Assembly. Required for all TM-associated part no's to be stand-mounted

Omni TM generator probes (not compatible with other Omni units)

Part no.	Description	Volume range
F2107-05	5 mm x 95 mm, flat bottom (fine to medium samples) (1)	200 µL - 5 mL
F2107	7 mm x 114 mm, flat bottom (fine to medium samples) (1)	200 µL - 5 mL
F2107-10	10 mm x 114 mm, flat bottom (fine to medium samples) (1)	1.5 mL - 100 mL



Omni Tip Generator Probes

Handheld rotor-stators and automated systems

Omni Tip™ plastic single-use probes (flat bottom with B-style drive connect)

Features Lexan® plastic outer tube & Ultem® plastic inner shaft

Note: Compatible as is with Omni THQ and automated systems. Requires separate purchase of Omni Tip adapters for use with Omni TH, micro, and GLH-850.



Part no.	Description	Volume range
30750 32750 34750 35750 3M750	7 mm soft tissue plastic probes (qty 25) 7 mm soft tissue plastic probes (qty 50) 7 mm soft tissue plastic probes (qty 100) 7 mm soft tissue plastic probes (qty 500) 7 mm soft tissue plastic probes (qty 1000)	250 µL - 30 mL
30750H 30750H-S 32750H 34750H 35750H 3M750H	7 mm hard tissue plastic probes (qty 25) 7 mm hard tissue plastic probes - sterile (qty 25) 7 mm hard tissue plastic probes (qty 50) 7 mm hard tissue plastic probes (qty 100) 7 mm hard tissue plastic probes (qty 500) 7 mm hard tissue plastic probes (qty 1000)	250 µL - 30 mL
301250 321250 341250 351250 3M1250	12 mm plastic probes (qty 25) 12 mm plastic probes (qty 50) 12 mm plastic probes (qty 100) 12 mm plastic probes (qty 500) 12 mm plastic probes (qty 1000)	>250 µL. Maximum recommended volume is application specific. Inquire.

Omni Tip hybrid probes (saw tooth B-style drive connect)

Features stainless steel outer tube with replaceable inner Ultem™ plastic shaft

Note: Compatible as is with Omni THQ and automated systems. Requires separate purchase of Omni Tip adapters for use with Omni TH, micro, and GLH-850.



Part no.	Description	Volume range
22-007 22-007-6	7 mm hybrid probe (qty 1), replacement shafts (2), brush (1), and cap (1) 7 mm hybrid probe (qty 6), replacement shafts (12), brushes and caps. Requires separate purchase of Omni Tip adapter for use with Omni TH.	250 µL - 40 mL
301725H	7 mm hybrid probe replacement plastic shafts (25)	250 µL - 40 mL
22-012 22-012-6	12 mm hybrid probe (qty 1), replacement shafts (2), brush (1), and cap (1) 12 mm hybrid probe (qty 6), replacement shafts (12), brushes and caps. Requires separate purchase of Omni Tip adapter for use with Omni TH.	250 µL - 40 mL
S-22-012	12 mm hybrid probe replacement plastic shafts (25)	250 µL - 40 mL

Industrial mixers and macros

Omni Mixer Homogenizer

- Powerful: 600-watt homogenizer, processing speeds up to 18,000 rpm
- Versatile: Wide range of sample volumes from 0.25 mL to 10 L. Compatible with 10 mm to 30 mm diameter stainless steel probes and sealed processing chambers with probe/blade assembly (glass, plastic, or stainless steel)



Units

Part no.	Description
HMG17105 (115V)	Includes stand assembly (1), generator probe adapter (1), and tool kit (1). Requires separate purchase of a generator probe or sealed chamber assembly to operate
HMG17106 (220V)	

Accessories (optional)

Part no.	Description
17204	Quick Connect adapter (1)
17210	Quick connect coupler (1)
17215	Quick connect kit (1 coupler, 3 adapters)

Omni Macro-ES Digital Programmable Homogenizer

- Powerful: 1800-watt homogenizer, processing speeds up to 18,000 rpm
- Programmable: digital display, programming, including cycling, ramping, and timer function
- Versatile: Wide range of sample volumes from 0.25 mL to 30 L. Compatible with 10 mm to 55 mm diameter stainless steel probes and sealed processing chambers with blade assembly (glass, plastic, or stainless steel)



Units and bundles

Part no.	Description
MES-115 (115V)	Includes stand assembly (1), generator probe adapter (1), Quick connect coupler (1), Quick connect adapters (2), and tool kit (1). Requires separate purchase of a generator probe or sealed chamber assembly to operate
MES-220 (220V)	
MESP-115 (115V)	Package additionally includes 3 stainless steel generator probes - 20 mm x 195 mm saw tooth (1), 30 mm x 195 mm flat bottom (1), 45 x 195 mm flat bottom (1)
MESP-220 (220V)	

Accessories

Part no.	Description
17204	Quick Connect adapter (1)
17210	Quick connect coupler (1)
17215	Quick connect kit (1 coupler, 3 adapters)



Omni Macro Homogenizer

- Powerful: 1800-watt homogenizer, processing speeds up to 20,000 rpm
- Versatile: Wide range of sample volumes from 0.25 mL to 30 L. Compatible with 10 mm to 55 mm diameter stainless steel probes and sealed processing chambers with blade included



Units

Part no.	Description
17505 (115V) 17506 (220V)	Includes stand assembly (1), generator probe adapter (1), and tool kit (1). Requires separate purchase of a generator probe or sealed chamber assembly to operate.

Accessories (optional)

Part no.	Description
17204	Quick Connect adapter (1)
17210	Quick connect coupler (1)
17215	Quick connect kit (1 coupler, 3 adapters)

Stainless steel generator probes for Omni Mixer, Macro, and Macro-ES

Part no.	Description	Volume range
15051 15051W 15010	10 mm x 95 mm, saw tooth (standard window/fine to medium samples) 10 mm x 95 mm, saw tooth (wide window/medium to coarse samples) 10 mm x 195 mm, saw tooth (standard window/fine to medium samples)	1.5 mL - 100 mL
15201 15401 15401W 15020 15020W	20 mm x 100 mm, saw tooth (standard window/fine to medium samples) 20 mm x 145 mm, saw tooth (standard window/fine to medium samples) 20 mm x 145 mm, saw tooth (wide window/medium to coarse samples) 20 mm x 195 mm, saw tooth (standard window/fine to medium samples) 20 mm x 195 mm, saw tooth (wide window/medium to coarse samples)	50 mL - 2 L
150-30NA-195 150-30WA-195	30 mm x 195 mm, flat bottom (standard window/fine to medium samples) 30 mm x 195 mm, flat bottom (wide window/medium to coarse samples)	75 mL - 10 L



Large volume stainless steel probes for Omni Macro and Macro-ES

Part no.	Description	Volume range
150-45NA-195	45 x 195 mm, flat bottom (standard window / fine to medium samples)	200 mL - 25 L
150-55NA-195	55 x 195 mm, flat bottom (wide window / medium to coarse samples)	300 mL - 35 L



Sealed chambers with stainless steel probe/blade assembly for Omni Mixer, Macro, and Macro-ES

Glass

Part No.	Description	Volume range
15080	20 mm probe (standard window), 473 mL chamber, particle size to 1 - 2 microns	150 mL - 375 mL
15081	20 mm probe (standard window), 946 mL chamber, particle size to 1 - 2 microns	200 mL - 750 mL
15081W	20 mm probe (wide window), 946 mL chamber, particle size to 1 - 2 microns	
17080	2" blade, 473 mL chamber, particle size to 10 - 15 microns	150 mL - 375 mL
17081	2" blade, 946 mL chamber, particle size to 10 - 15 microns	200 mL - 750 mL



Plastic/Polypropylene

Part No.	Description	Volume range
15065	20 mm probe (standard window), 473 mL chamber, particle size to 1 - 2 microns	180 mL - 375 mL
17065	2" blade, 473 mL chamber, particle size to 10 - 15 microns	150 mL - 375 mL
17098	2" blade, 1892 mL chamber, particle size to 10 - 15 microns	300 mL - 1.2 L



Stainless steel

Part No.	Description	Volume range
15050	10 mm probe (standard window), 50 mL chamber, particle size to 1 - 2 microns	10 mL - 50 mL
15400	20 mm probe (standard window), 400 mL chamber, particle size to 1 - 2 microns	90 mL - 350 mL
17077	1" blade, 50 mL chamber, particle size to 10 - 15 microns	10 mL - 50 mL



Other solutions from Revvity

Automated homogenizers and workstations



Omni Prep (Part No. 06-021)

- Benchtop unit
- up to 6 sample setup



Prep 96 (Part No. 51-02A-1)

- Walk-away homogenizer
- up to 96 sample setup



LH 96 (Inquire)

- Automated weighing, liquid handling, homogenization, and tube to well-plate reformatting
- Up to 96 sample setup

Single-probe sonicators



OM 50 (Part No. 29-060-050-xxx)

- 50 watt, 20 kHz
- Knob and switch controls



OM 100 (Part No. 29-060-100-xxx)

- 100 watt, 20 kHz
- Programmable interface



OM 500 (Part No. 29-060-450-xxx)

- 500 watt, 20 kHz
- Standard and microtips compatible

We also offer [high-speed, tabletop centrifuges](#) to ease workflow bottlenecks.



revvity

www.revvity.com

