

Omni TH

Tissue homogenizer

The Omni TH is a powerful tissue homogenizer that was designed to handle a wide range of samples from <200 μ L to 100 mL. It works quickly and efficiently to process a wide variety of sample types and is compatible with several types of generator probes for flexibility in application requirements.

Patented Omni Tips™ plastic homogenizing probes, in both soft tissue and hard tissue versions, are available for customers concerned about cross-contamination between samples. Stainless steel generator probes are recommended for larger, tougher sample types. Specially designed hybrid plastic/stainless steel probes are also available.

The Omni TH is ideal for a wide range of applications including tissue homogenization, emulsions and suspensions, cell disruption, protein extraction, and medical research. It is lightweight enough to be hand-held and can also be mounted to a motor stand (clamp included) for lengthier processing.

Specifications

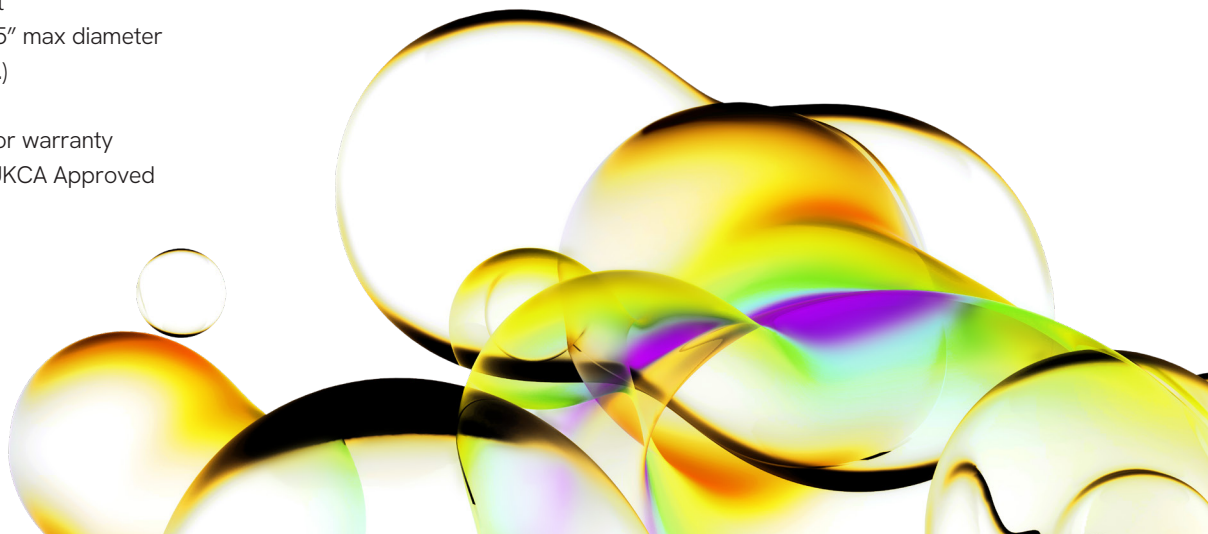
- **Processing range:** <200 μ L to 100 mL
- **Speed:** Variable from 500 to 35,000 rpm
- **Power rating:** 125 watt
- **Dimensions:** 9" H x 2.5" max diameter
- **Weight:** 466 g (16.4 oz.)
- **Sound level:** <72 db
- **Warranty:** 2 year motor warranty
- **Compliance:** CE and UKCA Approved






Part number: TH115 (120V) or TH220 (220V)
Bundle with stand: TH115-K or TH220-K

Key features

- Powerful 125 watt motor
- Light weight ergonomic design with optional post-mounting
- Quiet operation
- 1/4 turn quick connect coupling
- Works with a wide selection of stainless steel and Omni Tip™ plastic homogenizing probes



Stainless steel probes	Diameter	Processing range	Length	Type	Resulting particle size	Order no.
	5 mm	0.20 mL - 5 mL	75 mm	Flat bottom	Fine	G5-75
				Flat bottom - wide window	Coarse	G5-75W
	7 mm	0.25 mL - 30 mL	115 mm	Saw tooth	Fine	G7-115ST
				Saw tooth - wide window	Coarse	G7-115STW
	10 mm	1.5 mL - 100 mL	115 mm	Saw tooth	Fine	G10-115
				Saw tooth - wide window	Coarse	G10-115W

We offer a bundled package that includes 3 probes: 5 x 75 mm Stainless Steel Flat Bottom Probe, 7 x 115 mm Stainless Steel Saw Tooth Probe, and 10 x 115 mm Stainless Steel Saw Tooth Probe. Order Number THP115 (120V) or THP220 (220V)

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 significantly reduce risk of 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.

Features

- Economical: <1% of the cost of stainless steel probes
- Drastically reduces risk of cross contamination
- Ideal for PCR, RNA, and DNA sample preparation
- Maintenance free




“With Omni Tips I can be certain that all the tissue sample has been homogenized; there is no sample loss. I found that the Omni Tip hard tissue version worked as well as the stainless steel probe.”

Rob T., National Institutes of Health


“The Omni hard tissue tips have been indispensable to us in the analysis of heart, skeletal muscle and tumor tissue. Thank you for this product.”

Caryl A., Methodist Research Center

A selection of our best-selling plastic probes

	Omni Tips 7mm Soft Tissue Probes	Quantity	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

	Omni Tips 7 mm Hard Tissue Probes		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

	Omni Tips 12 mm Hard Tissue Probes		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

We offer additional cost-effective product bundles.

[Contact us](#) for more information or to get a quote.

