

Omni THQ

Digital tissue homogenizer

The Omni THQ with Whisper Drive Technology™ is a powerful tissue homogenizer that can handle a wide range of samples from <200µL to 120 mL. It provides reliable, repeatable performance with an ultra-quiet brushless motor. An integrated digital speed control stores the last set speed for ultimate repeatability. Tissue samples are usually processed in less than 30 seconds.

The Omni THQ operates with patented Omni Tip™ Plastic Homogenizing Probes, available in both soft tissue and hard tissue versions. Omni Tips are ideal for customers concerned about cross-contamination and carryover between samples. Stainless steel generator probes (5 mm to 10 mm diameter) are available for larger and tougher sample types. Specially designed hybrid steel/plastic probes are also available.

The Omni THQ is an ideal fit for a wide range of applications from tissue homogenization and dissociation, PD/PK studies, tablet disruption, emulsions and suspensions, cell disruption, food samples and much more. Lightweight enough to be hand-held it may also be stand mounted for lengthier processing.

Specifications

Processing Range: <200 µL to 120 mL

Speed: Variable from 8,000 to 32,000 rpm (increments of 500 rpm)

Power Rating: 40 watt max

Dimensions: 21.8 cm (8.6 in) height, 5.54 cm (2.18 in) max. diameter

Weight: 340 g (12.0 oz.)

Sound Level: <68 db

Warranty: 3 year motor warranty

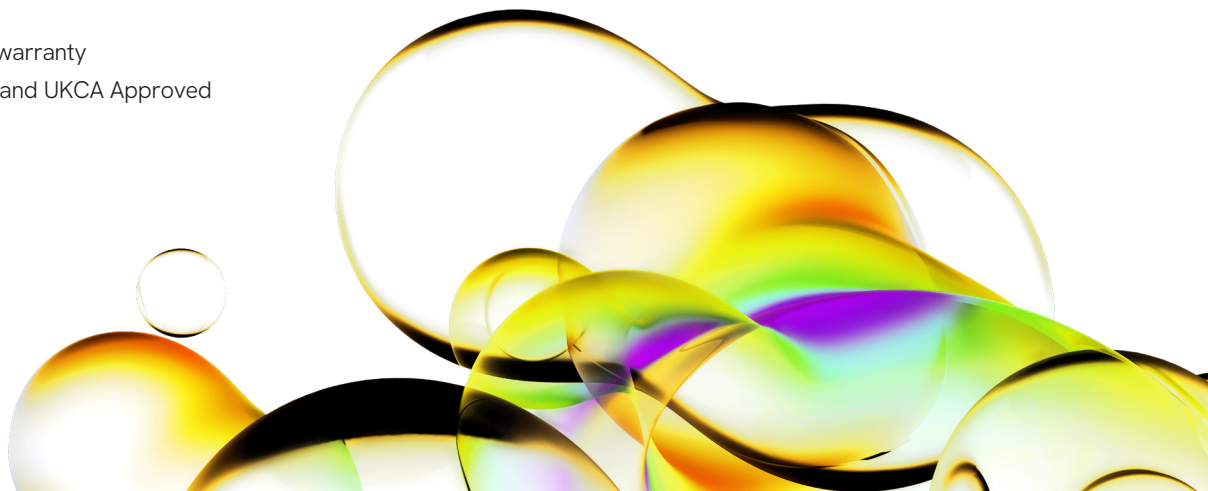
Standards Approval: CE and UKCA Approved



[Part Number: 12-500 \(stand sold separately\)](#)

Key features

- Three digit speed readout accurate to +/- <5 % at top speed
- Whisper Drive Technology™ provides ultra quiet operation
- On-board memory stores last set speed
- Bayonet probe mounting system for fast release
- Uses stainless steel probes or disposable Omni Tip plastic probes
- Motor stand mount included
- Light weight ergonomic design



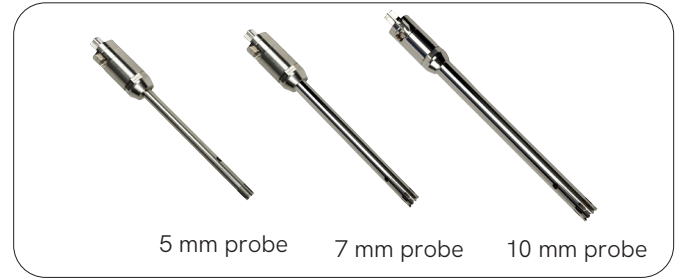
Omni International Stainless Steel Generator Probes

Our Stainless Steel Generator Probes come with PTFE upper and lower bearings. The entire probe can be autoclaved, or easily disassembled for cleaning. Optional Stainless steel upper bearings extend bearing life but are not autoclavable.

With 5 mm probe - [Order number B5-075](#)

With 7 mm probe - [Order number B7-110ST](#)

With 10 mm probe - [Order number B10-110ST](#)



Patented Omni Tip plastic disposable generator probes

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.

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

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

Omni THQ package includes: THQ motor, 3 year motor warranty, stand clamp, 6 Soft Tissue & 6 Hard Tissue Omni Tip Plastic Homogenizer Probes, tool kit, and instruction manual. Order number 12-500

