

Omni Tips

5, 7 and 12 mm disposable homogenizing probes

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

Proprietary Omni Tips probes are engineered from durable plastics and are economical enough to be disposed of after each use to reduce cross contamination during homogenization.

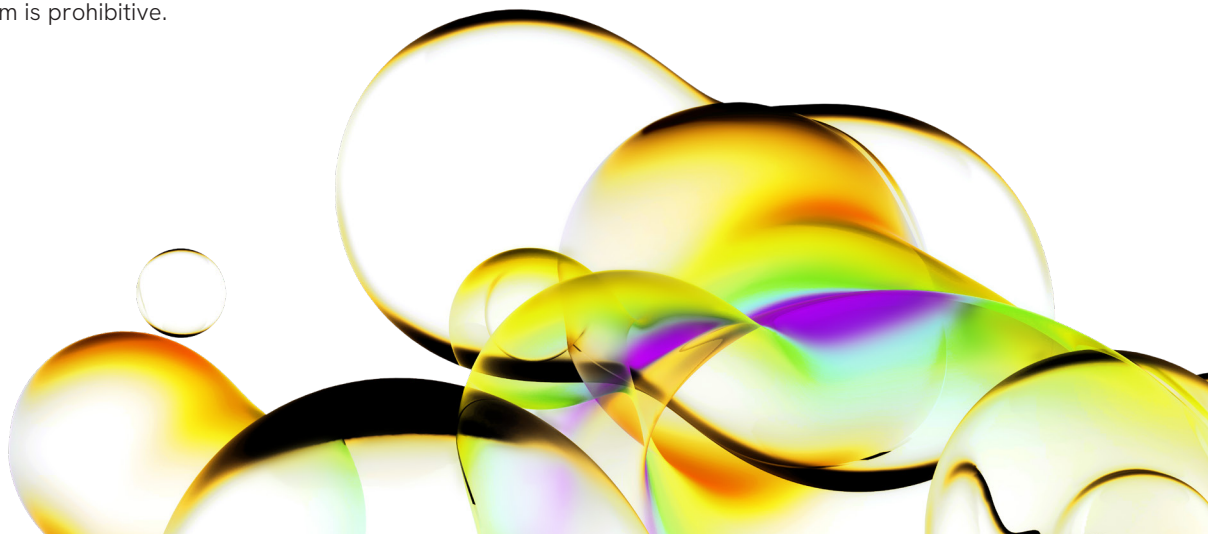
The outer tube is made of polycarbonate. The inner shafts are made of Ultem 1000, a polyetherimide. The patented design avoids the hassle of disassembly and maintenance of traditional stainless steel generators.

12 mm Omni Tips are suitable for homogenization where sample size or sample hardness dictates that processing with the 7 mm tip is prohibitive. 5 mm Omni Tips are suitable for homogenization where the sample size dictates that processing with the 7 mm is prohibitive.



Key features

- Economical: < 1% of the cost of stainless steel probes
- Mitigates cross contamination
- Ideal for PCR, RNA, and DNA sample preparation
- Maintenance free
- Pre-sterilized, electron beam treated
Omni Tips available



"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 Tips™ hard tissue version worked as well as the stainless steel probe."

Rob T., National Institutes of Health

A selection of our best-selling plastic probes

	Omni Tips 5 mm Soft Tissue Probes	Quantity	Order no.
	Omni Tips plastic homogenizing probes	25 pack	30550
	Omni Tips plastic homogenizing probes	50 pack	32550
	Omni Tips plastic homogenizing probes	100 pack	34550

	Omni Tips 7 mm 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	Quantity	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	Quantity	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

Note:

A1000SB adapter is required to use Omni Tips on TH motors with black nose ring, and GLH 850. 12-121 Omni Tip adapter is required use Omni Tips on TH motors with blue nose ring.



Scan or click
to shop now

revvity