

Omni Tip

Homogenizing Kits

Omni Tip™ homogenizing kits come complete with a redesigned TH motor, Omni Tip adapter, 12 hard or soft tissue Omni Tip plastic homogenizing probes, and tool kit. The kits can be used to process samples with a range of <math><200 \mu\text{L}</math> to 100 mL. The perfect solution for a small laboratory, this kit is cost effective and will handle a wide range of applications for multiple users. Kits are also available with optional stainless steel generator probes.

Proprietary Omni Tip probes are available in both soft tissue and hard tissue versions. Engineered from durable plastics, Omni Tip probes are economical enough to be disposed of after each use to eliminate cross contamination.



The outer tube of the Omni Tip probe is made of polycarbonate. The inner shafts are made of Ultem 1000, a polyetherimide. The Omni Tip probes' patented design eliminates the hassle of disassembly and maintenance of traditional stainless steel generators.

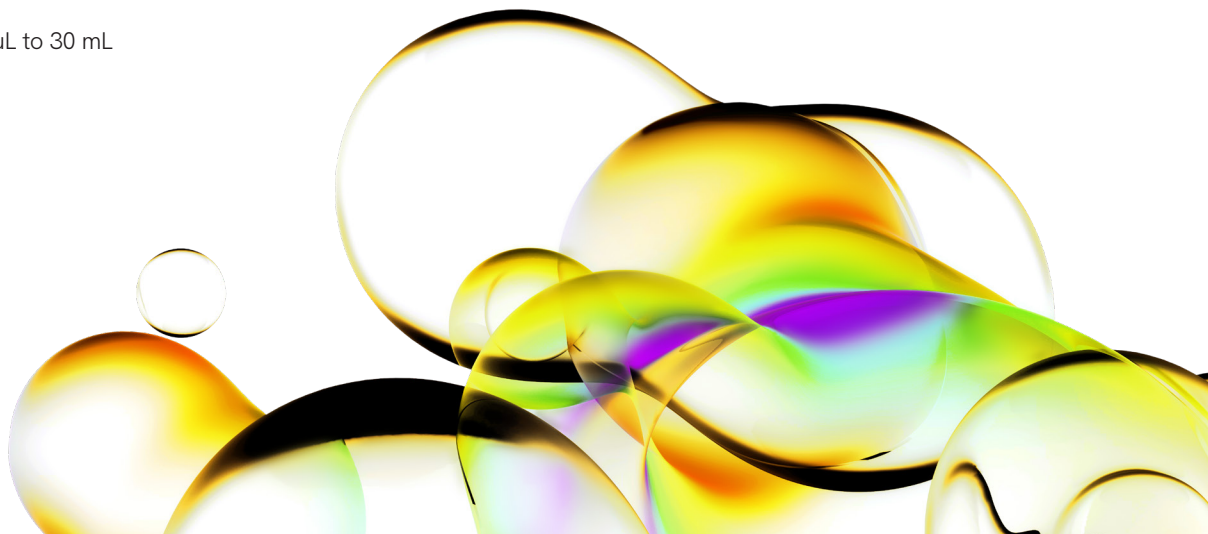
Treatable volume - 250 μL to 30 mL



Part Number: TH115-PCRD (120V) or TH220-PCRD (220V)

Key features

- Economical: <math><1\%</math> of the cost of stainless steel probes
- Drastically reduce cross contamination
- Ideal for PCR, RNA, and DNA sample preparation
- Maintenance Free





Part Number: TH115-PCR7D (120V) or TH220-PCR7D (220V)


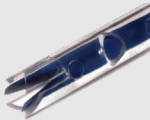
“The hard tissue tips have been indispensable to us in the analysis of heart, skeletal muscle and tumor tissue. Thank you for this product.”

Caryl A., Methodist Research Center

“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

Omni Tip homogenizing kits are also available with additional Stainless Steel Generator Probe for ultimate versatility

Omni Tip homogenizing kit selection guide		
	Soft Tissue Omni Tip Kits For processing liquids, soft tissues and samples that require gentle processing	Order no. (120V or 220V)
	Soft Tissue Omni Tip Kit (includes 7 mm Soft Tissue Omni Tips - quantity 12)	TH115-PCRD or TH220-PCRD
	With 5 x 75 mm stainless steel generator probe (G5-75)	TH115-PCR5D or TH220-PCR5D
	With 7 x 115 mm stainless steel generator probe (G7-115ST)	TH115-PCR7D or TH220-PCR7D
	With 10 x 115 mm stainless steel generator probe (G10-115)	TH115-PCR10D or TH220-PCR10D
	Hard Tissue Omni Tip Kits For processing hard or frozen tissue samples	Order no. (120V or 220V)
	Hard Tissue Omni Tip Kit (includes 7 mm Hard Tissue Omni Tips - quantity 12)	TH115-PCRH or TH220-PCRH
	With 5 x 75 mm stainless steel generator probe (G5-75)	TH115-PCR5H or TH220-PCR5H
	With 7 x 115 mm stainless steel generator probe (G7-115ST)	TH115-PCR7H or TH220-PCR7H
	With 10 x 115 mm stainless steel generator probe (G10-115)	TH115-PCR10H or TH220-PCR10H

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

