

Bioruptor tubes

Diagenode tubes	Volume	# rxns	Sterile	DNase-free	Temperature	Chemical resistance
Bioruptor Pico						
0,2 ml C30010020	20-100 µl	16	non-sterile	DNase-free	Stable from -80° to 120°C	Solvent-resistant
0,65 ml C30010011	100 µl	12	non-sterile	DNase-free	Stable from -80° to 120°C	Solvent-resistant
1,5 ml C30010016	100-300 µl	6	non-sterile	DNase-free	Stable from 4° to 60°C	Not compatible with solvents, strong acids and bases
15 ml C01020031	0,5-2 ml	6	sterile	DNase-free	Stable from 4° to 60°C	Not compatible with solvents, strong acids and bases
Bioruptor Plus						
0,5 ml C30010013	50-100 µl	12	non-sterile	DNase-free	Stable from -80° to 120°C	Solvent-resistant
1,5 ml C30010010	100-300 µl	6	non-sterile	DNase-free	Stable from -80° to 120°C	Solvent-resistant
10 ml C30010012	0,5-2 ml	6	non-sterile	DNase-free	Stable from -80° to 120°C	Solvent-resistant
15 ml C30010009	0,5-2 ml	6	sterile	DNase-free	Stable from -80° to 120°C	Solvent-resistant
Falcon™ 50mL Conical Centrifuge Tubes (Cat No Fisher Scientific 14-432-22)	15 ml	3			Stable from -80°C to ambient temperature	

⚠️ Autoclaving is not recommended for any sonication tubes since it can alter the properties of the plastics and affect the sonication efficiency.