

WHICH TUBES FOR WHICH BIORUPTOR®?

		APPLICATIONS	VALIDATED TUBES					
			0.1 ml	0.5/0.65 ml	1.5 ml	10 ml	15 ml	
BIORUPTOR®	BIORUPTOR® STANDARD/PLUS	Shearing	DNA Shearing	0.5 ml Bioruptor® Microtubes - C30010013				
			Chromatin Shearing	0.5 ml Bioruptor® Microtubes - C30010013	1.5 ml TPX Microtubes - C30010010	10 ml Tube - C30010012	15 ml TPX tubes - C30010009	
		Disruption Extraction	Cell lysis	1.5 ml TPX Microtubes - C30010010				
			Tissue disruption for Protein extraction	1.5 ml TPX Microtubes - C30010010 + Protein extraction beads C20000021				
			Tissue disruption for RNA extraction	1.5 ml TPX Microtubes - C30010010 + Protein extraction beads C20000021				
	Bacteria cell lysis	1.5 ml TPX Microtubes - C30010010	10 ml Tube - C30010012	15 ml TPX tubes - C30010009	15 ml TPX tubes - C30010009 + Protein extraction beads - C20000021			
	FFPE DNA Extraction	1.5 ml TPX Microtubes - C30010010						
	BIORUPTOR® PICO	Shearing	DNA Shearing	0.1 ml Bioruptor® Microtubes - C30010015	0.65 ml Bioruptor® Microtubes - C30010011			
			Chromatin Shearing	0.1 ml Bioruptor® Microtubes - C30010015	0.65 ml Bioruptor® Microtubes - C30010011	1.5 ml Bioruptor® Microtubes with Caps - C30010016	15 ml Bioruptor® Tubes & sonication beads - C01020031	
			RNA Shearing	0.65 ml Bioruptor® Microtubes - C30010011				
Disruption Extraction		Cell lysis	1.5 ml Bioruptor® Microtubes with Caps - C30010016			15 ml Bioruptor® Tubes - C30010017		
		Tissue disruption for Protein extraction	1.5 ml Bioruptor® Microtubes with Caps - C30010016 + Protein extraction beads C20000021			15 ml Bioruptor® Tubes - C30010017 + Protein extraction beads - C20000021		
		FFPE DNA Extraction	1.5 ml Bioruptor® Microtubes with Caps - C30010016					

Tubes and recommended sample volume

