

WHICH TUBES FOR YOUR BIORUPTOR® PICO?



**0.1 ML
BIORUPTOR
MICROTUBES**

**0.2 ML
BIORUPTOR
MICROTUBES**

**0.65 ML
BIORUPTOR
MICROTUBES**

**1.5 ML
BIORUPTOR
MICROTUBES
WITH CAPS**

**15 ML
BIORUPTOR
TUBES**

**15 ML
BIORUPTOR
TUBES AND
SONICATION
BEADS**

Catalog number	C30010015	C30010020	C30010011	C30010016	C30010017	C01020031
Recommended sample volume	5 µl - 50 µl	20 µl - 100 µl	100 µl	100 µl - 300 µl	500 µl - 2 ml	500 µl - 2 ml
APPLICATIONS	DNA shearing	✓	✓	✓	✓	
	Chromatin shearing	✓	✓	✓	✓	✓
	RNA shearing	✓	✓	✓		
	FFPE DNA Extraction				✓	
	Cell lysis		✓	✓	✓	✓
	Bacterial cell lysis		✓	✓	✓	
Tissue disruption for protein extraction				✓*	✓*	

✓ = validated tubes

* Protein extraction beads (C20000021) must be used



If you have any question, please contact the specialist at
<https://www.diagenode.com/en/pages/support>

WHICH TUBES FOR YOUR BIORUPTOR® PLUS?



**0.5 ML
BIORUPTOR
MICROTUBES**

**1.5 ML TPX
MICROTUBES**

**10 ML
BIORUPTOR
TUBES**

**15 ML
TPX TUBES**

**15 ML
TPX TUBES
AND RNA
EXTRACTION
BEADS**

**50ML
FALCON™
TUBES**

Catalog number	C30010013	C30010010	C30010012	C30010009	C20000010	14-432-22 <i>(Fisher Scientific)</i>
Recommended sample volume	50 µl - 100 µl	100 µl - 300 µl	500 µl - 2 ml	500 µl - 2 ml	500 µl - 2 ml	5 ml -15 ml

APPLICATIONS

DNA shearing	✓					
Chromatin shearing	✓	✓	✓	✓		
FFPE DNA Extraction		✓				
Cell lysis	✓	✓	✓	✓		✓
Bacterial cell lysis	✓	✓	✓			
Tissue disruption for protein extraction		✓*		✓*		
Tissue disruption for RNA extraction					✓	

✓ = validated tubes

* Protein extraction beads (C20000021) must be used



If you have any question, please contact the specialist at
<https://www.diagenode.com/en/pages/support>