

WHICH TUBES FOR BIORUPTOR® PICO?



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	C30010020	C30010011	C30010016	C30010017	C01020031
Recommended sample volume	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 Extraction			✓	
	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 (Fisher Scientific)
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>