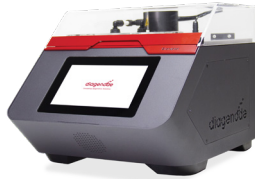


## WHICH TUBES FOR BIORUPTOR® PICO?



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>