

Recommended bead types by sample (2 mL)

Sample Type	Sample Toughness	Glass		Garnet		Ceramic			Metal
		0.1 mm	0.5 mm	0.15 mm	0.7 mm	0.1 mm	1.4 mm	2.8 mm	2.4 mm
Bacteria or yeast culture	Medium	19-621	19-622			19-632			
Plant leaf, root & stem	Medium-Hard							19-628	19-620
Soft soil, sediment, wastewater	Medium			19-623	19-624				
Stool	Medium					19-632			
GI tract with stool	Medium-Hard					19-636D		19-636D	
Soft tissues <small>Brain, liver, kidney</small>	Soft						19-627	19-628	19-620
Hard Tissues <small>Heart, muscle, skin</small>	Medium-Hard							19-628	19-620
Dense biological material <small>Hair, nails, bone, seeds</small>	Hard-Extra Hard						19-627	19-628	
Viruses	Soft					19-627	19-628		
Tissue containing bacteria <small>(tissue lysed, bacteria intact)</small>	Soft-Medium						19-628		
Fruit & vegetables	Soft-Medium						19-628		19-620

Each individual bead type is a potential option for selected sample type

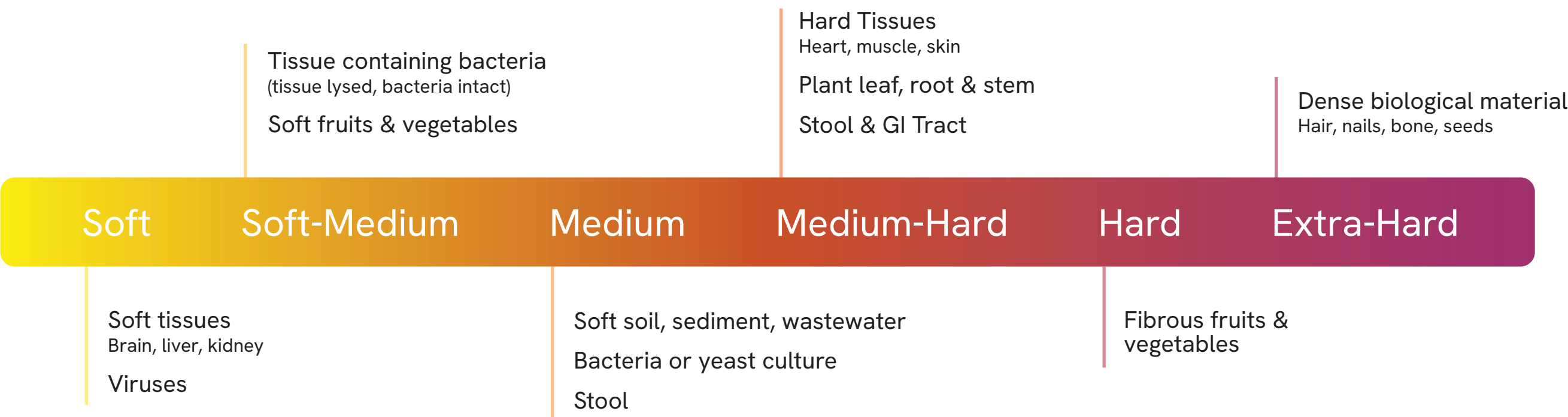
Combined bead types are recommended for use together per sample type

SoftSoft-MediumMediumMedium-HardHardExtra-Hard

Recommended tube size by sample volume

Tube Size	Volume Range							
	20 µL	100 µL	1 mL	2 mL	5 mL	7 mL	15 mL	30 mL
0.5 mL	20 µL - 300 µL							
1.5 mL		200 µL - 1.5 mL						
2 mL	20 µL - 800 µL							
7 mL			1 mL - 3 mL					
15 mL				2 mL - 10 mL				
30 mL					5 mL - 20 mL			
50 mL					5 mL - 35 mL			

Tissue toughness scale
How tough is it to ‘rupt’ your samples?



Which homogenizer is right for you?



Throughput	Less than 25 samples/week	BR 4	BR 12	BR 24	BR Elite	BR 96+	LH 96
	50-400 samples/week						
	500+ samples/week						
Tube size	0.5/2 mL Only						
	2/7 mL Only						
	2-50 mL Only						
	96 Well Plate						

Validate your choice with real data—
browse our app note database.

For application specific inquiries,
please email hmg-sales@revvity.com

