

PRODUCT NAME	
0.65 ml Bioruptor® Microtubes	
Cat. No: C30010011 (WA-005-0500)	Format: 500 tubes
Remark: Low Binding	

Description

The 0.65 ml Bioruptor® Microtubes for DNA and chromatin shearing strongly improve DNA and chromatin sonication efficiency. These tubes have been thoroughly validated for use on the Bioruptor® NGS (UCD-600) as well as on the Bioruptor® Pico (B01060001) and are strongly recommended for DNA and chromatin shearing applications using this instrument.

Characteristics

- Natural
- Graduated
- RNase and DNase free
- Non pyrogenic

Recommended shearing volume

Chromatin shearing: 100 µl

DNA shearing: 100 µl

Remark

Autoclaving is not required for standard use. 0.65 ml Bioruptor® Microtubes withstand 12,000 g.