

**PRODUCT NAME****0.5 ml Bioruptor® Microtubes****Cat. No:** C30010013 (WA-004-0500)**Format:** 500 tubes (2x250 tubes)**Remark:** Low Binding**Description**

The 0.5 ml Bioruptor® Microtubes for DNA shearing strongly improve DNA sonication efficiency. These tubes have been thoroughly validated for use on the Bioruptor® Standard (UCD-200) & Bioruptor® Plus (UCD-300) and are strongly recommended for DNA shearing applications using these instruments.

**Characteristics**

- Natural
- Low binding interior tube surface
- Flat caps surface for easier sample labeling
- Tight sealing caps that minimize evaporation
- Certified free of human DNA, RNase, DNase, and endotoxins
- Non pyrogenic

**Recommended shearing volume**

Chromatin shearing: not recommended

DNA shearing: 100 µl

**Remark**

Autoclaving is not required for standard use.