

PRODUCT NAME**1.5 ml Bioruptor® Microtubes****Cat. No:** C30010016**Format:** 300 tubes**Description**

The 1.5 ml Bioruptor® Microtubes for chromatin shearing strongly improve chromatin sonication efficiency. These tubes have been thoroughly validated for use on the Bioruptor® Pico (B01060001) and are strongly recommended for chromatin shearing applications using this instrument.

Characteristics

- Clear
- Non-steril

Recommended shearing volume: 100-300 µl

Remark

Autoclaving is not required for standard use. 1.5 ml Bioruptor® Microtubes withstand 3.000 g.