

PRODUCT NAME

0.1 ml Bioruptor® Microtubes

Cat. No: C30010015**Format:** 120 tubes**Description**

The 0.1 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® Pico (B01060001) and are strongly recommended for DNA and chromatin shearing applications using this instrument.

Characteristics

- Natural
- RNase and DNase free
- Non pyrogenic

Recommended shearing volume

Chromatin shearing: 10 µl

DNA shearing: 5, 10, or 50 µl

Remark

Autoclaving is not required for standard use).