

PRODUCT NAME	
15 ml Bioruptor® Tubes & sonication beads	
Cat. No: C01020031	Format: 50 rxns

Description

The 15 ml Bioruptor® Tubes & sonication beads for chromatin shearing strongly improve chromatin sonication efficiency. The tubes and beads have been thoroughly validated for use on the Bioruptor® Pico (B01060001) and are strongly recommended for chromatin shearing applications using this instrument.

Characteristics

- Clear
- Steril

Recommended shearing volume

Chromatin shearing: 500 µl - 2 ml (final sample volume incl. beads)

DNA shearing: not recommended

Remark

Autoclaving is not required for standard use.

Recommended amount of beads

500 µl final sample volume incl. beads: 250 mg; equivalent to first line of the 15 ml Bioruptor® Tubes graduation scale

1 ml final sample volume incl. beads: 500 mg; equivalent to second line of the 15 ml Bioruptor® Tubes graduation scale

2 ml final sample volume incl. beads: 800 mg; equivalent to third line of the 15 ml Bioruptor® Tubes graduation scale

Note

Prior to use, beads should be washed by vortexing in 3 volumes of PBS. All buffer should be removed before adding the samples.

