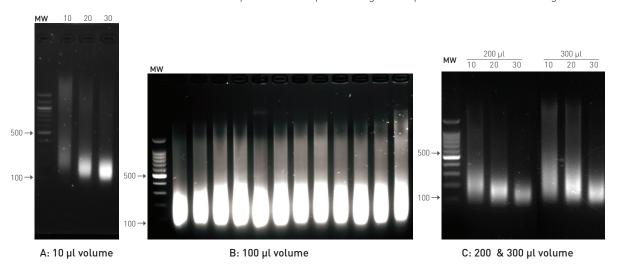


## Consistent and highly reproducible chromatin shearing with the Bioruptor® Pico 10 $\mu l$ - 300 $\mu l$

HeLa cells are fixed with 1% molecular grade formaldehyde (for 8 min at room temperature). Nucleus isolation is performed using buffers and reagents of Diagenode's Chromatin Shearing Optimization kit - Low SDS (Cat. No. AA-001-0100) 1x10e6 cells are then resuspended in 100 µl Shearing Buffer prior to chromatin shearing.



**Panel A**, 10  $\mu$ l volume: Chromatin samples are sheared for 10, 20 and 30 cycles of 30 sec ON/30 sec OFF with the Bioruptor® Pico using 0.1 ml Bioruptor® Microtubes (Cat. No. C30010015).

**Panel B**, 100 µl volume: Chromatin samples are sheared for 10 cycles of 30 sec ON/30 sec OFF with the Bioruptor® Pico using 0.65 ml Bioruptor® Microtubes (Cat. No. WA-005-0500).

**Panel C**, 200 & 300 µl volume: Chromatin samples are sheared for 10, 20 and 30 cycles of 30 sec ON/30 sec OFF with the Bioruptor® Pico using using 1.5 ml Bioruptor® Microtubes (Cat. No. C30010016).

Prior to de-crosslinking, samples are treated with RNase cocktail mixture at  $37^{\circ}$ C during 1 hour. The sheared chromatin is then de-crosslinked overnight and phenol/chloroform purified as described in the kit manual. 10  $\mu$ l of DNA (equivalent of 500, 000 cells) are analyzed on a 2% agarose gel (MW corresponds to the 100 bp DNA molecular weight marker).

(MW corresponds to the 100 bp DNA molecular weight marker)