

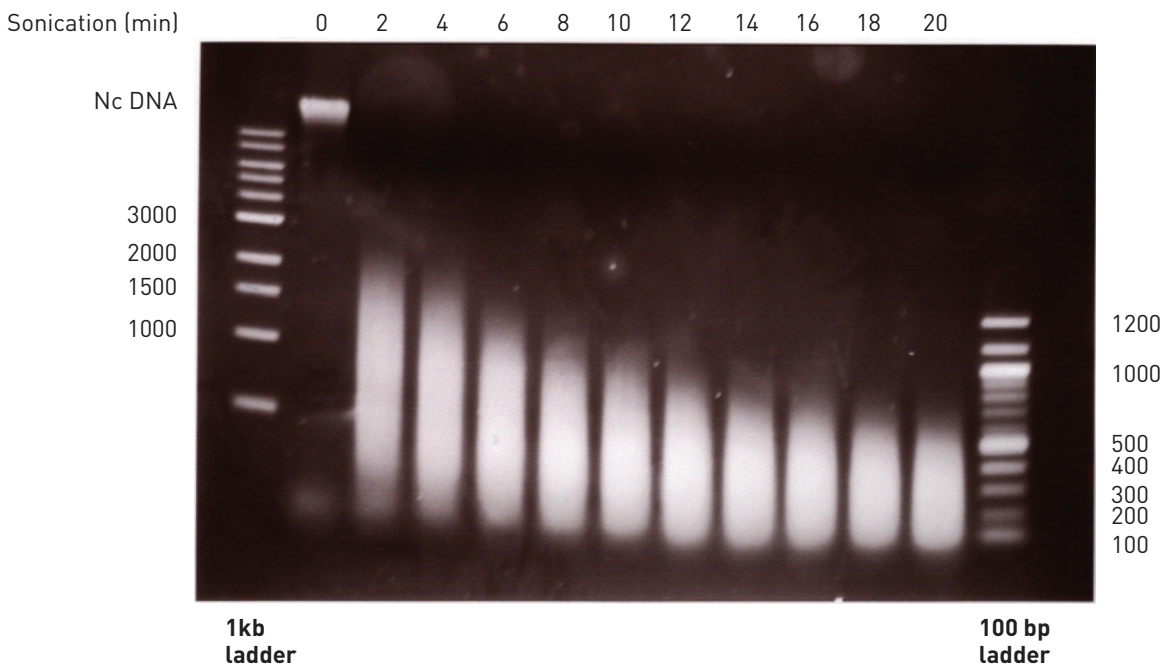
Next Generation SEQUENCING and Bioruptor

Best Solution for Shotgun libraries

Shearing of genomic DNA from *Neurospora*

Authors: Fei Yang, M.S., and Prof. Matthew S. Sachs, Ph.D., Texas A&M University, USA

“We observe a good correlation between sonication time and fragment-size ranges.”



Material: Purified genomic DNA from *Neurospora crassa*

Tubes: 1.5ml Diagenode TPX tubes

Sample: 15µg of purified genomic DNA in 300µl TE buffer (DNA concentration: 0.05µg/µl)

Bioruptor settings: Power: Low; Cycle: 15sec ON / 15sec OFF; Total time: 0-20min.

Cooling: Temperature in sonicator bath maintained stably at 5°C using recirculated chilled water (Lauda).

Agarose gel (1%TAE): 5µl (0.25 micrograms) of sheared DNA was removed at each timepoint for analysis.