

# Digestion with NEBNext dsDNA Fragmentase (M0348)

Protocols.io also provides an [interactive version of this protocol](#) where you can discover and share optimizations with the research community.

## Introduction

**Tip:** Adequate mixing of NEBNext dsDNA Fragmentase is important for the success of this reaction. NEBNext dsDNA Fragmentase should be vortexed for 3 seconds immediately prior to use.

For tough digestions, add 1  $\mu$ l of 200 mM MgCl<sub>2</sub> to the reaction. Additional MgCl<sub>2</sub> can be added if necessary.

The protocol listed below is for fragmentation of 5 ng–3  $\mu$ g of DNA.

## Protocol

- Vortex NEBNext dsDNA Fragmentase for 3 seconds, quick spin and place on ice.
- Combine the following components in a sterile PCR tube and vortex:

DNA (5 ng–3 $\mu$ g)	1–16 $\mu$ l
10X Fragmentase Reaction Buffer v2	2 $\mu$ l
Sterile Water	variable
Final Volume	18 $\mu$ l

- Add 2.0  $\mu$ l dsDNA Fragmentase and vortex the mixture for 3 seconds.  
**Note:** Fragmentase is very viscous and should be pipetted slowly. If the enzyme has been sitting for several minutes vortex it again before adding to the sample.
- Incubate at 37°C for the recommended times below to generate the desired fragment size. To determine the exact incubation time for a given sample type, a time course study should be performed.

Desired Fragment Size (bp)	Incubation Time (min)
50–200	25–35
200–1,000	15–25
1,000–2,000	10–15

\*If starting material is 100 ng or less, incubation times should be increased by 10 minutes.

- Add 5  $\mu$ l of 0.5 M EDTA to stop the reaction.
- Clean up the fragmented DNA with column purification or using SPRI beads. If using SPRI beads, it is recommended to dilute the sample 1:1 with sterile water for easier handling of the sample and faster collection of the beads to the magnet.

**Bioanalyzer:** Clean up the fragmented DNA prior to loading on a Bioanalyzer chip.

**End Repair:** Clean up the fragmented DNA then proceed with desired DNA end repair protocol.

**Polyacrylamide Gel Analysis:** Clean up the fragmented DNA prior to loading the samples on a PAGE gel.

**Long Term Storage:** Clean up the fragmented DNA prior to long term storage.

**Agarose Gel Size Selection/Analysis:** Samples can be loaded directly on to an agarose gel. It is not necessary to clean up the reactions prior to loading.