

## DNAFLUID+ KIT (E07020001)

### SOME MORE TIPS & TRICKS

## INTRODUCTION

The DNAFluid+ Kit (E07020001) is part of a pre-release program. The Diagenode pre-release program allows users to try out a pre-release of the kit and give feedback on their experience. We want to highlight user experience and improve the product accordingly.

The final version of the kit has not been released to the market by Diagenode yet, so please be aware that the protocol provides users with only very few guidelines, and the solution may not generate the expected results for your particular application. Therefore, we would suggest to test on few samples before using critical samples.

This document can constantly be improved thanks to your shared experience. We encourage you to visit our DNAFluid+ kit webpage to follow the new protocols and to share your experience with us. Feel free to take part in the improvements of the solution by giving us your comments, concerns, and suggestions.

## PROTOCOL

1. When the Circulomics Nanobind Kit and protocol is used to **extract (u)HMW DNA**:

- Use no more than 5 million cells/sample,
- Use 150 to 199  $\mu$ l as a final elution volume.

**Optional:** Incubate extracted DNA 10 min at 37°C.

### Remark

*Any viscous DNA extracted with other methods can be treated with the DNAFluid+ kit as long as its purity is acceptable (no cell debris and particle).*

2. Immediately **transfer** your samples to the Hydrotubes using wide bore 1 ml tips.
3. **Spin** your samples briefly (2 - 5 s) to make sure all the solution is in the bottom of the tube.
4. **Start** and **initialize** the Megaruptor.
5. **Go** on 'Protocols' and 'Go & Shear' to start the process.

6. **Set** parameters:

- Set the volume to a value between **150 and 199  $\mu$ l**,

**Remark**

*When the sample volume entered to the software is below the exact value of 200  $\mu$ l, 60 passes through the DNAFluid+ Hydropore will be applied during the homogenization process. It is however possible to treat a higher sample volume with the DNAFluid+ kit. In this case, the current software settings will apply 30 passes (see the manual of Megaruptor 3 for more details). Thus, to have 60 passes with a sample volume higher or equal to 200  $\mu$ l, the same run has to be applied twice.*

- Set the speed to **40**,

**Remark**

*This speed is a good starting point for most of the samples. Depending on the sample viscosity, other speeds between 35 and 59 can be used. Higher speeds are more optimal when the viscosity is higher and vice versa.*

- Set the concentration to any value below **50 ng/ $\mu$ l**.

7. **Plug** your samples to the Hydropore-Syringe.

8. **Insert** the samples in the cassette.

9. **Start** the protocol and check if the sample is passing through the hydropores during the first cycles.

10. Once the **protocol ended**, remove the assembly from the cassette and **unplug** carefully and slowly the Hydropore-Syringe from the Hydrotube. The samples are ready for any downstream step.

11. Don't forget to **share** your experience with us:

<https://www.diagenode.com/en/p/megaruptor-3-DNAFluid-kit>



For any assistance, please, contact us at [www.diagenode.com/en/pages/support](http://www.diagenode.com/en/pages/support) or scan:

