

Dear Customer,

Congratulations on getting early access to our DNAFluid+ Kit. Thank you for participating in the **pre-release program** and helping us to improve our solution.

The new DNAFluid+ Kit eliminates the challenge of handling **viscous Ultra High Molecular Weight DNA samples** for QC and any downstream applications. Using the **DNAFluid+ Kit** with the **Megaruptor 3** enables effective, automated homogenization of highly viscous DNA samples while keeping the integrity of the DNA as well as giving full control of the DNA size.

As part of the pre-release program, we encourage you to visit our DNAFluid+ kit webpage to follow the new upcoming protocols. Please share your experience with us. Feel free to provide your comments, concerns, and suggestions.

Sincerely,
The Diagenode Product Management Team.

PROTOCOL

Some guidelines

1. When the Circulomics Nanobind Kit and protocol is used to **extract (u)HMW DNA**:
 - o Use no more than 5 million cells/sample,
 - o Use 150 to 199 μl as a final elution volume.

Optional: Incubate extracted DNA 10 min at 37°C.
2. Immediately **transfer** your samples to the Hydrotubes using wide bore 1 ml tips.
3. **Spin** your samples briefly (2 - 5 s).
4. **Start** and **initialize** the Megaruptor.
5. **Go** on 'Protocols' and 'Go & Shear' to start the process.
6. **Set** parameters:
 - o Set the volume to a value between 150 and 199 μl ,
 - o Set the speed to 40 (as a good starting point),
 - o Set the concentration to any value below 50 ng/ μl .
7. **Plug** your samples to the Hydropore-Syringe.
8. **Insert** the samples in the cassette.
9. **Start** the protocol and **check** if the samples are passing through the Hydropores.
10. Once the **protocol ended**, 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>

Visit the
dedicated
webpage to
have more
tips and tricks
and share your
experience with us.

