

diagenode

Innovating Epigenetics Solutions

Megaruptor[®]: Your partner in long-read sequencing

Designed to provide simple and automated **DNA sample preparation**, the Megaruptor gives state-of-the-art shearing performance which is independent of the source, concentration, temperature, or salt content of a DNA sample, providing researcher with ultimate flexibility and ease of use. The shearing with the Megaruptor leads to optimal long-read sequencing using **PacBio[®]**, and **Oxford Nanopore[™]** technologies.

Recommended by  PACBIO[®]

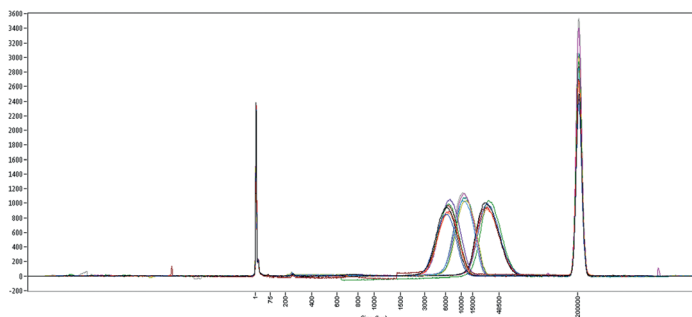


The NEW Megaruptor 3

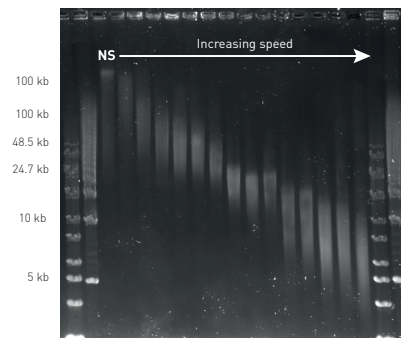
- **High throughput:** Processes **1 to 8 samples** simultaneously.
- **Automated:** Operates without user input and is easily tuneable to between **5 and 100 kb**.
- **Unmatched quality:** Provides the **tight fragment distribution** you need.
- **Safe from contamination and clogs:** Uses optimized consumables.
- **Ergonomic and space-efficient**

Excellent reproducibility and size distribution achieved with the Megaruptor 3

A. Reproducibility and size distribution



B. Versatile fragment range



A: Fragment Analyzer profiles of human genomic DNA samples (25 ng/μl; 200 μl/sample) sheared to 6, 10 and 30 kb. **B:** DNA samples sheared at different speed settings were analyzed by pulse-field gel electrophoresis in 1% agarose gel. (NS: No Sheared)

Megaruptor technical specifications



	Megaruptor 3 Cat. No. B06010003	Megaruptor 2 Cat. No. B06010002
Size range	5 - 100 kb	3 - 75 kb
Volume range	65 – 500 μl	50 – 400 μl
Concentration range	0 – 150 ng/μl	0 – 50 ng/μl
Throughput	1 to 8 samples in parallel	1 to 2 samples in series
Touch screen	✓	✓
Eliminate cross-contamination	with disposable elements in a completely closed-system	with disposable element and with washing sequences
Multiple-orifice disposables	✓	✓
Consumables	Megaruptor 3 Kit : Cat. No. E07010003	Hydropore - short: Cat. No. E07010001 Hydropore - long: Cat. No. E07010002 Hydro Tubes: Cat. No. C30010018

www.diagenode.com Please contact us for more information

Europe Diagenode sa / LIEGE SCIENCE PARK // Rue Bois Saint-Jean, 3 // 4102 Seraing (Ougrée) // Belgium // Phone: [+32] 4 364 20 50 // E-mail: info@diagenode.com
USA Diagenode Inc. / 400 Morris Avenue, Suite 101 // Denville, NJ 07834 // USA // Phone: [+1] 862 209-4680 // E-mail: info.na@diagenode.com