



diagenode

Innovating Epigenetic Solutions

# Bioruptor<sup>®</sup> Sonicator

The industry standard for DNA shearing: Essential for DNA methylation, chromatin immunoprecipitation and high-throughput sequencing studies.

ChIP-seq

ChIP-seq Input DNA

6000  
4000  
3000  
2000  
1500  
1000  
500.0  
200.0  
50.0  
L 1 2 3 4 5 6 7 8 9



# Bioruptor® Sonicator offers unsurpassed reproducibility & quality

## Bioruptor® Sonicator ensures success for a variety of applications

Diagenode's Bioruptor® Sonicator uses state-of-the-art ultrasound technology to disrupt, disperse, or shear a variety of sample types for biological, chemical, pharmaceutical, and industrial applications. The Bioruptor® Sonicator is widely used in biological settings and has proven success for efficient, reproducible sonication required for applications such as DNA/chromatin shearing, cell lysis, cell poration, and protein extraction. Researchers have confirmed the Bioruptor® Sonicator as an optimal shearing system, surpassing industry standards with high yields of superior quality material, as exemplified by over 200 publications.

## Features and benefits of the Bioruptor® Sonicator

The Bioruptor® Sonicator provides:

### ■ Unsurpassed quality, consistency, and efficiency

- Closed tube format prevents cross-contamination and aerosol formation
- Variable power range efficiently and evenly disrupts samples
- Unique cooling system retains integrity of biological complexes
- Gentle ultrasound method preserves samples
- Sample rotation of tubes in water bath ensures shearing and lysis consistency

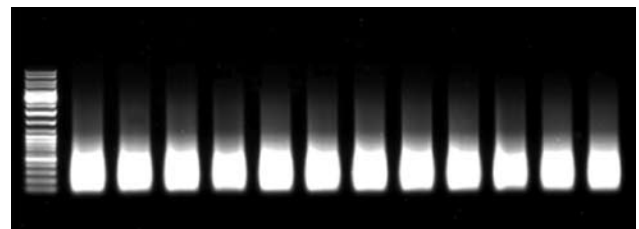
### ■ Compatibility with existing lab workflows: Uses standard microfuge and conical tubes

### ■ Ease of use: Easy set-up, operation, control parameters, and maintenance ensures success

### ■ High-throughput capability: Allows parallel processing of up to 48 samples

### ■ Scaling ability: Interchangeable sample holders allow for microliter to milliliter quantities.

### ■ Application flexibility: Efficient ultrasound technology enables high yields and reproducibility across biological, chemical, pharmaceutical, and industrial applications



**Figure 1. The Bioruptor® achieves highly consistent fragmentation** that ChIP requires in the shearing of chromatin DNA complexes.

### Customer Feedback

"Although many options exist, we routinely generate high quality libraries [using sonication]. The Bioruptor® is lower in cost, while achieving higher energy transfer efficiency and more reproducible performance than standard probe sonicators. In addition, multiple samples can be processed simultaneously in a uniform manner."

Hodges et al. [Nature Protocols 4, - 960 - 974, 2009] describing preparation of DNA for high-throughput sequencing

## Bioruptor® Sonicator shows high reproducibility for high-throughput sequencing library generation

### Application highlight: Bioruptor® in biology

Chromatin Immunoprecipitation (ChIP) and Methylated DNA Immunoprecipitation (MeDIP) Assays:

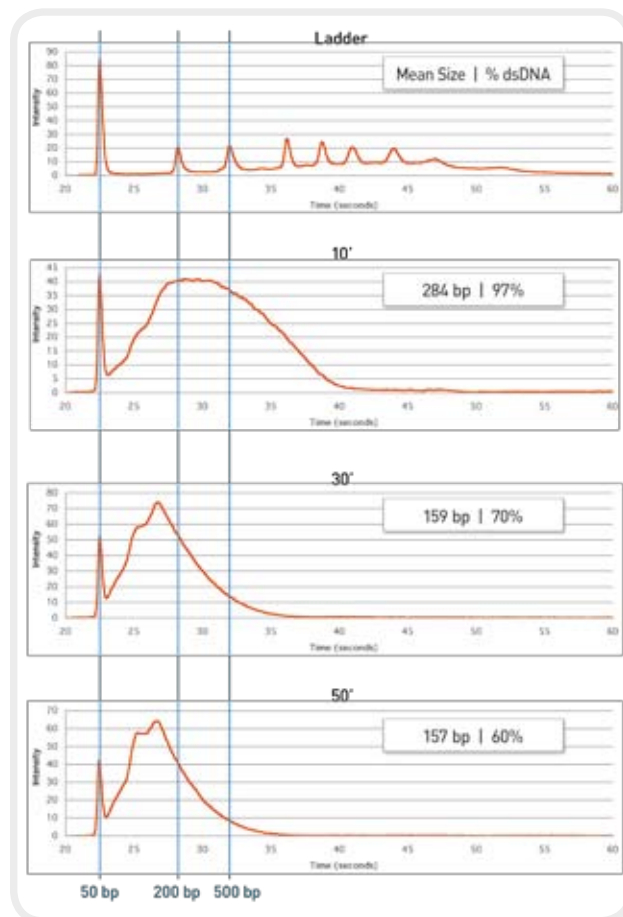
- Consistency and narrow size range of sheared DNA (between 300-600 bp) that ChIP and MeDIP demand
- Sample quality with intact chromatin complexes through the shearing process
- Multiplexing with standard lab tubes
- Compatibility with downstream ChIP and MeDIP applications (e.g. ChIP-qPCR, ChIP-chip, ChIP-seq, MeDIP-seq)
- Widespread success by several prominent chromatin research groups

For a complete list of customer publications and established protocols visit the Diagenode website at [www.diagenode.com](http://www.diagenode.com)

### High-throughput “next generation” sequencing

The Bioruptor® Sonicator is the DNA shearing device of choice for sequencing applications, providing optimal yields, lengths, and consistency as shown in Figure 2. Different fragment size ranges are frequently required for downstream applications (e.g. bridge amplification) for sequencing. The Bioruptor® can be easily programmed to modify duration of sonication for optimal fragmentation and produces:

- Desired narrow size distribution (e.g. 150–250 bp) crucial for sequencing accuracy
- High yields of double-stranded DNA needed for effective sequencing results
- Compatibility with downstream workflow steps (e.g. optimal linker additions, glycerol avoidance)



**Figure 2. Bioruptor® Sonicator achieves optimal size distributions and yields for high-throughput sequencing.** By manipulating the duration of sonication, the Bioruptor® ensures that most DNA fragments lie within the narrow size range required for applications such as ChIP or sequencing. Here, an automated shearing time of 50 minutes is best to maintain yield and complexity for optimal library construction for next generation sequencing, as shown by the narrow size distribution (compared to the DNA ladder) and high dsDNA yields (measured by Qubit; data not shown here).

#### Customer Feedback

“Our experiments indicate that the Bioruptor® provides a wider and more uniform size range compared to nebulization. Additionally, more of the DNA mass is sheared to the relevant size-range. This increases the flexibility of size selection and decreases the amount of DNA lost in the process. The Bioruptor® is also capable of processing up to 6 samples in parallel, while nebulization is performed one at a time, with a single sample per device.”

Shendure lab, University of Washington

# The Bioruptor<sup>®</sup> Sonicator is the instrument of choice

## Choosing the right Sonicator

The Bioruptor<sup>®</sup> Sonicator is the industry's most versatile sonicator. Tip/probe sonicators typically deliver focused energy, and the resulting heat generation may degrade the sample, leading to undesirable effects for ChIP (chromatin disaggregation) and high-throughput sequencing (low yields of double-stranded DNA). Other sonication instruments are not recommended for biological uses given the issues with low-throughput, high set-up costs, and variable sample quality.

	Bioruptor <sup>®</sup> Sonicator	Tip/Probe Sonicators
Wide range of applications	X	X
Enables random shearing vs. enzymatic methods	X	X
Multiplexing with standard lab tubes	X	
Scalable with holders for sample volume flexibility	X	
High-throughput processing ability	X	
Gentle ultrasound irradiation does not degrade sample	X	
Reproducible: Closed system prevents sample contamination	X	
High yields of dsDNA for high-throughput sequencing	X	

## Application versatility

The Bioruptor<sup>®</sup> can be used in industrial and pharmaceutical applications.

See [www.diagenode.com](http://www.diagenode.com) for a complete list of applications.

### Biological applications include:

- Shearing
- Cell disruption
- Lipid dispersion
- Protein solubilization
- Preparation of receptor specimens

### Selected chemical applications include:

- Chemical bond breaking
- Free radical formation
- Reaction catalysis
- Emulsification
- Degassing
- Polymerization

## Cavitation explained

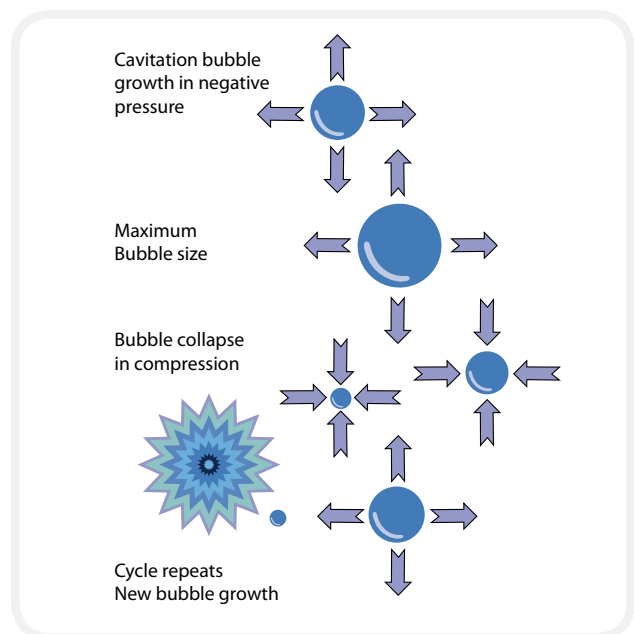
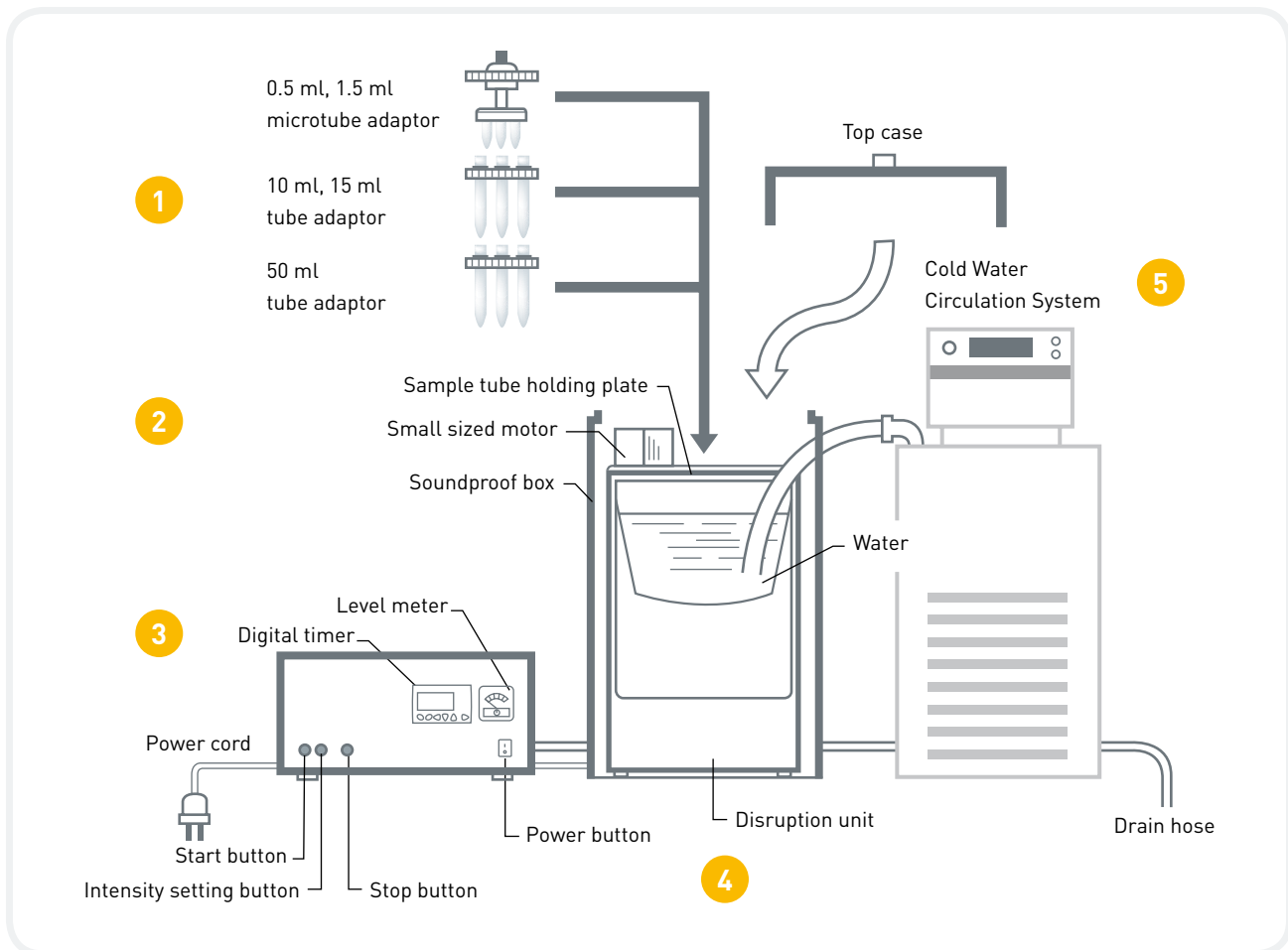


Figure 3. The Bioruptor<sup>®</sup> Sonicator uses ultrasound to create focused mechanical stress to lyse cells or shear DNA or chromatin. The ultrasound waves pass through the sample, expanding and contracting the liquid. During expansion, negative pressures pull the molecules away from one another and a form a cavity or bubble. The bubble continues to absorb energy until it can no longer sustain itself and then implodes, producing intense focused shearing forces, which disperses or breaks biomolecules.

## How the Diagenode Bioruptor® works

Some of the key design features of the Bioruptor® are the laboratory friendly format, ability to use many sample tube types in a water bath-based rotor, and flexible power controls. The walls of the waterbath reflect the ultrasound waves in a random but reproducible pattern. The samples in the adaptor are rotated through the ultrasound field to expose each sample to the same level and intensity of energy. This novel technology enables a wide range of applications for superior yields and quality.



- 1 Flexible sample Format**
- Fits into current workflow with standard tubes
  - Scales with flexible sample volume

- 2 Rotation**
- Prevents contamination with closed tubes
  - Continuous rotation through water bath guarantees equal distribution of energy

- 3 Flexible control**
- Easy to program
  - Power range effectively disrupts samples

- 4 Ultrasound**
- Gentle ultrasound method preserves sample

- 5 Cooling system**
- Cooling system maintains integrity of sensitive samples

# A flexible selection of Bioruptor® models suits your needs

## Attributes of our most popular Bioruptor® Sonicator models



STANDARD



NEXT-GEN



TWIN



XL

	<b>Bioruptor® UCD-200</b>	<b>Bioruptor® UCD-300</b>	<b>Bioruptor® UCD-400</b>	<b>Bioruptor® XL-2006</b>																																																
<b>Description summary</b>	Standard, basic model with reliable performance across applications	Upgrade of UCD-200 with improved displays and controls for greater timing accuracy and temperature control	Provides high-throughput ability and processes up to 24 samples simultaneously	Provides high-throughput ability and processes up to 48 samples simultaneously Ideal for cell lysis Not validated for all kind of DNA and chromatin samples																																																
<b>Throughput &amp; multiplexing</b>	<table border="1"> <thead> <tr> <th>Tube size</th> <th># of tubes processed</th> </tr> </thead> <tbody> <tr> <td>0.5ml</td> <td>12</td> </tr> <tr> <td>1.5ml</td> <td>6</td> </tr> <tr> <td>10ml</td> <td>6</td> </tr> <tr> <td>15ml</td> <td>6</td> </tr> <tr> <td>50ml</td> <td>3</td> </tr> </tbody> </table>	Tube size	# of tubes processed	0.5ml	12	1.5ml	6	10ml	6	15ml	6	50ml	3	<table border="1"> <thead> <tr> <th>Tube size</th> <th># of tubes processed</th> </tr> </thead> <tbody> <tr> <td>0.5ml</td> <td>12</td> </tr> <tr> <td>1.5ml</td> <td>6</td> </tr> <tr> <td>10ml</td> <td>6</td> </tr> <tr> <td>15ml</td> <td>6</td> </tr> <tr> <td>50ml</td> <td>3</td> </tr> </tbody> </table>	Tube size	# of tubes processed	0.5ml	12	1.5ml	6	10ml	6	15ml	6	50ml	3	<table border="1"> <thead> <tr> <th>Tube size</th> <th># of tubes processed</th> </tr> </thead> <tbody> <tr> <td>0.5ml</td> <td>24</td> </tr> <tr> <td>1.5ml</td> <td>12</td> </tr> <tr> <td>10ml</td> <td>12</td> </tr> <tr> <td>15ml</td> <td>12</td> </tr> <tr> <td>50ml</td> <td>6</td> </tr> </tbody> </table>	Tube size	# of tubes processed	0.5ml	24	1.5ml	12	10ml	12	15ml	12	50ml	6	<table border="1"> <thead> <tr> <th>Tube size</th> <th># of tubes processed</th> </tr> </thead> <tbody> <tr> <td>0.5ml</td> <td>48</td> </tr> <tr> <td>1.5ml</td> <td>24</td> </tr> <tr> <td>10ml</td> <td>24</td> </tr> <tr> <td>15ml</td> <td>24</td> </tr> <tr> <td>50ml</td> <td>12</td> </tr> </tbody> </table>	Tube size	# of tubes processed	0.5ml	48	1.5ml	24	10ml	24	15ml	24	50ml	12
Tube size	# of tubes processed																																																			
0.5ml	12																																																			
1.5ml	6																																																			
10ml	6																																																			
15ml	6																																																			
50ml	3																																																			
Tube size	# of tubes processed																																																			
0.5ml	12																																																			
1.5ml	6																																																			
10ml	6																																																			
15ml	6																																																			
50ml	3																																																			
Tube size	# of tubes processed																																																			
0.5ml	24																																																			
1.5ml	12																																																			
10ml	12																																																			
15ml	12																																																			
50ml	6																																																			
Tube size	# of tubes processed																																																			
0.5ml	48																																																			
1.5ml	24																																																			
10ml	24																																																			
15ml	24																																																			
50ml	12																																																			
<b>Timer &amp; controls</b>	Analog timer	Digital timer & controls	Digital timer & controls	Digital timer & controls																																																
<b>Sound-proofing</b>	Plastic soundproof box	Metal soundproof box	Metal soundproof box	Metal soundproof box																																																
<b>Monitoring ability &amp; control systems</b>		Control systems to automatically manage and monitor instrument components	Control systems to automatically manage and monitor instrument components																																																	

### Customer Feedback

"My general conclusion is that the Bioruptor® UCD-300 is much more appealing than the UCD-200 model. This is primarily due to the increased comfort of usability and the homogeneity of the sonication time programming. We find the UCD-300 model takes the robust platform of the UCD-200 and provides perfect sonication efficiency."

Jérémy Dufourt, Laboratory of Dr Chantal Vaury, University of Clermont-Ferrand, France

## Selected Bioruptor® references

The Bioruptor® has been cited in over 200 publications, and is trusted by leading researchers in epigenetics, ChIP, methylation studies, and for library preparation in next generation sequencing.

### ChIP

Cuthbert G. L., Daujat S., Snowden A. W., Erdjument-Bromage H., Hagiwara T., Yamada, M., Schneider R., Gregory P.D., Tempst P., Bannister A.J., Kouzarides T. Histone deimination antagonizes arginine methylation. *Cell*, **118** (5), p.545-553 (2004).

Acevedo, LG, Iniguez, AL, Holster, HL, Zhang, X., Green, R., Farnham, PJ. Genome-scale ChIP-chip analysis using 10,000 human cells. *Biotechniques*. **43**: 791-797 (2007).

Lopez-Serra L., Ballestar E., Fraga M.F, Alaminos M., Setien F. and Esteller M. A Profile of Methyl-CpG Binding Domain Protein Occupancy of Hypermethylated Promoter CpG Islands of Tumor Suppressor Genes in Human Cancer. *Cancer Res* **2006**; **66**: (17) (2006).

Sandmann T., Jakobsen J.S., Furlong, Eileen E.M. ChIP-on-chip protocol for genome-wide analysis of transcription factor binding in *Drosophila melanogaster* embryos. *Nature protocols*, **1** (6), p.2839-2855 (2006).

Navarro P., Page D.R., Avner P. and Rougeulle C. (2006). Tsix-mediated epigenetic switch of a CTCF-flanked region of the Xist promoter determines the Xist transcription program. *Genes & Development* **20**:2787-2792 (2006).

Unité de Génétique Moléculaire Murine, Institut Pasteur 75724, Paris Cedex 15, France.

Durant M. and Pugh B.F. NuA4-Directed Chromatin Transactions throughout the *Saccharomyces cerevisiae*. *Molecular and Cellular Biology*, p. 5327-5335 (2007).

Mikkelsen T.S., Ku M., Jaffe D.B., Issac B., Lieberman E., Giannoukos G., Alvarez P., Brockman W., Kim T.K., Koche R.P., Lee W., Mendenhall E., O'Donovan A., Presser A., Russ C., Xie X., Meissner A., Wernig M., Jaenisch R., Nusbaum C., Lander E.S., Bernstein B.E. Genome-wide maps of chromatin state in pluripotent and lineage-committed cells. *Nature*; **448**:553-560 (2007).

Rosenthal A.Z., Kim Y. and Gralla Y.D. Poising of *Escherichia coli* RNA Polymerase and Its Release from the  $\Delta 38$  C-Terminal Tail for *osmY* Transcription. *J. Mol. Biol.* **376**, 938-949 (2008).

### MeDIP

Yasui D.H., Peddada S., Bieda M.C., Vallero R.O., Hogart A., Nagarajan R.P., Thatcher K.N., Farnham P.J. and LaSalle J.M. Integrated epigenomic analyses of neuronal MeCP2 reveal a role for long-range interaction with active genes. *19416-19421 PNAS December 4, 2007 vol. 104 no. 49*.

### Library preparation for next generation sequencing

Lister R., O'Malley R.C., Tonti-Filippini J., Gregory B.D., Berry C.C., A. Harvey Millar C.A. and Ecker J.R. Highly Integrated Single-Base Resolution Maps of the Epigenome in Arabidopsis. *Cell* **133**, p523-536 (2008).

Gregory B.D., O'Malley R.C., Lister R., Urich M.A., Tonti-Filippini J., Chen H., Millar A.H. and Ecker J.R. A Link between RNA Metabolism and Silencing Affecting Arabidopsis Development. *Developmental Cell* **14**, 854-866 (2008).

Cokus S.J., Feng S., Zhang X., Chen Z., Merriman B., Haudenschild C.D., Pradhan S., Nelson S.F., Pellegrini M and Jacobsen S.E. Shotgun bisulphite sequencing of the Arabidopsis genome reveals DNA methylation patterning. *Nature* **452**, 215-219 (2008).

**Application Note:** Gregory B.D. Next generation sequencing library preparation following preparation of proper-sized DNA molecules using the Bioruptor® [complete application note [http://www.diagenode.com/pages/support\\_app\\_notes.html](http://www.diagenode.com/pages/support_app_notes.html)].

### Protein extraction

Bunk V., Kaehne T., Degtyarev D., Shcherbakova T. and Reiser G. Novel isoenzyme of 2-oxoglutarate dehydrogenase is identified in brain, but not in heart. *FEBS Journal* **275**, 4990-5006 (2008).

S. Lee, V. Horn, E. Julien, Y. Liu, J. Wysocka, B. Bowerman, M. O. Hengartner, W. Herr. Epigenetic Regulation of Histone H3 Serine 10 Phosphorylation Status by HCF-1 Proteins in *C. elegans* and Mammalian Cells. *PLoS ONE* **2**(11): e1213. doi:10.1371/journal.pone.0001213 (2007).

### Dissolution of aggregates

Diehl F., Li M., He Y., Kinzler K. W., Vogelstein B., Dressman D. BEAMing: single-molecule PCR on microparticles in water-in-oil emulsions. *Nature Methods*, **3** (7), p.551-559 (2006).

### Cell Dispersion

Sun B., Tanji Y., Unno H. Influences of iron and humic acid on the growth of the cyanobacterium *Anabaena circinalis*. *Biochemical Engineering Journal* **24**, 195-201 (2005).

### Bacterial Lysis

Md. Kabir S., Yamashita D., Koyama S., Oshima T., Kurokawa K., Maeda M., Tsunedomi R., Murat M., Wada C., Mori H. and Yamada M. Cell lysis directed by DE in early stationary phase and effect of induction of the *rpoE* gene on global gene expression in *Escherichia coli*. *Microbiology*, **151**, 2721-2735 (2005).

## Ordering information

Description	Reference
<b>Bioruptor® Models</b>	
Bioruptor® Standard	UCD-200 TM (1.5 mL) UCD-200 TO (1.5 mL & 15 mL)
Bioruptor® Standard (new)	UCD-300 TM (1.5 mL) UCD-300 TO (1.5 mL & 15 mL)
Bioruptor® Twin	UCD-400 TM (1.5 mL) UCD-400 TO (1.5 mL & 15 mL)
Bioruptor® XL	XL TM (1.5 mL) XL TO (1.5 mL & 15 mL)
<b>Sample Tube Adaptor Kits</b>	
0.5 mL (12) Microfuge tube accessory kit	UCD-pack 0.5
1.5 mL (6) Microfuge tube accessory kit	UCD-pack 1.5
10 mL (6) tube accessory kit	UCD-pack 10
15 mL (6) tube accessory kit	UCD-pack 15
50 mL (3) tube accessory kit	UCD-pack 50
<b>Cooling System</b>	
Water Cooler	BioAcc-cool
Connectors Kit for Water Cooler	CONN-7D1
Peristaltic pump (including Connectors kit)	TWI-pump
<b>Valve Kits</b>	
Valve kit for UCD-300	VB-100-0001
Valve kit for UCD-400	VB-101-0001



## Diagenode headquarters

### Diagenode s.a. BELGIUM | EUROPE

Avenue de l'hôpital, 1  
Tour GIGA, 3rd Floor  
4000 Liège - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

### Diagenode Inc. USA | NORTH AMERICA

376 Lafayette Road, Suite 202  
Sparta, NJ 07871 - USA  
Tel: +1 973-300-0976  
Fax: +1 973-300-1862  
orders.na@diagenode.com  
info.na@diagenode.com

## Regional offices

### | FRANCE

Centre d'affaire Genopole  
Rue Henri Desbruères, 5  
Bât. Genavenir 8  
91030 Evry cedex, France  
Tel: +33 1 60 87 14 82

### | UK

Wellington House, East Road, Room 101  
CB1 1BH Cambridge, UK  
Tel: +44 1223 911396  
Fax: +44 1223 451100

For a complete listing of Diagenode's international distributors visit:

<http://www.diagenode.com/pages/distributors.html>  
For the rest of the world, please contact Diagenode s.a.