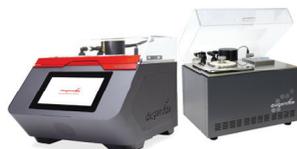


BIORUPTOR® FEATURES



Technical specification	Bioruptor® Pico	Bioruptor® Plus
Recommended sample volume	Excellent	Excellent
Fast processing*	+++	++
Shearing volume	(5 µl)** 20 µl - 2 ml	50 µl - 20 ml
Throughput	Up to 16 samples per run	Up to 12 samples per run
Low noise level	+++	+
Small footprint	+++	+

* Time required to simultaneously shear 12 genomic DNA (100ng/ul, 100ul samples) to 200 bp with Bioruptor Pico = 13 min. and Bioruptor Plus = 30min.

** -Pico 2013-2019



If you have any questions, please contact the specialist at
<https://www.diagenode.com/en/pages/support>

BIORUPTOR® APPLICATION VERSALITY



Applications	Bioruptor® Pico	Bioruptor® Plus
DNA shearing	Excellent	Good
Chromatin shearing	Excellent	Excellent
Cell and tissue disruption (mass spectrometry or other applications)	Excellent	Excellent
RNA shearing	Excellent	Less suitable
DNA extraction from FFPE samples	Excellent	Excellent
Cell and tissue disruption for RNA extraction	Not suitable	Excellent
Bacterial cell lysis	Excellent but limited to small volumes 20x16-300x6 µl	Excellent volumes 50 µl*12-15 mlx3
Compound solubilization	Less suitable	Excellent
Liposome size reduction	Less suitable	Excellent
Protein aggregation (seeding standardization)	Excellent	Excellent



If you have any questions, please contact the specialist at <https://www.diagenode.com/en/pages/support>

BIORUPTOR® COMPETITIVE LANDSCAPE

diagenode

SONICATION

QIMAGEN

Covaris

Features/benefits	Bioruptor®	Probe / water bath sonicators	Probe / cup horn sonicators	Focused ultra-sonicators
Wide range of applications	✓	✓	✓	✓
Enables random shearing	✓	✓	✓	✓
Closed system prevents sample contamination	✓		✓	✓
High-throughput processing	✓		✓	✓
Reproducibility between samples	✓			✓
Scalability and sample volume flexibility	✓		✓	
Low consumable cost	✓		✓	
Gentle ultrasound and temperature control prevent sample degradation	✓			
High yields of chromatin and dsDNA for high throughput sequencing	✓			



If you have any questions, please contact the specialist at <https://www.diagenode.com/en/pages/support>