

EpiShear™ Sonicator

consistent preparation of sheared chromatin samples ensures reproducible ChIP results

Active Motif's EpiShear™ Sonicator is ideal for preparing sheared chromatin from small or large sample sizes for use in chromatin immunoprecipitation (ChIP), as well as for standard cell disruption, DNA/RNA shearing and other homogenization applications.

The EpiShear is a microprocessor-based, ultrasonic processor that offers both programmable and manual operation. Features include pulse mode and a digital display of both wattage and joules. The unit is supplied with an ultrasonic electric generator, a piezoelectric converter, a 1/8" tip probe, and a comprehensive manual. It is backed by a two-year warranty.

A variety of accessories are also available, including different sized microtip probes, a support stand / convert clamp for hands-free use and a sound enclosure to reduce noise. For more complete details, please visit www.activemotif.com/sonicator.



The EpiShear generator, converter and 1/8" tip probe.

EpiShear Features

- **Programmable operation**
Set the time and amplitude so you can perform hands-free operation
- **Pulse mode**
Prevents the buildup of heat in temperature-sensitive samples
- **Digital amplitude / intensity control**
Output intensity can be set from 20-100%
- **Elapsed time indicator**
Displays duration of sonication
- **Display of wattage and joules**
Real-time energy monitoring
- **Overload protection**
Prevents damage to circuitry in the event of a fault occurring
- **Compact design**
Takes up less space than other units

Technical Specifications

Sample size: 200 µl - 50 ml
(500 µl - 15 ml with the included 1/8" probe)

Power Rating: 120 watts

Frequency: 20 kHz

Programmable Timer: 10 hours

Adjustable Pulse: 1 second to 59 seconds

Dimensions: 8" W x 13.75" L x 5.75" H
(203 mm x 349 mm x 146 mm)

Voltage: 110V @ 50/60 Hz or
230V @ 50/60 Hz

EpiShear Includes

- Generator
- Converter
- 1/8" tip probe
- Power cable
- Converter cable
- Wrench set

Product	Voltage	Catalog No.
EpiShear™ Sonicator	110V	53051
	230V	53052