

Quick Guide: PIXUL™ Multi-Sample Sonicator

This Quick Guide provides instructions for how to operate the PIXUL Multi-Sample Sonicator (cat. no. 53130).

IMPORTANT: PIXUL sonication requires using PIXUL 96-well round bottom plates (cat. no. 53139). Using any other plates may result in inefficient sonication, may damage the instrument transducers, and will void the instrument warranty.

How to Turn On

1. Turn on the large switch on the back of the instrument to the ON position. The light around the center power button on the front of the instrument should illuminate.
2. Press the center power button. The touchscreen will begin initializing and load the home screen.

How to Get Ready for Sonication

1. Once PIXUL is on, the first step is to check the Coupling Fluid level (cat. no. 53136).
2. Coupling Fluid should be just above the circular cut-out in the plastic column within reservoir. The reservoir capacity is approximately 500 mL, and the fluid level should be topped off as needed.

How to Chill the PIXUL Prior to Sonication

1. Place the plate to be sonicated on the grid in the PIXUL instrument.
NOTE: Using PIXUL 96-well round bottom plates (cat. no. 53139) is required and all wells without samples in the columns containing samples being sonicated must be filled with liquid (water, buffer, etc.) prior to sonicating.
2. Secure plate by closing lid and fastening tension rods.
3. Once secured, press the “circulate” button and allow the instrument to cool to at least 17 °C. This should take 10-15 minutes.

How to Program the Instrument

1. Use the touchscreen to select the wells you want to sonicate.
2. We recommend using the following parameter specifications as a starting point, and adjusting Process Time to optimize for your specific sample and application:

Sonication Parameter	Setting
Pulse [N]	50
PRF [kHz]	1.00
Process Time [min]	36:00
Burst Rate [Hz]	20.00

3. Once programmed and the plate is secure, make sure the coupling fluid is circulating and press “start” on the touchscreen.
4. The time to completion will be on the top left corner of the display.
5. Once sonication is complete, open the lid and allow coupling fluid to drain from under the plate (this will happen automatically, and will take 5-10 seconds).
6. Remove plate from instrument and close lid but do not fasten the tension rods.
7. Turn off instrument in reverse order as it was turned on.

Technical Support

If you need assistance at any time, please contact Active Motif Technical Support at tech_service@activemotif.com.