

PIXUL™ Chromatin Shearing Kit Manual

Catalog No. 53132

(version A1)

Active Motif North America

1914 Palomar Oaks Way, Suite 150
Carlsbad, California 92008, USA
Toll free: 877 222 9543
Telephone: 760 431 1263
Fax: 760 431 1351

Active Motif Europe

Waterloo Atrium
Drève Richelle 167 – boîte 4
BE-1410 Waterloo, Belgium
UK Free Phone: 0800 169 31 47
France Free Phone: 0800 90 99 79
Germany Free Phone: 0800 181 99 10
Telephone: +32 (0)2 653 0001
Fax: +32 (0)2 653 0050

Active Motif Japan

Azuma Bldg, 7th Floor
2-21 Ageba-Cho, Shinjuku-Ku
Tokyo, 162-0824, Japan
Telephone: +81 3 5225 3638
Fax: +81 3 5261 8733

Active Motif China

787 Kangqiao Road
Building 10, Suite 202, Pudong District
Shanghai, 201315, China
Telephone: (86)-21-20926090
Hotline: 400-018-8123

Copyright 2020 Active Motif, Inc.

TABLE OF CONTENTS	Page
Overview	1
Kit Components and Storage	2
Reagents for Chromatin Shearing	2
Additional Materials Required	2
Protocol	3
Technical Services	5

Information in this manual is subject to change without notice and does not constitute a commitment on the part of Active Motif, Inc. It is supplied on an “as is” basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time.

This documentation may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Active Motif, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties.

PIXUL™ is sold under an exclusive license to patents owned by Matchstick Technologies Inc. and University of Washington, specifically Matula, T.J.; Bomsztyk, K.; Darlington, G.P., Maxwell, A.D.; MacConaghy, B.E., Reed, J. “Ultrasound system for shearing cellular material”. US Patent US10809166. European Patent EP3169451.

The manufacturer of this documentation is Active Motif, Inc.

© 2020 Active Motif, Inc., 1914 Palomar Oaks Way, Suite 150; Carlsbad, CA 92008. All rights reserved.

All trademarks, trade names, service marks or logos referenced herein belong to their respective companies.

For Research Use Only. Not for use in diagnostic procedures.

Overview

Highly consistent shearing of chromatin is critical for successful downstream chromatin immunoprecipitation assays. Manual shearing with a probe sonicator or shearing with an inconsistent high-throughput sonicator can lead to sample-to-sample variability.

The PIXUL Chromatin Shearing Kit delivers highly consistent average fragment lengths across entire plates time after time. This kit is compatible with a wide range of fixed cells (10,000 - 5,000,000) with very low standard deviations in sonication efficiency.

PIXUL Advantages

- Process 1 - 96 samples in parallel
- Extremely consistent sonication
- Up to 12 different sonication conditions per run

product	format	catalog no.
PIXUL™ Chromatin Shearing Kit	1 X 96 reactions	53132

Kit Components and Storage

The kit contains sufficient reagents to sonicate 96 wells in one run. The reagents in this kit have multiple storage temperatures. Please store components according to the storage conditions below. All reagents are guaranteed stable for 6 months from date of receipt when stored properly.

Reagents	Quantity	Storage
Fixation Buffer	75 μ L	4°C
Stop Solution	560 μ L	RT
10X PBS	6 mL	4°C
Protease Inhibitor Cocktail	2 X 100 μ L	-20°C
Cell Shearing Buffer	11 mL	RT
PIXUL™ 96-well plate with sealer	1 unit	RT

Additional Materials Required

- 100% Ethanol
- Nuclease-free water
- 16% methanol-free formaldehyde
- 1X PBS
- PIXUL Multi-Sample Sonicator (Cat. No. 53130)
- Microcentrifuge tubes
- PCR tube strips (300 μ L well volume capacity)
- Centrifuge to spin 96-well plate
- Thermal cycler or other 37°C incubator
- 96-well plate magnet or bar magnet
- Single channel and multi-channel pipettors and tips
- Hemocytometer and light microscope

PIXUL Chromatin Shearing Kit Protocol

IMPORTANT: PIXUL sonication requires the use of PIXUL 96-well round bottom plates. These plates are available as Active Motif Catalog No. 53139, and these are the same plate as Corning Catalog No. 3799. Using any other plates will result in inefficient sonication, may damage the instrument, and will void the instrument warranty.

Buffer Preparation

Prepare the following buffers before starting the protocol.

1X PBS

- 4.5 mL 10 X PBS
- 40.5 mL nuclease-free water

Fixation Solution (Stock Solution)

- 293 μ L nuclease-free water
- 90 μ L Fixation Buffer
- 867 μ L 16% methanol-free formaldehyde

Fixation Solution in PBS (Working Solution)

- 1 mL Fixation Solution (Stock Solution)
- 10 mL 1X PBS

1X Stop Solution

- 0.6 mL Stop Solution
- 11.4 mL 1X PBS

Cell Shearing Buffer with Inhibitors

- 10 mL Cell Shearing Buffer
- 100 μ L Protease Inhibitor Cocktail

Cell Fixation

1. Culture cells directly in the PIXUL 96-well round-bottom plate or cell culture dish or flask of your choice.
2. If you are using the PIXUL 96-well plates for cell culture, carefully aspirate media and add 100 μ L Fixation Solution (Working Solution).
3. If you are not using the PIXUL 96-well plate for cell culture, collect cells, and resuspend the appropriate number of cells in 100 μ L Fixation Solution (Working Solution) and transfer to the PIXUL 96-well plate.
4. Incubate for 10 minutes at room temperature. If you are working with suspension cells, collect cells by centrifugation.
5. Aspirate Fixation Solution and add 100 μ L 1X Stop Solution.
5. Incubate for 5 minutes at room temperature. If you are working with suspension cells, collect cells by centrifugation.
6. Aspirate 1X Stop Solution and wash the cells twice with 100 μ L 1X PBS. If you are working with suspension cells, collect cells by centrifugation before aspiration.

PIXUL Sonication of Fixed Cells

1. Aspirate PBS and add 100 μ L Cell Shearing Buffer with Inhibitors to the fixed cells and seal the plate with a PIXUL plate seal.

Note: All wells lacking sample in the columns being sonicated **MUST** be filled with liquid (water, coupling buffer, etc.) prior to starting the run.

2. Turn on and run the instrument as specified in the PIXUL Multi-Sample Sonicator User Manual.
3. We recommend using the following parameter specifications as a starting point, and adjusting Process Time to optimize for your specific sample and application:

Labile marks, like phosphorylated epitopes, may be preserved better by discontinuous sonication (for example, rather than 30 minutes, 4 rounds of 4 - 6 minutes each, where the PIXUL Coupling Fluid is allowed to circulate and cool between runs).

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

Note: The instrument software will not allow settings that could damage the transducers. The upper limits are a maximum Pulse of 50 or a maximum PRF of 1.00. We recommend a minimum Burst Rate of 20.00. Varying the Process Time is typically the only parameter that needs adjusting, with longer Time needed for difficult samples such as adipose tissue.

Analysis of Sonicated Chromatin

Remove 10 μ L of the sonicated chromatin and use the PIXUL Chromatin Input Kit (Cat. No. 53134) to determine yield and sonication efficiency. We recommend shearing chromatin to 200 - 500 bp in length for ChIP assays.

Using Chromatin Sheared Using PIXUL in ChIP Assays

Chromatin sheared with PIXUL is compatible with many ChIP protocols, but for best results we recommend performing ChIP assays using the ChIP-IT High Sensitivity® Kit (Cat. No. 53040) or High-Throughput ChIP-IT® Kit (Cat. No. 53146).

Technical Services

If you need assistance at any time, please call Active Motif Technical Service at one of the numbers listed below.

Active Motif North America

1914 Palomar Oaks Way, Suite 150

Carlsbad, CA 92008, USA

E-mail: tech_service@activemotif.com

Toll Free: 877 222 9543

Direct: 760 431 1263

Fax: 760 431 1351

Active Motif Europe

Waterloo Atrium

Drève Richelle 167 – boîte 4

BE-1410 Waterloo, Belgium

E-mail: eurotech@activemotif.com

Direct: +32 (0)2 653 0001

Germany Free Phone: 0800 181 99 10

France Free Phone: 0800 90 99 79

UK Free Phone: 0800 169 31 47

Fax: +32 (0)2 653 0050

Active Motif Japan

Azuma Bldg, 7th Floor

2-21 Ageba-Cho, Shinjuku-Ku

Tokyo, 162-0824, Japan

Direct: +81 3 5225 3638

Fax: +81 3 5261 8733

E-mail: japantech@activemotif.com

Active Motif China

787 Kangqiao Road

Building 10, Suite 202

Pudong District

Shanghai, 201315, China

Direct: (86)-21-20926090

Hotline: 400-018-8123

E-mail: techchina@activemotif.com

Visit Active Motif online at <http://www.activemotif.com>

At this site:

- Read about who we are, where we are, and what we do
- Review data supporting our products and the latest updates
- Share your ideas and results with us
- View our job opportunities

Don't forget to bookmark our site for easy reference!