

PIXUL™ Chromatin Input Prep Kit Manual

Catalog No. 53134

(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 Input Prep	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

This kit is designed to analyze yield and sonication efficiency of the chromatin sheared with the PIXUL™ Chromatin Shearing Kit (Cat. No. 53132) using the PIXUL™ Multi-Sample Sonicator (Cat. No. 53130).

The kit contains sufficient reagents to prepare ChIP-ready chromatin for 96 IP reactions.

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 Input Prep Kit	1 X 96 reactions	53134

Kit Components and Storage

The kit contains sufficient reagents to prepare analyze ChIP-ready chromatin for 96 IP reactions. 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
SPRI Beads	3.98 mL	4°C
Elution Buffer AM4	1.1 mL	RT
Proteinase K	56 µL	-20°C
RNase A	2 X 45 µL	-20°C
Low EDTA TE	2 X 1.2 mL	RT

Additional Materials Required

- 100% Ethanol
- Nuclease-free water
- 1X PBS
- Microcentrifuge tubes
- PCR tube strips (300 µL well volume capacity)
- Centrifuge to spin 96-well plate
- Thermal cycler
- 96-well plate magnet or bar magnet
- Single channel and multi-channel pipettors and tips
- Hemocytometer and light microscope

PIXUL Chromatin Input Prep Kit Protocol

Buffer Preparation

1. Prepare TE-RNase by combining 95 μL Low EDTA TE and 5 μL RNase A (10 $\mu\text{g}/\mu\text{L}$).
2. Prepare Elution Buffer with Proteinase K (EB-PK) by combining 95 μL Elution Buffer AM4 with 5 μL Proteinase K (10 $\mu\text{g}/\mu\text{L}$).
3. Prepare 80% ethanol (4.8 mL 100% ethanol + 1.2 mL nuclease-free water).

Reverse Crosslinking

1. Remove 10 μL of the chromatin that has been sonicated with the PIXUL Chromatin Shearing Kit (cat. no. 53132) and transfer to a PCR tube, strip-well, or 96-well plate
2. Add 10 μL TE-RNase to each sample and incubate for 30 minutes at 37°C.
3. Add 10 μL EB-PK to each sample and incubate for 30 minutes at 55°C, followed by 90 minutes at 65°C.

DNA Purification using SPRI Beads

Take SPRI Beads out of 4°C and warm to room temperature by simply placing the vial of SPRI beads on the benchtop.

1. Add 0.5 μL RNase A to each well and incubate for 30 minutes at 37°C.
2. Add 5 μL Proteinase K to each well and incubate for 30 minutes at 55°C.
3. Add 120 μL (1.2 volumes) of SPRI Beads to each well. Pipette the bead-lysate mixture 10 times to mix, and incubate for 5 minute at room temperature.
4. Place plate on a magnet for 5 minutes to collect beads, and then aspirate supernatant.
5. Keeping the plate on the magnet, add 180 μL of 80% ethanol and incubate for 30 seconds at room temperature.
6. Aspirate the supernatant and perform a second wash with 180 μL of 80% ethanol.
7. Completely aspirate the supernatant and air dry the beads until they are no longer shiny (2-5 minutes).
8. Remove the plate from the magnet and elute DNA by adding 50 μL of Low EDTA TE.
9. Pipette the beads 10 times and incubate for 3 minutes at room temperature.
10. Place tube or plate back on the magnet for 5 minutes to collect beads.
11. Carefully remove eluted DNA in the supernatant and transfer to a new tube or plate.
12. Analyze gDNA yield using a spectrophotometric (Nanodrop) or fluorometric (Qubit) method of your choice and analyze fragmentation efficiency using either agarose gel electrophoresis, Agilent TapeStation, or Agilent Fragment Analyzer. For most applications, we recommend shearing DNA to 200 - 600 bp.

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!