

# Chromatin IP DNA Purification Kit

(version A2)

Catalog No. 58002

## **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**

Avenue Reine Astrid, 92

B-1310 La Hulpe, 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

Copyright 2014 Active Motif, Inc.

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.

The manufacturer of this documentation is Active Motif, Inc.

© 2014 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.

<b>TABLE OF CONTENTS</b>	<b>Page</b>
<b>Overview</b> .....	1
<b>Kit Components and Storage</b> .....	2
Additional Materials Required. ....	2
<b>Chromatin IP DNA Purification Kit Experimental Design</b> .....	3
<b>Protocols</b>	
A. Purification of DNA from Chromatin IP Samples .....	4
<b>Appendix</b>	
Section A. Supplementary Chromatin IP Data .....	5
Section B. DNA Recovery .....	6
Section C. Troubleshooting Guide .....	8
Section D. Frequently Asked Questions (FAQ) .....	8
Section E. Related Products .....	9
<b>Technical Services</b> .....	12

## Overview

---

Chromatin Immunoprecipitation (ChIP) is a powerful, well established technique for studying interactions between chromatin-associated proteins and specific regions of the genome. The use of ChIP in combination with genome-wide analysis techniques can yield a tremendous amount of information regarding the distribution of transcription factors and histone modifications. But, these techniques require DNA that has been purified away from the components and contaminants present in an eluted ChIP sample. Active Motif's Chromatin IP DNA Purification Kit enables you to quickly clean up your ChIP DNA samples and get them ready for analysis, without the need for messy, labor intensive and time consuming phenol/chloroform extraction.

The Chromatin IP DNA Purification Kit is designed to be used to purify DNA following chromatin immunoprecipitation. It contains all necessary reagents to purify DNA from 50 ChIP reactions and can be used with material derived from any of the Active Motif line of ChIP-IT® Kits, including our ChIP-IT® Protein G Magnetic Beads (Catalog No. 53014).

<b>product</b>	<b>format</b>	<b>catalog no.</b>
Chromatin IP DNA Purification Kit	50 rxns	58002

## Kit Components and Storage

---

Please store each component at the temperature indicated in the table below.

Reagents	Quantity	Storage / Stability
DNA Purification Binding Buffer	50 ml	Room temperature
DNA Purification Wash Buffer*	50 ml	Room temperature
DNA Purification Elution Buffer	5 ml	Room temperature
3M Sodium Acetate**	500 µl	Room temperature
DNA purification columns	50	Room temperature
DNA column collection tubes	50	Room temperature

Columns and buffers (with the exception of the 3M Sodium Acetate) can be stored at 4°C, but all components must be warmed to room temperature before use. Some of the salts present in the buffers may precipitate out of solution during storage and will need to be resuspended prior to use.

\* DNA Purification Wash Buffer must be reconstituted to a final concentration of 80% ethanol prior to use.

\*\* **Do not** store the 3M Sodium Acetate at 4°C.

### Additional materials required

- Ethanol, 100%
- Microcentrifuge
- 1.5 or 2 ml microcentrifuge tubes

## Chromatin IP DNA Purification Kit Experimental Design

---

### PLEASE READ THE ENTIRE PROTOCOL BEFORE STARTING!

Prior to using the Chromatin IP DNA Purification Kit for the first time, the DNA Purification Wash Buffer must be reconstituted to a final concentration of 80% ethanol using 100% ethanol supplied by the user. This only needs to be performed once, prior to first using the kit. After reconstitution, store the buffer at room temperature. See step A1 in on page 4.

The Chromatin IP DNA Purification Kit is designed for use with a microcentrifuge for sample processing but can also be used with a vacuum manifold. Please consult the manufacturer's instructions for adaptation of the kit for vacuum manifold use.

**Compatibility:** Once your ChIP experiments are complete, the DNA purification procedure can be started immediately. The entire procedure should take no more than five to ten minutes, depending upon the number of samples to purify. The Chromatin IP DNA Purification Kit is compatible with samples from all Active Motif ChIP-IT Kits, or from any standard chromatin IP kit or procedure, using either mechanical or enzymatic shearing of chromatin, agarose or paramagnetic beads. Before purification, DNA samples should be eluted from the beads, cross-link reversed, and, if the ChIP procedure calls for it, Proteinase K and/or RNase digested.

**Binding of DNA to Column:** One critical aspect of the procedure is the pH of the sample when it is added to the column. Binding of DNA to the silica matrix of the DNA purification column will only occur at or below pH 7.5. A pH indicator dye is present in the DNA Purification Binding Buffer so that you can easily monitor that your samples are at the correct pH. To ensure successful purification of your ChIP DNA, sodium acetate (pH 5.4) is supplied in case the pH of your sample needs to be adjusted lower.

**Washing:** DNA bound to the column is extensively washed after binding.

**Elution:** After washing, the DNA is eluted using the included DNA Purification Elution Buffer. Elution can be carried out using highly pure water (pH 7) if desired.

**Purity:** Use of the Chromatin IP DNA Purification Kit results in recovery of purified DNA of high quality suitable for a number of downstream analysis techniques, including PCR (endpoint or quantitative), Southern blotting, microarray, standard or Next-Gen sequencing or cloning.

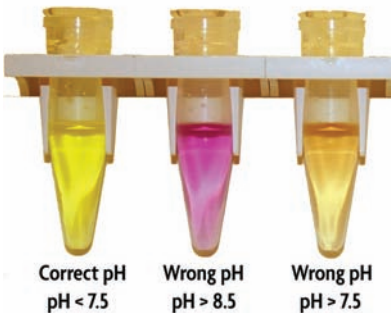
**Recovery:** DNA fragments below 50 base pairs in length are not recovered efficiently.

**Yield:** Recovery of DNA will be 85-100%. Depending upon the number of cell equivalents of chromatin used in each ChIP reaction, this typically will be from 100 ng to 1 µg. DNA can be successfully recovered from ChIP experiments starting with as few as 10,000 cells.

## Protocols

### A. Purification of DNA from Chromatin IP Samples

1. **ADD ETHANOL BEFORE THE FIRST USE ONLY!** The final concentration of ethanol in the DNA Purification Wash Buffer must be 80%. Add 40 ml of fresh 100% ethanol to the DNA Purification Wash Buffer bottle. Invert repeatedly. The buffer can be stored at room temperature after the addition of ethanol.
2. Your eluted ChIP samples should be in an Eppendorf tube. To each sample, add 5 volumes DNA Purification Binding Buffer for every volume of DNA sample. For example, if your ChIP sample is 100  $\mu$ l, add 500  $\mu$ l Binding Buffer.
3. Add 5  $\mu$ l 3M Sodium Acetate and mix. Check that the color of the DNA sample / DNA Purification Binding Buffer mixture is bright yellow, not light orange or violet. If the color is light orange or violet, this indicates that its pH is too high. Add more 3M Sodium Acetate 5  $\mu$ l at a time, mixing after addition, until the color is bright yellow (see Figure 1). As the elution buffers used in many ChIP protocols (including those in the ChIP-IT Kits) have a basic pH, **this step is crucial to the successful binding and purification of your ChIP DNA sample.**



**Figure 1: Solution color as a function of pH.**

The DNA Purification Binding Buffer has a pH indicator dye, so that the pH of the solution can be easily determined.

Only apply the sample to the column when the solution is bright yellow (left), indicating a pH under 7.5. DNA will not bind to the column if the pH is higher than 7.5.

**Note:** A full-color PDF of this manual can be downloaded from the Active Motif website, which will enable you to see this figure in color.

4. For each sample, place a DNA purification column in the collection tube and add each sample to its own column. Close the cap on each column, place them with the collection tubes in a microcentrifuge and spin them at 14,000 rpm for 1 minute. Alternatively, a compatible vacuum manifold may be used for sample processing.
5. Remove the column from the collection tube, then remove and discard the flow through from the collection tube. Return the column to the collection tube.
6. Add 750  $\mu$ l of Wash Buffer to the column and cap the column.
7. Spin at 14,000 rpm for 1 minute in a microcentrifuge.
8. Remove the column from the collection tube, then remove and discard the flow through from the collection tube. Return the column to the collection tube.
9. With the column cap open, spin at 14,000 rpm for 2 minutes in a microcentrifuge to remove any residual Wash Buffer from the column.

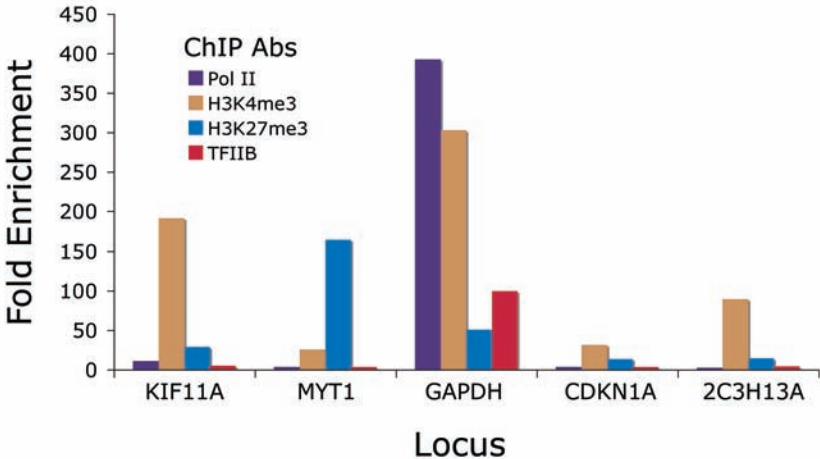
10. Transfer the column to a clean Eppendorf tube.
11. Add 50  $\mu$ l of DNA Purification Elution Buffer (or purified water with a pH equal to or greater than 7) to the center of the column matrix and wait for 1 minute.
 

**Note:** The amount of Elution Buffer used can be reduced to less than 50  $\mu$ l in order to attain a more concentrated DNA sample, if required. Reducing the elution volume will result in an overall reduction in the recovery efficiency of the purification. For more information, please see Appendix, Section B, Figure 5.
12. Spin at 14,000 rpm for 1 minute in a microcentrifuge.
13. Discard column. The DNA eluted into the Eppendorf tube is purified and ready to use in your application.

## Appendix

### Section A. Supplementary Chromatin IP Data

The Chromatin IP DNA Purification Kit is designed for purification of DNA after performing a ChIP experiment. Below are examples of quantitative PCR (qPCR) data from DNA purified using the Chromatin IP DNA Purification Kit after ChIP.



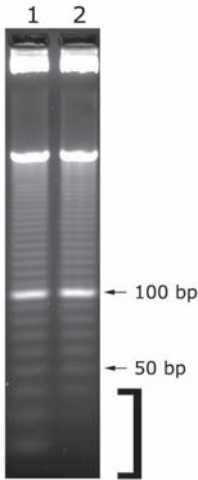
**Figure 2: Quantitative PCR performed on ChIP DNA purified with the Chromatin IP DNA Purification Kit.**

Quantitative PCR was performed using DNA purified with the Chromatin IP DNA Purification Kit after ChIP using the indicated antibodies (ChIP Abs). Chromatin IP experiments were performed using the ChIP-IT<sup>®</sup> Express Kit (Catalog No. 53008) and Ready-to-ChIP HeLa Chromatin (Catalog No. 53015,  $7.5 \times 10^5$  cell equivalents per ChIP). Quantitative PCR was carried out using primers specific for the indicated gene (Locus) and normalized data represent fold enrichment over ChIP experiments carried out with control IgG.



## Section B. DNA Recovery

The Chromatin IP DNA Purification Kit enables recovery of DNA fragments as small as 50 base pairs (Figure 3). As nucleosomal DNA is on the order of 150 base pairs, all of your chromatin IP DNA should be recovered efficiently.

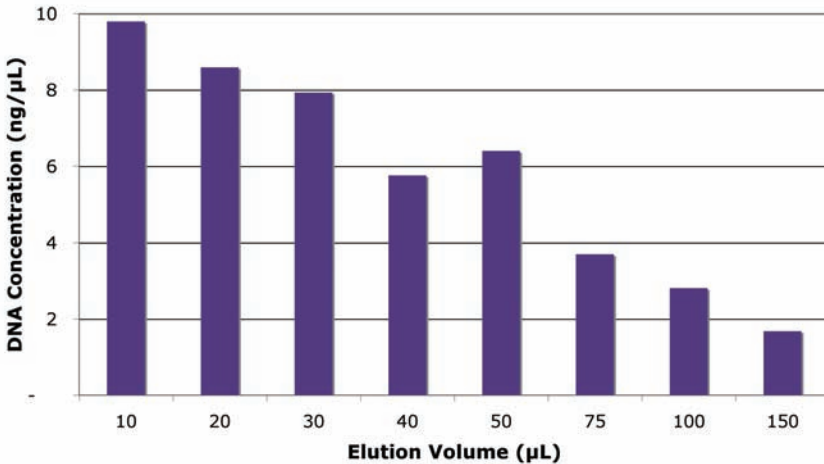


**Figure 3: DNA recovery as a function of fragment size.**

The Chromatin IP DNA Purification Kit enables efficient recovery of DNA fragments as small as 50 base pairs. DNA molecular weight markers (10 base pair ladder) were purified with the Chromatin IP DNA Purification Kit and run on a 5% agarose gel and stained with ethidium bromide to visualize the recovered DNA fragments.

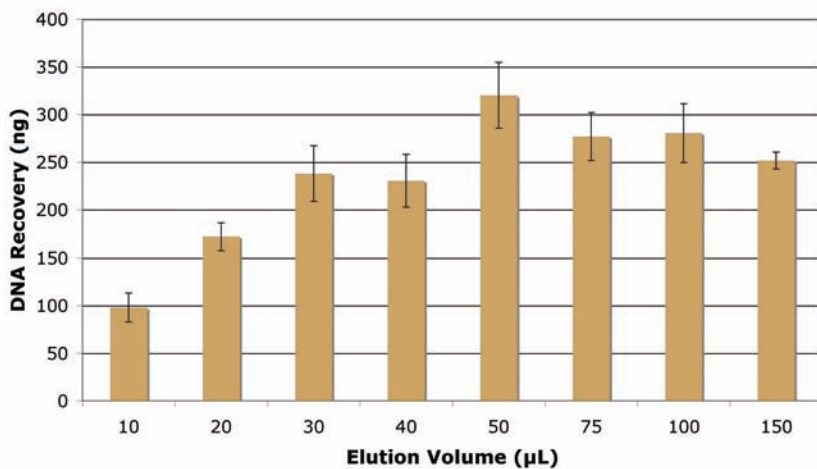
Lane 1: DNA not purified.  
Lane 2: DNA after purification.

As the figure illustrates, DNA larger than 50 base pairs is represented equally in the purified and non-purified DNA samples. DNA fragments smaller than 50 base pairs (bracket) are not efficiently purified, and thus are not represented in the purified DNA sample (Lane 2, bracket).



**Figure 4: Concentration of DNA recovered as a function of elution volume.**

Human genomic DNA was purified using the Chromatin IP DNA Purification Kit and eluted using the indicated volume of Elution Buffer and the concentration (ng/μl) of the recovered DNA was determined.



**Figure 5: Yield of DNA recovered as a function of elution volume.**

Human genomic DNA was purified using the Chromatin IP DNA Purification Kit and eluted using the indicated volume of Elution Buffer and the quantity (ng) of the recovered DNA was determined.

## Section C. Troubleshooting Guide

Problem/question	Recommendation
No amplification by PCR of ChIP DNA	<p>If you are experiencing problems with downstream analysis of ChIP DNA, it is possible that the purification process was compromised and your DNA samples were not recovered. This could be due to one of several factors.</p> <ol style="list-style-type: none"><li><b>The pH of the Binding Buffer was too high.</b> Refer to Figure 1 on page 4 to make sure that the pH of your sample is below 7.5 to ensure proper binding of ChIP DNA to the column. Your sample should be bright yellow (not orange, peach or purple). Add 3M Sodium Acetate in aliquots of 5 <math>\mu</math>l until the pH reaches the proper point and the sample is bright yellow. (A full-color PDF of this manual can be downloaded from the Active Motif website, which will enable you to see what is meant by "bright yellow".)</li><li><b>No ethanol in the DNA Purification Wash Buffer.</b> Make sure that ethanol has been added to the DNA Purification Wash Buffer prior to use. See step A1 on page 4.</li><li><b>Residual DNA Purification Wash Buffer in the column prior to elution.</b> Make certain to spin the column one extra time after washing to remove any residual Wash Buffer, as it will interfere with elution of the DNA from the column. Open the cap of the column to enable proper removal of residual Wash Buffer.</li><li><b>Salt in the Elution Buffer.</b> DNA will not be eluted from the silica matrix of the DNA purification column if the buffer used for elution contains high salt. Use the included DNA Purification Elution Buffer or purified water (pH 7) for elution.</li></ol>

## Section D. Frequently Asked Questions (FAQ)

Can the Chromatin IP DNA Purification Kit be...

used after 5-hydroxymethylcytidine MeDIP? Yes.

used after 5-methylcytidine MeDIP? Yes.

**Note:** The MeDIP procedure using the 5-Methylcytidine antibody (Catalog No. 39649) requires the DNA to be made single stranded. Single-stranded DNA does not bind as efficiently to the silica matrix of the column so recovery may not be 100%, but our kit has been shown to work as well as the leading competitor's kit.

used with other Active Motif kits, such as the hMeDIP Kit and the MethylCollector™ family? Yes.

used to purify DNA samples from other manufacturer's ChIP kits? Yes.

used to purify DNA samples from non-kit based ChIP protocols? Yes.

used with a vacuum manifold? Yes.

used to purify RNA? Not recommended. The buffers are not guaranteed to be RNase-free.

used to purify PCR products? Yes.

used to clean up restriction enzyme reactions? Yes.

## Section E. Related Products

Chromatin Immunoprecipitation	Format	Catalog No.
ChIP-IT® Express	25 rxns	53008
ChIP-IT® Express Enzymatic	25 rxns	53009
ChIP-IT® Express Shearing Kit	10 rxns	53032
ChIP-IT® Express Enzymatic Shearing Kit	10 rxns	53035
ChIP-IT® High Sensitivity	16 rxns	53040
ChIP-IT® qPCR Analysis Kit	10 rxns	53029
ChIP-IT® ChIP-Seq	10 libraries	53041
ChIP-IT® FFPE	16 rxns	53045
ChIP-IT® FFPE Chromatin Preparation Kit	5 rxns	53030
ChIP-IT® Express HT	96 rxns	53018
Re-ChIP-IT®	25 rxns	53016
RNA ChIP-IT®	25 rxns	53024
Chromatin IP DNA Purification Kit	50 rxns	58002
EpiShear™ Probe Sonicator	110 V	53051
EpiShear™ Cooled Sonication Platform, 1.5 ml	1 platform	53080
ChIP-IT® Protein G Magnetic Beads	25 rxns	53014
Protein G Agarose Columns	30 rxns	53039
Siliconized Tubes, 1.7 ml	25 tubes	53036
ChIP-IT® Control qPCR Kit – Human	5 rxns	53026
ChIP-IT® Control qPCR Kit – Mouse	5 rxns	53027
ChIP-IT® Control qPCR Kit – Rat	5 rxns	53028
ChIP-IT® Control Kit – Human	5 rxns	53010
ChIP-IT® Control Kit – Mouse	5 rxns	53011
ChIP-IT® Control Kit – Rat	5 rxns	53012
Ready-to-ChIP HeLa Chromatin	10 rxns	53015
Ready-to-ChIP Hep G2 Chromatin	10 rxns	53019
Ready-to-ChIP K-562 Chromatin	10 rxns	53020
Ready-to-ChIP NIH/3T3 Chromatin	10 rxns	53021
Bridging Antibody for Mouse IgG	500 µg	53017
Dounce Homogenizer	1 ml	40401
Dounce Homogenizer	15 ml	40415

### ChIP-validated Antibodies

For an up-to-date list of over 135 ChIP-validated antibodies, please visit [www.activemotif.com/chipabs](http://www.activemotif.com/chipabs).

Whole Genome Amplification	Format	Catalog No.
GenoMatrix™ Whole Genome Amplification Kit	1 kit	58001

Modified Histones Array	Format	Catalog No.
MODified™ Histone Peptide Array	1 array	13001

### Histone Antibodies

For an up-to-date list of over 140 antibodies against histones and modified histones, please visit [www.activemotif.com/histoneabs](http://www.activemotif.com/histoneabs).

<b>Histone ELISAs</b>	<b>Format</b>	<b>Catalog No.</b>
Total Histone H3 ELISA	1 x 96 rxns	53110
Histone H3 monomethyl Lys4 ELISA	1 x 96 rxns	53101
Histone H3 dimethyl Lys4 ELISA	1 x 96 rxns	53112
Histone H3 trimethyl Lys4 ELISA	1 x 96 rxns	53113
Histone H3 acetyl Lys9 ELISA	1 x 96 rxns	53114
Histone H3 dimethyl Lys9 ELISA	1 x 96 rxns	53108
Histone H3 trimethyl Lys9 ELISA	1 x 96 rxns	53109
Histone H3 phospho Ser10 ELISA	1 x 96 rxns	53111
Histone H3 acetyl Lys14 ELISA	1 x 96 rxns	53115
Histone H3 monomethyl Lys27 ELISA	1 x 96 rxns	53104
Histone H3 trimethyl Lys27 ELISA	1 x 96 rxns	53106
Histone H3 phospho Ser28 ELISA	1 x 96 rxns	53100

<b>Histone Purification &amp; Chromatin Assembly</b>	<b>Format</b>	<b>Catalog No.</b>
Histone Purification Kit	10 rxns	40025
Histone Purification Mini Kit	10 rxns	40026
Chromatin Assembly Kit	10 rxns	53500
HeLa Core Histones	36 µg	53501

### **Recombinant Methylated, Acetylated and Phosphorylated Histone Proteins**

For an up-to-date list of Recombinant Histone Proteins, please visit [www.activemotif.com/recombhis](http://www.activemotif.com/recombhis).

<b>Histone Acetyltransferase and Deacetylase Activity</b>	<b>Format</b>	<b>Catalog No.</b>
HAT Assay Kit (Fluorescent)	1 x 96 rxns	56100
Recombinant p300 protein, catalytic domain	5 µg	31205
Recombinant GCN5 protein, active	5 µg	31204
HDAC Assay Kit (Fluorescent)	1 x 96 rxns	56200
HDAC Assay Kit (Colorimetric)	1 x 96 rxns	56210

<b>Histone Demethylase Activity</b>	<b>Format</b>	<b>Catalog No.</b>
Histone Demethylase Assay (Fluorescent)	48 rxns	53200

<b>DNA Methylation</b>	<b>Format</b>	<b>Catalog No.</b>
hMeDIP	10 rxns	55010
MeDIP	10 rxns	55009
MethylDetector™	50 rxns	55001
MethylCollector™ Ultra	30 rxns	55005
UnMethylCollector™	30 rxns	55004
DNMT Activity / Inhibition Assay	96 rxns	55006
Methylated DNA Standard Kit	3 x 2.5 µg	55008
Fully Methylated Jurkat DNA	10 µg	55003
Jurkat genomic DNA	10 µg	55007

### **DNA Methylation Antibodies**

For an up-to-date list of over 25 DNA Methylation antibodies, visit [www.activemotif.com/dnamethabs](http://www.activemotif.com/dnamethabs).

<b>Co-Immunoprecipitation</b>	<b>Format</b>	<b>Catalog No.</b>
Nuclear Complex Co-IP Kit	50 rxns	54001
Universal Magnetic Co-IP Kit	25 rxns	54002

<b>SUMOylation</b>	<b>Format</b>	<b>Catalog No.</b>
SUMOlink™ SUMO-1 Kit	20 rxns	40120
SUMOlink™ SUMO-2/3 Kit	20 rxns	40220

<b>In-cell Phospho-specific ELISAs</b>	<b>Format</b>	<b>Catalog No.</b>	<b>Catalog No.</b>
FACE™ AKT	1 x 96 rxns	48120	48220
FACE™ ATF-2	1 x 96 rxns	48115	48215
FACE™ Bad	1 x 96 rxns	48165	48265
FACE™ c-Jun (S63)	1 x 96 rxns	48125	48225
FACE™ c-Jun (S73)	1 x 96 rxns	48135	48235
FACE™ c-Src	1 x 96 rxns	48155	48255
FACE™ EGFR (Y845)	1 x 96 rxns	48340	48440
FACE™ EGFR (Y992)	1 x 96 rxns	48150	48250
FACE™ EGFR (Y1173)	1 x 96 rxns	48190	48290
FACE™ ErbB-2 (Y877)	1 x 96 rxns	48130	48230
FACE™ ErbB-2 (Y1248)	1 x 96 rxns	48105	48205
FACE™ ERK1/2	1 x 96 rxns	48140	48240
FACE™ FAK	1 x 96 rxns	48145	48245
FACE™ FKHR (FOXO1)	1 x 96 rxns	48160	48260
FACE™ HSP27	1 x 96 rxns	48350	48450
FACE™ JAK1	1 x 96 rxns	48185	48285
FACE™ JNK	1 x 96 rxns	48110	48210
FACE™ MEK1/2	1 x 96 rxns	48180	48280
FACE™ NFκB p65 Profiler	3 x 96 rxns	48300	48400
FACE™ p38	1 x 96 rxns	48100	48200
FACE™ PI3 Kinase p85	1 x 96 rxns	48175	48275
FACE™ STAT2	1 x 96 rxns	48310	48410
FACE™ STAT4	1 x 96 rxns	48320	48420
FACE™ STAT6	1 x 96 rxns	48330	48430
FACE™ Maker	1 x 96 rxns	48000	48050
Suspension Cell FACE™	2 x 96 rxns	48305	48405

For a complete, up-to-date list of available FACE™ Kits, please visit [www.activemotif.com/face](http://www.activemotif.com/face)

## 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

Toll Free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

E-mail: [tech\\_service@activemotif.com](mailto:tech_service@activemotif.com)

### Active Motif Europe

Avenue Reine Astrid, 92

B-1310 La Hulpe, 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

E-mail: [eurotech@activemotif.com](mailto:eurotech@activemotif.com)

### 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

E-mail: [japantech@activemotif.com](mailto:japantech@activemotif.com)

Visit Active Motif on the worldwide web 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
- Enter your name into our mailing list to receive our catalog, *MotifVations* newsletter and notification of our upcoming products
- Share your ideas and results with us
- View our job opportunities

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