

# FunctionELISA™

IκBα

For the detection and analysis of IκBα phosphorylation

(version B1)

Catalog Nos. 48005 & 48505

## 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

104 Avenue Franklin Roosevelt

B-1330 Rixensart, 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

**Background**

Overview .....	1
Advantages .....	2
Example Experiment .....	2

**Kit Components and Storage**

Additional Materials Required .....	3
-------------------------------------	---

**Preparation of Reagents** ..... 4**Preparation of Whole-Cell Extract** ..... 4**I $\kappa$ B $\alpha$  Assay**

Flow Chart of Process .....	5
-----------------------------	---

## Protocol

Step 1: Preparation of Phospho-I $\kappa$ B $\alpha$ Standard Curve .....	6
Step 2: Binding of I $\kappa$ B $\alpha$ .....	8
Step 3: Binding of I $\kappa$ B $\alpha$ -Detecting Antibody .....	8
Step 4: Binding of Secondary Antibody .....	8
Step 5: Chemiluminescent Detection .....	8

**References** ..... 9**Appendix**

Section A. Troubleshooting .....	10
Section B. Related Products .....	11

**Technical Services** ..... 12

## Background

The transcription factor NF $\kappa$ B is implicated in the regulation of many genes that code for mediators of the immune, acute phase and inflammatory responses<sup>1</sup>. NF $\kappa$ B is composed of homo- and heterodimeric complexes of members of the Rel (NF $\kappa$ B) family. There are five subunits of the NF $\kappa$ B family in mammals: p50, p65 (RelA), c-Rel, p52 and RelB<sup>2</sup>. These proteins share a conserved 300 amino acid sequence in the N-terminal region, known as the Rel homology domain, that mediates DNA binding, protein dimerization and nuclear localization. This domain is also a target of the I $\kappa$ B inhibitors, which include I $\kappa$ B $\alpha$ , I $\kappa$ B $\beta$ , I $\kappa$ B $\gamma$ , Bcl-3, p105 and p100<sup>3</sup>. In the majority of cells, NF $\kappa$ B exists in an inactive form in the cytoplasm, bound to the inhibitory I $\kappa$ B proteins<sup>4</sup>. I $\kappa$ B proteins are phosphorylated by the I $\kappa$ B kinase complex, which consists of at least three proteins: IKK1/ $\alpha$ , IKK2/ $\beta$  and IKK3/ $\gamma$ <sup>4-7</sup>. External stimuli, such as tumor necrosis factor (TNF) or other cytokines, initiate a signal transduction cascade that leads to the activation of the I $\kappa$ B kinase complex. This complex specifically phosphorylates I $\kappa$ B $\alpha$  on Serine-32 and Serine-36, which leads to ubiquitination of I $\kappa$ B $\alpha$  and its subsequent degradation by the 26S proteasome. Degradation of I $\kappa$ B $\alpha$  results in unmasking of the nuclear localization signal of NF $\kappa$ B dimers, which subsequently translocate to the nucleus and activate target genes<sup>8,9</sup>. Thus, analysis of the phosphorylation state of I $\kappa$ B $\alpha$  provides a correlation to the activity of the IKK complex as well as the activation and nuclear localization of NF $\kappa$ B.

## Overview

FunctionELISA™ Kits provide a useful tool to capture and quantifiably measure the amount of a specific protein in a sample. They utilize the “Sandwich ELISA” method for detecting a protein in a sample, which uses two antibodies that each recognize a distinct epitope on the protein. An ELISA plate is coated with the first antibody, called the Capture Antibody, which is used to capture the protein from the sample. The second antibody, called the Detecting Antibody, is used

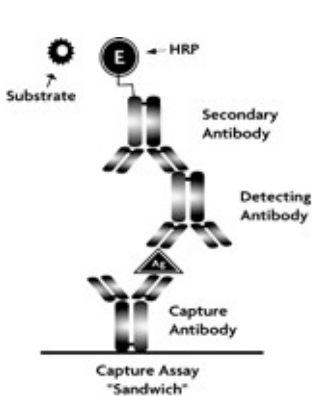


Figure 1: Sandwich ELISA schematic.

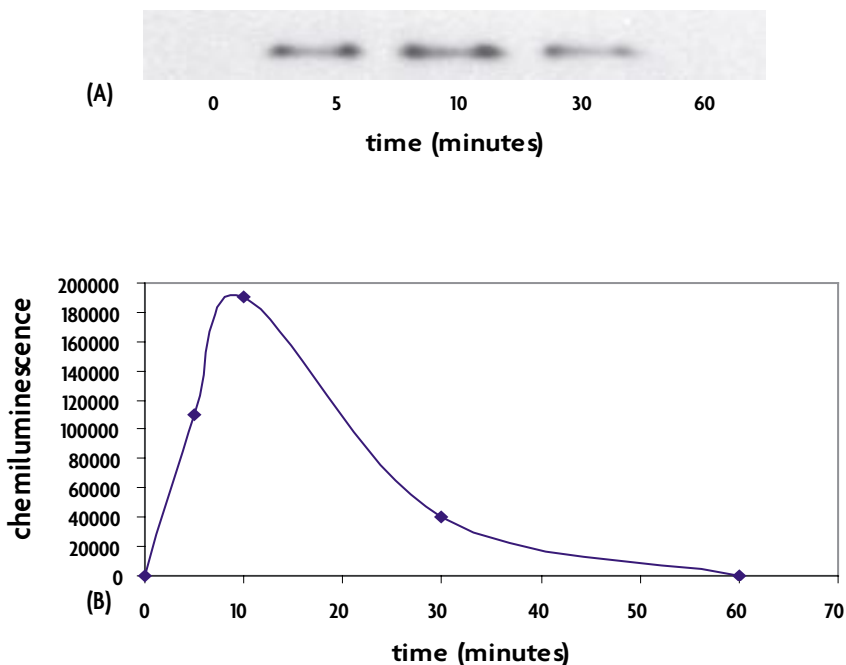
to detect the protein bound by the Capture Antibody. An HRP-conjugated Secondary Antibody is then used to quantitate the amount of bound Detecting Antibody (Figure 1). By using Capture Antibody that recognizes a phosphorylation site on a protein, one can specifically detect phosphorylated protein. Because many proteins involved in signaling cascades are regulated by phosphorylation, this assay proves useful in monitoring and analyzing the phosphorylation state of a given protein. Compared to Western blots, FunctionELISA provides a quicker, more quantifiable way to monitor a protein's phosphorylation state.

## Advantages

- Multiple samples can be analyzed in low-volume, high-throughput experiments
- Analysis is complete in just hours
- Quantitative nature of the assay enables direct measurement of changes in phosphorylation

## Example Experiment

To monitor the phosphorylation state of I $\kappa$ B $\alpha$ , Jurkat cells were grown to  $2 \times 10^6$  cells/ml and treated with 1 nM TNF- $\alpha$ . The cells were harvested at various time points, lysed and Western blot and the FunctionELISA I $\kappa$ B $\alpha$  Kit were used to monitor and quantify I $\kappa$ B $\alpha$  phosphorylation using 100  $\mu$ g of lysate per well. TNF- $\alpha$ -induced phosphorylation of I $\kappa$ B $\alpha$  was very rapid, peaking at around 10 minutes and disappearing by 30 minutes (see Figure 2).



**Figure 2: Induction of I $\kappa$ B $\alpha$  phosphorylation with TNF- $\alpha$  in Jurkat cells.** Jurkat cells, grown to a density of  $2 \times 10^6$  cells/ml, were treated with 1 nM TNF- $\alpha$ . The cells were harvested at various indicated time points, and cell lysates were tested in Western analysis (Fig. 2A) using an anti-Phospho-I $\kappa$ B $\alpha$  antibody (Active Motif Cat. No. 40904) and in a FunctionELISA I $\kappa$ B $\alpha$  assay (Fig. 2B).

## Kit Components and Storage

Reagents	Quantity 1 plate / 5 plates	Composition	Storage
Detecting Antibody (200 µg/ml)	24 µl / 120 µl	PBS, 0.1% sodium azide 0.2% gelatin	4°C for 1 year
HRP-conjugated Secondary Antibody	12 µl / 60 µl	PBS, 0.1% sodium azide	4°C for 1 year
Lysis Buffer	10 ml / 50 ml	20 mM Hepes, pH 7.5 400 mM NaCl 20% glycerol 1% Triton X-100 0.5% NP-40 0.1 mM EDTA 10 mM NaF 10 µM Na <sub>2</sub> MoO <sub>4</sub> 1 mM NaVO <sub>3</sub> 10 mM PNPP 10 mM β-glycerophosphate	4°C for 6 months
100X Protease Inhibitor	100 µl / 500 µl	Proprietary Mix	-20°C for 1 year
1M DTT	100 µl / 500 µl	solution in H <sub>2</sub> O	-20°C for 1 year
BSA powder	0.3 g / 1.5 g		4°C for 1 year
20X Wash Buffer	25 ml / 125 ml	20X PBS, pH 7.4 1% Tween-20 4 mM Sodium Orthovanadate	4°C for 6 months
HRP Substrate Solution A	2.5 ml / 12.5 ml		4°C for 6 months
HRP Substrate Solution B	2.5 ml / 12.5 ml		4°C for 6 months
FunctionELISA 1κBα plate	1 plate / 5 plates	12 x 8-microwell strips w/ holder	4°C for 1 year
Plate sealer	1 roll / 5 rolls		
Phospho-1κBα Control (20 ng/ml)	1 ml / 5 ml	Blocking Buffer	-20°C for 6 months

**Note:** The included buffers and reagents have been optimized for use with this kit. Substitution with other reagents may not give optimal results.

### Additional Materials Required

- Multi-channel pipettor
- Multi-channel pipettor reservoirs
- Orbital shaker
- Microplate luminometer
- Distilled water (dH<sub>2</sub>O)
- Phosphate buffered saline (PBS, pH 7.4)

## Preparation of Reagents

---

### Complete Lysis Buffer

An excess of Lysis Buffer is provided for preparing cell extracts for the assay. Prepare the amount of Complete Lysis Buffer required for the assay by adding 1  $\mu$ l DTT and 10  $\mu$ l Protease Inhibitor Cocktail per ml of Lysis Buffer.

**Note:** Some protease inhibitors lose activity within 24 hours of dilution. Therefore, we recommend using the Complete Lysis Buffer immediately for cell lysis. The remaining amount should be discarded if not used on the same day.

### Wash Buffer

1X Wash Buffer may be stored at 4°C. If you do not plan on using the entire ELISA plate at once, we recommend preparing only enough Wash Buffer to meet the needs of the experiment.

- Dilute 20X Wash Buffer to 1X with dH<sub>2</sub>O. The following chart can be used as a guide.

1X Wash Buffer	20X Wash Buffer	dH <sub>2</sub> O	Sufficient for
25 ml	1.25 ml	23.75 ml	5 rxns
50 ml	2.5 ml	47.5 ml	10 rxns
100 ml	5 ml	95 ml	20 rxns
250 ml	12.5 ml	237.5 ml	48 rxns
500 ml	25 ml	475 ml	96 rxns

### Blocking Buffer

Prepare Blocking Buffer by dissolving 0.3 g BSA in 30 ml 1X Wash Buffer in a sterile bottle.

**Note:** Sodium azide has not been included as it inhibits Horseradish Peroxidase. To prevent microbial growth, store Blocking Buffer at 4°C and reduce exposure to contaminants. Dissolve the BSA in 1X Wash Buffer only when ready to use.

## Preparation of Whole-Cell Extract

---

This procedure can be used for a confluent cell layer of 25 cm<sup>2</sup> (60 mm dish). The yield is approximately 0.3 mg of total cellular protein.

1. Treat the cells as required for activation.
2. Wash the cells 2X with 8 ml ice-cold PBS (10 mM phosphate buffer, pH 7.4, 150 mM NaCl).
3. Collect the cells in 3 ml ice-cold PBS. Use a cell scraper for adherent cells.
4. Centrifuge the cells for 10 min at 1,000 rpm at 4°C.

5. Discard the supernatant and add a volume of Complete Lysis Buffer equivalent to the volume of the cell pellet. Pipette gently to mix.
 

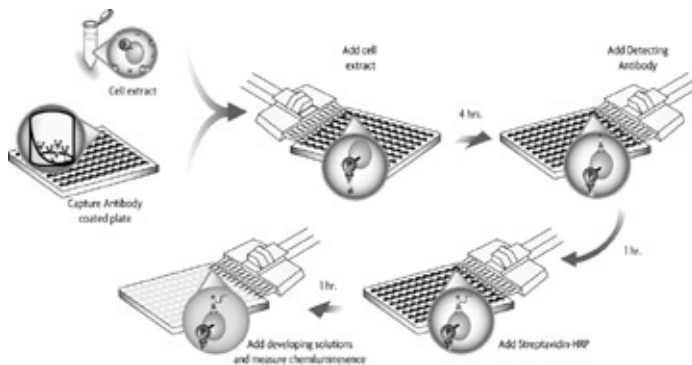
**Note:** Preparing such a concentrated lysate will aid in the capture of phosphorylated proteins, whose concentration is very low.
6. Transfer into a 1.5 ml microcentrifuge tube.
7. Incubate on ice for 30 minutes.
8. Centrifuge for 20 min at 14,000 x g (full speed in a microcentrifuge) at 4°C.
9. Collect the supernatant at 4°C, aliquot and store at -80°C. Avoid freeze/thaw cycles.
10. Measure the protein content by a Bradford-based assay.

## I $\kappa$ B $\alpha$ Assay

---

I $\kappa$ B $\alpha$  protein is phosphorylated in response to stimulation of the NF $\kappa$ B pathway, and then broken down very rapidly. Thus, capture and detection of this transient form of I $\kappa$ B $\alpha$  is very difficult, as is the long-term storage of extracts. Because of these limitations, this kit's positive control is a synthetic antigen generated by coupling a phospho-peptide to full-length I $\kappa$ B $\alpha$  protein. The conjugated phospho-peptide is the same one used to generate the monoclonal Capture Antibody. The Phospho-I $\kappa$ B $\alpha$  Control is stable and can be used to generate a standard curve.

Determine the appropriate number of microwell strips required for testing your samples, controls and blanks in duplicate. If less than 8 wells in a strip are to be used, cover the unused wells with a portion of the plate sealer before you perform the assay. The contents of these wells are stable at room temperature and, if kept dry, can be used later for a separate assay. Store the unused strips in the aluminum pouch at 4°C. Use the strip holder for the assay. Prepare the Washing and Blocking Buffers as described in the section **Preparation of Reagents**. Multi-channel pipettor reservoirs may be used for dispensing the various buffers and Substrate Solutions into the wells being used.



**Figure 3:** Flow Chart of the FunctionELISA™ I $\kappa$ B $\alpha$  Process.

## Protocol

**PLEASE READ THE ENTIRE PROTOCOL BEFORE STARTING!**

### Step 1: Preparation of Phospho-I $\kappa$ B $\alpha$ Standard Curve

1. Set up a Phospho-I $\kappa$ B $\alpha$  Standard Curve in duplicate using the following concentrations: 10, 5, 2.5, 1.25, 0.63, 0.32, 0.15 and 0.0 ng/ml. (See table below for suggested layout.)

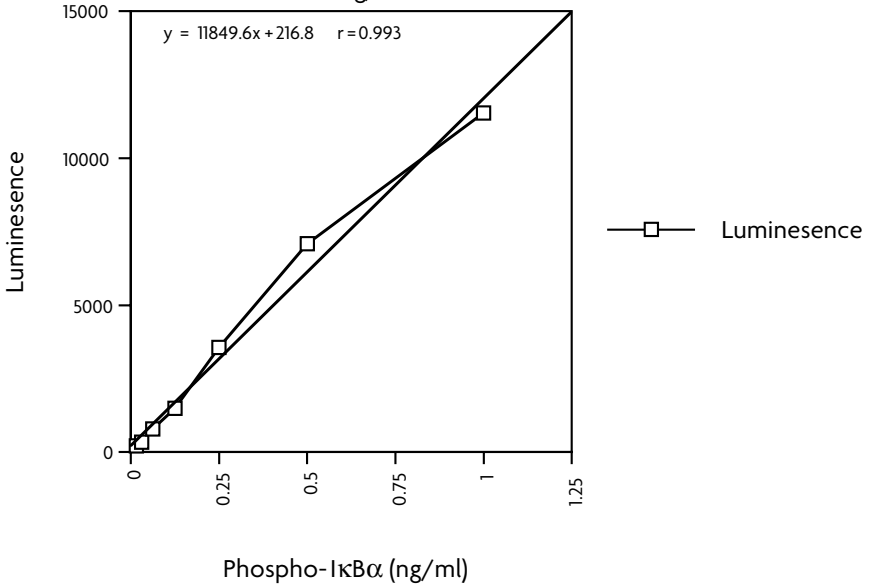
**Note:** For higher sensitivity, set up one or both Phospho-I $\kappa$ B $\alpha$  Standard Curve(s) using a 10-fold dilution of the 20 ng/ml Phospho-I $\kappa$ B $\alpha$  Control in Blocking Buffer to 2 ng/ml (final standard curve concentrations of 1, 0.5, 0.25, 0.125, 0.063, 0.032, 0.015 and 0.0 ng/ml).

2. Add 100  $\mu$ l of Blocking Buffer to wells A1 through H1 and A2 through H2.
3. Pipette 100  $\mu$ l Phospho-I $\kappa$ B $\alpha$  Control solution (20 ng/ml) into wells A1 and A2.
4. Mix wells A1 and A2 by pipetting.
5. Transfer 100  $\mu$ l from well A1 to B1 and A2 to B2.
6. Mix wells B1 and B2 by pipetting.
7. Transfer 100  $\mu$ l from well B1 to C1 and B2 to C2.
8. Continue this procedure to wells G1 and G2. After mixing, discard 100  $\mu$ l of solution from wells G1 and G2.
9. Wells H1 and H2 are blanks and should contain only 100  $\mu$ l of Blocking Buffer.

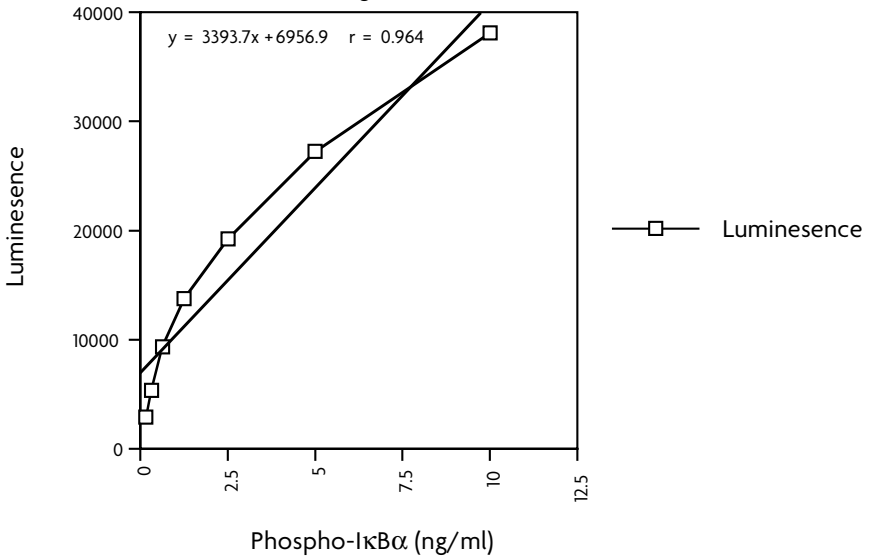
	1	2	3	4	5	6	7	8	9	10	11	12
A	10 (1) ng/ml	10 (1) ng/ml	-	-	-	-	-	-	-	-	-	-
B	5(0.5) ng/ml	5(0.5) ng/ml	-	-	-	-	-	-	-	-	-	-
C	2.5 (0.25) ng/ml	2.5 (0.25) ng/ml	-	-	-	-	-	-	-	-	-	-
D	1.25 (0.125) ng/ml	1.25 (0.125) ng/ml	-	-	-	-	-	-	-	-	-	-
E	0.63 (0.063) ng/ml	0.63 (0.063) ng/ml	-	-	-	-	-	-	-	-	-	-
F	0.32 (0.032) ng/ml	0.32 (0.032) ng/ml	-	-	-	-	-	-	-	-	-	-
G	0.15 (0.015) ng/ml	0.15 (0.015) ng/ml	-	-	-	-	-	-	-	-	-	-
H	Blank	Blank	-	-	-	-	-	-	-	-	-	-



Phospho-IκBα Standard Curve  
1 - 0.015 ng/ml



Phospho-IκBα Standard Curve  
10 - 0.15 ng/ml



### **Step 2: Binding of I $\kappa$ B $\alpha$**

1. **Sample wells:** Dilute concentrated cell lysates to 1 mg/ml in Blocking Buffer. Pipette 100  $\mu$ l lysate into each well.
2. Cover the plate with plate sealer and incubate at 4°C for 4 hours.
3. Remove cell lysate from wells and wash four times (4X) with 250  $\mu$ l Wash Buffer. After final wash, tap plate upside down several times to remove any residual Wash Buffer.

### **Step 3: Binding of I $\kappa$ B $\alpha$ -Detecting Antibody**

1. Dilute Detecting Antibody 1:250 in Blocking Buffer. For one entire plate, use 20  $\mu$ l Detecting Antibody in 5 ml Blocking Buffer.
2. Add 50  $\mu$ l diluted Detecting Antibody to each well.
3. Cover the plate and incubate at room temperature for 1 hour.
4. Remove Detecting Antibody and wash wells four times (4X) with 250  $\mu$ l Wash Buffer. After final wash, tap plate upside down several times to remove any residual Wash Buffer.

### **Step 4: Binding of Secondary Antibody**

1. Dilute Secondary Antibody 1:1000 in Blocking Buffer. For one entire plate, use 5  $\mu$ l Secondary Antibody in 5 ml Blocking Buffer.
2. Add 50  $\mu$ l diluted Secondary Antibody to each well.
3. Cover the plate and incubate at room temperature for 1 hour.
4. While Secondary Antibody is reacting, mix equal amounts of HRP Substrate Solution A and HRP Substrate Solution B. For one entire plate, combine 2.5 ml of each solution. Mix and allow temperature to equilibrate to room temperature.
5. Remove Secondary Antibody solution and wash thoroughly five times (5X) with 250  $\mu$ l Wash Buffer to remove any unbound conjugate. Let plate sit for 1 minute and then tap plate upside down several times to remove any residual Wash Buffer.

### **Step 5: Chemiluminescent Detection**

1. Add 50  $\mu$ l Substrate Solution mixture (See Step 4, No.4) per well.
2. Use a 96-well luminometer to measure each well and take readings for 2-second and for 20-second intervals.

## References

---

1. Bauerle P.A. and Baichwal V.R. (1997) *Adv. Immunol.* 65: 111-137.
2. Gilmore T.D. and Morin P.J. (1993) *Trends Genetics* 9: 427-433.
3. Bauerle P.A. and Baltimore D. (1988) *Cell* 53: 211-217.
4. DiDonato, J.A., Hayakawa, M., Rothwarf, D.M. Zandi, E., and Karin, M. (1997) *Nature* 388: 548.
5. Regnier, C.H., Ho, Y.S., Gao, X., Goeddel, D.V., Cao, Z., and Rothe, M. (1997) *Cell* 90: 373.
6. Mercurio, F., Zhu, H., Murray, B.W., Shevchenko, A., Bennett, B.L., Li, J., Young, D.B., Barbosa, M. Mann, M., Manning, A., and Rao, A. (1997) *Science* 278: 860.
7. Yamaoka, S., Courtois, G., Bessia, C., Whiteside, S.T., Weil, R., Agou, F., Kirk, H.E., Kay, R.J., and Israel, A. (1998) *Cell* 93: 1231-1240.
8. Ghosh, S., May, M.J., and Kopp, E.B. (1998) *Annu Rev. Immunol.* 16: 225-260.
9. Baldwin, A.S. (1996) *Annu. Rev. Immunol.* 14: 694-681.

## Appendix

### Section A. Troubleshooting

PROBLEM	POSSIBLE CAUSE	RECOMMENDATION
No signal or weak signal in all wells	Omission of key reagent	Check that all reagents have been added in the correct order
	Substrate or conjugate is no longer active	Test conjugate and substrate for activity
	Enzyme inhibitor present	Sodium azide will inhibit the peroxidase reaction, follow our recommendations when preparing buffers
	Plate reader settings not optimal	Verify the wavelength and filter settings of the plate reader
	Incorrect assay temperature	Bring substrate to room temperature
	Inadequate volume of Substrate Solution	Check to make sure that correct volume is delivered by pipette
High background in all wells	Concentration of antibodies too high	Increase antibody dilutions
	Inadequate washing	Ensure all wells are filled with Wash Buffer and follow washing recommendations
Uneven signal development	Incomplete washing of wells	Ensure all wells are filled with Wash Buffer and follow washing recommendations
	Well cross-contamination	Follow washing recommendations
High background in sample wells	Too much cell extract per well	Decrease amount of cell extract to 50 µg/well
	Concentration of antibodies too high	Perform antibody titration to determine optimal working concentration. Start using 1:1000 for Detecting Antibody and 1:2000 for the Secondary Antibody. The sensitivity of the assay will be decreased
No signal or weak signal in sample wells only	Not enough cell extract per well	Increase amount of cell extract to 150 µg/well
	IκBα is poorly activated or is inactivated	Perform a time course for IκBα activation in the studied cell line
	Cell extracts are not from rat, mouse or human origin	Perform study with a rat, mouse or human model

## Section B. Related Products

NFκB PRODUCTS	Application	Format	Catalog No.
<b>TransAM™</b>			
NFκB p65	Transcription Factor Assay	96-well plate	40096
NFκB p50	Transcription Factor Assay	96-well plate	41096
<b>Cell extracts</b>			
Unstimulated HeLa whole-cell extract		125 µg	40050
IL-1α-stimulated HeLa whole-cell extract		125 µg	40100
TNF-α-stimulated HeLa whole-cell extract		125 µg	40200
TNF-α-stimulated HeLa nuclear extract		125 µg	40210
Nuclear Extract Kit		100 rxns	40010
<b>Antibodies</b>			
A20 mAb	WB, IP	100 µg	40901
Phosphorylated-Akt1 mAb	WB, IP	100 µg	40902
IκBα mAb	WB, IP	100 µg	40903
Phosphorylated-IκBα mAb	WB, IP	100 µg	40904
IKKα mAb	WB	100 µg	40905
IKKβ mAb (Clone 10AG2)	WB	100 µg	40906
IKKβ mAb (Clone 10A9B6)	IP	100 µg	40907
IKKγ mAb	WB	100 µg	40908
IKKi/IKKε mAb	WB	100 µg	40950
IKKi/IKKε pAb	WB	100 µg	40957
TBK1 mAb	WB	100 µg	40909
IRAK-1 pAb	WB, IP	100 µg	40910
IRAK-2 pAb	WB	100 µg	40912
MyD88 pAb	WB	100 µg	40914
NFκB (p65) mAb	WB	100 µg	40916
RANK mAb	WB	100 µg	40917
TAB1 mAb	WB	100 µg	40918
TRAF2 mAb	WB	100 µg	40919
TRAF5 mAb	WB	100 µg	40920
TRAF6 mAb	WB	100 µg	40921
TLR2 pAb	WB	100 µg	40981
TLR3 mAb	WB	100 µg	40952
TLR6 mAb	WB, FC	100 µg	40954
TLR8 mAb	WB, FC	100 µg	40955
TLR9 mAb	WB, FC	100 µg	40956
TRAIL mAb	WB, FC, E	100 µg	40966

## 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

104 Avenue Franklin Roosevelt

B-1330 Rixensart, 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, *Motivations* 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!**