

FACE™

HSP27 ELISA Kits

(version A)

Catalog Nos. 48350 & 48850
Catalog Nos. 48450 & 48950 (Chemi)

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

TABLE OF CONTENTS	Page
Overview	1
Flow Chart of Process	2
Introduction	
HSP27	3
Traditional Kinase Assays	4
FACE HSP27	5
Kit Performance and Benefits	6
FACE HSP27 Experimental Design	7
Kit Components and Storage - Colorimetric Assay	8
Kit Components and Storage - Chemiluminescent Assay	9
Protocols - Colorimetric Assay	
Buffer Preparation and Recommendations	10
Quick Chart for Preparing Buffers	12
Adherent Cell Protocol	13
Non-Adherent Cell Protocol	15
Protocols - Chemiluminescent Assay	
Buffer Preparation and Recommendations	16
Quick Chart for Preparing Buffers	18
Adherent Cell Protocol	19
Non-Adherent Cell Protocol	21
References	21
Appendix	
Section A. Troubleshooting Guide	22
Section B. Related Products	24
Technical Services	25

Overview

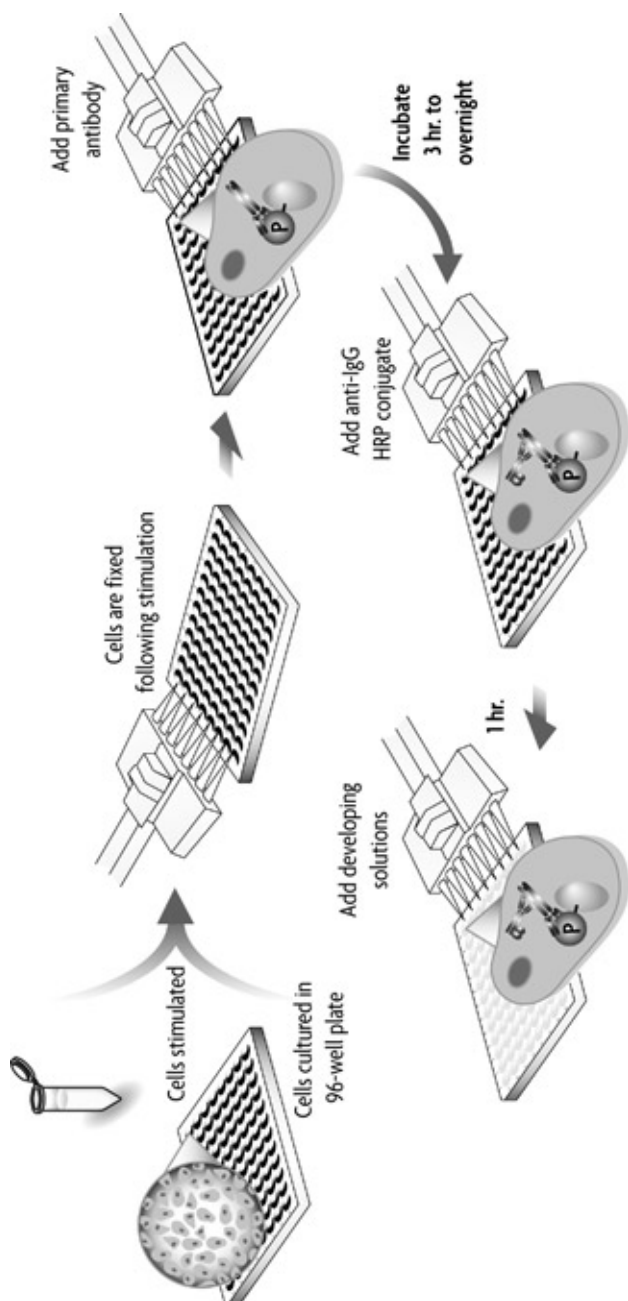
Heat Shock Protein 27 (HSP27) is a 27 kDa protein whose expression and function are stimulated by heat shock and other stress stimuli. It has been implicated in various biological processes including the response to heat shock, oxidative stress and cytokine treatment. Like other small heat shock proteins, HSP27 is regulated at both the transcriptional and post-translational levels¹. A major function of these proteins is to serve as chaperones that bind to and stabilize the active conformation of other proteins. HSP27, along with other members of the small HSP group, possesses a C-terminal α -crystalline homology domain². HSP27 was postulated to inhibit apoptosis through its activity as a molecular chaperone³.

Fast Activated Cell-based ELISA (FACE™)* Kits provide a simple, efficient, cell-based method to monitor proteins activated by phosphorylation.⁴ FACE HSP27 Kits are designed specifically to quantify activated (phosphorylated) HSP27 and/or total HSP27. In the FACE method, cells are cultured in 96-well plates and stimulated to induce the pathway of interest. Following stimulation, the cells are rapidly fixed to preserve activation-specific protein modifications. Each well is then incubated with a primary antibody that recognizes either phosphorylated HSP27 or total HSP27. Subsequent incubation with secondary HRP-conjugated antibody and developing solution provides an easily quantified colorimetric or chemiluminescent readout. The relative number of cells in each well is then determined using the provided Crystal Violet solution. The 96-well plate format is suitable for high-throughput screening applications. FACE HSP27 Kits are available in both colorimetric and chemiluminescent formats for studying the Ser 82 phosphorylation site of HSP27:

product	format	catalog no.
FACE HSP27 (S82)	1 x 96 rxns	48350
	5 x 96 rxns	48850
FACE HSP27 (S82) Chemi	1 x 96 rxns	48450
	5 x 96 rxns	45950

* Developed in collaboration with Dr. M. Peppelenbosch and Dr. H. Versteeg.

Flow Chart of Process



Introduction

HSP27

In response to stress, the expression level of HSP27 increases several fold to confer cellular resistance to the adverse environmental change. HSP27 is localized to the cytoplasm of unstressed cells but can redistribute to the nucleus in response to stress, where it may function to stabilize DNA and/or the nuclear membrane⁵. Cytoplasmic HSP27 exists in multiple complexes. One complex consists of HSP27, AKT (PKB), MAPKAPK-2 and p38 MAPK⁶. The presence of HSP27 in this complex is required for AKT activation by stress stimuli. Another complex consists of HSP27 and the IKK complex⁷. It is also an actin capping protein that binds to the barbed (growing) ends of actin filaments, thereby inhibiting filament extension^{8,9}. It has been demonstrated that phosphorylation and increased concentration of HSP27 modulates actin polymerization and reorganization^{9,10}.

HSP27 is phosphorylated at serines 15, 78 and 82 by MAPKAPK-2 as a result of the activation of the p38 MAPK pathway^{11,12}. AKT only phosphorylates HSP27 on Ser82, but not on Ser15 and Ser78 both *in vitro* and in intact cells³. Phosphorylation of HSP27 causes a change in its tertiary structure, which shifts from large homotypic multimers to dimers and monomers. Thus, phosphorylated HSPs have significantly decreased ability to act as molecular chaperones suppressing thermal denaturation and facilitating refolding of citrate synthase *in vitro*¹³. Phosphorylation of HSP27 on Ser82 by MAPKAPK-2 or AKT leads to HSP dissociation from the AKT/MAPKAPK2/p38 complex, which may promote independent survival signals³. The phosphorylation of HSP27 also enhances its association with IKK β to result in decreased IKK activity⁷.

Traditional Kinase Assays

To date, two methods are widely used to perform kinase assays:

1. One method used is the in-gel kinase assay, which is an activity staining technique used to study protein kinases¹⁴. A given protein substrate is immobilized on a gel and phosphorylated by protein kinases, which are separated by SDS-PAGE. The bands of incorporated [³²P]phosphate are then visualized by autoradiography. While this method is sensitive, it is also cumbersome and is not suitable for high-throughput applications. In-gel kinase assays also require special precautions and equipment for handling radioactivity.
2. Another method used is Western blot analysis. Western blots are performed using antibodies that recognize only the phosphorylated version of the protein of interest. Although less tedious than in-gel kinase assays, Western blotting, like in-gel kinase, requires the preparation of nuclear or whole-cell extract and separation by SDS-PAGE. Furthermore, this process is expensive due to the large quantity of phospho-specific antibody required.

FACE HSP27

To date efforts to study HSP27 signaling have been hampered by the lack of convenient and high-throughput assays suitable for quantifying HSP27 activation (phosphorylation).

To overcome this, Active Motif has introduced FACE HSP27 to its line of FACE™ (Fast Activated Cell-based ELISA) Kits. These are highly sensitive 96-well assays designed for detecting activated proteins within mammalian cells. Unlike Western blot, FACE assays do not require cell extracts, electrophoresis or membrane blotting. And, unlike typical kinase assays, FACE assays are non-radioactive and simple to perform. FACE HSP27 Kits contains two 96-well plates and two primary antibodies. The FACE phospho-HSP27 (Ser82) antibody was raised in rabbits against a synthetic phospho-peptide corresponding to residues surrounding phosphorylated Ser 82 of human HSP27. The phospho-HSP27 (Ser82) antibody recognizes HSP27 only when phosphorylated at Ser82. The total-HSP27 antibody supplied in the FACE HSP27 Kits recognizes HSP27 protein regardless of its phosphorylation state.

FACE HSP27 Kits can be used to study phosphorylated HSP27 relative to cell number. In this application, cells are cultured in the wells of one of the provided 96-well plates, treated as desired and then assayed using the FACE protocol with only the phospho-HSP27 antibody. The relative number of cells in each well is then determined through use of the Crystal Violet reagent. In this application, the second 96-well plate can be kept on reserve in case of culturing problems or two 48-well assays can be performed.

FACE HSP27 Kits can also be used to determine HSP27 phosphorylation relative to the total HSP27 protein found in the cells. In this application, the two 96-well plates are cultured as replicates, with the wells within each plate treated with reagents that may affect the phosphorylation state of HSP27. After the cells are fixed, one plate is studied with the phospho-HSP27 antibody, while the other plate is studied with the total-HSP27 antibody. The relative number of cells in each well is then determined through use of the Crystal Violet reagent. Once the phospho-HSP27 and total-HSP27 signals have been normalized for cell number, a comparison of the ratio of phosphorylated HSP27 to total HSP27 for each of the cell growth conditions can be made.

In the FACE HSP27 assay, the provided total-HSP27 antibody can be used as a positive control to demonstrate that the cells contain HSP27, the kit reagents are functional and that the protocol is performed correctly. Also, because fixed cells are stable for several weeks, you can prepare many plates simultaneously and then perform the FACE assay when desired. Fixed cells should be stored refrigerated in a zip-lock or heat-sealed bag with the formaldehyde solution in the wells.

Kit Performance and Benefits

FACE HSP27 Kits are for research use only. Not for use in diagnostic procedures.

Antibody specificity: The FACE phospho-HSP27 (S82) antibody was raised in rabbits against a synthetic phospho-peptide corresponding to residues surrounding phosphorylated Ser82 of human HSP27. The phospho-HSP27 (S82) antibody recognizes HSP27 only when phosphorylated at Ser82. The total-HSP27 antibody supplied in the FACE HSP27 Kits recognizes HSP27 protein regardless of its phosphorylation state.

Cross-reactivity: The phospho-HSP27 (S82) antibody detects human phosphorylated HSP27. The total-HSP27 antibody detects HSP27 from human and mouse origin.

Assay time: < 3 hours of hands-on time.

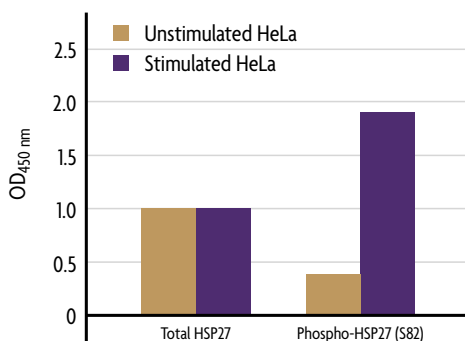


Figure 1: Measurement of phosphorylated HSP27 (S82) and total HSP27.

Human HeLa cells were cultured in 96-well plates and serum-starved (0.5% FBS) for 16 hours. Cells were then stimulated with 25 µg/ml of Anisomycin for 30 minutes and fixed. Total HSP27 and HSP27 phosphorylated on Ser82 were each assayed in duplicate using the phospho-HSP27 (S82) and total-HSP27 antibodies included in the FACE HSP27 Kit. Data were plotted after correction for cell number (performed through use of Crystal Violet).

Note on data interpretation

The phospho-HSP27 and total-HSP27 antibodies can be used on equivalent cell cultures to determine the effects of various cell treatments on the ratio of phosphorylated HSP27 to total HSP27. However, if the signals obtained with the phospho-HSP27 antibody and the total-HSP27 antibody are identical, one cannot conclude that the treatment resulted in phosphorylation of 100% of the HSP27.

FACE HSP27 Experimental Design

The FACE HSP27 assay is a high-throughput method for quantifying cellular levels of HSP27 and phosphorylated HSP27. It should be used with cell types that have been shown to contain readily detectable levels of HSP27 and, under appropriate induction conditions, phosphorylated HSP27.

Before starting a FACE assay, it is necessary to determine the experimental conditions for each well of the 96-well plate to maximize the information obtained.

Points to consider:

1. Are you working with adherent or non-adherent cells? Protocol modifications for use of non-adherent cells are given after the protocol for adherent cells.
2. Do you want to compare phosphorylated HSP27 to total HSP27? If so, replicate wells must be cultured so that the two different antibodies can be used on equivalently grown cells.
3. Which wells will be used as positive controls (*i.e.* incubated with the total-HSP27 antibody) and which will be used as negative controls (*i.e.* incubated with secondary antibody alone)?
4. Each experimental condition should be performed in duplicate or in triplicate to control for possible errors.
5. FACE assays are most easily performed when all 96 wells of the assay plate are used. This makes it possible to perform washing steps by “flicking” liquid from the plate into a sink. The inverted plate is then tapped gently onto several layers of paper towel to remove the remaining liquid. See “Kit Components” section if you need additional 96-well plates.
6. Fixed cells are stable for several weeks, so you can prepare many plates simultaneously and then perform the FACE assay when desired. Fixed cells should be stored with the formaldehyde solution in the wells and then sealed in a zip-lock bag or, preferably, a heat-sealed bag and refrigerated.

After planning the experiment, determine the amount of each buffer/reagent required and prepare according to the Quick Chart for Preparing Buffers. Multi-channel pipettors and pipettor reservoirs should be used when appropriate. The volumes given are appropriate for multi-channel pipetting if the assay is performed on 48 wells or more. Volumes may need to be adjusted if the assay is performed on less than 48 wells.

Kit Components and Storage - Colorimetric Assay

FACE Colorimetric Kit components can be stored at -20°C prior to first use. Then, we recommend storing each component at the temperature indicated in the table below.

Reagents	Quantity	Storage / Stability
	1 x 96 rxns / 5 x 96 rxns	
Phospho-HSP27 (S82) antibody	9 µl / 45 µl	-20°C for 6 months
Total-HSP27 antibody	9 µl / 45 µl	-20°C for 6 months
Anti-rabbit HRP-conjugated IgG	11 µl / 55 µl	4°C for 6 months
1X Antibody Blocking Buffer	22 ml / 110 ml	-20°C for 6 months
1X Antibody Dilution Buffer	30 ml / 150 ml	-20°C for 6 months
10X PBS (pH 7.4)	120 ml / 600 ml	Room temperature for 6 months
10% Triton X-100	9 ml / 45 ml	Room temperature for 6 months
Crystal Violet Solution	22 ml / 110 ml	4°C for 6 months
Developing Solution	22 ml / 110 ml	4°C for 6 months
Stop Solution	22 ml / 110 ml	4°C for 6 months
1% SDS Solution	22 ml / 110 ml	Room temperature for 6 months
96-well tissue culture plate*	2 / 10	
Plate sealing tape	2 / 10	

* Suitable tissue culture plates are Greiner part no. 655180 and Corning Costar part no. 3596.

Additional materials required

- Multi-channel pipettor
- Multi-channel pipettor reservoirs
- Rocking platform
- Parafilm
- Microplate spectrophotometer capable of reading at 595 nm and at 450 nm (655 as optional reference wavelength)
- Fresh 10% hydrogen peroxide (H₂O₂) in dH₂O (3 ml are required)
- 10 mg/ml poly-L-Lysine (if using non-adherent cells)
- 10% Sodium Azide (NaN₃) in dH₂O (250 µl are required)
- 37% Formaldehyde (2.5 ml are required for adherent cells; 5.0 ml required for non-adherent cells)

WARNING: Sodium Azide and Formaldehyde are highly toxic chemicals. Appropriate safety precautions (gloves and eye protection) should be used. In addition, formaldehyde is highly toxic by inhalation and should be used only in a ventilated hood.

Kit Components and Storage - Chemiluminescent Assay

FACE Chemi Kit components can be stored at -20°C prior to first use. Then, we recommend storing each component at the temperature indicated in the table below.

Reagents	Quantity	Storage / Stability
	1 x 96 rxns / 5 x 96 rxns	
Phospho-HSP27 (S82) antibody	9 µl / 45 µl	-20°C for 6 months
Total-HSP27 antibody	9 µl / 45 µl	-20°C for 6 months
Anti-rabbit HRP-conjugated IgG	11 µl / 55 µl	4°C for 6 months
1X Antibody Blocking Buffer	22 ml / 110 ml	-20°C for 6 months
1X Antibody Dilution Buffer	30 ml / 150 ml	-20°C for 6 months
10X PBS (pH 7.4)	120 ml / 600 ml	Room temperature for 6 months
10% Triton X-100	9 ml / 45 ml	Room temperature for 6 months
Crystal Violet Solution	22 ml / 110 ml	4°C for 6 months
Chemiluminescent Reagent	4 ml / 20 ml	4°C for 6 months
Reaction Buffer	8 ml / 40 ml	4°C for 6 months
1% SDS Solution	22 ml / 110 ml	Room temperature for 6 months
96-well tissue culture plate*	2 / 10	
Plate sealing tape	2 / 10	

* Suitable tissue culture plates are Greiner part no. 655098.

Additional materials required

- Multi-channel pipettor
- Multi-channel pipettor reservoirs
- Rocking platform
- Parafilm
- Microplate spectrophotometer capable of reading at 595 nm for Crystal Violet staining
- Microplate luminometer or CCD camera-coupled imaging system for chemiluminescent detection
- Fresh 10% hydrogen peroxide (H₂O₂) in dH₂O (3 ml are required)
- 10 mg/ml poly-L-Lysine (if using non-adherent cells)
- 10% Sodium Azide (NaN₃) in dH₂O (250 µl are required)
- 37% Formaldehyde (2.5 ml are required for adherent cells; 5.0 ml required for non-adherent cells)

WARNING: Sodium Azide and Formaldehyde are highly toxic chemicals. Appropriate safety precautions (gloves and eye protection) should be used. In addition, formaldehyde is highly toxic by inhalation and should be used only in a ventilated hood.

Protocols - Colorimetric Assay

Buffer Preparation and Recommendations

We provide an excess of buffer components in order to perform one 96-well FACE assay with the phospho-HSP27 antibody and one 96-well FACE assay with the total-HSP27 antibody. Required reagents that are not supplied are listed on the previous page. Please review the Quick Chart for Preparing Buffers in this section prior to preparing the assay buffers.

Preparation of 1X PBS

1X PBS is the basis of several buffers used in the FACE protocol. 1X PBS is also used in several of the wash steps in the protocol (see the Quick Chart for Preparing Buffers). It is prepared by adding 1 volume of 10X PBS (pH 7.4) to 9 volumes of dH₂O and mixing thoroughly.

Preparation of Fixing Buffer (4% or 8% Formaldehyde in PBS)

Fixing Buffer is used to fix cells after cell culturing. It is prepared by adding formaldehyde to 1X PBS and mixing well. 4% formaldehyde is used with adherent cells, 8% formaldehyde is used with non-adherent cells. The recipe in the Quick Chart for Preparing Buffers is written for use with a stock solution of 37% formaldehyde.

Preparation of Wash Buffer (0.1% Triton X-100 in PBS)

Wash Buffer is used throughout the FACE protocol and is prepared by adding the provided 10% Triton X-100 solution to 1X PBS and mixing thoroughly.

Quenching Buffer (Wash Buffer containing 1% H₂O₂ and 0.1% Azide)

Quenching Buffer is used to inactivate the cells' endogenous peroxidase activity. It is prepared by adding fresh Sodium Azide and fresh hydrogen peroxide to the Wash Buffer.

Blocking Buffer

This is supplied ready-to-use. A small amount of white precipitate may form if thawed in a warm water bath. This does not interfere with buffer function.

Antibody Dilution Buffer

This is supplied ready-to-use. A small amount of white precipitate may form if thawed in a warm water bath. This does not interfere with buffer function.

Diluted phospho-HSP27 antibody

The phospho-HSP27 (S82) antibody recognizes only the Ser82 phosphorylated form of the HSP27 protein. The phospho-antibody will be diluted 1/500 in Antibody Dilution Buffer (see the Quick Chart for Preparing Buffers in this section).

Diluted total-HSP27 antibody

The total-HSP27 antibody recognizes both the non-phosphorylated and the phosphorylated forms of HSP27 protein. The supplied antibody will be diluted 1/500 in Antibody Dilution Buffer (see the Quick Chart for Preparing Buffers in this section).

Diluted HRP-conjugated secondary antibody

HRP-conjugated anti-rabbit IgG is used as the secondary antibody to detect bound primary antibodies. The supplied antibody will be diluted 1/2000 in Antibody Dilution Buffer (see the Quick Chart for Preparing Buffers in this section).

1% SDS Solution

1% SDS Solution is used in the Crystal Violet counting procedure to solubilize cells and release the dye for subsequent quantification at 595 nm. This buffer is supplied ready-to-use.

Crystal Violet Solution

This is supplied ready-to-use. Crystal Violet is used to determine the relative number of cells in each well. This stain binds to cell nuclei and gives an OD₅₉₅ reading that is proportional to cell number.

Developing Solution

The Developing Solution must be warmed to room temperature before use. This solution is light sensitive, therefore, we recommend avoiding direct exposure to intense light during storage. The Developing Solution may develop a yellow hue over time. This does not affect product performance. A blue color present in the solution indicates that it has been contaminated and must be discarded. Prior to use, transfer the amount of Developing Solution required for the assay into a secondary container (see the Quick Chart for Preparing Buffers in this section), avoid direct exposure to intense light and leave at room temperature for at least 1 hour. After use, discard any remaining solution that was transferred into the secondary container.

Stop Solution

Prior to use, transfer the amount of Stop Solution required for the assay into a secondary container (see the Quick Chart for Preparing Buffers in this section). After use, discard any remaining Stop Solution that was transferred into the secondary container.

WARNING: The Stop Solution is corrosive. Wear personal protective equipment when handling, *i.e.* labcoat, gloves and eye protection.

Quick Chart for Preparing Buffers - Colorimetric Assay

Reagents to prepare	Components	1 well	48 wells	96 wells	192 wells
Fixing Buffer for adherent cells	1X PBS	98 µl	4.7 ml	9.41 ml	18.82 ml
	37% Formaldehyde	12 µl	576 µl	1.15 ml	2.30 ml
	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Fixing Buffer for non-adherent cells	1X PBS	86.0 µl	4.13 ml	8.26 ml	16.51 ml
	37% Formaldehyde	24.0 µl	1.15 ml	2.30 ml	4.61 ml
	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Wash Buffer	1X PBS	3.376 ml	162 ml	310 ml	620 ml
	10% Triton X-100	34.1 µl	1.64 ml	3.13 ml	6.26 ml
	TOTAL REQUIRED	3.41 ml	163.7 ml	313 ml	626 ml
Quenching Buffer	Wash Buffer	97.9 µl	4.7 ml	9.40 ml	18.8 ml
	10% H ₂ O ₂	11 µl	528 µl	1.06 ml	2.11 ml
	10% Azide	1.1 µl	52.8 µl	106 µl	211 µl
	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Blocking Buffer	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Diluted total-HSP27 antibody	Antibody Dilution Buffer	45 µl	2160 µl	4320 µl	-
	Total-HSP27 antibody	0.09 µl	4.32 µl	8.64 µl	-
	TOTAL REQUIRED	45.09 µl	2164.32 µl	4328.64 µl	-
Diluted phospho-HSP27 antibody	Antibody Dilution Buffer	45 µl	2160 µl	4320 µl	-
	Phospho-HSP27 antibody	0.09 µl	4.32 µl	8.64 µl	-
	TOTAL REQUIRED	45.09 µl	2164.32 µl	4328.64 µl	-
Diluted HRP-conjugated secondary antibody	Antibody Dilution Buffer	110 µl	5.28 ml	10.56 ml	21.12 ml
	HRP-conjugated secondary ab	0.055 µl	2.64 µl	5.28 µl	10.56 µl
	TOTAL REQUIRED	110.055 µl	5.282 ml	10.56 ml	21.13 ml
1X PBS (for wash steps)	10X PBS	154 µl	7.39 ml	14.11 ml	28.22 ml
	dH ₂ O	1.39 ml	66.53 ml	127.01 ml	254.02 ml
	TOTAL REQUIRED	1.54 ml	73.92 ml	141.12 ml	282.24 ml
1% SDS Solution	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Developing Solution	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Stop Solution	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Crystal Violet Solution	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml

Adherent Cell Protocol - Colorimetric Assay

PLEASE READ THE ENTIRE PROTOCOL BEFORE STARTING

Step 1: Culture, fix and block cells

1. Seed cells in the 96-well plate so that they will be approximately 80% confluent at the time of fixing, after they have been treated as desired. The growth area in each well of the 96-well plate is 0.32 cm². The provided plates are sterile and treated for tissue culture.
2. Grow and treat cells as desired.
3. Fix cells by replacing the growth medium with 100 µl of 4% formaldehyde in PBS. To minimize the escape of formaldehyde vapors, place a 10 cm x 17 cm piece of parafilm over the plate and then cover the plate with the lid. The covered plate can also be placed in a zip-lock bag. Incubate for 20 minutes at room temperature.

WARNING: Formaldehyde is highly toxic. Confine vapors to a chemical hood and wear appropriate gloves and eye protection when using this chemical.

4. Remove formaldehyde solution and wash cells 3 times with 200 µl Wash Buffer. Each wash step should be performed for 5 minutes with gentle shaking.
5. Remove Wash Buffer, add 100 µl Quenching Buffer and incubate for 20 minutes at room temperature.
6. Remove Quenching Buffer and wash cells 2 times for 5 minutes each with 200 µl Wash Buffer.
7. Remove Wash Buffer, add 100 µl Antibody Blocking Buffer and incubate 1 hour at room temperature.

Step 2: Binding of primary and secondary antibodies

NOTE: Depending on experiment design, some wells may be incubated with diluted phospho-HSP27 antibody, some with total-HSP27 antibody and some with secondary antibody alone (negative controls). For negative control wells, incubate with 40 µl Antibody Dilution Buffer during primary antibody incubation step.

1. Remove Antibody Blocking Buffer and wash cells 2 times with 200 µl Wash Buffer.
2. Remove Wash Buffer, add 40 µl of diluted primary antibody (or Antibody Dilution Buffer for negative control wells) and seal plate with sealing tape. Place a 10 cm x 17 cm piece of parafilm over the plate, cover with lid and incubate overnight at 4°C. Be sure that the plate is level and that each well is tightly sealed with the sealing tape to prevent evaporation.

NOTE: In cells known to generate high amounts of phosphorylated-HSP27, a three hour primary antibody incubation is sufficient. For maximum sensitivity an overnight incubation is recommended.

3. Remove primary antibody, wash cells 3 times for 5 minutes each with 200 µl Wash Buffer.

4. Remove Wash Buffer, add 100 μ l diluted secondary antibody, cover plate with tissue culture plate lid or sealing tape, and incubate 1 hour at room temperature.
5. During this incubation, transfer the amount of Developing Solution required for the assay into a secondary container and leave at room temperature for at least an hour (avoid light).

Step 3: Colorimetric reaction

1. Remove secondary antibody, wash cells 3 times for 5 minutes with 200 μ l Wash Buffer and then 2 times for 5 minutes with 200 μ l 1X PBS.
2. Transfer the amount of Developing Solution required for the assay into a secondary container. Remove PBS from plate wells and add 100 μ l Developing Solution to each well.
3. Incubate 2-20 minutes at room temperature protected from direct light. Monitor the blue color development until the darkest-staining wells are medium- to dark-blue. Do not over-develop.
4. Add 100 μ l Stop Solution. This acidic solution turns the blue color to yellow. Take care with pipetting to ensure that each well is developed for the same amount of time.

WARNING: The Stop Solution is corrosive. Wear personal protective equipment when handling, *i.e.* labcoat, gloves and eye protection.

5. Read absorbance on a spectrophotometer within 5 minutes at 450 nm with an optional reference wavelength of 655 nm.

OPTIONAL - Crystal Violet cell staining

Crystal Violet is an intense stain that binds to the cell nuclei and gives an OD₅₉₅ reading that is proportional to cell number. If you wish to normalize your readings from above simply follow the steps below.

1. After reading at 450 nm is complete, wash wells twice with 200 μ l Wash Buffer and 2 times with 200 μ l 1X PBS. Tap plates onto paper towels to remove excess liquid from wells and air-dry at room temperature for 5 minutes.
2. Add 100 μ l Crystal Violet solution to each well and incubate 30 minutes at room temperature.

WARNING: Crystal Violet is an intense stain. Avoid contact with skin and clothing.

3. Wash wells 3 times with 200 μ l 1X PBS for 5 minutes each.
4. Add 100 μ l of 1% SDS Solution to each well and incubate on shaker for 1 hour at room temperature.
5. Read absorbance on a spectrophotometer at 595 nm. If the signals obtained are greater than the range of your spectrophotometer, the signal can be reduced by removing some (*e.g.* 50 μ l) of the liquid from each well and replacing with an equivalent volume of dH₂O.
6. The measured OD₄₅₀ readings are corrected for cell number by dividing the OD₄₅₀ reading for a given well by the OD₅₉₅ reading for that well.

Non-Adherent Cell Protocol - Colorimetric Assay

The protocol given above can be modified for use with non-adherent cells by culturing and fixing the cells as follows:

1. Treat the 96-well culture plate with 10 µg/ml poly-L-Lysine for 30 minutes at 37°C. Wash twice for 5 minutes with PBS.
2. Seed 17,000 cells/well, or whatever amount is appropriate for your particular cell line.
3. Grow and treat cells as desired.
4. Fix cells by replacing the growth medium with 100 µl of 8% formaldehyde in PBS. Incubate 20 minutes at room temperature.
5. Continue with Step 1, No. 4 of the Adherent Cell Protocol above.

Protocols - Chemiluminescent Assay

Buffer Preparation and Recommendations

We provide an excess of buffer components in order to perform one 96-well FACE assay with the phospho-HSP27 antibody and one 96-well FACE assay with the total-HSP27 antibody. Required reagents that are not supplied are listed on the previous page. Please review the Quick Chart for Preparing Buffers in this section prior to preparing the assay buffers.

Preparation of 1X PBS

1X PBS is the basis of several buffers used in the FACE protocol. 1X PBS is also used in several of the wash steps in the protocol (see the Quick Chart for Preparing Buffers). It is prepared by adding 1 volume of 10X PBS (pH 7.4) to 9 volumes of dH₂O and mixing thoroughly.

Preparation of Fixing Buffer (4% or 8% Formaldehyde in PBS)

Fixing Buffer is used to fix cells after cell culturing. It is prepared by adding formaldehyde to 1X PBS and mixing well. 4% formaldehyde is used with adherent cells, 8% formaldehyde is used with non-adherent cells. The recipe in the Quick Chart for Preparing Buffers is written for use with a stock solution of 37% formaldehyde.

Preparation of Wash Buffer (0.1% Triton X-100 in PBS)

Wash Buffer is used throughout the FACE protocol and is prepared by adding the provided 10% Triton X-100 solution to 1X PBS and mixing thoroughly.

Quenching Buffer (Wash Buffer containing 1% H₂O₂ and 0.1% Azide)

Quenching Buffer is used to inactivate the cells' endogenous peroxidase activity. It is prepared by adding fresh Sodium Azide and fresh hydrogen peroxide to the Wash Buffer.

Blocking Buffer

This is supplied ready-to-use. A small amount of white precipitate may form if thawed in a warm water bath. This does not interfere with buffer function.

Antibody Dilution Buffer

This is supplied ready-to-use. A small amount of white precipitate may form if thawed in a warm water bath. This does not interfere with buffer function.

Diluted phospho-HSP27 antibody

The phospho-HSP27 (S82) antibody recognizes only the Ser82 phosphorylated form of the HSP27 protein. The phospho-antibody will be diluted 1/500 in Antibody Dilution Buffer (see the Quick Chart for Preparing Buffers in this section).

Diluted total-HSP27 antibody

The total-HSP27 antibody recognizes both the non-phosphorylated and the phosphorylated forms of HSP27 protein. The supplied antibody will be diluted 1/500 in Antibody Dilution Buffer (see the Quick Chart for Preparing Buffers in this section).

Diluted HRP-conjugated secondary antibody

HRP-conjugated anti-rabbit IgG is used as the secondary antibody to detect bound primary antibodies. The supplied antibody will be diluted 1/2000 in Antibody Dilution Buffer (see the Quick Chart for Preparing Buffers in this section).

Preparation of Chemiluminescent Working Solution

The Chemiluminescent Reagent and Reaction Buffer should be warmed to room temperature before use. These components are light sensitive, therefore, we recommend avoiding direct exposure to intense light during storage. Prior to use, place the Chemiluminescent Reagent and Reaction Buffer at room temperature for at least 1 hour. In a separate container, mix 1 volume of Chemiluminescent Reagent with 2 volumes of Reaction Buffer to prepare the Chemiluminescent Working Solution (see the Quick Chart for Preparing Buffers in this section). The Chemiluminescent Working Solution is stable for several hours. After the Chemiluminescent Working Solution is aliquoted into the wells, discard the remaining solution.

1% SDS Solution

1% SDS Solution is used in the Crystal Violet counting procedure to solubilize cells and release the dye for subsequent quantification at 595 nm. This buffer is supplied ready-to-use.

Crystal Violet Solution

This is supplied ready-to-use. Crystal Violet is used to estimate the relative number of cells in each well. This stain binds to cell nuclei and gives an OD₅₉₅ reading that is proportional to cell number.

Quick Chart for Preparing Buffers - Chemiluminescent Assay

Reagents to prepare	Components	1 well	48 wells	96 wells	192 wells
Fixing Buffer for adherent cells	1X PBS	98 µl	4.7 ml	9.41 ml	18.82 ml
	37% Formaldehyde	12 µl	576 µl	1.15 ml	2.30 ml
	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Fixing Buffer for non-adherent cells	1X PBS	86.0 µl	4.13 ml	8.26 ml	16.51 ml
	37% Formaldehyde	24.0 µl	1.15 ml	2.30 ml	4.61 ml
	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Wash Buffer	1X PBS	3.376 ml	162 ml	310 ml	620 ml
	10% Triton X-100	34.1 µl	1.64 ml	3.13 ml	6.26 ml
	TOTAL REQUIRED	3.41 ml	163.7 ml	313 ml	626 ml
Quenching Buffer	Wash Buffer	97.9 µl	4.7 ml	9.40 ml	18.8 ml
	10% H ₂ O ₂	11 µl	528 µl	1.06 ml	2.11 ml
	10% Azide	1.1 µl	52.8 µl	106 µl	211 µl
	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Blocking Buffer	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Diluted total-HSP27 antibody	Antibody Dilution Buffer	45 µl	2160 µl	4320 µl	-
	Total-HSP27 antibody	0.09 µl	4.32 µl	8.64 µl	-
	TOTAL REQUIRED	45.09 µl	2164.32 µl	4328.64 µl	-
Diluted phospho-HSP27 antibody	Antibody Dilution Buffer	45 µl	2160 µl	4320 µl	-
	Phospho-HSP27 antibody	0.09 µl	4.32 µl	8.64 µl	-
	TOTAL REQUIRED	45.09 µl	2164.32 µl	4328.64 µl	-
Diluted HRP-conjugated secondary antibody	Antibody Dilution Buffer	110 µl	5.28 ml	10.56 ml	21.12 ml
	HRP-conjugated secondary ab	0.055 µl	2.64 µl	5.28 µl	10.56 µl
	TOTAL REQUIRED	110.055 µl	5.282 ml	10.56 ml	21.13 ml
1X PBS (for wash steps)	10X PBS	154 µl	7.39 ml	14.11 ml	28.22 ml
	dH ₂ O	1.39 ml	66.53 ml	127.01 ml	254.02 ml
	TOTAL REQUIRED	1.54 ml	73.92 ml	141.12 ml	282.24 ml
Chemiluminescent Working Solution	Chemiluminescent Reagent	18 µl	864 µl	1.728 ml	3.46 ml
	Reaction Buffer	36 µl	1.728 ml	3.456 ml	6.91 ml
	TOTAL REQUIRED	54 µl	2.592 ml	5.184 ml	10.37 ml
1% SDS Solution	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml
Crystal Violet Solution	TOTAL REQUIRED	110 µl	5.28 ml	10.56 ml	21.12 ml

Adherent Cell Protocol - Chemiluminescent Assay

PLEASE READ THE ENTIRE PROTOCOL BEFORE STARTING

Step 1: Culture, fix and block cells

1. Seed cells in the 96-well plate so that they will be approximately 80% confluent at the time of fixing, after they have been treated as desired. The growth area in each well of the 96-well plate is 0.32 cm². The provided plates are sterile and treated for tissue culture.
2. Grow and treat cells as desired.
3. Fix cells by replacing the growth medium with 100 µl of 4% formaldehyde in PBS. To minimize the escape of formaldehyde vapors, place a 10 cm x 17 cm piece of parafilm over the plate and then cover the plate with the lid. The covered plate can also be placed in a zip-lock bag. Incubate for 20 minutes at room temperature.

WARNING: Formaldehyde is highly toxic. Confine vapors to a chemical hood and wear appropriate gloves and eye protection when using this chemical.

4. Remove formaldehyde solution and wash cells 3 times with 200 µl Wash Buffer. Each wash step should be performed for 5 minutes with gentle shaking.
5. Remove Wash Buffer, add 100 µl Quenching Buffer and incubate for 20 minutes at room temperature.
6. Remove Quenching Buffer and wash cells 2 times for 5 minutes each with 200 µl Wash Buffer.
7. Remove Wash Buffer, add 100 µl Antibody Blocking Buffer and incubate 1 hour at room temperature.

Step 2: Binding of primary and secondary antibodies

NOTE: Depending on experiment design, some wells may be incubated with diluted phospho-HSP27 antibody, some with total-HSP27 antibody and some with secondary antibody alone (negative controls). For negative control wells, incubate with 40 µl Antibody Dilution Buffer during primary antibody incubation step.

1. Remove Antibody Blocking Buffer and wash cells 2 times with 200 µl Wash Buffer.
2. Remove Wash Buffer, add 40 µl of diluted primary antibody (or Antibody Dilution Buffer for negative control wells) and seal plate with sealing tape. Place a 10 cm x 17 cm piece of parafilm over the plate, cover with lid and incubate overnight at 4°C. Be sure that the plate is level and that each well is tightly sealed with the sealing tape to prevent evaporation.

NOTE: In cells known to generate high amounts of phosphorylated-HSP27, a three hour primary antibody incubation is sufficient. For maximum sensitivity an overnight incubation is recommended.

3. Remove primary antibody, wash cells 3 times for 5 minutes each with 200 µl Wash Buffer.
4. Remove Wash Buffer, add 100 µl diluted secondary antibody, cover plate with tissue culture

plate lid or sealing tape, and incubate 1 hour at room temperature.

5. During this incubation, place the Chemiluminescent Reagent and Reaction Buffer at room temperature.

Step 3: Chemiluminescent detection

1. Remove secondary antibody, wash cells 3 times for 5 minutes with 200 μ l Wash Buffer and then 2 times for 5 minutes with 200 μ l 1X PBS.
2. Remove PBS from plate wells and add 50 μ l room temperature Chemiluminescent Working Solution to each well.
3. Read chemiluminescence using a luminometer or CCD camera system. Readings should be taken within 10 minutes to minimize changes in signal intensity.

OPTIONAL - Crystal Violet cell staining

Crystal Violet is an intense stain that binds to the cell nuclei and gives an OD₅₉₅ reading that is proportional to cell number. If you wish to normalize your readings from above simply follow the steps below.

1. After reading chemiluminescence, wash wells twice with 200 μ l Wash Buffer and 2 times with 200 μ l 1X PBS. Tap plates onto paper towels to remove excess liquid from wells and air-dry at room temperature for 5 minutes.
2. Add 100 μ l Crystal Violet solution to each well and incubate 30 minutes at room temperature.
WARNING: Crystal Violet is an intense stain. Avoid contact with skin and clothing.
3. Wash wells 3 times with 200 μ l 1X PBS for 5 minutes each.
4. Add 100 μ l of 1% SDS Solution to each well and incubate on shaker for 1 hour at room temperature.
5. Read absorbance on a spectrophotometer at 595 nm. If the signals obtained are greater than the range of your spectrophotometer, the signal can be reduced by removing some (e.g. 50 μ l) of the liquid from each well and replacing with an equivalent volume of dH₂O.
6. The measured OD₅₉₅ readings indicate the relative number of cells in each well. This relative cell number is then used to normalize each reading from Step 3.

Non-Adherent Cell Protocol - Chemiluminescent Assay

The protocol given above is suitable for use with non-adherent cells if the cells are cultured and fixed as follows:

1. Treat the 96-well culture plate with 10 µg/ml poly-L-Lysine for 30 minutes at 37°C. Wash twice for 5 minutes with PBS.
2. Seed 17,000 cells/well, or whatever amount is appropriate for your particular cell line.
3. Grow and treat cells as desired.
4. Fix cells by replacing the growth medium with 100 µl of 8% formaldehyde in PBS. Incubate 20 minutes at room temperature.
5. Continue with Step 1, No. 4 of the Adherent Cell Protocol above.

References

1. Arrigo, A.P. and Landry, J. (1994) Cold Spring Harbor Laboratory Press, NY, 335-373.
2. De Jong, W.W., *et al.* (1988) *J. Biol. Chem.* 263: 5141-5149.
3. Rane, M.J., *et al.* (2003) *J. Biol. Chem.* 278: 27828-27835.
4. Versteeg H.H. *et al.* (2000) *Biochem J.* 350(Pt3): 717-22.
5. Pantos, C., *et al.* (2003) *Eur. J. Pharmacol.* 478: 53-60.
6. Rane, M.J., *et al.* (2001) *J. Biol. Chem.* 276: 3517-3523.
7. Park, K.J., *et al.* (2003) *J. Biol. Chem.* 278: 35272-35278.
8. Lavoie, J.N., *et al.* (1993) *J. Biol. Chem.* 268: 3420-3429.
9. Lavoie, J.N., *et al.* (1993) *J. Biol. Chem.* 268: 24210-24214.
10. Rousseau, S., *et al.* (1997) *Oncogene* 15: 2169-2177.
11. Landry, J., *et al.* (1992) *J. Biol. Chem.* 267: 794-803.
12. Rouse, J., *et al.* (1994) *Cell* 78: 1027-1037.
13. Rogalla, T., *et al.* (1999) *J. Biol. Chem.* 274: 18947-18956.
14. Kameshita I., *et al.* (1989) *Analytical Biochem.* 183: 139-143.

* The FACE method was developed in the laboratory of Dr. Maikel P. Peppelenbosch, Laboratory for Experimental Internal Medicine, Academic Medical Centre, Amsterdam, The Netherlands. We thank Dr. Henri H. Versteeg and Dr. Peppelenbosch for their assistance in developing the FACE Kits.

Appendix

Section A. Troubleshooting Guide

PROBLEM	POSSIBLE CAUSE	RECOMMENDATION
No signal or weak signal in wells incubated with either phospho-HSP27 antibody or total-HSP27 antibody	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 to prepare buffers
	Plate reader or CCD camera settings not optimal	Verify the wavelength (measurement mode) and filter settings in the plate reader
	Developing Solution was cold	Bring Developing Solution to room temperature
	Inadequate volume of Developing Solution	Check to make sure that correct volume is delivered by pipette
	Cells do not contain detectable levels of phospho HSP27 and/or total HSP27	Use Western blotting to confirm that cells contain detectable levels of protein(s) of interest. If you do not require all of the included antibodies for FACE assays, they can be used in Colorimetric Western blotting at a 1:500 dilution for the total and a 1:100 for the phospho antibody. Use the secondary antibody provided in the FACE kit at 1 in 2000 for Western blotting. Alternatively, the customer can use their own secondary antibody.
	Insufficient number of cells were plated	Plate cells so that they are 80% confluent at time of fixing
	Cells did not adhere correctly to plate	Follow protocol for use of non-adherent cells
	Cells are not from correct origin	The phospho HSP27 (S82) antibody detects human HSP27 (S82) and the total HSP27 antibody detects human and mouse HSP27. .
	Excessive washing	Wash steps should be 5 minutes each
Incubation of secondary antibody was too long	Incubate secondary antibody for 1 hour	
High background in all wells	Developing time too long (Colorimetric Assay)	Stop enzymatic reaction as soon as the positive wells turn medium-dark blue
	Measurement time too long (Chemiluminescent Assay)	Reduce integration time or exposure time on luminometer or CCD camera
	Concentration of antibodies too high	Perform antibody titration to determine optimal working concentration. Use primary antibody only, or secondary antibody only to confirm that the high background came from the high antibody concentration. Start with 1 in 1000 for the primary antibody (total and phospho), while keeping the secondary at 1 in 2000. If the background is still high, then try the secondary antibody at 1 in 4000, while keeping the primary at 1 in 1000. Different cell lines react differently with the antibodies. The sensitivity of the assay may be decreased

High background in all wells (cont.)	Inadequate washing	Ensure all wells are filled with Wash Buffer and follow washing recommendations
	Inadequate quenching or blocking	Ensure that quenching and blocking steps were performed according to the protocol
Uneven color development	Incomplete washing of wells	Ensure all wells are filled with Wash Buffer and follow washing recommendations
	Well cross-contamination	Follow washing recommendations
No signal or weak signal in wells incubated with phospho-HSP27 antibody	Cell culture conditions did not induce phosphorylation of HSP27	Perform Western blot with phospho-HSP27 antibody to confirm that cells contain detectable levels of phosphorylated HSP27
Antibody solution evaporates from well during overnight incubation with primary antibody	Sealing tape was incorrectly applied	Ensure that each well is sealed when sealing tape is applied and ensure that the parafilm sheet covers the plate completely before the lid is placed on the plate. The plate can also be placed in a zip-lock or heat-sealed bag
Insufficient sensitivity	Antibody concentration incorrect	If the cells studied have very low levels of the protein of interest, the sensitivity of detection may be improved by increasing the concentration of primary antibody used and by minimizing the incubation volume. It is possible to perform the overnight incubation in as little as 25 μ l, however, this will make multichannel pipetting difficult and requires the plate be carefully sealed and incubated on a level surface. Alternatively, if the cells have easily detectable levels of the phosphorylated protein and the detection of small changes in phosphorylation is desired, sensitivity of the assay may be improved by decreasing the concentration of the phospho antibody used
Poor precision	Cross-well read through	The 96-well plates provided are designed to minimize signal cross-well contamination. If possible, do not use the phospho and total antibodies in adjoining wells. If this is not possible, use the total antibody at a higher dilution

Section B. Related Products

TransAM™ Kits	Unit	Catalog No.
TransAM™ CREB	1 x 96 rxns	42096
	5 x 96 rxns	42596
TransAM™ pCREB	1 x 96 rxns	43096
	5 x 96 rxns	43596
TransAM™ c-Myc	1 x 96 rxns	43396
	5 x 96 rxns	43896
TransAM™ Elk-1	1 x 96 rxns	44396
	5 x 96 rxns	44896
TransAM™ MAPK Family	2 x 96 rxns	47296
TransAM™ STAT Family	2 x 96 rxns	42296
TransAM™ STAT3	1 x 96 rxns	45196
	5 x 96 rxns	45696
TransAM™ IRF Family	2 x 96 rxns	45296
Cell-based ELISAs	Unit	Catalog No.
FACE™ AKT	1 x 96 rxns	48120
	5 x 96 rxns	48620
FACE™ AKT Chemi	1 x 96 rxns	48220
	5 x 96 rxns	48720
FACE™ ErbB-2 (Y877)	1 x 96 rxns	48130
	5 x 96 rxns	48630
FACE™ ErbB-2 (Y877) Chemi	1 x 96 rxns	48230
	5 x 96 rxns	48730
FACE™ ErbB-2 (Y1248)	1 x 96 rxns	48105
	5 x 96 rxns	48605
FACE™ ErbB-2 (Y1248) Chemi	1 x 96 rxns	48205
	5 x 96 rxns	48705
FACE™ ERK1/2	1 x 96 rxns	48140
	5 x 96 rxns	48640
FACE™ ERK1/2 Chemi	1 x 96 rxns	48240
	5 x 96 rxns	48740
FACE™ JNK	1 x 96 rxns	48110
	5 x 96 rxns	48610
FACE™ JNK Chemi	1 x 96 rxns	48210
	5 x 96 rxns	48710
FACE™ p38	1 x 96 rxns	48100
	5 x 96 rxns	48600
FACE™ p38 Chemi	1 x 96 rxns	48200
	5 x 96 rxns	48700
Sandwich ELISAs	Unit	Catalog No.
FunctionELISA™ IκBα	1 x 96 rxns	48005
	5 x 96 rxns	48505
FunctionELISA™ TRAIL	1 x 96 rxns	48010
	5 x 96 rxns	48510
FunctionELISA™ Cytochrome c	1 x 96 rxns	48006
	5 x 96 rxns	48506

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

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

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

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!