

Product Information

p38 MAPK Activity Assay Kit

Product Number **CS0250**

Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

The mitogen-activated protein kinase (MAPK) p38 pathway is a signaling cascade activated by pro-inflammatory stimuli and cellular stresses, playing a critical role in the translational regulation of pro-inflammatory cytokine synthesis. p38 MAPK regulates gene expression in response to various extracellular stimuli including growth factors, hormones, ligands of G protein coupled receptors, inflammatory cytokines (IL-1, IL-8, TNF α) and stress (UV, osmotic shock, LPS). p38 MAPK is activated by dual specificity MAPK kinase, including MAPK Kinase 3 (MKK3), MAPK Kinase 6 (MKK6) and Jun N-terminal Kinase Kinase 1 (JNKK1) that phosphorylate it on Thr¹⁸⁰ and Tyr¹⁸². p38 MAPK stimulates the activity of protein kinases: MAPK Activated Protein Kinase-2 and 3 (MAPKAP2, MAPKAP3) and MAPK-interacting Kinase 1 (MNK1), which in turn activate several transcription factors including Activating Transcription Factor 2 (ATF2), C/EBP Homologous Protein (CHOP) and Myocyte Enhancer Factor 2C (MEF2C).^{1,2}

To date four members of the p38 MAPK family have been cloned: p38 α , β , γ and δ . The α and β isoforms are characterized by their sensitivity to the pyridinyl imidazole inhibitors such as SB 203580.²

p38 MAPK is of great interest in both basic and therapeutic research due to its critical role in inflammation and stress response.²

The p38 MAPK Assay Kit provides an easy method to assay p38 MAPK activity and to explore new p38 MAPK stimuli, inhibitors and activators.^{3,4} The kit provides all the reagents required for the straightforward detection and measurement of p38 MAPK activity in cell lysates, tissue homogenates, column fractions or the purified enzyme. In addition, the kit provides a specific p38 MAPK inhibitor to enable the verification of the specificity of the kinase activity observed. The kit assay is based on immunoprecipitation of the active form of the p38 α kinase, and detection of its phosphorylation

activity on the substrate ATF2 by immunoblotting, without the need for a secondary antibody. An alternative protocol for radioactive measurement of the p38 MAPK activity is also provided.

Reagents

The kit is sufficient for 50 reactions.

- Anti-p38 MAP Kinase antibody. 0.1ml, Product Number M 0800
- Detection antibody, 0.5 mg, Product Number A 4728 - Monoclonal anti-phospho-ATF2 (pThr^{69,71}), Peroxidase conjugate
- Assay Buffer For Kinase Activity, 1 ml, Product Number A 4603 - 75 mM β -glycerophosphate (pH 7.3), 3.75mM EGTA, 30mM MgCl₂, 4.5 mM DTT, 0.15 mM Na-Vanadate, 0.3 mM ATP
- Wash Buffer 10X, 8 ml, Product Number W 3264 - 500 mM β -glycerophosphate (pH 7.3), 15mM EGTA, 10mM MgCl₂, 10 mM DTT, 1 mM Na-Vanadate
- EZviewTM Red Protein A Affinity Gel, 1ml, Product Number P 6486
- ATF2, 0.5 mg, Product Number A 2353
- p38 Inhibitor (SB 203580), 0.05 ml, Product Number P 2496

Reagents and Equipment required but not provided

- Microfuge centrifuge e.g. Eppendorf[®] microcentrifuge 5415 Series (Product Number Z60,406-2) or equivalent.
- Dulbecco's Phosphate Buffered Saline (PBS, Product Number D 8537)
- Sample buffer 4X for western blot (preparation instructions are detailed in the Appendix below).
- Materials and equipment required for immunoblot analysis procedure including the ProteoQwestTM, chemiluminescent Western blotting kit (Product Code PQ0201) or equivalent.
- Materials and equipment required for the radioactive assay (see below).

Precautions and Disclaimer

The kit is for R&D use only, not for drug, household or other uses. Please refer to the Material Safety Data Sheet (MSDS) for information regarding hazards and safe handling practices.

Preparation Instructions

1. ATF2 Substrate

Reconstitute the ATF2 substrate by adding 0.9 ml of 17 MOhm water to the ATF2 substrate bottle, to obtain about 0.6-0.7 mg/ml ATF2. Mix well by pipetting. Aliquot the suspended ATF2 substrate and store at -20°C.

Note: ATF2 protein is a Maltose binding protein (MBP) fusion protein of 53 kDa with an apparent MW of 60 kDa.

2. Detection antibody

Reconstitute the detection antibody, monoclonal anti-phospho-ATF2, by addition of 100 µl 17 MOhm water to the vial. Mix well by pipetting. Aliquot the suspended antibodies and store at -20°C.

3. Wash buffer 1x

Dilute the Wash buffer 10X ten fold with 17 MOhm water. Dilute 160 µl of Wash buffer 10x per each sample. For each set of test samples calculate accordingly.

Storage/Stability

The kit is shipped on dry ice and stored at -20 °C. Upon initial thawing, reconstitute the ATF2 substrate and the detection antibodies according to the procedure instructions and freeze them in working aliquots to avoid multiple freeze-thaw cycles.

Procedure

I. Procedure for detection by Immunoblotting

Principle of the assay

p38 MAPK is immunoprecipitated from the sample with anti-p38 antibodies and EZview Red Protein A Affinity Gel. Then, the immunoprecipitated p38 MAPK phosphorylates ATF2 and the level of ATF2 phosphorylation is detected by immunoblotting using anti-phospho-ATF2 (pThr^{69,71}), Peroxidase conjugate.

The kit can be used for detection of p38 MAPK activity in lysates of induced (e.g. UV, anisomycin etc.) versus non-induced cells (as a control), and for the analysis of p38 MAPK inhibitors using purified activated p38 MAPK.

General comments

- For each test set two negative control reactions:
 - Without the addition of the anti-p38 MAP Kinase antibody to detect non-specific protein binding to the agarose.
 - Without the addition of the ATF2 substrate to determine the existence of endogenous phosphorylated product.
- For ascertaining that the phosphorylation of the ATF2 is performed by p38 MAPK and not by any other kinase, set a p38 MAPK inhibition reaction: Set an additional reaction tube to which the p38 MAPK specific inhibitor [p38 Inhibitor (SB 203580) Product Number P 2496] will be added.

All preparations should be performed on ice unless otherwise stated.

Reaction scheme (example)

	Immunoprecipitation			Phosphorylation			Termination
	Cell lysate	Anti-p38	EZview Protein A 50% slurry	Assay buffer 2X	ATF2	Inhibitor SB203580	Sample buffer 4X
Control 1	250-1000 µl	-----	30 µl	15 µl	15 µl		12 µl
Control 2	250-1000 µl	2 µl	30 µl	15 µl	-----		12 µl
Reaction 1	250-1000 µl	2 µl	30 µl	15 µl	15 µl		12 µl
Reaction + inhibitor	250-1000 µl	2 µl	30 µl	15 µl	15 µl	2 µl	12 µl

The procedure described is for one test. Adjust the procedure according to your experiment.

Immunoprecipitation

1. Transfer up to 1 ml cell lysate (200-500 µg protein) into an Eppendorf tube. The optimal cell lysate volume is around 250 µl.
2. Add 2 µl of anti-p38 MAP Kinase antibody. Mix gently and set the tube on ice while equilibrating the EZview Red Protein A Affinity Gel beads.
3. For equilibrating the EZview Red Protein A Affinity Gel beads:
 - a. Carefully mix the gel beads until uniformly suspended. Aliquot 30 µl of the 50 % slurry into clean 1.5 ml microtube. For beads dispensing, use a wide orifice pipette tip or cut about 1 mm off the tip to enlarge the opening and allow unrestricted flow of the bead suspension.
 - b. Wash/equilibrate beads with PBS: Add 750 µl PBS to the tube, vortex and centrifuge in a microcentrifuge for 30 seconds at around 8,000xg. Carefully remove the supernatant with a micropipette (or carefully aspirate the supernatant).
 - c. Repeat the wash step as indicated above. After removing the supernatant, set the washed bead pellet on ice.
4. Briefly centrifuge the tube containing the cell lysate and the antibody (from steps 1-2) for several seconds at around 8000xg to collect all the liquid to the bottom part of the microcentrifuge tube. Carefully remove all the lysate with a 1 ml micropipette and transfer into the tube containing the washed EZview Red Protein A Affinity Gel beads from step 3c.
5. Vortex briefly and incubate with thorough, gentle rocking for 4 hours at 2 - 8 °C to allow the antibody-antigen complexes to bind the Protein A on the EZview Red Protein A Affinity Gel beads.
6. Centrifuge the tube in a microcentrifuge for 30 seconds at around 8,000xg. Set the tube on ice.
7. Aspirate the supernatant carefully (or remove with a micropipette) and set the tube (containing the bead pellet) on ice.

Note: Because of the enhanced visibility of the red affinity resin beads, it is easy to see if beads were accidentally removed during the washing steps. If

this happens, transfer the washed supernatant back into the tube and repeat the centrifugation step to pellet the resin once again.

8. Wash the bead pellet by adding 400 µl Wash buffer 1 X. Vortex briefly and incubate with a thorough and gentle mixing at 2 - 8°C, for 1 minute.
9. Centrifuge the tube in a microcentrifuge for 30 seconds at around 8,000xg. Aspirate the supernatant carefully (or remove with a micropipette) and set the tube with the bead pellet on ice.

Perform two-three additional washes as indicated in step 8-9.

ATF2 phosphorylation

1. Add to the bead pellet 15 µl of Assay buffer and 15 ml µl of reconstituted ATF2 substrate. Suspend the pellet by gentle pipetting.
2. Incubate for 30 minutes at 30°C.
3. Terminate the reaction with 12µl of SDS sample buffer 4X. Mix well.
4. Boil the sample for 5 minutes, and then spin down for 30 seconds. Save the supernatant.

Immunoblot analysis

1. Load the sample (15-20 µl) on SDS-PAGE gel (10-12%), run the gel and then transfer the proteins to nitrocellulose. Block the nitrocellulose paper using an appropriate blocking solution.
2. To detect the 53 kDa (~60 kDa apparent size) phospho-ATF2 work according to the protocol of the ProteoQwest™, chemiluminescent Western blotting kit (or equivalent). Incubate the membrane with a 1:2000 dilution in blocking buffer of the reconstituted Detection antibody [anti-phospho-ATF2 (pThr^{69,71}), Peroxidase conjugate] . During this incubation step the blot should be gently agitated at room temperature for 2 hours (or alternatively, overnight at 4°C).
3. Expose the membrane to an x-ray film. An initial 10-second exposure will indicate the need for a different exposure time.

Note: A dot blot procedure may be performed in order to get a sneak preview of the results. Refer to the appendix for procedure.

II. Procedure for radioactive detection

Perform steps 1-9 as described in the procedure for immunoprecipitation section. Continue with the following:

1. Add 1 μl [$\gamma^{32}\text{P}$]-ATP with a specific activity of 10mCi/ml to 100 μl of Assay buffer.
2. Suspend the immunoprecipitated pellet by gentle pipetting in 15 μl radioactive Assay buffer and 15 μl reconstituted ATF2 substrate.
Note: For performing p38 inhibition reaction only, add at this stage 2 μl of the p38 inhibitor (SB203580) into the "inhibition reaction" tube. We recommend to first suspend the pellet with 15 μl radioactive Assay buffer, then add the p38 inhibitor and finally to add the ATF2 substrate.
3. Incubate the tube for 30 minutes at 30°C. From this step on, work at room temperature.
4. Terminate the reaction by spotting 25 μl of the liquid phase of the sample on 2 cm x 2 cm phosphocellulose P81 squares.
5. Soak the phosphocellulose squares in 0.5% phosphoric acid.
6. Wash the phosphocellulose squares 4 times with 0.5% phosphoric acid. For each wash, agitate gently for 5-6 min.
7. Wash once with ethanol for 1 min.
8. Wash once with acetone for 1 min.
9. Dry the phosphocellulose squares at room temperature or under a heat lamp and count the radioactivity incorporated using Cerenkov mode (i.e. count the emission without scintillation liquid, using tritium channel).

Troubleshooting

Problem	Possible cause	Solution
The signal is very poor or no signal is observed	The amount of activated p38 MAPK in the sample is very low	<ul style="list-style-type: none"> • Add more than 2 μl of anti-p38 MAP Kinase antibody to the cell lysate in step 2 of the immunoprecipitation procedure. • Increase the sample volume - increasing the sample volume up to 1 ml usually has a minor effect on the interaction between the antibody and the activated p38 MAPK. • Use a lower dilution of the anti-phospho-ATF2 antibody, Peroxidase conjugate. • Increase the reaction incubation time from 30 min up to 90 min (Step 2 in the ATF2 phosphorylation section or step 3 in the radioactive procedure). • If possible, load on the gel a higher volume of reaction sample.
	There is no activated p38 MAPK in the sample	<ul style="list-style-type: none"> • Prepare a fresh lysate. • Add the appropriate phosphatase inhibitors to the sample (Product Numbers P 2850 and P 5726) or increase their concentration to prevent dephosphorylation of activated p38 MAPK. • Add the appropriate protease inhibitors to the sample (Product Number P 8340) or increase their concentration to prevent degradation of activated p38 MAPK. • Make sure the extraction buffer is not interfering with the kinase activity. Some extraction buffers that could be used for immunoprecipitation are not suitable for activity assays. We highly recommend using the CellLytic™ Cell Lysis Reagent (Product Number C 2978) for cell extract preparation. • Verify that the sample is appropriate or that the induction procedure for p38 MAPK activation is appropriate. Determine the presence of activated p38 MAPK in the sample by immunoblotting of the sample using anti activated p38 MAP Kinase (Product Number M 8177) prior to the performance of the immunoprecipitation.
	Incubation time is inadequate	<ul style="list-style-type: none"> • Prolong the incubation duration of the anti p38 antibody with the EZview Red Protein A (from several hours to overnight).
	Interfering substance present in sample	<ul style="list-style-type: none"> • Excessive detergent concentration may interfere with the interaction between the antibody and the activated p38 MAPK.
Signal is too strong or is spread out of the lane	There is an excess of ATF2 in the reaction	<ul style="list-style-type: none"> • Dilute ATF2 substrate with 17 MOhm water before adding it to the reaction tube (dilution from 1:2 to 1:10 is allowable). • Dilute the terminated reaction with more sample buffer or load a smaller volume of reaction sample.
Background is too high or additional bands are visible	Proteins bind non-specifically to Protein A, the resin beads or the microcentrifuge tube	<ul style="list-style-type: none"> • Pre-clear the sample once or several times by pre-incubation with EZview red protein A affinity gel (without the anti-p38 MAP Kinase antibody) to remove proteins that may bind non-specifically. • During the final wash (step 9 in the immunoprecipitation procedure), after suspending the resin, transfer the entire sample to a clean microcentrifuge tube before centrifuging the sample.

	Insufficient washes in the immunoprecipitation step	<ul style="list-style-type: none"> • Increase the number of washes • Prolong the duration of the washes to at least 15 minutes incubation. • Centrifuge at a lower speed to avoid non-specific trapping of lysate proteins during the initial centrifugation of Protein A/antigen complexes.
	Detection antibody is in excess	<ul style="list-style-type: none"> • Perform a higher dilution of the HRP-conjugated anti-phospho-ATF2 antibody.

Appendix

I. Preparation of 10 ml Sample buffer 4X

To 4 ml of 0.5M Tris-HCl pH-6.8 (Trizma base, Product Number T 1503 titrated to pH-6.8) add:

- 1.2 g SDS (Product Number L 3771)
- 616 mg DTT (Product Number D 0632)
- 10 mg Bromophenol blue (Product Number B 0126)
- 4 g Glycerol (Product Number G 9012).

Complete the volume to 10 ml with water. Mix well.

II. Dot Blot

The dot blot procedure may be performed in order to get a sneak preview of the results. This procedure does not replace the requirement for a western blot.

1. Cut nitrocellulose membrane to oblong strips of 5 cm x 2 cm (total of 10 cm²). This is sufficient for 10 samples. Each additional sample (spot) requires an additional membrane of 1 cm² area.
2. Spot 1 µl of the sample that is ready for loading on the gel (from step 4 in the ATF2 phosphorylation section) on the nitrocellulose membrane. Let it dry for a few minutes.
3. Rinse briefly with PBS or equivalent.
4. Incubate the membrane in Blocking Solution for 15 minutes.
5. Incubate the membrane with a 1:2000 dilution in blocking buffer of the reconstituted Detection antibody [anti-phospho-ATF2, Peroxidase conjugate]. Incubate with gentle agitation for 20 minutes at room temperature.

6. Wash the membrane five times with PBS containing 0.2% Tween 20 (or an equivalent buffer), for 4 minutes each.
7. Briefly rinse in a wash buffer (with no detergent).
8. Prepare Chemiluminescent Peroxidase Substrate solution, enough to cover the membrane (based on 0.125 ml/cm² membrane). Incubate the membrane in the reagent, with gentle agitation for 1 minute at room temperature.
9. Drain the membrane of excess reagent solution, wrap in Saran Wrap and expose to x-ray film. An initial 10 seconds exposure will indicate the need for different exposure times.

References:

1. Johnson, G.L, and Lapadat, R., Mitogen-activated protein kinase pathways mediated by ERK, JNK, and p38 protein kinases. *Science*, **298**, 1911-1912 (2002).
2. Clark, A.R., et al., Post-transcriptional regulation of gene expression by mitogen-activated protein kinase p38. *FEBS Lett.*, **546**, 37-44 (2003).
3. Eliopoulos, A.G, et al., Activation of the p38 Mitogen-activated Protein Kinase Pathway by Epstein-Barr Virus-encoded Latent Membrane Protein 1 co-regulates Interleukin-6 and Interleukin-8 production. *J. Biol. Chem.*, **274**, 16085-16096 (1999).
4. Torcia, M., et al., Nerve Growth Factor Inhibits Apoptosis in Memory B Lymphocytes via Inactivation of p38 MAPK, prevention of Bcl-2 phosphorylation, and Cytochrome c release. *J. Biol. Chem.*, **276**, 39027-39036 (2001).

EB/JK 06/2004

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.