

ab126459 – STAT3 (pY705) + total STAT3 ELISA Kit

Instructions for Use

For the semi-quantitative measurement of Human, Mouse and Rat phosphorylated STAT3 (Tyr705) and total STAT3 concentrations in cell lysates.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Introduction	2
2. Assay Summary	3
3. Kit Contents	4
4. Storage and Handling	5
5. Additional Materials Required	6
6. Preparation of Samples	7
7. Preparation of Reagents	8
8. Assay Method	10
9. Data Analysis	13
10. Troubleshooting	15

1. Introduction

ab126459 is a very rapid, convenient and sensitive assay kit that can monitor the activation or function of important biological pathways in cell lysates. By determining phosphorylated STAT3 protein in your experimental model system, you can verify pathway activation in your cell lysates. You can simultaneously measure numerous different cell lysates without spending excess time and effort in performing a Western Blot analysis.

This Sandwich ELISA kit is an in vitro enzyme-linked immunosorbent assay for the measurement of phospho-STAT3 (Tyr705) and total STAT3 in human, mouse and rat cell lysates (help normalize the results of phospho-STAT3 from different cell lysates being compared). A pan STAT3 antibody has been coated onto a 96-well plate. Samples are pipetted into the wells and STAT3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and anti-phospho-STAT3 (Tyr705) is used to detect phosphorylated STAT3 or biotinylated anti-total-STAT3 is used to detect total STAT3. After washing away unbound antibody, HRP-conjugated anti-rabbit IgG or HRP- streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of STAT3 (Tyr705) or total STAT3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

2. Assay Summary

Prepare all reagents, samples and standards as instructed.



Add 100 μ l sample or positive control to each well. Incubate 2.5 hours at room temperature or overnight at 4°C with gentle shaking.



Add 100 μ l prepared primary antibody to each well. Incubate 1 hour at room temperature with gentle shaking.



Add 100 μ l prepared 1X HRP-conjugated secondary antibody solution to corresponding wells. Incubate 1 hour at room temperature with gentle shaking.



Add 100 μ l TMB One-Step Substrate Reagent to each well. Incubate 30 minutes at room temperature.



Add 50 μ l Stop Solution to each well. Read at 450 nm immediately.

3. Kit Contents

- STAT3 Microplate (Item A): 96 wells (12 strips x 8 wells) coated with anti-pan STAT3 antibody.
- Wash Buffer Concentrate (20x) (Item B): 25 ml of 20x concentrated solution.
- Assay Diluent (Item E): 15 ml of 5x concentrated buffer. For diluting cell lysate sample, detection antibody (Item C-1 and Item C-2) and secondary antibody (Item D-1 and Item G) concentrates.
- Detection Antibody STAT3 (Tyr705) (Item C-1): 1 vial of anti-phospho-STAT3 (Tyr705) (each vial is enough to assay half microplate).
- Detection Antibody STAT3 (Item C-2): 1 vial of biotinylated anti-total-STAT3 (1 vial is enough to assay half microplate).
- HRP-conjugated Anti-rabbit IgG (Item D-1): 25 μ l of 2,000x HRP-conjugated anti-rabbit IgG concentrate.
- HRP-streptavidin concentrate (Item G): 200 μ l of 80 fold HRP-Streptavidin concentrate.
- TMB One-Step Substrate Reagent (Item H): 12 ml of 3,3',5,5'-tetramethylbenzidine (TMB) in buffered solution.

- Stop Solution (Item I): 8 ml of 0.2 M sulfuric acid.
- Cell Lysis Buffer (Item J): 5 ml 2x cell lysis buffer (not including protease and phosphatase inhibitors).
- Positive Control (Item K): 1 vial of lyophilized powder from A431 cell lysate.

4. Storage and Handling

Upon receipt, the kit should be stored at -20°C and used within 6 months of shipment date.

Return unused wells to the pouch containing desiccant pack, reseal along entire edge and return to -20°C .

After dilution, Wash Buffer Concentrate (Item B), Assay Diluent (Item E), TMB One-Step Substrate Reagent (Item H), Stop Solution (Item I) and Cell Lysis Buffer (Item J) can be stored at 4°C for up to 1 month.

Reconstituted Positive Control (Item K) should be stored at -80°C for up to 1 week.

After preparation, the detection Antibody STAT3 (Tyr705) (Item C-1) concentrate and Detection Antibody STAT3 (Item C-2) concentrate can be stored at 4°C for 5 days or at -80°C for one month.

After dilution, HRP-conjugated Anti-rabbit IgG (Item D-1) and HRP-streptavidin concentrate (Item G) working solutions should be used immediately and not re-used.

5. Additional Materials Required

- Microplate reader capable of measuring absorbance at 450 nm.
- Protease and Phosphatase inhibitors.
- Shaker.
- Precision pipettes to deliver 2 μ l to 1 ml volumes.
- Adjustable 1-25 ml pipettes for reagent preparation.
- 100 ml and 1 L graduated cylinders.
- Distilled or deionized water.
- Tubes to prepare sample dilutions.

6. Preparation of Samples

Cell lysates - Rinse cells with PBS, making sure to remove any remaining PBS before adding the Cell Lysis Buffer. Solubilize cells at 4×10^7 cells/ml in 1x Cell Lysis Buffer (we recommend adding protease and phosphatase inhibitors to Cell Lysis Buffer prior to sample preparation). Pipette up and down to resuspend and incubate the lysates with shaking at 2-8°C for 30 minutes. Microcentrifuge at 13,000 rpm for 10 minutes at 2-8°C, and transfer the supernatants into a clean test tube. Lysates should be used immediately or aliquoted and stored at -80°C. Avoid repeated freeze-thaw cycles. Thawed lysates should be kept on ice prior to use.

For the initial experiment, we recommend to do a serial dilution testing such as 5-fold and 50-fold dilution for your cell lysates with Assay Diluent (Item E) before use.

Note: The fold dilution of sample used depends on the abundance of phosphorylated proteins and should be determined empirically. More of the sample can be used if signals are too weak. If signals are too strong, the sample can be diluted further.

7. Preparation of Reagents

1. Bring all reagents and samples to room temperature (18-25°C) before use.
2. Assay Diluent (Item E) should be diluted 5-fold with deionized or distilled water before use.
3. Cell Lysis Buffer (Item J) should be diluted 2-fold with deionized or distilled water (for cell lysate and tissue lysate). We also recommend the addition of protease and phosphatase inhibitors (not included) to the Lysis Buffer prior to use.
4. Preparation of Positive Control (Item K): Briefly spin the Positive Control Vial. Add 400 μ L of prepared 1X Assay Diluent into Positive Control. Gently mix the powder to allow it to dissolve thoroughly. If a precipitate is seen in the solution after mixing, this can be removed by a quick centrifuge of the positive control vial and then pipetting the supernatant only for the assay.
5. If the Wash Concentrate (20x) (Item B) contains visible crystals, warm to room temperature and mix gently until dissolved. Dilute 20 ml of Wash Buffer Concentrate into deionized or distilled water to yield 400 ml of 1x Wash Buffer.

6. Preparation of Detection Antibodies:

- a. Preparation of Detection Antibody STAT3 (Tyr705) (Item C-1): Briefly spin the vial of rabbit anti-phospho-Stat3 (Tyr705). Add 100 μ L of 1X Assay Diluent into the vial to prepare a phospho detection antibody concentrate. Pipette up and down to mix gently (the concentrate can be stored at 4°C for 5 days or at -80°C for one month). The concentrate should then be diluted 55-fold with 1X Assay Diluent and used in step 5 of the Assay Procedure.
- b. Preparation of Detection Antibody STAT3 (Item C-2): Briefly spin the vial of biotinylated anti-pan-Stat3. Add 100 μ L of 1X Assay Diluent into the vial to prepare a detection antibody concentrate. Pipette up and down to mix gently (the concentrate can be stored at 4°C for 5 days or at -80°C for one month). The concentrate should then be diluted 55-fold with 1X Assay Diluent.

7. Preparation of HRP-conjugated concentrates:

- a. Preparation of HRP-conjugated Anti-rabbit IgG (Item D-1): Briefly spin the vial of HRP-conjugated anti-rabbit IgG (Item D-1)

before use. Prepare a 1:2000 working solution by diluting the HRP-conjugated anti-rabbit IgG concentrate in 1X Assay Diluent.

For example, add 5 μ L of HRP-conjugated anti rabbit IgG concentrate to 10 mL of 1 \times Assay Diluent and mix gently

- b. Preparation of HRP-streptavidin concentrate (Item G): Briefly spin the vial of HRP-Streptavidin concentrate before use. Prepare a 1:80 working solution by diluting HRP Streptavidin concentrate in 1X Assay Diluent

For example, add 100 μ L of HRP-conjugated anti rabbit IgG concentrate to 8 mL of 1 \times Assay Diluent and mix gently

8. Assay Method

1. Bring all reagents and samples to room temperature (18 - 25°C) before use. It is strongly recommended to run all positive controls and samples in at least duplicate. It is also recommended to run the positive controls in singlet for each of the pan and phospho-specific antibodies.

2. Label removable 8-well strips as appropriate for your experiment
3. Add 100 μL of positive control (Item K) (see Reagent Preparation step 4) or sample into appropriate wells. Cover the wells and incubate for 2.5 hours at room temperature or overnight at 4°C with gentle shaking.
4. Discard the solution and wash 4 times with 1X Wash Solution (Item B). Wash by filling each well with Wash Buffer (300 μL) using a multi-channel Pipette or autowasher. Complete removal of liquid at each step is essential for good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
5. Add 100 μL of prepared 1X Detection Antibody STAT3 (Tyr705) (Item C-1) (see Reagent Preparation step 6a) into the wells designated to detect phosphorylated protein. Add 100 μL of prepared 1X Detection Antibody STAT3 (Item C-2) (see Reagent Preparation step 6b) to the remaining wells to detect pan protein. Incubate for 1 hour at room temperature with gentle shaking.
6. Discard the solution. Repeat the wash as in step 4.
7. Add 100 μL of prepared HRP-conjugated Anti-rabbit IgG (Item D-1) (see Reagent Preparation step 7a) to the wells corresponding with Detection Antibody STAT3 (Tyr705) (Item

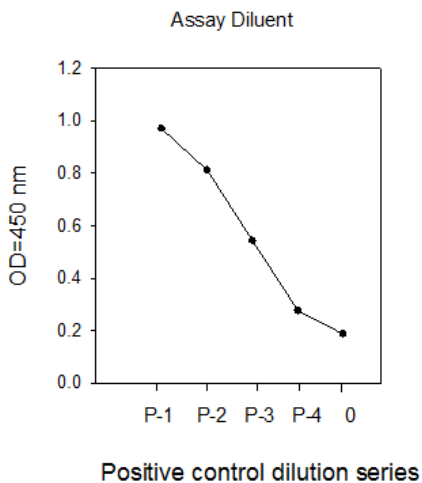
- C-1) in order to detect phosphorylated protein. To the remaining wells (corresponding with biotinylated anti-pan-Stat3), add 100 μ L of prepared HRP-Streptavidin solution (see Reagent Preparation step 7b) in order to detect pan protein. Incubate for 1 hour at room temperature with gentle shaking.
8. Discard the solution. Repeat the wash as in step 4.
 9. Add 100 μ L of TMB One-Step Substrate Reagent (Item H) to each well. Incubate for 30 minutes at room temperature in the dark with gentle shaking.
 10. Add 50 μ L of Stop Solution (Item I) to each well. Read at 450 nm immediately.

9. Data Analysis

ELISA data analysis: Calculate the mean absorbance for each sample. Then, subtract the average zero (blank) optical density from each sample mean and set of singlet positive controls.

Positive Control

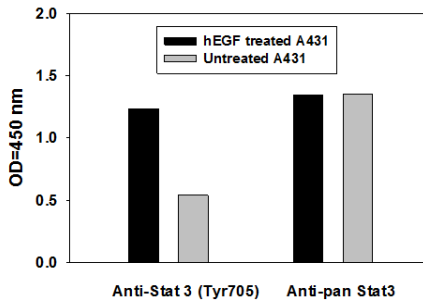
A431 cells were treated with recombinant human EGF at 37°C for 20 min. Solubilize cells at 4×10^7 cells/ml in Cell Lysis Buffer. Serial dilutions of lysates were analyzed in this ELISA.



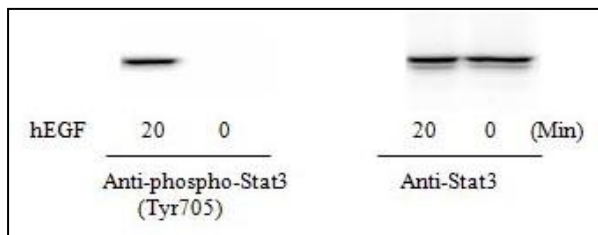
i. Recombinant Human EGF Stimulation of A431 Cell Lines

A431 cells were treated or untreated with 100 ng/ml recombinant human EGF for 20 min. Cell lysates were analyzed using this phosphoELISA and Western Blot.

a) ELISA



b) Western Blot Analysis



10. Troubleshooting

Problem	Cause	Solution
Sample signals	Too low: Sample concentration is too low.	Increasing sample concentration.
	Too high: Sample concentration is too high.	Reduce sample concentration.
Large CV	Inaccurate pipetting	Check pipettes
High background	Plate is insufficiently washed.	Review the manual for proper washing. If using an automated plate washer, check that all ports are unobstructed.
	Contaminated wash buffer.	Make fresh wash buffer.

Positive Control: Low signal	Improper storage of the ELISA kit.	Upon receipt, the kit should be stored at -20°C. Store the positive control at -70°C after reconstitution.
	Stop solution	Stop solution should be added to each well before measurement and read OD immediately.
	Improper primary or secondary antibody dilution.	Ensure correct dilution.

Technical Support

Copyright © 2026 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)