



ab126458 - STAT3 (pY705) ELISA Kit

Instructions for Use

For the quantitative measurement of human, mouse and rat phosphorylated STAT3 (Tyr705) 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	12
10.	Troubleshooting	14

1. Introduction

ab126458 is a very rapid, convenient and sensitive assay kit that can monitor the activation or function of important biological pathways in human, mouse and rat 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 Blotting analysis.

This Sandwich ELISA kit is an in vitro enzyme-linked immunosorbent assay for the measurement of human, mouse and rat phospho-STAT3 (Tyr705). An anti-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-STAT3 (Tyr705) antibody is used to detect phosphorylated STAT3 (Tyr705). After washing away unbound antibody, HRP-conjugated anti-rabbit IgG 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) 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 anti-rabbit IgG solution.

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 HRP-conjugated anti-rabbit IgG Concentrate (Item G).
- Detection Antibody STAT3 (Tyr705) (Item C-1): 2 vials of rabbit anti-STAT3 (Tyr705) (each 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.
- 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 .

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 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) working solution 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 $2 - 8^{\circ}\text{C}$ for 30 minutes. Microcentrifuge at 13,000 rpm for 10 minutes at $2 - 8^{\circ}\text{C}$, and transfer the supernatants into a clean test tube. Lysates should be used immediately or aliquoted and stored at -70°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 1x 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 Cell Lysis Buffer (Item J) 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 Antibody STAT3 (Tyr705) (Item C-1): Briefly spin the vial of Detection Antibody STAT3 (Tyr705) (Item C-1). Add 100 μ L of 1X Assay Diluent (Item E) 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.
7. 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 1 $\times$ 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

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.
2. Label removable 8-well strips as appropriate for your experiment.
3. Add 100 μ L 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. Wash by filling each well with Wash Buffer (300 μ L) 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 μ L of prepared 1X Detection Antibody STAT3 (Tyr705) (Item C-1) (see Reagent Preparation step 6) to each well. Incubate for 1 hour at room temperature with gentle shaking.
6. Discard the solution. Repeat the wash as in step 4.

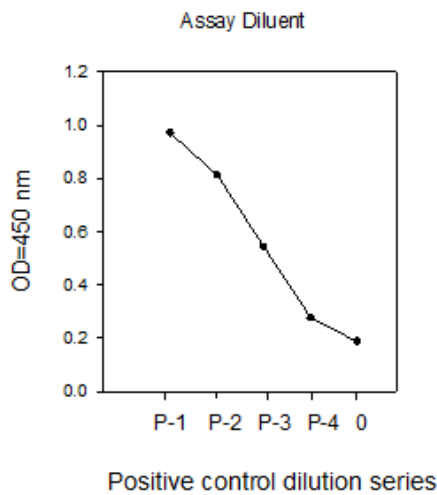
7. Add 100 μ L of prepared HRP-conjugated Anti-rabbit IgG (Item D-1) (see Reagent Preparation step 7) to each well. Incubate for 1 hour at room temperature with gentle shaking.
8. Discard the solution. Repeat the wash as in step 4.
9. Add 100 μ L of TMB One-Step Substrate Reagent (Item H) to each well. Incubate for 30 minutes at room temperature in the dark with 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: Average the duplicate readings for each sample or positive, then subtract the average zero (blank) optical density.

i. **Positive Control**

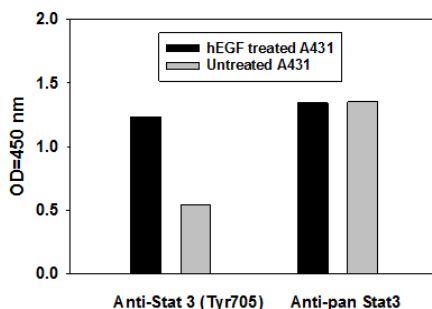
A431 cells were treated with recombinant human EGF at 37°C for 20 min. Cells were solubilized at 4×10^7 cells/ml in Cell Lysis Buffer. Serial dilutions of lysates were analyzed in this ELISA.



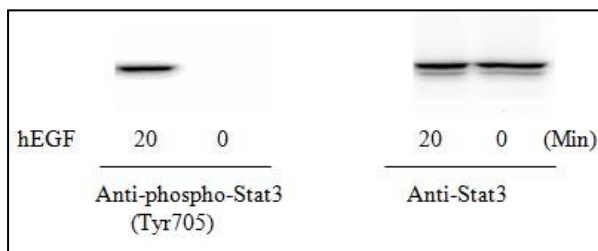
ii. Recombinant Human EGF Stimulation of A431 Cell Lines

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)