



EGFR (pY1086) (Human) ELISA Kit

Catalog Number KA2159

96 assays

Version: 01

Intended for research use only

www.abnova.com

I. INTRODUCTION

EGFR (pY1086) (Human) ELISA (Enzyme-Linked Immunosorbent Assay) kit 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 EGFR 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 human phospho-EGFR (Tyr 1086). An anti-EGFR antibody has been coated onto a 96-well plate. Samples are pipetted into the wells and phosphorylated and unphosphorylated EGFR present in a sample is bound to the wells by the immobilized antibody. The wells are washed and anti-phospho-EGFR (Tyr 1086) antibodies are used to detect phosphorylated EGFR. After washing away unbound antibody, HRP-conjugated anti-Rabbit IgG (secondary antibody) 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 EGFR bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

II. REAGENTS

1. EGFR Microplate (Item A): 96 wells (12 strips x 8 wells) coated with monoclonal anti-EGFR.
2. Wash Buffer Concentrate (20x) (Item B): 25 ml of 20x concentrated solution
3. Anti-phospho-EGFR (Tyr 1086) (Item C): 11µl rabbit anti-EGFR (Tyr 1086).
4. HRP-conjugated Anti-rabbit IgG (Item D), 25 µl HRPconjugated Anti-rabbit IgG.
5. Assay Diluent (Item E): 15 ml of 5x concentrated buffer. For diluting cell lysate sample and reagent (Item C and Item D).
6. TMB One-Step Substrate Reagent (Item H): 12 ml of 3,3',5,5'-tetramethylbenzidine (TMB) in buffered solution.
7. Stop Solution (Item I): 8 ml of 2 M sulfuric acid.
8. Cell Lysate Buffer (Item J): 5 ml 2x Cell Lysate Buffer (not including protease and phosphatase inhibitors).
9. Positive Control A431S001-1 (Item K): 1 vial of lyophilized powder from 100 µl cell lysate.

III. STORAGE

Upon receipt, the kit should be stored at -20°C. Please use within 6 months from the date of shipment. After initial use, Wash Buffer Concentrate (Item B), HRP-conjugated Anti-rabbit IgG (Item D), Assay Diluent (Item E), TMB One-Step Substrate Reagent (Item H), Stop Solution (Item I) and Cell Lysate Buffer (Item J) should be stored at 4°C to avoid repeated freeze-thaw cycles. Anti-phospho-EGFR (Tyr 1086) (Item C) should be stored at -20°C. Return unused wells to the pouch containing desiccant pack, reseal along entire edge and store at -20°C. Reconstituted Positive Control A4S001 (Item K) should be stored at -70°C.

IV. ADDITIONAL MATERIALS REQUIRED

1. Microplate reader capable of measuring absorbance at 450 nm.
2. Protease and Phosphatase inhibitors.
3. Shaker.
4. Precision pipettes to deliver 2 µl to 1 ml volumes.
5. Adjustable 1-25 ml pipettes for reagent preparation.
6. 100 ml and 1 liter graduated cylinders.
7. Distilled or deionized water.
8. Tubes to prepare sample dilutions.

V. SAMPLE PREPARATION

Cell lysates - Rinse cells with PBS, making sure to remove any remaining PBS before adding the Cell Lysate Buffer. Solubilize cells at 4×10^7 cells/ml in 1x Cell Lysate Buffer (we recommend adding protease and phosphatase inhibitors to Cell Lysate 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 supernates 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 100-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.

Cell Lysate Buffer should be diluted 2-fold with deionized or distilled water before use (recommend to add protease and phosphatase inhibitors).

VI. REAGENT PREPARATION

1. Bring all reagents and samples to room temperature (18 - 25 °C) before use.
2. Item E, Assay Diluent should be diluted 5-fold with deionized or distilled water before use.
3. Preparation of Positive Control: Briefly spin the Positive Control vial of Item K. Add 500 µl 1x Assay Diluent (Item E, Assay Diluent should be diluted 5-fold with deionized or distilled water before use) into Item K vial to prepare a Positive Control Stock Solution. Dissolve the powder thoroughly by a gentle mix. Add 50 µl prepared Positive Control Stock Solution from the vial of Item K, into a tube with 350 µl 1x Assay Diluent Buffer to prepare P-1 (See i. Positive control of part IX. TYPICAL DATA for a typical result). Pipette 240 µl 1x Assay Diluent into each tube. Use the Positive Control (1) to produce a dilution series (shown below). Mix each tube thoroughly before the next transfer. 1x Assay Diluent serves as the background.

50µl Positive Control

Stock Solution + 350 µl

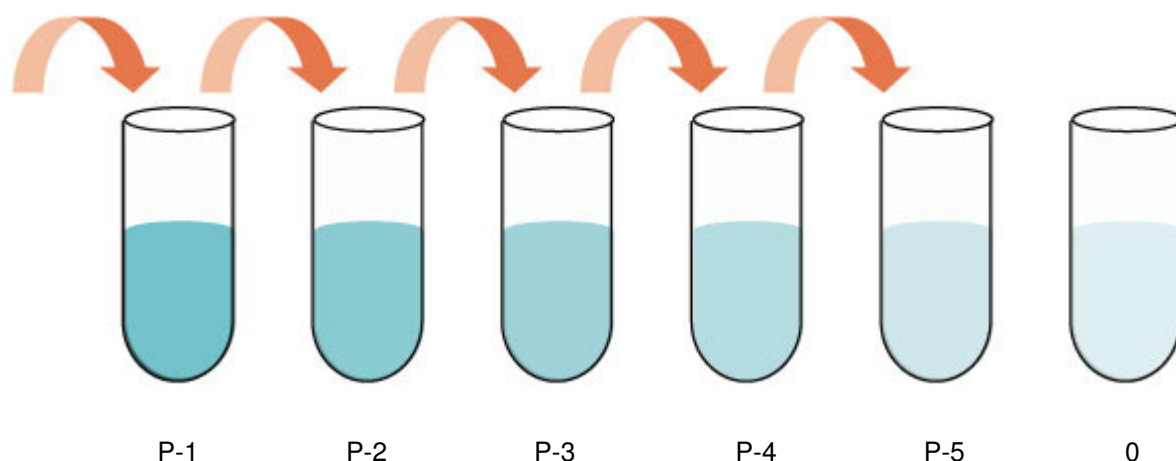
1x Assay Diluent

60 µl

60 µl

60 µl

60 µl








4. 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.
5. Briefly spin the anti-phospho-EGFR (Tyr 1086) (Item C) before use. Pipette up and down to mix gently. The antiphospho-EGFR (Tyr 1086) should be diluted 1,000-fold with 1x Assay Diuent. For example, add 5 µl anti-phospho-EGFR (Tyr 1086) into a tube with 5.0 ml 1x Assay Diluent to prepare a 1,000-fold diluted antibody.
6. Briefly spin the HRP-conjugated anti-rabbit IgG (Item D) before use. Pipette up and down to mix gently. HRP-conjugated antirabbit IgG concentrate should be diluted 1,000-fold with 1x Assay Diuent. For example: Briefly spin the vial (ItemD) and pipette up and down to mix gently. Add 5 µl of HRP-conjugated antirabbit IgG concentrate into a tube with 5.0 ml 1x Assay Diluent to prepare a 1,000-fold diluted HRP-conjugated anti-rabbit IgG solution.

7. Cell Lysate Buffer should be diluted 2-folds with deionized or distilled water before use (recommend to add protease and phosphatase inhibitors).

VII. ASSAY PROCEDURE

1. Bring all reagents to room temperature (18 - 25°C) before use. It is recommended that all samples or Positive Control should be run at least in duplicate.
2. Add 100 µl of each sample or positive control into appropriate wells. Cover well with plate holder and incubate for 2.5 hours at room temperature or over night at 4 °C with shaking.
3. Discard the solution and wash 4 times with 1x Wash Solution. 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 to 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.
4. Add 100 µl of prepared 1,000 fold diluted anti-phospho-EGFR (Tyr 1086) (Reagent Preparation step 6) to each well. Incubate for 1.5 hour at room temperature with shaking.
5. Discard the solution. Repeat the wash as in step 3.
6. Add 100 µl of prepared 1,000 fold diluted HRP-conjugated anti-rabbit IgG (see Reagent Preparation step 7) to each well. Incubate for 1 hour at room temperature with shaking.
7. Discard the solution. Repeat the wash as in step 3.
8. Add 100 µl of TMB One-Step Substrate Reagent (Item H) to each well. Incubate for 30 minutes at room temperature in the dark with shaking.
9. Add 50 µl of Stop Solution (Item I) to each well. Read at 450 nm immediately.

VIII. ASSAY PROCEDURE SUMMARY

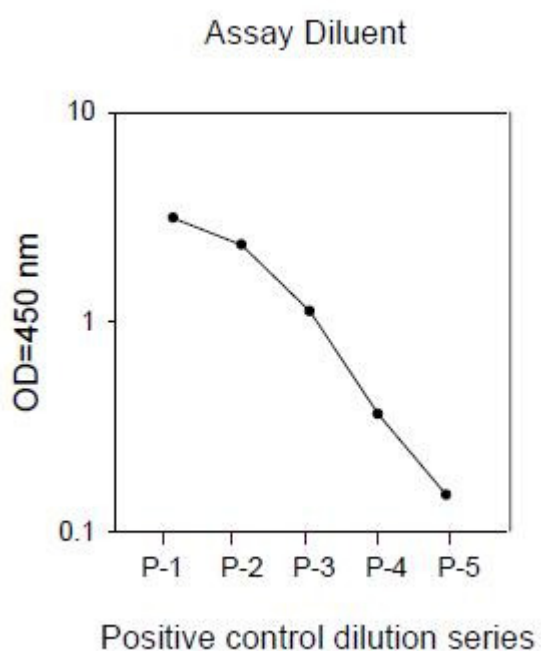
1. Prepare all reagents, samples and standards as instructed.

2. Add 100 µl sample or positive control to each well. Incubate 2.5 hours at room temperature or over night at 4 °C.

3. Add 100 µl prepared primary antibody to each well. Incubate 1.5 hours at room temperature.

4. Add 100 µl prepared secondary antibody solution. Incubate 1 hour at room temperature.

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

6. Add 50 µl Stop Solution to each well. Read at 450 nm immediately.

IX. TYPICAL DATA

ELISA data analysis: Average the duplicate readings for each sample or positive control then subtract the average blank optical density.

A. POSITIVE CONTROL

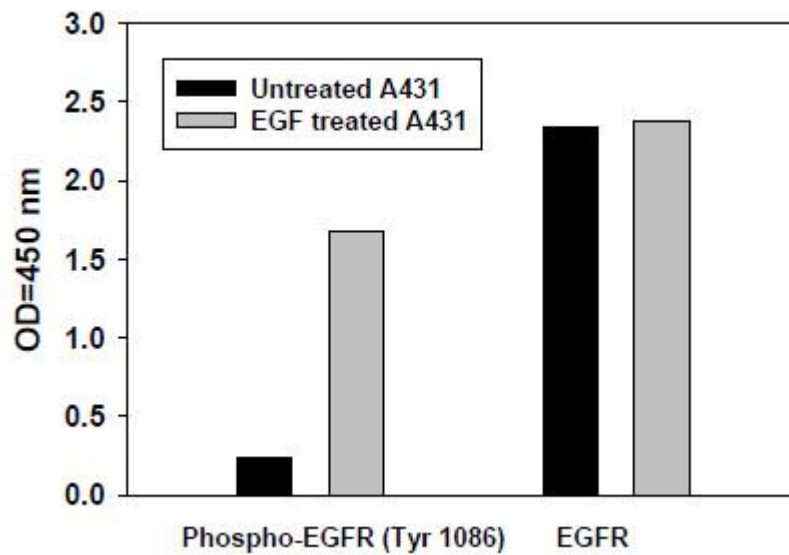
A431 cells were treated with recombinant human EGF at 37°C for 20 min. Solubilize cells at 4×10^7 cells/ml in Cell Lysate Buffer. Serial dilutions of lysates were analyzed in this ELISA. Please see step 3 of Part VI Reagent Preparation for detail.



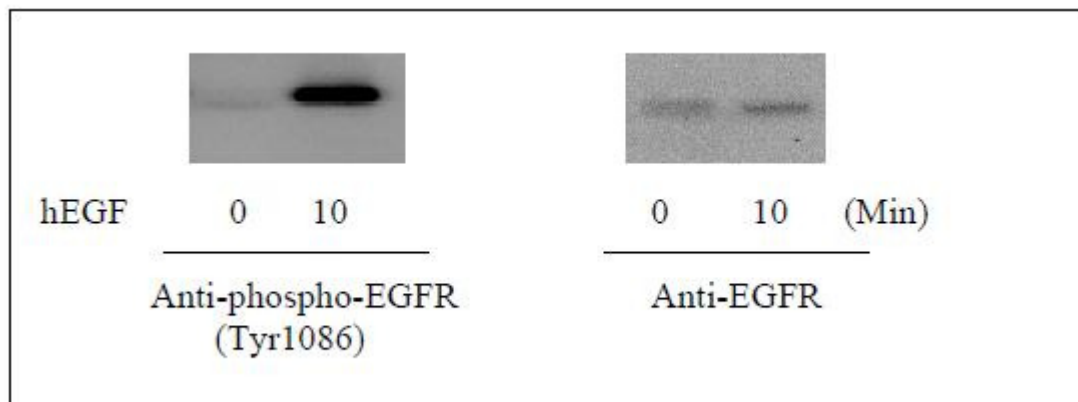
B. Recombinant Human EGF Stimulation of A431 Cell Lines

A431 cells were treated or untreated with 100 ng/ml recombinant human EGF for 10 min. Cell lysates were analyzed using this phosphoELISA and Western Blot.

a. ELISA



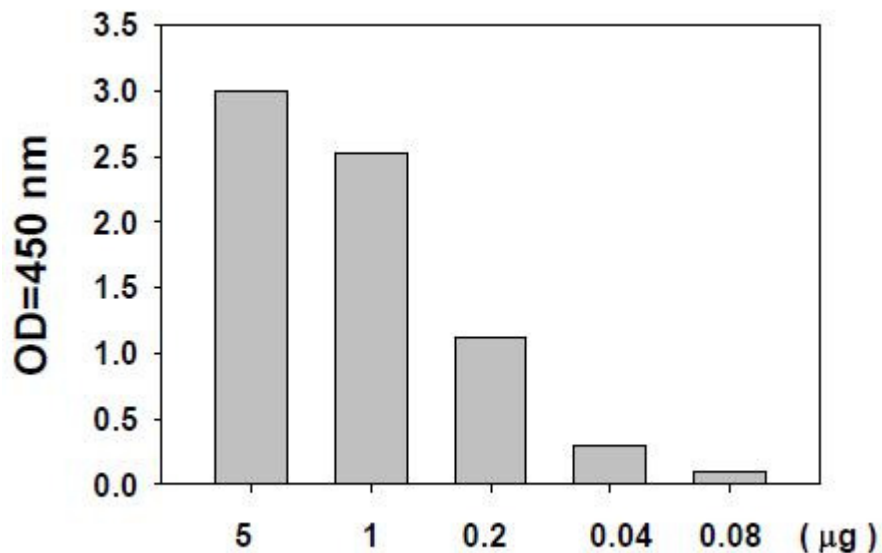
b. Western-Blot



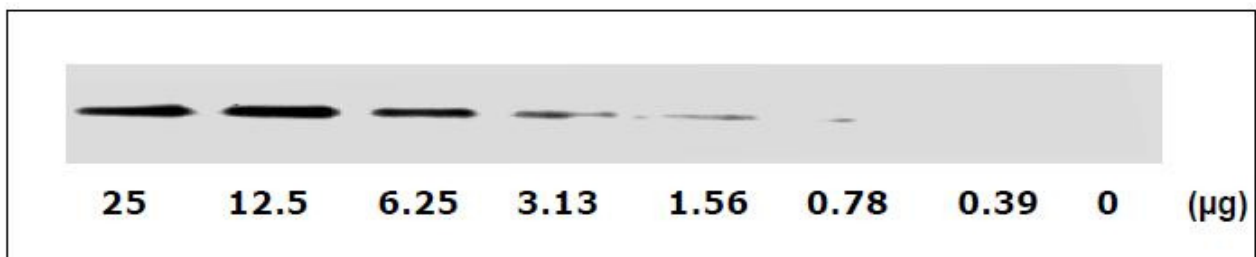
C. SENSITIVITY

The A431 cells were treated with 100 ng/mL recombinant human EGF for 20 minutes to induce phosphorylation of EGF R. Serial dilutions of lysates were analyzed in this ELISA and by Western blot. Immunoblots were incubated with anti-phospho-EGFR (Tyr 1086).

a. ELISA



b. Western-Blot



D. REFERENCES:

1. Hackel, P.O. et al. (1999) Curr. Opin. Cell Biol. 11, 184-189.
2. Alroy, I. and Y. Yarden (1997) FEBS Lett. 410:83.
3. Cooper, J.A. and Howell, B. (1993) Cell 73, 1051-1054.
4. Riedemann, J. et al. (2007) Biochem. Biophys. Res. Commun. **355**:707.

X. TROUBLESHOOTING GUIDE

Problem	Cause	Solution
1. Sample signals: a. Too low b. Too high	a. Sample concentration is too low b. Sample concentration is too high	a. Increasing sample concentration b. Reducing sample concentration
2. Large CV	a. Inaccurate pipetting	a. Check pipettes
3. High background	a. Plate is insufficiently washed b. Contaminated wash buffer	a. Review the manual for proper wash. If using an automated plate washer, check that all ports are unobstructed. b. Make fresh wash buffer
4. Positive Control: Low signal	a. Improper storage of the ELISA kit b. Stop solution c. Improper primary or secondary antibody dilution	a. Upon receipt, the kit should be stored at -20°C. Store the positive control at -70°C after reconstitution. b. Stop solution should be added to each well before measurement and read OD immediately. c. Ensure correct dilution