

Version 12b Last updated 16 August 2022

ERK1 (phospho T202 + Y204) + ERK2 (phospho T185 + Y187) + Total ELISA Kit

For the semi-quantitative measurement of ERK1/2 (pT202/Y204) and Total ERK1/2 in Human and mouse cell lysates.

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

On the 11th of December 2014, we removed the Antibody diluent (2 x 3 mL), Capture (phospho and total) antibody (2 x 300 µL), Detector (phospho and total) antibody (2 x 300 µL) and (1 vial) Positive control lysate and replaced with phospho and total capture antibody (2 x 1.5 mL) and phospho and total detector antibody (2 x 1.5 mL) prepared in 2 mL tubes.

Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	5
8. Technical Hints	6
9. Reagent Preparation	7
10. Standard Preparation	9
11. Sample Preparation	11
12. Plate preparation	14
13. Assay Procedure	15
14. Calculations	17
15. Typical Data	18
16. Typical Sample Values	20
17. Assay Specificity	23
18. Species Reactivity	23
19. Troubleshooting	24
20. Notes	25
Technical Support	30

1. Overview

ERK1 (phospho T202 + Y204) + ERK2 (phospho T185 + Y187) + Total ELISA kit is designed for the semi-quantitative measurement of ERK1/2 (pT202/Y204) and Total ERK1/2 protein in Human and mouse cells.

The SimpleStep ELISA™ employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. TMB substrate is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

ERK1 and ERK2, belong to a conserved family of serine/threonine protein kinases. ERK1/2 form part of a classical MAPK signaling cascade, consisting of a MAP kinase kinase kinase (MAPKKK), a MAP kinase kinase (MAPKK), and a MAP kinase (MAPK).

ERK1/2 is involved in cellular signaling events associated with a range of stimuli, including GPCRs, growth factors, and cytokines. The kinase activity of ERK proteins is regulated by dual phosphorylation at Thr202/Tyr204 in ERK1, and Threonine 185/Tyrosine 187 in ERK2. MEK1 and MEK2 are the primary upstream kinases responsible for ERK1/2 phosphorylation in this pathway. Several targets of ERK1/2 have been identified, including downstream kinases and transcription factors.

2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add standard or sample to appropriate wells.



Add Antibody Cocktail to all wells. Incubate at room temperature.



Aspirate and wash each well.



Add TMB Substrate to each well and incubate.



Add Stop Solution at a defined endpoint.

Alternatively, record color development kinetically after TMB substrate addition.

3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

4. Storage and Stability

Store kit at +4°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

Item	Quantity	Storage Condition
ERK1/2 (pT202/Y204) Capture Antibody	1.5 mL	+4°C
ERK1/2 (pT202/Y204) Detector Antibody	1.5 mL	+4°C
ERK1/2 (Total) Capture Antibody	1.5 mL	+4°C
ERK1/2 (Total) Detector Antibody	1.5 mL	+4°C
Lyophilized ERK1/2 Control Lysate	1 Vial	+4°C
10X Wash Buffer PT	15 mL	+4°C
5X Cell Extraction Buffer PTR	10 mL	+4°C
50X Cell Extraction Enhancer Solution	1 mL	+4°C
TMB Substrate	12 mL	+4°C

Stop Solution	12 mL	+4°C
Pre-Coated 96 Well Microplate (12 x 8 well strips)	96 Wells	+4°C
Plate Seal	1	+4°C

Control Lysate:

Prepared from A431 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 200ng/mL EGF for 10min.

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at 450 or 600 nm.
- Method for determining protein concentration (BCA assay recommended).
- Deionized water.
- PBS (1.4 mM KH_2PO_4 , 8 mM Na_2HPO_4 , 140 mM NaCl , 2.7 mM KCl , pH 7.4).
- Multi- and single-channel pipettes.
- Tubes for standard dilution.
- Plate shaker for all incubation steps.

8. Technical Hints

- This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.
- The provided 50X Cell Extraction Enhancer Solution may precipitate when stored at + 4°C. To dissolve, warm briefly at + 37°C and mix gently. The 50X Cell Extraction Enhancer Solution can be stored at room temperature to avoid precipitation.
- The provided 5X Cell Extraction Buffer contains phosphatase inhibitors. Protease inhibitors can be added if required.
- Selected components in this kit are supplied in surplus amount to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Make sure all buffers and solutions are at room temperature before starting the experiment.
- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Make sure you have the right type of plate for your detection method of choice.
- Make sure the heat block/water bath and microplate reader are switched on before starting the experiment.
- Complete removal of all solutions and buffers during wash steps is necessary to minimize background.

9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. **The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.**
- Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the formulations provided.

9.1 1X Cell Extraction Buffer PTR

Prepare 1X Cell Extraction Buffer PTR by diluting 5X Cell Extraction Buffer PTR and 50X Cell Extraction Enhancer Solution to 1X with deionized water. To make 10 mL 1X Cell Extraction Buffer PTR combine 7.8 mL deionized water, 2 mL 5X Cell Extraction Buffer PTR and 200 μ L 50X Cell Extraction Enhancer Solution Mix thoroughly and gently. If required protease inhibitors can be added.

Alternative – Enhancer may be added to Cell Extraction Buffer after extraction of cells or tissue. Refer to note in Section 19.

9.2 5X Cell Extraction/Enhancer Buffer PTR

Prepare 5X Cell Extraction/Enhancer Buffer PTR by adding 1/10th volume of 50X Cell Extraction Enhancer Solution. To prepare 1 mL, add 100 μ L of 50X Cell Extraction Enhancer Solution to 900 μ L of 5X Cell Extraction Buffer PTR. This concentrated mix is used for lysing cells directly in cell culture medium.

9.3 1X Wash Buffer

Prepare 1X Wash Buffer PT by diluting 10X Wash Buffer PT with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL 10X Wash Buffer PT with 45 mL deionized water. Mix thoroughly and gently.

9.4 Antibody Cocktail

Prepare Antibody Cocktails for phospho and total ERK1/2 separately, by combining an appropriate volume of the capture and detector antibodies immediately prior to assay. To make 3 mL of the Antibody Cocktail combine 1.5 mL Capture Antibody with 1.5 mL Detector Antibody. Mix thoroughly and gently.

10. Control Lysate Preparation

Kit Control lysates are provided at a concentration that give consistent signal between different lots. Lysates are produced and formulated by signal intensity to be consistent to within 30% of the previous lot. As such, Control lysates are not provided with a protein concentration.

- Always prepare a fresh set of positive controls for every use.
- Prepare serially diluted control lysates immediately prior to use.
- Discard working lysate dilutions after use as they do not store well.
- The following section describes the preparation of a lysate dilution series for duplicate measurements (recommended).

10.1 Reconstitute the Lyophilized ERK1/2 Control Lysate* by adding 250 μ L water. Mix thoroughly and gently. Hold at room temperature for 1 minute and mix gently. This is the 100% **Stock Lysate** Solution. Remaining stock material should be aliquoted and stored at -80 °C.

10.2 Label eight tubes, Controls 1– 8.

10.3 Add 300 μ L of 1X Cell Extraction Buffer PTR into tube number 1 and 100 μ L of 1X Cell Extraction Buffer PTR into numbers 2-8.

10.4 Use the Stock Lysate to prepare the following dilution series. Control #8 contains no protein and is the Blank control:

*Control lysates are supplied as a control reagent - not an absolute quantitation measure. A 3 - 4 point lysate dilution series is sufficient for this purpose.

Note: the extent of the dilution series appropriate for your samples needs to be determined empirically. Further dilutions than those shown below may be required.

Control #	Volume to dilute (μL)	Volume Diluent N (μL)	Lyophilized ERK1/2 (pT202/Y204) (%)
1	100 μL Stock Lysate	300	25
2	100 μL Control #1	100	12.5
3	100 μL Control #2	100	6.25
4	100 μL Control #3	100	3.125
5	100 μL Control #4	100	1.563
6	100 μL Control #5	100	0.781
7	100 μL Control #6	100	0.391
8 (Blank)	N/A	100	0

11. Sample Preparation

A cell density that yields 10,000 – 40,000 cells/well is suitable for the analysis of many cell lines. The lysis buffer volume should be adjusted so that lysates are in the range of 100-500 µg/mL of protein.

11.1 Preparation of extracts from adherent cells:

- 11.1.1 Remove growth media and rinse adherent cells 2 times in PBS.
- 11.1.2 Solubilize the cells by addition of chilled 1X Cell Extraction Buffer PTR directly to the plate (for cells cultured in 96-well microplates, lyse the cells with 100 µL* of 1X Cell Extraction Buffer PTR).
*Lysis volume should be adjusted depending on the desired lysate concentration. Lysates in the range of 100 - 500 µg/mL protein are usually sufficient.

11.2 Preparation of extracts from non-adherent cells:

- 11.2.1 Collect non-adherent cells by centrifugation and resuspend at an appropriate density in RPMI containing 10% FBS. Typical centrifugation conditions for cells are 500 x g for 5 minutes at RT.
- 11.2.2 Return cells to a 37°C incubator for 1 - 2 hours. For certain pathways, this can allow handling-mediated pathway activation to subside. This step is optional, and depends on the activation status of your cells following re-suspension.
- 11.2.3 At the completion of the cell treatment, harvest cells by centrifugation and lyse with 1X Cell Extraction Buffer PTR.*
- 11.2.4 Alternatively, in the absence of centrifugation cells may be lysed directly with a 20% volume of 5X Cell Extraction/Enhancer Buffer PTR (e.g. for 80 µL of cells, use 20 µL of 5X Cell Extraction/Enhancer Buffer PTR).
- 11.2.5 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.
- 11.2.6 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

*For best results, we recommend centrifugation and lysis of cells with 1X Cell Extraction Buffer PTR (11.2.3). Matrix effects may be observed in different cell media using the direct lysis approach (11.2.4). FBS may contain levels of ERK1/2 (Total) protein.

11.3 Preparations of extracts from tissue homogenates:

- 11.3.1 Tissue lysates are typically prepared by homogenization of tissue that is first minced and thoroughly rinsed in PBS to remove blood (dounce homogenizer recommended).
- 11.3.2 Homogenize 100 to 200 mg of wet tissue in 500 μ L - 1 mL of chilled 1X Cell Extraction Buffer PTR. For lower amounts of tissue adjust volumes accordingly.
- 11.3.3 Incubate on ice for 20 minutes.
- 11.3.4 Centrifuge at 18,000 x g for 20 minutes at 4°C.
- 11.3.5 Transfer the supernatants into clean tubes and discard the pellets.
- 11.3.6 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.
- 11.3.7 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

Refer to Dilution Guidelines for further instruction.

Guidelines for Dilutions of 100-fold or Greater <i>(for reference only; please follow the insert for specific dilution suggested)</i>	
100x	10000x
<p>4 μl sample + 396 μl buffer (100X) = 100-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 400 μl</i></p>	<p>A) 4 μl sample + 396 μl buffer (100X) B) 4 μl of A + 396 μl buffer (100X) = 10000-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 400 μl</i></p>
1000x	100000x
<p>A) 4 μl sample + 396 μl buffer (100X) B) 24 μl of A + 216 μl buffer (10X) = 1000-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 240 μl</i></p>	<p>A) 4 μl sample + 396 μl buffer (100X) B) 4 μl of A + 396 μl buffer (100X) C) 24 μl of A + 216 μl buffer (10X) = 100000-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 240 μl</i></p>

12. Plate preparation

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.
- For each assay performed, a minimum of two wells must be used as the zero control.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Differences in well absorbance or "edge effects" have not been observed with this assay.

13. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
 - We recommend that you assay all standards, controls and samples in duplicate.
 - To assay phospho and total ERK1/2, please use separate wells.
- 13.1** Prepare all reagents, working standards, and samples as directed in the previous sections.
 - 13.2** Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
 - 13.3** Add 50 µL of all samples and standards to appropriate wells.
 - 13.4** Add 50 µL of the corresponding Antibody Cocktail (phospho or total ERK1/2) to each well.
 - 13.5** Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
 - 13.6** Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and blot it against clean paper towels to remove excess liquid.
 - 13.7** Add 100 µL of TMB Substrate to each well and incubate for 15 minutes in the dark on a plate shaker set to 400 rpm.
 - 13.8** Add 100 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.

Alternative to 13.7 – 13.8: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

Mode:	Kinetic
Wavelength:	600 nm
Time:	up to 15 min
Interval:	20 sec - 1 min
Shaking:	Shake between readings

ΔNote: *An endpoint reading can also be recorded at the completion of the kinetic read by adding 100 μL Stop Solution to each well and recording the OD at 450nm.*

13.9 Analyze the data as described below.

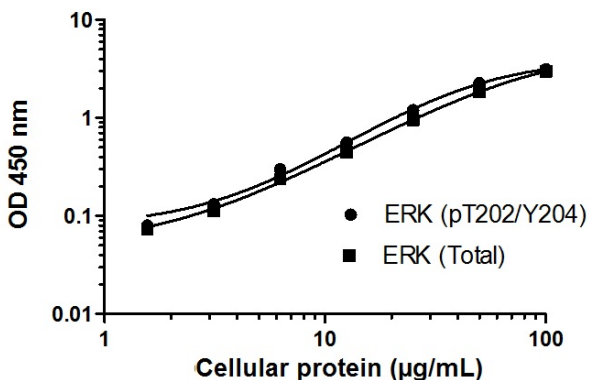
14. Calculations

Subtract average zero lysate control from all readings. Average the duplicate readings of the positive control dilutions.* If required, draw the best smooth curve through these points to fit a curve. Most plate reader software or graphing software can plot these values and curve fit. A four parameter algorithm (4PL) usually provides the best fit, though other equations can be examined to see which provides the most accurate (e.g. linear, semi-log, log/log, 4 parameter logistic).

*Control Lysates are supplied as a control reagent - not an absolute quantitation measure.

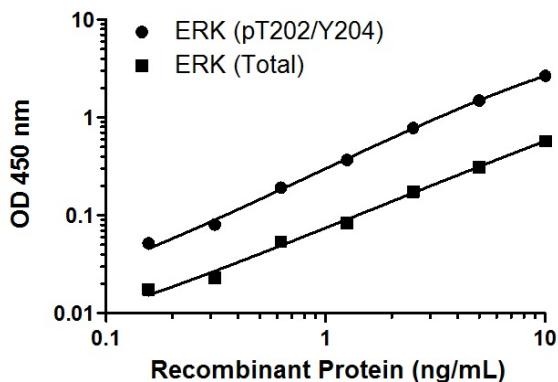
15. Typical Data

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Lysate Dilution Series Measurements						
Conc. (µg/mL)	ERK1/2 (pT202/Y204)			ERK1/2 (Total)		
	O.D. 450 nm		Mean O.D.	O.D. 450 nm		Mean O.D.
	1	2		1	2	
0.000	0.043	0.044	0.044	0.072	0.068	0.070
1.563	0.124	0.123	0.124	0.146	0.142	0.144
3.125	0.175	0.175	0.175	0.183	0.182	0.183
6.250	0.348	0.338	0.343	0.308	0.309	0.309
12.50	0.604	0.605	0.605	0.522	0.516	0.519
25.00	1.266	1.221	1.244	1.013	1.034	1.024
50.00	2.326	2.267	2.297	1.894	1.931	1.913
100.0	3.162	3.183	3.173	3.054	3.062	3.058

Figure 1. Example of a typical ERK1/2 (pT202/Y204) and ERK1/2 (Total) cell lysate dilution series. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Recombinant Protein Standard Curve Measurements						
Conc. (ng/mL)	ERK1/2 (pT202/Y204)			ERK1/2 (Total)		
	O.D. 450 nm		Mean O.D.	O.D. 450 nm		Mean O.D.
	1	2		1	2	
0.000	0.047	0.046	0.047	0.072	0.072	0.072
0.156	0.102	0.095	0.099	0.090	0.089	0.090
0.313	0.126	0.128	0.127	0.094	0.096	0.095
0.625	0.244	0.233	0.239	0.127	0.125	0.126
1.250	0.413	0.415	0.414	0.163	0.149	0.156
2.500	0.844	0.811	0.828	0.264	0.231	0.248
5.000	1.506	1.573	1.540	0.375	0.394	0.385
10.00	2.727	2.707	2.717	0.626	0.668	0.647

Figure 2. Example of a typical ERK1/2 (pT202/Y204) and ERK1/2 (Total) recombinant protein standard curve. The proportion of total protein that is phosphorylated is unknown - data is indicative only. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed

16. Typical Sample Values

SENSITIVITY –

The calculated minimal detectable (MDD) dose of ERK1/2 (pT202/Y204) and ERK1/2 (Total) recombinant protein is 100 pg/mL and 600 pg/mL respectively. The MDD was determined by calculating the mean of zero standard replicates (n=16) and adding 2 standard deviations then extrapolating the corresponding concentrations. Data for ERK1/2 (pT202/Y204) is indicative only – the proportion of total protein that is phosphorylated is unknown.

The minimal detectable (MDD) dose of ERK1/2 (pT202/Y204) and ERK1/2 (Total) is estimated in cellular extracts. As a guide, ERK1/2 (pT202/Y204) and ERK1/2 (Total) are detectable in activated A431 extracts with a total cellular protein concentration < 1.0 µg/mL (see Figure 1).

PRECISION –

Mean coefficient of variations of interpolated values from 3 concentrations of A431 extracts within the working range of the assay.

Target		Intra-Assay	Inter-Assay
ERK1/2 (pT202/Y204)	n=	6	3
	CV (%)	3.3	1.3
ERK1/2 (Total)	n=	6	3
	CV (%)	3.0	4.9

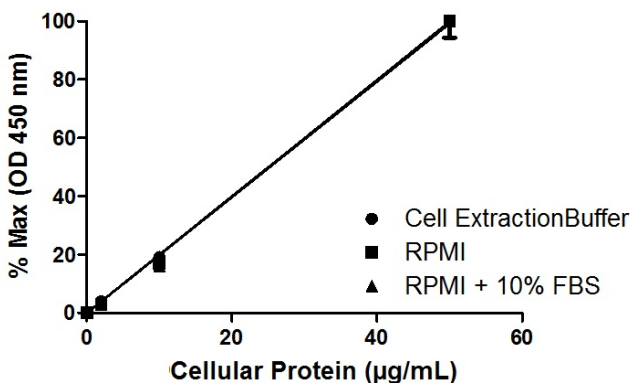


Figure 3: Linearity of dilution in representative sample matrices. Cellular lysates were prepared at 3 concentrations in common media containing 1X Cell Extraction Buffer PTR. Data from duplicate measurements of ERK1/2 (pT202/Y204) are normalized and plotted.

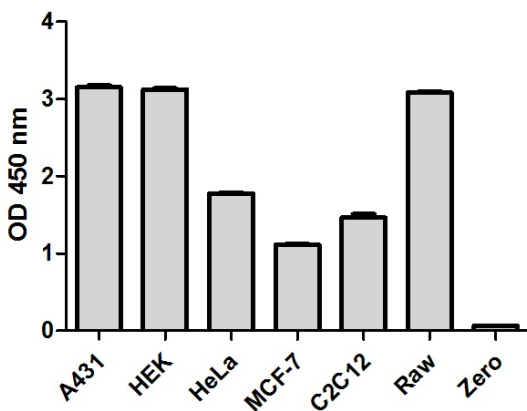


Figure 4: Cell line analysis for Total ERK1/2 from 100 µg/mL preparations of cell extracts. Data from triplicate measurements (mean +/- SD) are plotted and compared to 1X Cell Extraction Buffer PTR (zero).

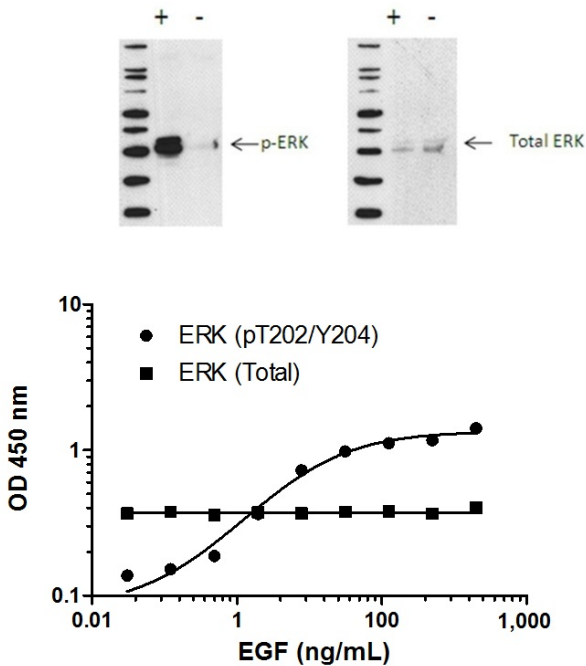


Figure 5: Induction of ERK1/2 (pT202/Y204) phosphorylation in MCF-7 cells in response to EGF treatment. MCF-7 cells were cultured in 96-well tissue culture plates, serum starved and treated (10 min) with a dose-range of EGF before cell lysis. Data from quadruplicate measurements of ERK1/2 (pT202/Y204) are plotted and compared against Total ERK1/2 protein levels. Comparative ERK1/2 (pT202/Y204) and ERK1/2 (Total) data also shown by Western Blot.

17. Assay Specificity

The ERK1/2 (pT202/Y204) assay detects endogenous levels of ERK1/2 (GenBank Accessions NP 002737.2 [ERK1], NP 620407 [ERK2]) in cellular lysates, only when phosphorylated at Thr202/Tyr204.

The ERK1/2 (Total) assay detects endogenous levels of ERK1/2 (GenBank Accessions NP 002737.2 [ERK1], NP 620407 [ERK2]) in cellular lysates. Total ERK1/2 assay kits detect ERK1/2, irrespective of phosphorylation status.

18. Species Reactivity

This kit detects ERK1/2 (pT202/Y204) and ERK1/2 (Total) in Human and mouse cell culture extracts. Detection in rat samples is also expected. Other species should be tested on a case-by-case basis.

Serum and plasma samples have not been tested with this kit.

Please contact our Technical Support team for more information.

19. Troubleshooting

Problem	Cause	Solution
Difficulty pipetting lysate; viscous lysate.	Genomic DNA solubilized	Prepare 1X Cell Extraction Buffer PTR (without enhancer). Add enhancer to lysate after extraction.
Poor standard curve	Inaccurate pipetting	Check pipettes
	Improper standard dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low signal	Incubation times too brief	Ensure sufficient incubation times; increase to 2 or 3 hour standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
	Incubation times with TMB too brief	Ensure sufficient incubation time until blue color develops prior addition of Stop solution
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions.
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the ELISA kit	Store your reconstituted standards at -80°C , all other assay components 4°C . Keep TMB substrate solution protected from light.

20. Notes

Technical Support

Copyright © 2022 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:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)