


NF kappaB p65 ELISA Kit ab176648

SimpleStep ELISA

22 References 6 Images

Overview

| Product name       | NF kappaB p65 ELISA Kit                                                                                                                                                                                                                                                                                                                                                 |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
|--------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|------|----|------|--------|---|------|----|-----|--------------|---|--|--|------|--------|---|------|----|-----|--------------|---|--|--|------|
| Detection method   | Colorimetric                                                                                                                                                                                                                                                                                                                                                            |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Precision          | <div>Intra-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>HeLa extract</td><td>6</td><td></td><td></td><td>2.8%</td></tr></table> <div>Inter-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>HeLa extract</td><td>3</td><td></td><td></td><td>8.7%</td></tr></table> |      |    |      | Sample | n | Mean | SD | CV% | HeLa extract | 6 |  |  | 2.8% | Sample | n | Mean | SD | CV% | HeLa extract | 3 |  |  | 8.7% |
| Sample             | n                                                                                                                                                                                                                                                                                                                                                                       | Mean | SD | CV%  |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| HeLa extract       | 6                                                                                                                                                                                                                                                                                                                                                                       |      |    | 2.8% |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Sample             | n                                                                                                                                                                                                                                                                                                                                                                       | Mean | SD | CV%  |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| HeLa extract       | 3                                                                                                                                                                                                                                                                                                                                                                       |      |    | 8.7% |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Sample type        | Cell Lysate, Tissue Homogenate                                                                                                                                                                                                                                                                                                                                          |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Assay type         | Semi-quantitative                                                                                                                                                                                                                                                                                                                                                       |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Sensitivity        | 5 µg/ml                                                                                                                                                                                                                                                                                                                                                                 |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Range              | 6 µg/ml - 600 µg/ml                                                                                                                                                                                                                                                                                                                                                     |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Assay time         | 1h 30m                                                                                                                                                                                                                                                                                                                                                                  |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Assay duration     | One step assay                                                                                                                                                                                                                                                                                                                                                          |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |
| Species reactivity | <b>Reacts with:</b> Mouse, Human<br><b>Predicted to work with:</b> Rat                                                                                                                                                                                                               |      |    |      |        |   |      |    |     |              |   |  |  |      |        |   |      |    |     |              |   |  |  |      |

**Product overview**

Abcam's NFκB p65 *in vitro* SimpleStep ELISA™ (Enzyme-Linked Immunosorbent Assay) kit is designed for the semi-quantitative measurement of NFκB p65 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.

**As of October 2019, this kit was reformulated with new antibodies to maintain continued long term supply.**

## Notes

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

## Platform

Microplate

## Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

| Components                                          | 1 x 96 tests | 1 x 96 tests |
|-----------------------------------------------------|--------------|--------------|
| 10X Wash Buffer PT                                  | 1 x 15ml     | 1 x 15ml     |
| 50X Cell Extraction Enhancer Solution               | 1 x 1ml      | 1 x 1ml      |
| 5X Cell Extraction Buffer PTR                       | 1 x 12ml     | 1 x 12ml     |
| Lyophilized NF kappaB p65 Control Lysate            | 1 vial       | 1 vial       |
| NF kappaB p65 (Total) Capture Antibody              | 1 x 3ml      | 1 x 3ml      |
| NF kappaB p65 (Total) Detector Antibody             | 1 x 3ml      | 1 x 3ml      |
| Plate Seal                                          | 1 unit       | 1 unit       |
| SimpleStep Pre-Coated 96-Well Microplate (ab206978) | 1 unit       | 1 unit       |
| Stop Solution                                       | 1 x 12ml     | 1 x 12ml     |
| TMB Substrate                                       | 1 x 12ml     | 1 x 12ml     |

## Function

NF-kappa-B is a pleiotropic transcription factor which is present in almost all cell types and is involved in many biological processes such as inflammation, immunity, differentiation, cell growth, tumorigenesis and apoptosis. NF-kappa-B is a homo- or heterodimeric complex formed by the Rel-like domain-containing proteins RELA/p65, RELB, NFkB1/p105, NFkB1/p50, REL and NFkB2/p52 and the heterodimeric p65-p50 complex appears to be most abundant one. The dimers bind at kappa-B sites in the DNA of their target genes and the individual dimers have distinct preferences for different kappa-B sites that they can bind with distinguishable affinity and specificity. Different dimer combinations act as transcriptional activators or repressors, respectively. NF-kappa-B is controlled by various mechanisms of post-translational modification and subcellular compartmentalization as well as by interactions with other cofactors or corepressors. NF-kappa-B complexes are held in the cytoplasm in an inactive state complexed with members of the NF-kappa-B inhibitor (I-kappa-B) family. In a conventional activation pathway,

I-kappa-B is phosphorylated by I-kappa-B kinases (IKKs) in response to different activators, subsequently degraded thus liberating the active NF-kappa-B complex which translocates to the nucleus. NF-kappa-B heterodimeric p65-p50 and p65-c-Rel complexes are transcriptional activators. The NF-kappa-B p65-p65 complex appears to be involved in invasin-mediated activation of IL-8 expression. The inhibitory effect of I-kappa-B upon NF-kappa-B the cytoplasm is exerted primarily through the interaction with p65. p65 shows a weak DNA-binding site which could contribute directly to DNA binding in the NF-kappa-B complex. Associates with chromatin at the NF-kappa-B promoter region via association with DDX1.

### Sequence similarities

Contains 1 RHD (Rel-like) domain.

### Domain

the 9aaTAD motif is a transactivation domain present in a large number of yeast and animal transcription factors.

### Post-translational modifications

Ubiquitinated, leading to its proteasomal degradation. Degradation is required for termination of NF-kappa-B response.

Monomethylated at Lys-310 by SETD6. Monomethylation at Lys-310 is recognized by the ANK repeats of EHMT1 and promotes the formation of repressed chromatin at target genes, leading to down-regulation of NF-kappa-B transcription factor activity. Phosphorylation at Ser-311 disrupts the interaction with EHMT1 without preventing monomethylation at Lys-310 and relieves the repression of target genes.

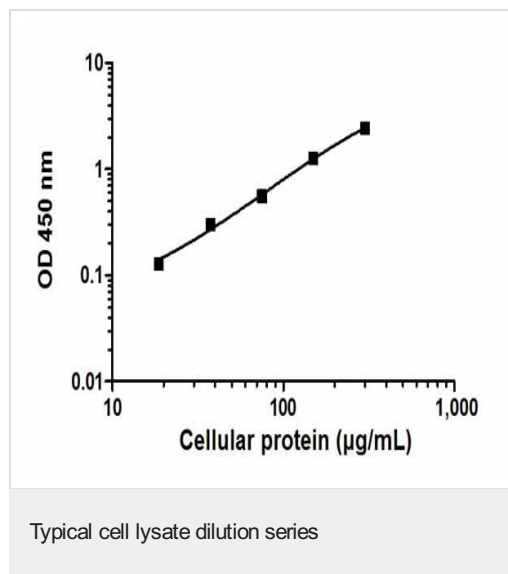
Phosphorylation at Ser-311 disrupts the interaction with EHMT1 and promotes transcription factor activity (By similarity). Phosphorylation on Ser-536 stimulates acetylation on Lys-310 and interaction with CBP; the phosphorylated and acetylated forms show enhanced transcriptional activity.

Reversibly acetylated; the acetylation seems to be mediated by CBP, the deacetylation by HDAC3. Acetylation at Lys-122 enhances DNA binding and impairs association with NFKBIA. Acetylation at Lys-310 is required for full transcriptional activity in the absence of effects on DNA binding and NFKBIA association. Acetylation can also lower DNA-binding and results in nuclear export. Interaction with BRMS1 promotes deacetylation of 'Lys-310'.

### Cellular localization

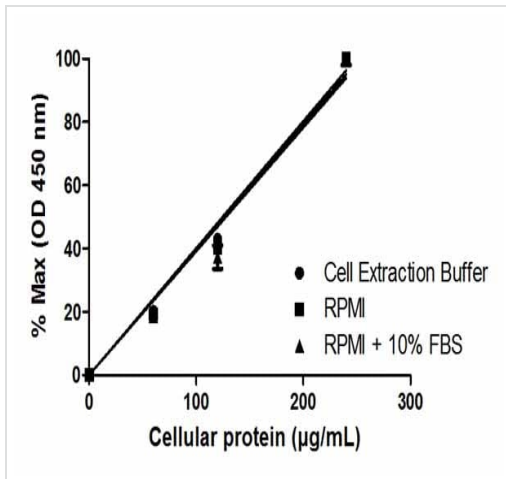
Nucleus. Cytoplasm. Nuclear, but also found in the cytoplasm in an inactive form complexed to an inhibitor (I-kappa-B). Colocalized with RELA in the nucleus upon TNF-alpha induction.

### Images



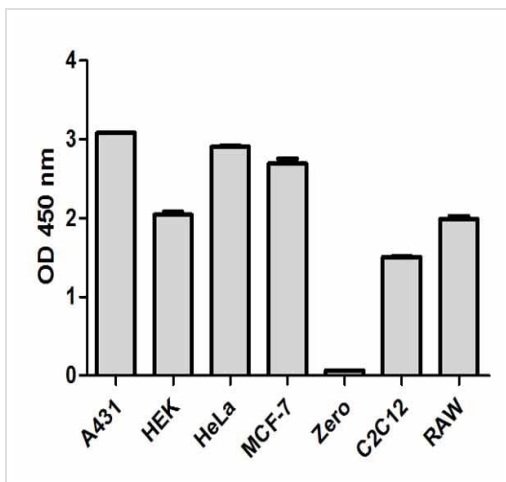
Example of a typical NFkB p65 cell lysate dilution series.

Background-subtracted data values (mean +/- SD) are graphed.



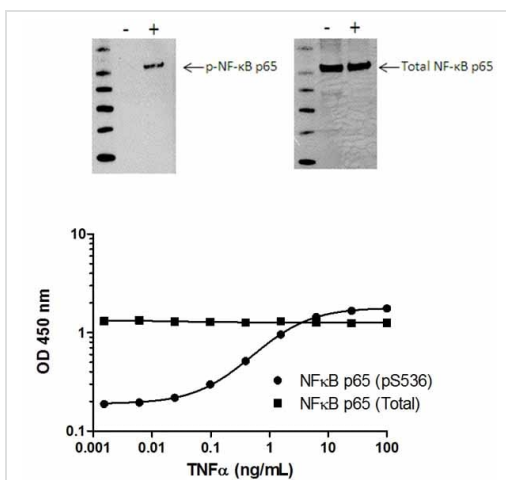
Linearity of dilution

Linearity of dilution in representative sample matrices. Cellular lysates were prepared at 3 concentrations in common media containing 1 x Cell Extraction Buffer PTR. Data from duplicate measurements of NFκB p65 (Total) are normalized and plotted.



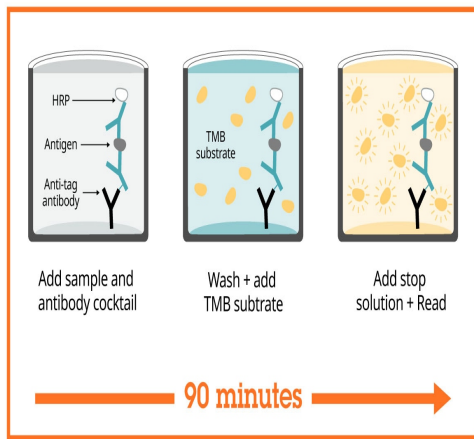
Comparison of NFκB p65 expression in different cell lines

Cell line analysis for Total NFκB p65 from 200 μg/mL preparations of cell extracts. Data from triplicate measurements (mean +/- SD) are plotted and compared to 1X Cell Extraction Buffer PTR (zero).



NFκB p65 phosphorylation in response to TNFα treatment

Induction of NFκB p65 (pS536) phosphorylation in MCF-7 cells in response to TNFα treatment. HeLa cells were cultured in 96-well tissue culture plates, and treated (10 min) with a dose-range of TNFα before cell lysis. Data from quadruplicate measurements of NFκB p65 (pS536) are plotted and compared against total NFκB p65 protein levels. Comparative NFκB p65 (pS536) and NFκB p65 (Total) data also shown by Western Blot.



Sandwich ELISA - NF kappaB p65 ELISA Kit  
(ab176648)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

#### Get more done with SimpleStep ELISA



**Easy to use**  
Single-wash 90-minute protocol



**Flexible**  
Matched antibody pairs available



**Precision antibodies**  
High sensitivity, specificity and reproducibility



**Scalable**  
Now in 10-pack and 384-well formats

Sandwich ELISA - NF kappaB p65 ELISA Kit  
(ab176648)

To learn more about the advantages of SimpleStep ELISA® kits see [here](#).

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

#### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors