



NFkB p50 ELISA Kit (Colorimetric)

Catalog Number TE-0002

(For Research Use Only)

Introduction

NF-kappaB (NFkB) proteins comprise a family of eukaryotic transcription factors that are involved in the control of a large number of cellular and organismal processes. In addition, these transcription factors are associated with many diseases including cancer and arthritis. NFkB commonly refers specifically to a p50-RelA(p65) heterodimer, which is the major Rel/NF-kB complex in most cells. P65-p65 and p50-p50 heterodimers have been demonstrated to bind on DNA as well. NF-kB is present as a latent, inactive, IkB-bound complex in the cytoplasm. When a cell receives any of a multitude of extracellular signals, NF-kB rapidly enters the nucleus and activates gene expression. Signosis developed the NFkB p50 ELISA kits for sensitive and specific analysis of the activities of NFkB in a high throughput way. The kit can be used for human, mouse and rat samples.

Principle of the assay

NFkB p50 ELISA kit is high sensitive and specific assay with a simple and optimized procedure. The 96-well (8X12 strip) clear plate is pre-immobilized with the NFkB consensus sequencing oligo. The activated NFkB in nuclear extract or the whole cell lysate is added in the well and binds to the oligo. The activated NFkB is detected with a specific antibody against p50 subunit and a HRP conjugated secondary antibody. The assay utilizes colorimetric detection method, which can be easily measured by spectrophotometry.

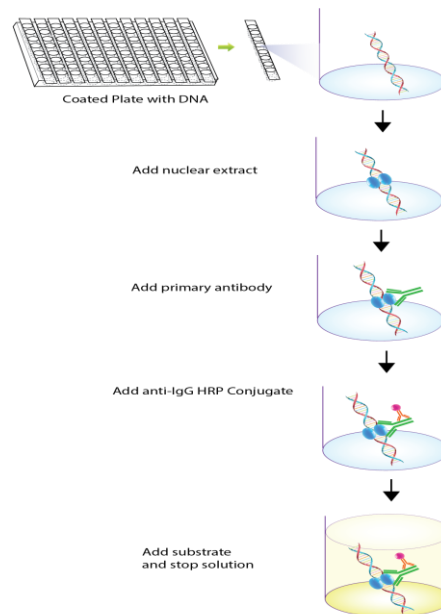


Diagram of TF ELISA

Materials provided with the kit

- 8x12 96-well microplate coated with NFkB consensus oligo (4°C).
- Antibody against NFkB p50 (4°C).
- HRP conjugate secondary antibody (4°C).
- 2X TF binding buffer (-20°C).
- 1X Nuclear extract dilution buffer (RT).
- NFkB p50 positive control (-20°C).
- 1X Diluent buffer (4°C).
- 5X Assay wash buffer (RT).
- Substrate (4°C).
- Stop Solution (4°C).

Material required but not provided

- Microplate reader capable of measuring absorbance at 450 nm
- Deionized or distilled water.

Reagent preparation before starting experiment

- Dilute the 5x Assay wash buffer to 1x buffer
40ml 5x Assay wash buffer
160ml ddH₂O
- Dilute 100 times of antibody against NFκB p50 with 1X Diluent buffer before use.
- Dilute 1000 times of HRP conjugate secondary antibody with 1X Diluent buffer before use.

Assay procedure

1. Calculate the number of samples to decide how many strips need to be used.
2. Make TF binding mix

30ul 2X TF binding buffer
X Nuclear extract (2-10ug)
X ul Nuclear extract dilution buffer
Total 60ul

-
- For the positive control, add 2ul positive control, and 28ul nuclear extract dilution buffer. Add the mix on a well and incubate with gentle shaking at room temperature for 1-2 hours.
3. Discard the contents and wash by adding 200μl of 1X Assay wash buffer. Repeat the process for a total of three washes. Complete removal of liquid at each wash. After the last wash, remove any remaining liquid by inverting the plate against clean paper towels.
 4. Add 60μl of diluted antibody against NFκB to each well and incubate for 2 hours at room temperature with gentle shaking, or 4°C overnight without shaking.
 5. Repeat the aspiration/wash as in step 4.
 6. Add 60 μl of diluted anti-mouse IgG HRP conjugate secondary antibody to each well and incubate for 45 minutes at room temperature with gentle shaking (DO NOT incubate longer than 45 mins to prevent from high background)
 7. Repeat the aspiration/wash as in step 4.
 8. Add 60μl of substrate to each well and incubate for 5-15 minutes or until positive wells turn blue.
 9. Add 30μl of stop solution to each well. The color in the wells should change from blue to yellow.
 10. Determine the optical density of each well with a microplate reader at 450 nm immediately.