



Human EGF ELISA

Catalog Number EA-0403

(For Research Use Only)

Introduction

Epidermal growth factor (EGF) is a strong mitogen for many cells, such as fibroblasts. EGF also stimulates the proliferation of embryonic cells. The proliferation of some cell lines has been shown to be inhibited by EGF. EGF acts as a differentiation factor for some cell types. To a limited extent EGF also augments angiogenesis because it is mitogenic for endothelial cells. The mitogenic activity of EGF for endothelial cells can be potentiated by thrombin. EGF is a strong chemoattractant for fibroblasts and epithelial cells. EGF alone and also in combination with other cytokines is an important factor mediating wound healing processes.

Principle of the assay

EGF ELISA is based on the principle of a solid phase enzyme-linked immunosorbent assay. The assay utilizes a mouse anti-human EGF antibody for immobilization on the microtiter wells and goat anti-human EGF antibody along with streptavidin conjugated to horseradish peroxidase (HRP) for detection. The test sample is allowed to react simultaneously with the two antibodies, resulting in the EGF molecules being sandwiched between the solid phase and enzyme-linked antibodies. After incubation, the wells are washed to remove unbound-labeled antibodies. A HRP substrate, TMB, is added to result in the development of a blue color. The color development is then stopped with the addition of Stop Solution changing the color to yellow. The concentration of EGF is directly proportional to the color intensity of the test sample. Absorbance is measured spectrophotometrically at 450 nm.

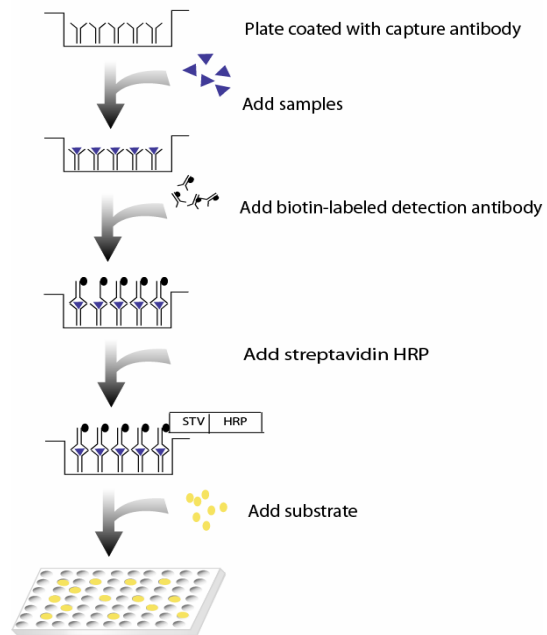


Diagram of ELISA

Materials provided with the kit

- 96 well microplate coated with a mouse anti-human EGF antibody (4°C).
- Biotin labeled goat anti-human EGF antibodies (9µg/ml) (-20°C).
- Streptavidin-HRP conjugate (4°C).
- Recombinant EGF standard (20ng/ml) (-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 200 times of human recombinant EGF (20ng/ml) with 1X Diluent buffer to 100pg/ml and then 2-fold serial dilutions.
- Dilute 400 times of biotin labeled goat anti-human EGF antibody with 1X Diluent buffer before use.
- Dilute 200 times of streptavidin-HRP with 1X Diluent buffer before use.

Assay procedure

1. Cut the sealing film over the plate and remove it from the desired number of well strips. Make sure the rest of wells are well sealed.
2. Add 100 μ l of Standard, control, or sample per well and incubate for 1 hour at room temperature with gentle shaking.
3. Aspirate each well and wash by adding 200 μ l of 1X Assay wash buffer. Repeat the process three times 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 100 μ l of diluted biotin-labeled goat anti-human EGF antibody to each well and incubate for 1 hour at room temperature with gentle shaking.
5. Repeat the aspiration/wash as in step 3.
6. Add 100 μ l of diluted streptavidin-HRP conjugate to each well and incubate for 45 min at room temperature with gentle shaking.
7. Repeat the aspiration/wash as in step 3.
8. Add 100 μ l substrate to each well and incubate for 5-30 minutes.
9. Add 50 μ l 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 within 30 minutes.

Example of standard curve

