

Fibronectin (human) ELISA Kit

rev. 6/14

(Catalog # K3631-100, 100 assays; Store at -20°C)

I. Introduction:

Fibronectin is a 430 KDa dimeric glycoprotein that exists in 2 forms - cellular and plasma fibronectin. BioVision's human Fibronectin ELISA Kit is based on the standard principle of a sandwich enzyme-linked immunosorbent assay. A rabbit polyclonal antibody specific for human Fibronectin is coated on a 96-well plate. Standards and test samples are added to the wells and Fibronectin present in a sample is bound by the immobilized antibody. A biotinylated polyclonal antibody from goat specific for Fibronectin is added subsequently. After washing away the unbound biotinylated antibody with PBS or TBS buffer, avidin-Biotin-Peroxidase Complex is added to the wells. The wells are again washed with PBS or TBS buffer to remove the unbound conjugates. HRP substrate TMB is used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue product that changes into yellow after adding acidic stop solution. The density of yellow color is proportional to the human Fibronectin captured onto the plate. This ELISA kit shows no cross-reactivity with other relevant proteins. Detection Range: 156 pg/ml – 10,000 pg/ml. Sensitivity: < 10 pg/ml.

II. Application:

Quantitative protein detection, establishing normal range etc.

III. Specificity:

Native and recombinant human Fibronectin

IV. Sample Type:

- Serum & plasma (heparin, EDTA, Citrate)
- Cell culture supernatants

V. Kit Contents:

Components	K3631-100	Part No.
Plate Coated with Fibronectin Ab	12 strips x 8 wells	K3631-100-1
Human Fibronectin Standard	10 ng x 2	K3631-100-2
Biotinylated anti-human Fibronectin Ab	130 µl	K3631-100-3
Avidin-Biotin-Peroxidase Complex	130 µl	K3631-100-4
Sample Diluent Buffer	30 ml	K3631-100-5
Antibody Diluent Buffer	12 ml	K3631-100-6
ABC Diluent Buffer	12 ml	K3631-100-7
TMB	10 ml	K3631-100-8
TMB Stop Solution	10 ml	K3631-100-9

VI. User Supplied Reagents and Equipment:

- Microplate reader capable of measuring absorbance at 450 nm.
- Absorbent paper.
- Adjustable pipettes and pipette tips. Multichannel pipettes are recommended for large number of samples.
- Washing buffer (neutral PBS or TBS).
 - Preparation of 0.01M TBS: Add 1.2 g Tris, 8.5 g NaCl; 450 µl of purified acetic acid or 700 µl of concentrated hydrochloric acid to 1000 ml H₂O and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.
 - Preparation of 0.01 M PBS: Add 8.5 g sodium chloride, 1.4 g Na₂HPO₄ and 0.2 g NaH₂PO₄ to 1000 ml distilled water and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

VII. Storage Conditions and Reagent Preparation:

Store kit at 4°C for 6 months or at -20°C for 12 months. Avoid repeated freeze-thaw cycles. Centrifuge tubes briefly to spin down all components to the bottom before opening.

- **Reconstitution of the human Fibronectin standard:** Two vials of Fibronectin standard (10 ng per vial) are included in each kit. Use one vial for each experiment. Prepare 10,000 pg/ml of human Fibronectin standard solution by adding 1ml of sample diluent buffer into one of the vials. Keep the tube at room temperature for 10 min. and mix thoroughly. Label 6 Eppendorf tubes with 5000 pg/ml, 2500 pg/ml, 1250 pg/ml, 625 pg/ml, 312 pg/ml & 156 pg/ml respectively. Aliquot 0.3 ml of the sample diluent buffer into each tube. Add 0.3 ml of the 10,000 pg/ml Fibronectin standard solution into 1st tube and mix. Transfer 0.3 ml from 1st tube to 2nd tube and mix. Transfer 0.3 ml from 2nd tube to 3rd tube and mix, and so on.

Note: The standard solutions are best used within 2 hrs. The 10 ng/ml standard solution should be stored at 4°C for up to 12 hrs, or at -20°C for up to 48 hrs. Avoid repeated freeze-thaw cycles.

- **Preparation of biotinylated anti-human Fibronectin antibody working solution:** Dilute 1:100 with the antibody diluent buffer and mix thoroughly. Prepare 0.1 ml of Fibronectin antibody working solution for each well. Solution should be prepared no more than 2 hrs. prior to the experiment.
- **Preparation of Avidin-Biotin-Peroxidase Complex (ABC) working solution:** Dilute 1:100 with the ABC dilution buffer and mix thoroughly. Prepare 0.1 ml of ABC working solution for each well. Solution should be prepared no more than 1 hr. prior to the experiment.

VIII. Sample Preparation and Storage:

Centrifuge cell culture supernatants to remove particulates, assay immediately or aliquot and store at -20°C. Allow the serum to clot in a serum separator tube (about 4 hrs) at room temperature. Centrifuge at approximately 1000 x g for 15 min. Analyze the serum immediately

or aliquot and store frozen at -20°C. Collect plasma using anticoagulants such as heparin, EDTA or Citrate. Centrifuge for 15 min. at 1000 x g within 30 min. of collection. Analyze immediately or aliquot and store frozen at -20°C.

Notes:

- Store samples to be assayed within 24 hrs. at 2-8°C. For long-term storage, aliquot and freeze samples at -20°C. Avoid repeated freeze-thaw cycles.
- Sample dilution guidelines: The user needs to estimate the concentration of the target protein in the sample and select a proper dilution factor so that the diluted target protein concentration falls near the middle of the linear regime of the standard curve. Dilute the sample using the provided diluent buffer. The sample must be well mixed with the diluent buffer. The following is a guideline for sample dilution. Several trials may be necessary to optimize sample dilution. For high target protein concentration (100-1000 ng/ml): dilute 1:100. For medium target protein concentration (10-100 ng/ml): dilute 1:10. For low target protein concentration (156-10,000 pg/ml): dilute 1:2. For very low target protein concentration (≤ 156 pg/ml). No dilution necessary or dilute 1:2.

IX. Assay Protocol:

The ABC working solution and TMB color developing agent must be kept warm at 37°C for 30 min. before use. When diluting samples and reagents, they must be mixed completely and evenly. The 96-well plate should not be dry at any time, as drying will inactivate the active components on the plate.

- Aliquot 0.1ml per well of the 10,000 pg/ml, 5000 pg/ml, 2500 pg/ml, 1250 pg/ml, 625 pg/ml, 312 pg/ml and 156 pg/ml human Fibronectin standard solutions into the precoated 96-well plate. Add 0.1ml of the sample diluent buffer into the control well (Zero well). Add 0.1ml each of the properly diluted samples of human cell culture supernates, serum or plasma (heparin, EDTA, citrate) to each empty well. See "Sample Dilution Guideline" for details.

Notes:

- We recommend that each human Fibronectin standard solution and each sample is measured in duplicate.
 - We recommend doing a pilot experiment using standards and a small number of samples to validate the assay procedure with the specific samples and optimize the appropriate sample dilution.
- Seal the plate with the cover and incubate at 37°C for 90 min. Remove the cover, discard plate contents, and blot the plate onto paper towels or other absorbent material. Do not let the wells completely dry at any time.
 - Add 0.1ml of biotinylated anti-human Fibronectin antibody working solution into each well and incubate the plate at 37°C for 60 min. Wash plate 3 times with 0.01M TBS or 0.01M PBS, and each time leave the washing buffer in the wells for 1 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material. (Plate Washing Method: Discard the solution in the plate without touching the side walls. Blot the plate onto paper towels or other absorbent material. Soak each well with at least 0.3 ml PBS or TBS buffer for 1~2 min. Repeat this process two additional times for a total of three washes. Note: For automated washing, aspirate all wells and wash three times with PBS or TBS buffer, overfilling wells with buffer each time. Blot the plate onto paper towels or other absorbent material.)
 - Add 0.1ml of prepared ABC working solution into each well and incubate the plate at 37°C for 30 min. Wash plate 5 times with 0.01M TBS or 0.01 M PBS, and each time leave washing buffer in the wells for 1-2 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material. (See Step 3 for plate washing method).
 - Add 90 μ l of prepared TMB color developing agent into each well and incubate plate at 37°C in dark for 20-25 min.
Note: These guidelines are for reference only; the optimal incubation time should be determined by end user. The shades of blue can be seen in the wells with the four most concentrated human Fibronectin standard solutions; the other wells show no obvious color.
 - Add 0.1ml of prepared TMB stop solution into each well. The color changes into yellow immediately.
 - Read absorbance at 450 nm in a microplate reader within 30 min. after adding the stop solution.
 - Calculation: Relative O.D.₄₅₀ = O.D.₄₅₀ of each well – O.D.₄₅₀ of Zero well. The standard curve can be plotted as the relative O.D.₄₅₀ of each standard solution (Y) vs. the respective concentration of the standard solution (X). The human Fibronectin concentration of the samples can be interpolated from the standard curve. **Note:** if the samples were diluted, multiply the dilution factor to the concentrations from interpolation to obtain the original concentration in the sample.

Typical Data Obtained from Human Fibronectin
(TMB reaction incubated at 37°C for 20 min.)

Concentration(pg/ml)	0	156	312	625	1250	2500	5000	10,000
O.D	0.056	0.134	0.193	0.336	0.593	0.904	1.596	2.042

Human Fibronectin ELISA Kit

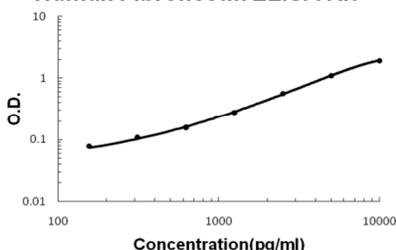


Figure: Standard Curve: This standard curves is for demonstration only. A standard curve must be run with each assay.

X. RELATED PRODUCTS:

FOR RESEARCH USE ONLY!



Fibronectin Antibody (3630-100)

Fibronectin (mouse) ELISA Kit (K3632)

Fibronectin (rat) ELISA Kit (K3633)

FOR RESEARCH USE ONLY! Not to be used on humans.