

Human β IG-H3 ELISA Kit

(Catalog # K926-100; 100 assays; Store at -20°C)

I. Introduction: BioVision's Human β IG-H3 (also known as Kerato-epithelin, RGDCAP) ELISA (Enzyme-Linked Immunosorbent Assay) kit is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of human β IG-H3 in serum, plasma, cell culture supernatants and urine. This assay employs an antibody specific for human β IG-H3 coated on a 96-well plate. Standards and samples are pipetted into the wells and β IG-H3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human β IG-H3 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of β IG-H3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm. The minimum detectable dose of β IG-H3 is typically less than 25 pg/ml. This ELISA kit shows no cross-reactivity with the $\text{m}\beta$ IGH3. The inter assay reproducibility CV is $<12\%$ and the intra assay CV is $<10\%$.

II. Sample Type: Serum, Plasma, Cell culture supernatants and urine.

III. Kit Contents:

Components	K926-100	Part No.
β IG-H3 Microplate (Item A): 96 wells coated with anti-human β IG-H3	12 stripsx8 wells	K926-100-1
Wash Buffer Concentrate (Item B) (20x)	25 ml	K926-100-2
Standards (Item C): recombinant human β IG-H3	2 vials	K926-100-3
Assay Diluent A (Item D): for Standard/Sample (serum/plasma) diluent. 0.09% sodium azide as preservative	30 ml	K926-100-4
Assay Diluent B [5X] (Item E): for Standard/Sample (cell culture medium/urine) diluent	15 ml	K926-100-5
Detection Antibody β IG-H3 (Item F): biotinylated anti-human β IG-H3 (each vial enough for half plate)	2 vial	K926-100-6
HRP-Streptavidin Concentrate [300X] (Item G)	200 μ l	K926-100-7
TMB One-Step Substrate Reagent (Item H): 3,3',5,5'-tetramethylbenzidine (TMB) in buffer solution	12 ml	K926-100-8
Stop Solution (Item I): 0.2 M sulfuric acid	8 ml	K926-100-9

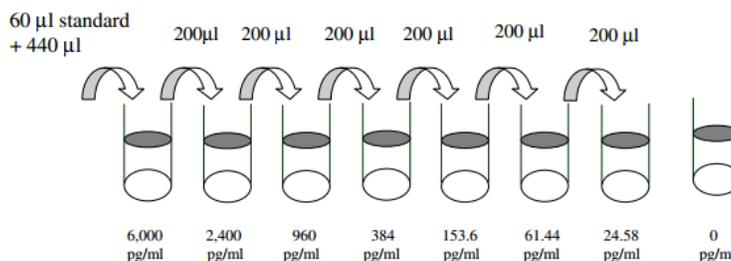
IV. User Supplied Reagents and Equipment:

- Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 2 μ l to 1 ml volumes.
- Adjustable 1-25 ml pipettes for reagent preparation.
- 100 ml and 1 liter graduated cylinders.
- Absorbent paper.
- Distilled or deionized water.
- Log-log graph paper or computer and software for ELISA data analysis.
- Tubes to prepare standard or sample dilutions.

V. Storage and Handling: May be stored for up to 6 months at 2° to 8°C from the date of shipment. Standard (recombinant protein) should be stored at -20°C or -80°C (recommended at -80°C) after reconstitution. Opened Microplate Wells or reagents may be stored for up to 1 month at 2° to 8°C . Return unused wells to the pouch containing desiccant pack, reseal along entire edge. Note: the kit can be used within one year if the whole kit is stored at -20°C . Avoid repeated freeze-thaw cycles.

VI. Reagent Preparation and Storage Conditions:

- Bring all reagents and samples to room temperature ($18 - 25^{\circ}\text{C}$) before use.
- Sample dilution: If your samples need to be diluted, Assay Diluent A (Item D) should be used for dilution of serum/plasma samples. 1x Assay Diluent B (Item E) should be used for dilution of culture supernatants and urine. Suggested dilution for normal serum/plasma: 400-4,000 fold. Please note that levels of the target protein may vary between different specimens. Optimal dilution factors for each sample must be determined by the investigator.
- Assay Diluent B should be diluted 5-fold with deionized or distilled water before use.
- Preparation of standard: Briefly spin the vial of Item C and then add 400 μ l Assay Diluent A (for serum/plasma samples) or 1x Assay Diluent B (for cell culture supernates/urine) into Item C vial to prepare a 50 ng/ml standard. Dissolve the powder thoroughly by a gentle mix. Add 60 μ l β IG-H3 standard (50 ng/ml) from the vial of Item C, into a tube with 440 μ l Assay Diluent A or 1x Assay Diluent B to prepare a 6,000 pg/ml standard solution. Pipette 300 μ l Assay Diluent A or 1x Assay Diluent B into each tube. Use the 6,000 pg/ml standard solution to produce a dilution series (shown below). Mix each tube thoroughly before the next transfer. Assay Diluent A or 1x Assay Diluent B serves as the zero standard (0 pg/ml).

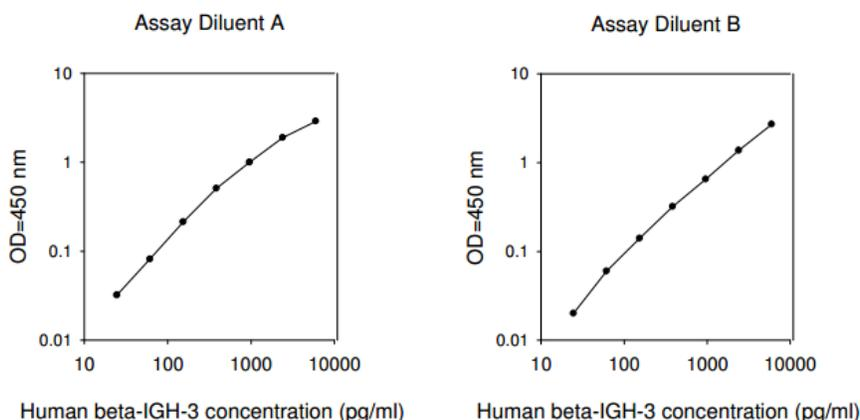


- If the Wash Concentrate (20x) (Item B) contains visible crystals, warm to room temperature and mix gently until dissolved. Dilute 20 ml of Wash Buffer Concentrate into deionized or distilled water to yield 400 ml of 1x Wash Buffer.
- Briefly spin the Detection Antibody vial (Item F) before use. Add 100 μ l of 1x Assay Diluent B into the vial to prepare a detection antibody concentrate. Pipette up and down to mix gently (the concentrate can be stored at 4°C for 5 days). The detection antibody concentrate should be diluted 80-fold with 1x Assay Diluent B and used in step 4 of Part VI Assay Procedure.

- Briefly spin the HRP-Streptavidin concentrate vial (Item G) and pipette up and down to mix gently before use. HRP-Streptavidin concentrate should be diluted 300-fold with 1x Assay Diluent B. For example: Briefly spin the vial (Item G) and pipette up and down to mix gently. Add 50 μ l of HRP-Streptavidin concentrate into a tube with 15 ml 1x Assay Diluent B to prepare a 300-fold diluted HRP-Streptavidin solution (don't store the diluted solution for next day use). Mix well.

VII. Assay Protocol:

- Bring all reagents and samples to room temperature (18 - 25°C) before use. It is recommended that all standards and samples be run at least in duplicate.
- Add 100 μ l of each standard (see Reagent Preparation step 2) and sample into appropriate wells. Cover well and incubate for 2.5 hours at room temperature or O/N at 4°C with gentle shaking.
- Discard the solution and wash 4 times with 1x Wash Solution. Wash by filling each well with Wash Buffer (300 μ l) using a multi-channel Pipette or autowasher. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
- Add 100 μ l of 1x prepared biotinylated antibody (Reagent Preparation step 6) to each well. Incubate for 1 hour at room temperature with gentle shaking. Discard the solution. Repeat the wash as in step 3.
- Add 100 μ l of prepared Streptavidin solution (see Reagent Preparation step 7) to each well. Incubate for 45 minutes at room temperature with gentle shaking. Discard the solution. Repeat the wash as in step 3.
- Add 100 μ l of TMB One-Step Substrate Reagent (Item H) to each well. Incubate for 30 minutes at room temperature in the dark with gentle shaking. Add 50 μ l of Stop Solution (Item I) to each well. Read at 450 nm immediately.
- Calculate the mean absorbance for each set of duplicate standards, controls and samples, and subtract the average zero standard optical density. Plot the standard curve on log-log graph paper or using Sigma plot software, with standard concentration on the x-axis and absorbance on the y-axis. Draw the best-fit straight line through the standard points.



- VIII. Recovery:** Recovery was determined by spiking various levels of human β IG-H3 into human serum, plasma and cell culture media. Mean recoveries are as in the table:

Sample type	Average % recovery	Range %
Serum	92.44	76-113
Plasma	86.91	69-104
Cell culture media	104.8	96-114

IX. Linearity:

Sample type	Serum	Plasma	Cell culture media
1:2 Average % of expected Range (%)	107.9 100-116	104.9 97-113	109.1 101-117
1:4 Average % of expected Range (%)	116.2 108-124	113.0 95-121	78.96 70-87

X. RELATED PRODUCTS:

TGF- β 1 Antibody
 Recombinant proteins
 Apoptosis Assay Kits and Reagents
 HDAC and HAT Assay Kits
 siRNA Expression Vectors
 Polyclonal and monoclonal antibodies

TGF- β 1 blocking peptide
 Antibodies and ELISA kits Metabolism Assay kits and Reagents
 Cellular Fractionation Kits
 DNA Damage, SOD Quantification Kits
 Recombinant Growth Factors and Cytokines

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