

Progranulin (Rat) ELISA Kit

(Catalog # K4735-100; 100 assays; Store at 4°C)

- I. Introduction:** Progranulin (PGRN) also called epithelin precursor, proepithelin (PEPI), PC cell-derived growth factor (PCDGF), acrogranin, or paraganulin is a 593aa cysteine-rich protein of 68.5kDa, that is typically secreted in a highly glycosylated 88kDa form. BioVision's Progranulin (rat) ELISA kit is a sandwich Enzyme Linked-Immunoassay (ELISA) for quantitative determination of rat progranulin in biological fluids. A polyclonal antibody specific for progranulin has been precoated onto the 96-well microtiter plate. Standards and samples are pipetted into the wells for binding to the coated antibody. After extensive washing to remove unbound compounds, progranulin is recognized by the addition of a biotinylated polyclonal antibody specific for progranulin (Detection Antibody). After removal of excess biotinylated antibody, HRP labeled streptavidin (Detector) is added. Following a final washing, peroxidase activity is quantified using the substrate 3,3',5,5'-tetramethylbenzidine (TMB). The intensity of the color reaction is measured at 450 nm after acidification and is directly proportional to the concentration of progranulin in the samples. The lowest level of progranulin that can be detected by this assay is 40 pg/ml. The Assay range is 0.063 ng/ml – 4 ng/ml. This ELISA is specific for the measurement of natural and recombinant rat progranulin. It does not cross-react with human progranulin, human granulin C, human granulin F, rat RBP4, rat adiponectin, rat Nampt, rat ANGPTL4, rat lipocalin-2, rat resistin, mouse RBP4, mouse adiponectin, human RBP4, human adiponectin.
- II. Sample Type:** Serum and Cell culture supernatants.
- III. Kit Contents:**

Components	K4735-100	Part No.
Plate coated with rat Progranulin antibody	6 stripsx16 wells	K4735-100-1
Wash Buffer (10x)	2x30 ml	K4735-100-2
Diluent (10X)	2x30 ml	K4735-100-3
Detection Antibody	20 µl	K4735-100-4
Detector (HRP Labeled Streptavidin) (lyophilized)	1 vial	K4735-100-5
rat progranulin Standard (lyophilized)	8 ng	K4735-100-6
TMB Substrate solution	12 ml	K4735-100-7
plate sealers (plastic film)	2	K4735-100-8
Stop Solution	12 ml	K4735-100-9

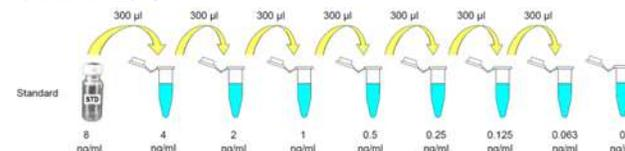
IV. User Supplied Reagents and Equipment:

- Microtiterplate reader at 450 nm, with the correction wavelength set at 540 nm or 570 nm
- Calibrated precision single and multi-channel pipettes. Disposable pipette tips
- Deionized water
- Microtubes or equivalent for preparing dilutions
- Disposable plastic containers for preparing working buffers
- Plate washer: automated or manual
- Glass or plastic tubes for diluting and aliquoting standard

- V. Storage and Handling:** Reagent must be stored at 2-8°C when not in use. Plate and reagents should be at room temperature before use. Do not expose reagents to temperatures greater than 25°C.

VI. Reagent Preparation and Storage Conditions:

- Bring all reagents and samples to room temperature (18 - 25°C) before use.
- Wash Buffer 10X has to be diluted with deionized water 1:10 before use (e.g. 50 ml Wash Buffer 10X + 450 ml water) to obtain Wash Buffer 1X.
- Diluent 10X has to be diluted with deionized water 1:10 before use (e.g. 20 ml Diluent 10X + 180 ml water) to obtain Diluent 1X.
- Detection Antibody has to be diluted 1:1000 in 1X Diluent (e.g.10 µl Detection Antibody + 9990 µl 1X Diluent). The diluted antibody cannot be stored.
- Detector (HRP Labeled Streptavidin) has to be reconstituted with 100 µl 1X Diluent. NOTE: After reconstitution, prepare aliquots and store at -20°C. Avoid freeze/thaw. Dilute the reconstituted Detector to the working concentration by adding 50 µl in 10 ml 1X Diluent (1:200 dilution). The diluted Detector is to be used within one hour of preparation.
- Rat Progranulin Standard (STD) has to be reconstituted with 1 ml of deionized water. This reconstitution produces a stock solution of 8 ng/ml. Mix the standard to ensure complete reconstitution and allow the standard to sit for a minimum of 15 minutes. Mix well prior to making dilutions. NOTE: The reconstituted standard is aliquoted and stored at -20°C
 - Dilute the standard protein concentrate (STD) (8 ng/ml) in Diluent 1X. A seven-point standard curve using 2-fold serial dilutions in Diluent 1X is recommended.
 - Suggested standard points are: 4, 2, 1, 0.5, 0.25, 0.125, 0.063 and 0 ng/ml.



VII. Assay Protocol:

- Determine the number of 16-well strips needed for the assay and insert them in the frame for current use. The extra strips should be resealed in the foil pouch bag and stored at 4°C. NOTE: Remaining 16-well strips coated with progranulin antibody when opened can be stored at 4°C for up to 1 month.
- Add 100 µl of the different standards into the appropriate wells in duplicate! At the same time, add 100 µl of appropriately diluted serum, plasma, urine, cell culture supernatant or CSF samples in duplicate to the wells.

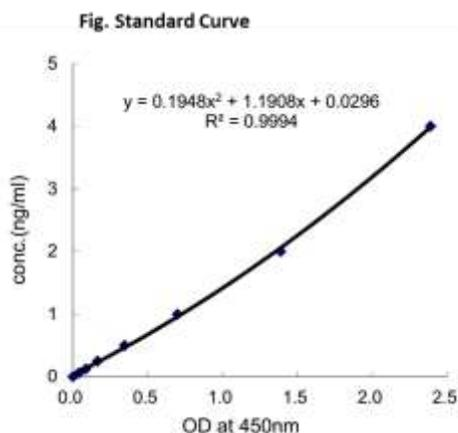
- Cover the plate with plate sealer and incubate for 1 hour at 37°C.
- Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of three washes. After the last wash, complete removal of liquid is essential for good performance.
- Add 100 µl to each well of the Detection Antibody.
- Cover the plate with plate sealer and incubate for 1 hour at 37°C.
- Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of three washes. After the last wash, complete removal of liquid is essential for good performance.
- Add 100 µl to each well of the diluted Detector.
- Cover the plate with plate sealer and incubate for 1 hour at 37°C.
- Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of five washes. After the last wash, complete removal of liquid is essential for good performance.
- Add 100 µl to each well of TMB Substrate Solution.
- Allow the color reaction to develop at room temperature (RT°C) in the dark for 10 minutes.
- Stop the reaction by adding 100 µl of Stop Solution. Tap the plate gently to ensure thorough mixing. The substrate reaction yields a blue solution that turns yellow when Stop Solution is added.
- Measure the OD at 450 nm in an ELISA reader within 30 minutes.

VIII. Calculation of Results:

- Average the duplicate readings for each standard, control and sample and subtract the average blank value (obtained with the 0 ng/ml point).
- Generate the standard curve by plotting the average absorbance obtained for each standard concentration on the horizontal (X) axis vs. the corresponding progranulin concentration (ng/ml) on the vertical (Y) axis (see Typical Data).
- Calculate the progranulin concentrations of samples by interpolation of the regression curve formula as shown above in a form of a quadratic equation.
- If the test samples were diluted, multiply the interpolated values by the dilution factor to calculate the concentration of human progranulin in the samples.

IX. Typical Data:

The following data are obtained using the different concentrations of standard as described in this protocol:



rProgranulin (ng/ml)	O.D.
4	2.389
2	1.388
1	0.699
0.5	0.346
0.25	0.165
0.125	0.087
0.063	0.042
0	0

X. RELATED PRODUCTS:

Progranulin, Human recombinant
 Progranulin, Mouse recombinant
 ELISA kits
 Apoptosis Assay Kits and Reagents
 Recombinant Growth Factors and Cytokines

Progranulin, Rat recombinant
 Recombinant proteins and enzymes
 Metabolism Assay kits and Reagents
 Cellular Fractionation Kits
 Polyclonal and monoclonal antibodies

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