



# B2M (Human) ELISA Kit

Catalog Number KA1045

96 assays

Version: 09

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Background .....	3
Principle of the Assay .....	3
<b>General Information .....</b>	<b>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	5
<b>Assay Protocol .....</b>	<b>6</b>
Reagent Preparation .....	6
Sample Preparation .....	6
Assay Procedure .....	7
<b>Data Analysis .....</b>	<b>9</b>
Calculation of Results .....	9
Performance Characteristics .....	10
<b>Resources .....</b>	<b>11</b>
References .....	11
Plate Layout .....	12

## **Introduction**

### **Background**

Beta-2-Microglobulin ( $\beta$ 2M) is a small serum protein that constitutes the light chain of the major histocompatibility class I human leukocyte antigen (HLA class I), an integral membrane protein involved in the immune response. The protein is 99 amino acid residues in length and has molecular mass of 12 kDa (1 - 4).  $\beta$ 2M is released from the cell surface of HLA class I into the serum and carried to the kidneys for degradation and secretion (5). In chronic renal failure,  $\beta$ 2M accumulates as insoluble amyloid aggregates and causes arthralgias, destructive osteoarthropathies, carpal tunnel syndrome and dialysis-related amyloidosis (6 - 8). Elevated serum  $\beta$ 2M levels are associated with poor prognosis in multiple myeloma and lymphoma (9 - 12).

### **Principle of the Assay**

The B2M (Human) ELISA Kit is designed for detection of human  $\beta$ 2M in plasma, serum, milk, saliva, urine, CSF and cell culture samples. This assay employs a quantitative sandwich enzyme immunoassay technique that measures  $\beta$ 2M in less than 4 hours. A polyclonal antibody specific for  $\beta$ 2M has been pre-coated onto a 96-well microplate with removable strips.  $\beta$ 2M in standards and samples is sandwiched by the immobilized antibody and biotinylated polyclonal antibody specific for  $\beta$ 2M, which is recognized by a streptavidin-peroxidase conjugate. All unbound material is then washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

## General Information

### Materials Supplied

List of component

Component	Amount
Human beta-2-Microglobulin Microplate: A 96-well polystyrene microplate coated with a polyclonal antibody against human $\beta$ 2M.	96 (8 x 12) wells
Sealing Tapes: Pre-cut, pressure-sensitive sealing tapes, which can be cut to fit the format of the individual assay.	3 slices
Human beta-2-Microglobulin Standard: Human $\beta$ 2M in a buffered protein base (lyophilized).	60 ng
Biotinylated Human beta-2-Microglobulin Antibody (50x): A 50-fold biotinylated polyclonal antibody against $\beta$ 2M.	140 $\mu$ L
MIX Diluent Concentrate (10x): A 10-fold concentrated buffered protein base.	30 mL
Wash Buffer Concentrate (20x): A 20-fold concentrated buffered surfactant.	30 mL x 2
Streptavidin-Peroxidase Conjugate (SP Conjugate): A 100-fold concentrate.	80 $\mu$ L
Chromogen Substrate: A ready-to-use stabilized peroxidase chromogen substrate tetramethylbenzidine.	8 mL
Stop Solution: A 0.5 N hydrochloric acid to stop the chromogen substrate reaction.	12 mL

### Storage Instruction

- ✓ Upon arrival, immediately store components of the kit at recommended temperatures up to the expiration date.
- ✓ Store SP Conjugate and Biotinylated Antibody at  $-20^{\circ}\text{C}$ .
- ✓ Store Microplate, Diluent Concentrate (10x), Wash Buffer, Stop Solution, and Chromogen Substrate at  $2-8^{\circ}\text{C}$ .
- ✓ Unused microplate wells may be returned to the foil pouch with the desiccant packs and resealed. May be stored for up to 30 days in a vacuum desiccator.
- ✓ Diluent (1x) may be stored for up to 30 days at  $2-8^{\circ}\text{C}$ .
- ✓ Store Standard at  $2-8^{\circ}\text{C}$  before reconstituting with Diluent and at  $-20^{\circ}\text{C}$  after reconstituting with diluent.

### Materials Required but Not Supplied

- ✓ Microplate reader capable of measuring absorbance at 450 nm.
- ✓ Pipettes (1-20  $\mu$ L, 20-200  $\mu$ L, 200-1000  $\mu$ L and multiple channel).
- ✓ Deionized or distilled reagent grade water.

**Precautions for Use**

- ✓ Prepare all reagents (working diluent buffer, wash buffer, standard, biotinylated antibody, and SP conjugate) as instructed, prior to running the assay.
- ✓ Prepare all samples prior to running the assay. The dilution factors for the samples are suggested in this protocol. However, the user should determine the optimal dilution factor.
- ✓ Spin down the SP conjugate vial and the biotinylated antibody vial before opening and using contents.
- ✓ This kit is for research use only.
- ✓ The kit should not be used beyond the expiration date.
- ✓ The Stop Solution is an acidic solution.

## Assay Protocol

### Reagent Preparation

- ✓ Freshly dilute all reagents and bring all reagents to room temperature before use.
- ✓ MIX Diluent Concentrate (10x): If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the MIX Diluent Concentrate 1:10 with reagent grade water. Store for up to 30 days at 2-8°C.
- ✓ Human beta-2-Microglobulin Standard: Reconstitute the 60 ng (4301 mU) of human beta-2-Microglobulin Standard with 1.2 mL of MIX Diluent to generate a 50 ng/mL (3584 mU/mL) standard solution. Allow the standard to sit for 10 minutes with gentle agitation prior to making dilutions. Prepare duplicate or triplicate standard points by serially diluting the Standard solution (50 ng/mL) 1:4 with MIX Diluent to produce 12.5, 3.125, 0.781, 0.195 and 0.049 ng/mL solutions. MIX Diluent serves as the zero standard (0 ng/mL). Any remaining solution should be frozen at -20°C and used within 30 days.

Standard Point	Dilution	[β2M] (ng/mL)	[β2M] (mU/mL)
P1	Standard (50 ng/mL)	50.00	3584
P2	1 part P1 + 3 part MIX Diluent	12.50	896
P3	1 part P2 + 3 part MIX Diluent	3.125	224
P4	1 part P3 + 3 part MIX Diluent	0.781	56
P5	1 part P4 + 3 part MIX Diluent	0.195	14
P6	1 part P5 + 3 part MIX Diluent	0.049	3.5
P7	MIX Diluent	0.000	0.0

- ✓ Biotinylated Human beta-2-Microglobulin Antibody (50x): Spin down the biotinylated antibody briefly and dilute the desired amount of the antibody 1:50 with MIX Diluent. Any remaining solution should be frozen at -20°C.
- ✓ Wash Buffer Concentrate (20x): If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the Wash Buffer Concentrate 1:20 with reagent grade water.
- ✓ SP Conjugate (100x): Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 1:100 with MIX Diluent. Any remaining solution should be frozen at -20°C.

### Sample Preparation

- ✓ Plasma: Collect plasma using one-tenth volume of 0.1 M sodium citrate as an anticoagulant. Centrifuge samples at 3000 x g for 10 minutes. Dilute samples 1:1000 into MIX Diluent and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles. (EDTA or Heparin can also be used as anticoagulant.)
- ✓ Serum: Samples should be collected into a serum separator tube. After clot formation, centrifuge

samples at 3000 x g for 10 minutes and remove serum. Dilute samples 1:1000 into MIX Diluent and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

- ✓ Cell Culture Supernatants: Centrifuge cell culture media at 3000 x g for 10 minutes to remove debris. Collect supernatants and assay. Store samples at -20°C or below. Avoid repeated freeze-thaw cycles.
- ✓ Urine: Collect urine using sample pot. Centrifuge samples at 800 x g for 10 minutes. Dilute samples 1:100 into MIX Diluent and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ Saliva: Collect saliva using sample tube. Centrifuge samples at 800 x g for 10 minutes. Dilute saliva 1:200 into MIX Diluent and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ Milk: Collect milk using sample tube. Centrifuge samples at 800 x g for 10 minutes. Dilute milk 1:4000 into MIX Diluent and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ CSF: Collect cerebrospinal fluid (CSF) using sample tube. Centrifuge samples at 3000 x g for 10 minutes. Dilute samples 1:1000 into MIX Diluent and assay. The undiluted samples can be stored at -80°C for up to 3 months. Avoid repeated freeze-thaw cycles.

### **Assay Procedure**

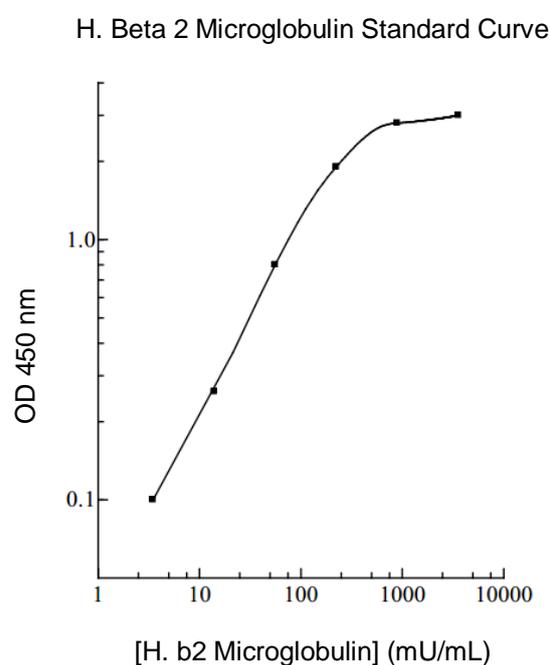
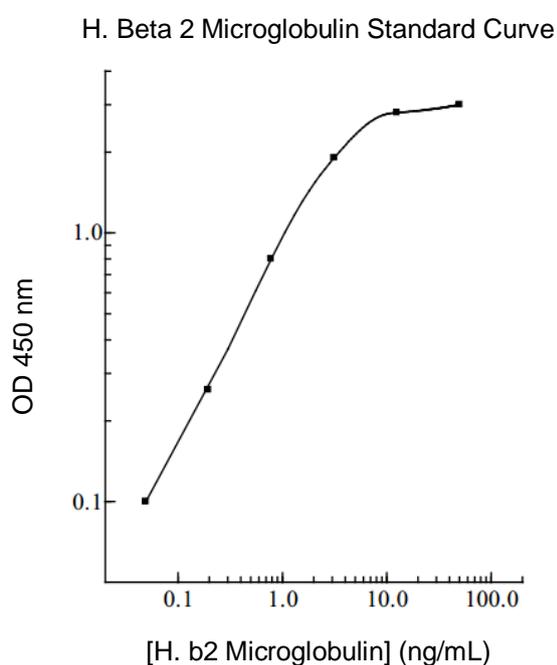
1. Prepare all reagents, standard solutions, and samples as instructed. Bring all reagents to room temperature before use. The assay is performed at room temperature (20-25°C).
2. Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccant inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
3. Add 50 µL of Human beta-2-Microglobulin Standard or sample per well. Cover wells with a sealing tape and incubate for 2 hour. Start the timer after the last sample addition.
4. Wash five times with 200 µL of Wash Buffer manually. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a machine wash six times with 300 µL of Wash Buffer and then invert the plate, decant the contents; hit 4-5 times on absorbent paper towel to completely remove the liquid.
5. Add 50 µL of Biotinylated Human beta-2-Microglobulin Antibody to each well and incubate for 1 hour.
6. Wash the microplate as described above.
7. Add 50 µL of Streptavidin-Peroxidase Conjugate per well and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
8. Wash the microplate as described above.
9. Add 50 µL of Chromogen Substrate per well and incubate for about 10 minutes or till the optimal blue color density develops. Gently tap plate to ensure thorough mixing and break the bubbles in the well with pipette tip.

10. Add 50  $\mu\text{L}$  of Stop Solution to each well. The color will change from blue to yellow.
11. Read the absorbance on a microplate reader at a wavelength of 450 nm immediately. If wavelength correction is available, subtract readings at 570 nm from those at 450 nm to correct optical imperfections. Otherwise, read the plate at 450 nm only. Please note that some unstable black particles may be generated at high concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

## Data Analysis

### Calculation of Results

- ✓ Calculate the mean value of the triplicate readings for each standard and sample.
- ✓ To generate a Standard Curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance on the y-axis. The best-fit line can be determined by regression analysis using log-log or four-parameter logistic curve-fit.
- ✓ Determine the unknown sample concentration from the Standard Curve and multiply the value by the dilution factor.



The curve is provided for illustration only. A standard curve should be generated each time the assay is performed.

### Performance Characteristics

- ✓ The minimum detectable dose of  $\beta$ 2M is typically ~ 0.04 ng/mL.
- ✓ Intra-assay and inter-assay coefficients of variation were 4.9 % and 7.2% respectively.
- ✓ Kit standard has been calibrated against WHO International Standard.

#### ✓ Linearity

Sample Dilution	Average Percentage of Expected Value		
	Plasma	Serum	Milk
1:1000	95%	97%	-
1:2000	99%	101%	94%
1:4000	101%	104%	98%
1:8000	-	-	104%

Sample Dilution	Urine	Saliva
1:50	97%	-
1:100	99%	85%
1:200	93%	95%
1:400	-	105%

#### ✓ Recovery

Standard Added Value	0.195 – 12.5 ng/mL
Recovery %	84-111 %
Average Recovery %	97 %

#### ✓ Cross-Reactivity

Species	% Cross Reactivity
Canine	None
Monkey	50%
Mouse	None
Rat	None
Swine	None
Bovine	None
Rabbit	None
Human	100%

#### ✓ Reference Value

Normal human  $\beta$ 2M plasma levels are < 2.7 ug/mL.

## Resources

### References

1. Gussow D *et al.* (1987) *J. Immunol.* 139:3132-3138
2. Krangel MS *et al.* (1979) *Cell* 18:979-991
3. Cattoni-Celli S *et al.* (1992) *Cancer Res.*1:1201-1204
4. Saper MA *et al.* (1991) *J. Mol. Biol.* 219:277-319
5. Trinh CH *et al.* (2002) *Proc. Natl. Acad. Sci. USA* 99:9771-9776
6. Dorthe BC *et al.* (2005) *Clinical Chemistry* 51:1177-1184
7. Eakin CM *et al.* (2006) *Nat. Struct. Mol. Biol.* 13:202-208
8. Drüeke TB (2000) *Nephrol. Dial. Transplant.* 15:17-24
9. Spati B. *et al.* (1980) *Acta Haematol.* 64:79-86
10. Bataille R *et al.* (1982) *Eur. J. Cancer Clin. Oncol.* 18:59-66
11. Norfolk D *et al.* (1979) *Br. J. Cancer* 39:510-515
12. Kantarjian HM *et al.* (1992) *Am. J. Med.* 93:599-604

**Plate Layout**

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H