

# ab137973

## Human Factor B

### ELISA Kit

#### Instructions for Use

For the quantitative measurement of Human Factor B in plasma, serum, saliva, milk, CSF, cell culture supernatants, cell lysates and tissue samples.

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

## INTRODUCTION

1. BACKGROUND	2
2. ASSAY SUMMARY	3

## GENERAL INFORMATION

3. PRECAUTIONS	4
4. STORAGE AND STABILITY	4
5. MATERIALS SUPPLIED	4
6. MATERIALS REQUIRED, NOT SUPPLIED	5
7. LIMITATIONS	5
8. TECHNICAL HINTS	6

## ASSAY PREPARATION

9. REAGENT PREPARATION	7
10. STANDARD PREPARATIONS	10
11. SAMPLE PREPARATION	13
12. PLATE PREPARATION	15

## ASSAY PROCEDURE

13. ASSAY PROCEDURE	16
---------------------	----

## DATA ANALYSIS

14. CALCULATIONS	18
15. TYPICAL DATA	18
16. TYPICAL SAMPLE VALUES	19
17. ASSAY SPECIFICITY	20

## RESOURCES

18. TROUBLESHOOTING	21
---------------------	----

## 1. BACKGROUND

Abcam's Factor B Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Human Factor B in plasma, serum, saliva, milk, CSF, cell culture supernatants, cell lysates and tissue samples.

A Factor B specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor B specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Factor B captured in plate.

Factor B (FB) is a component of the alternative pathway of complement activation. The zymogen circulates in the blood as a 93 kDa single chain glycoprotein with 739 amino acids. In the presence of C3b, it is cleaved by factor D into a 30 kDa N terminal noncatalytic Ba fragment and a 63 kDa C terminal catalytic Bb fragment. The active subunit Bb associates with C3b to form the alternative pathway C3 convertase. Human Factor B plays a major role in the initiation of the alternative pathway and in amplification of C3 cleavage. The polymorphism of Factor B influences C3 convertase formation, and is associated with age-related macular degeneration and polypoidal choroidal vasculopathy.

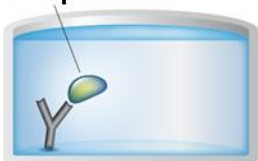
## 2. ASSAY SUMMARY

### Primary capture antibody



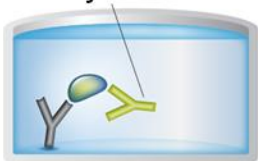
Prepare all reagents, samples and standards as instructed.

### Sample



Add standard or sample to each well used. Incubate at room temperature.

### Primary detector antibody



Wash and add prepared biotin antibody to each well. Incubate at room temperature.

### Streptavidin Label



Wash and add prepared Streptavidin-Peroxidase Conjugate. Incubate at room temperature.

### Substrate **Colored product**



Add Chromogen Substrate to each well. Incubate at room temperature. Add Stop Solution to each well. Read immediately.

## 3. PRECAUTIONS

Please read these instructions carefully prior to beginning the assay.

Modifications to the kit components or procedures may result in loss of performance.

## 4. STORAGE AND STABILITY

Store kit at 4°C immediately upon receipt, apart from the SP Conjugate & Biotinylated Antibody, which should be stored at -20°C.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in sections 9 & 10.

## 5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
Factor B Microplate (12 x 8 well strips)	96 wells	4°C
Factor B Standard	1 vial	4°C
10X Diluent M Concentrate	30 mL	4°C
Biotinylated Human Factor B Antibody	1 vial	-20°C
100X Streptavidin-Peroxidase Conjugate (SP Conjugate)	80 µL	-20°C
Chromogen Substrate	7 mL	4°C
Stop Solution	11 mL	4°C
20X Wash Buffer Concentrate	2 x 30 mL	4°C
Sealing Tapes	3	N/A

### 6. MATERIALS REQUIRED, NOT SUPPLIED

These materials are not included in the kit, but will be required to successfully utilize this assay:

- 1 Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 1  $\mu$ 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.
- 8 tubes to prepare standard or sample dilutions.

### 7. LIMITATIONS

- Do not mix or substitute reagents or materials from other kit lots or vendors.

### 8. TECHNICAL HINTS

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps.
- This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.

## 9. REAGENT PREPARATION

Equilibrate all reagents to room temperature (18-25°C) prior to use. Prepare fresh reagents immediately prior to use. If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved.

### 9.1 1X Diluent M

Dilute the 10X Diluent M Concentrate 10-fold with reagent grade water. Mix gently and thoroughly. *Store for up to 1 month at 4°C.*

### 9.2 1X Wash Buffer

Dilute the 20X Wash Buffer Concentrate 20-fold with reagent grade water. Mix gently and thoroughly.

### 9.3 1X Biotinylated Factor B Detector Antibody

9.3.1 The stock Biotinylated Factor B Antibody must be diluted with 1X Diluent M according to the label concentration to prepare 1X Biotinylated Factor B Antibody for use in the assay procedure. Observe the label for the "X" concentration on the vial of Biotinylated Factor B Antibody.

9.3.2 Calculate the necessary amount of 1X Diluent M to dilute the Biotinylated Factor B Antibody to prepare a 1X Biotinylated Factor B Antibody solution for use in the assay procedure according to how many wells you wish to use and the following calculation:

Number of Wells Strips	Number of Wells	(V <sub>T</sub> ) Total Volume of 1X Biotinylated Detector Antibody (μL)
4	32	1,760
6	48	2,640
8	64	3,520
10	80	4,400
12	96	5,280



*The undiluted antibody should be stored at -20°C.*

Where:

$C_S$  = Starting concentration (X) of stock Biotinylated Factor B Antibody (variable)

$C_F$  = Final concentration (always = 1X) of 1X Biotinylated Factor B Antibody solution for the assay procedure

$V_T$  = Total required volume of 1X Biotinylated Factor B Antibody solution for the assay procedure

$V_A$  = Total volume of (X) stock Biotinylated Factor B Antibody

$V_D$  = Total volume of 1X Diluent M required to dilute (X) stock Biotinylated Factor B Antibody to prepare 1X Biotinylated Antibody solution for assay procedures

Calculate the volume of (X) stock Biotinylated Antibody required for the given number of desired wells:

$$(C_F / C_S) \times V_T = V_A$$

Calculate the final volume of 1X Diluent M required to prepare the 1X Biotinylated Factor B Antibody:

$$V_T - V_A = V_D$$

Example:

**NOTE: This example is for demonstration purposes only. Please remember to check your antibody vial for the actual concentration of antibody provided.**

$C_S$  = 50X Biotinylated Factor B Antibody stock

$C_F$  = 1X Biotinylated Factor B Antibody solution for use in the assay procedure

$V_T$  = 3,520  $\mu$ L (8 well strips or 64 wells)

$$(1X/50X) \times 3,520 \mu\text{L} = 70.4 \mu\text{L}$$

$$3,520 \mu\text{L} - 70.4 \mu\text{L} = 3,449.6 \mu\text{L}$$

$V_A$  = 70.4  $\mu$ L total volume of (X) stock Biotinylated Factor B Antibody required

$V_D = 3,449.6 \mu\text{L}$  total volume of 1X Diluent M required to dilute the 50X stock Biotinylated Antibody to prepare 1X Biotinylated Factor B Antibody solution for assay procedures

9.3.3 First spin the Biotinylated Factor B Antibody vial to collect the contents at the bottom.

9.3.4 Add calculated amount  $V_A$  of stock Biotinylated Factor B Antibody to the calculated amount  $V_D$  of 1X Diluent M. Mix gently and thoroughly.

### 9.4 1X SP Conjugate

Spin down the 100X Streptavidin-Peroxidase Conjugate (SP Conjugate) briefly and dilute the desired amount of the conjugate 100-fold with 1X Diluent M.

*The undiluted conjugate should be stored at  $-20^\circ\text{C}$*

## 10. STANDARD PREPARATIONS

- Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use.
- Any remaining standard should be stored at -20°C after reconstitution and used within 5 days.
- This procedure prepares sufficient standard dilutions for duplicate wells.

10.1 Reconstitution of the Factor B Standard vial to prepare a 140 ng/mL Factor B **Standard #1**:

10.1.1 First consult the Factor B Standard vial to determine the mass of protein in the vial.

10.1.2 Calculate the appropriate volume of 1X Diluent M to add when resuspending the Factor B Standard vial to produce a 140 ng/mL Factor B **Standard #1** by using the following equation:

$C_S$  = Starting mass of Factor B Standard (see vial label)  
(ng)

$C_F$  = 140 ng/mL Factor B **Standard #1** final required  
concentration

$V_D$  = Required volume of 1X Diluent M for reconstitution  
( $\mu$ L)

Calculate total required volume 1X Diluent M for resuspension:

$$(C_S / C_F) \times 1,000 = V_D$$

Example:

**NOTE:** This example is for demonstration purposes only. Please remember to check your standard vial for the actual amount of standard provided.

$C_S$  = 280 ng of Factor B Standard in vial

$C_F$  = 140 ng/mL Factor B **Standard #1** final concentration

$V_D$  = Required volume of 1X Diluent M for reconstitution

$$(280 \text{ ng} / 140 \text{ ng/mL}) \times 1,000 = 2,000 \text{ }\mu\text{L}$$

- 10.1.3 Reconstitute the Factor B Standard vial by adding the appropriate calculated amount  $V_D$  of 1X Diluent M to the vial to generate the 140 ng/mL Factor B **Standard #1**. Mix gently and thoroughly.
- 10.2 Allow the reconstituted 140 ng/mL Factor B **Standard #1** to sit for 10 minutes with gentle agitation prior to making subsequent dilutions
- 10.3 Label eight tubes #2-8.
- 10.4 Add 120  $\mu\text{L}$  of 1X Diluent M to tubes #2 – 8.
- 10.5 To prepare **Standard #2**, add 120  $\mu\text{L}$  of the **Standard #1** into tube #2 and mix gently.
- 10.6 To prepare **Standard #3**, add 120  $\mu\text{L}$  of the **Standard #2** into tube #3 and mix gently.
- 10.7 Using the table below as a guide, prepare subsequent serial dilutions.
- 10.8 1X Diluent M serves as the zero standard, 0 ng/mL (tube #8).

## ASSAY PREPARATION

Standard Dilution Preparation Table

Standard #	Volume to Dilute (μL)	Volume Diluent M (μL)	Total Volume (μL)	Starting Conc. (ng/mL)	Final Conc. (ng/mL)
1	Step 10.4				140
2	120	120	240	140	70.00
3	120	120	240	70.00	35.00
4	120	120	240	35.00	17.50
5	120	120	240	17.50	8.750
6	120	120	240	8.750	4.375
7	120	120	240	4.375	2.188
8	-	120	120	-	0

## 11. SAMPLE PREPARATION

### 11.1 **Saliva**

Collect saliva using sample tube. Centrifuge samples at  $800 \times g$  for 10 minutes. Using an undiluted sample is suggested; however, the user should determine the optimal dilution factor into 1X Diluent M depending on the application needs, in the range 1:2-1:20. The undiluted samples can be stored at  $-20^{\circ}\text{C}$  or below for up to 3 months. Avoid repeated freeze-thaw cycles.

### 11.2 **Milk**

Collect milk using sample tube. Centrifuge samples at  $800 \times g$  for 10 minutes. A 50-fold sample dilution is suggested into 1X Diluent M; however, the user should determine optimal dilution factor depending on application needs in the range 1:5-1:500. The undiluted samples can be stored at  $-20^{\circ}\text{C}$  or below for up to 3 months. Avoid repeated freeze-thaw cycles.

### 11.3 **Plasma**

Collect plasma using one-tenth volume of 0.1 M sodium citrate as an anticoagulant. Centrifuge samples at  $3000 \times g$  for 10 minutes and collect plasma. A 4000-fold sample dilution is suggested into 1X Diluent M and assay, however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at  $-20^{\circ}\text{C}$  or below for up to 3 months. Avoid repeated freeze-thaw cycles (EDTA or heparin can also be used as an anticoagulant).

### 11.4 **Serum**

Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at  $3,000 \times g$  for 10 minutes and remove serum. A 16000-fold sample dilution is suggested into 1X Diluent M; however, user should determine optimal dilution factor depending on application needs. The undiluted

samples should be aliquoted to limit repeated freeze-thaw cycles and stored at  $-80^{\circ}\text{C}$  for up to 3 months. When needed, the frozen sample should be thawed rapidly in a water bath at  $37^{\circ}\text{C}$  and immediately placed on ice until use to prevent complement activation.

### 11.5 Cell Culture Supernatants

Centrifuge cell culture media at  $3,000 \times g$  for 10 minutes at  $4^{\circ}\text{C}$  to remove debris and collect supernatants. If necessary, dilute samples into 1X Diluent M; user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at  $-80^{\circ}\text{C}$ . Avoid repeated freeze-thaw cycles.

### 11.6 Cell lysate

Rinse cell with cold PBS and then scrape the cell into a tube with 5 ml of cold PBS and 0.5 M EDTA. Centrifuge suspension at 1500 rpm for 10 minutes at  $4^{\circ}\text{C}$  and aspirate supernatant. Resuspend pellet in ice-cold Lysis Buffer (PBS, 1% Triton X-100, protease inhibitor cocktail). For every  $1 \times 10^6$  cells, add approximately 100  $\mu\text{l}$  of ice-cold Lysis Buffer. Incubate on ice for 60 minutes. Centrifuge at 13000 rpm for 30 minutes at  $4^{\circ}\text{C}$  and collect supernatant. If necessary, dilute samples into EIA Diluent; user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at  $-80^{\circ}\text{C}$ . Avoid repeated freeze-thaw cycles.

### 11.7 Tissue

Extract tissue samples with 0.1 M phosphate-buffered saline (pH 7.4) containing 1% Triton X-100 and centrifuge at  $14,000 \times g$  for 20 minutes. Collect the supernatant and measure the protein concentration. If



necessary, dilute samples into EIA Diluent; user should determine optimal dilution factor depending on application needs. Store remaining extract at -80°C. Avoid repeated freeze-thaw cycles.

### 11.8 CSF

Collect cerebrospinal fluid (CSF) using sample pot. Centrifuge samples at 3000 x g for 10 minutes. A 2-fold sample dilution is suggested into 1X Diluent M; however, user should determine optimal dilution factor depending on application needs in the range 1:1-1:40. The undiluted samples can be stored at -80°C for up to 3 months. Avoid repeated freeze-thaw cycles.

## 12. PLATE PREPARATION

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused well plate strips should be returned to the plate packet and stored at 4°C.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section.

### 13. ASSAY PROCEDURE

- Equilibrate all materials and prepared reagents to room temperature (18 - 25°C) prior to use.
  - It is recommended to assay all standards, controls and samples in duplicate.
- 13.1 Prepare all reagents, working standards and samples as instructed. Equilibrate reagents to room temperature before use. The assay is performed at room temperature (18-25°C).
  - 13.2 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.
  - 13.3 Add 50 µL of Factor B Standard or sample per well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for two hours. Start the timer after the last addition.
  - 13.4 Wash five times with 200 µL of 1X Wash Buffer manually. Invert the plate each time and decant the contents; tap it 4-5 times on absorbent paper towel to completely remove the liquid. If using a machine wash six times with 300 µL of 1X Wash Buffer and then invert the plate, decant the contents; tap it 4-5 times on absorbent paper towel to completely remove the liquid.
  - 13.5 Add 50 µL of 1X Biotinylated Factor B Antibody to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed.

Cover wells with a sealing tape and incubate for 1 hour.

- 13.6 Wash microplate as described above.
- 13.7 Add 50  $\mu\text{L}$  of 1X SP Conjugate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
- 13.8 Wash microplate as described above.
- 13.9 Add 50  $\mu\text{L}$  of Chromogen Substrate per well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Incubate in ambient for 25 minutes or until the optimal blue color density develops.
- 13.10 Add 50  $\mu\text{L}$  of Stop Solution to each well. The color will change from blue to yellow. Gently tap plate to ensure thorough mixing. Break any bubbles that may have formed.
- 13.11 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.

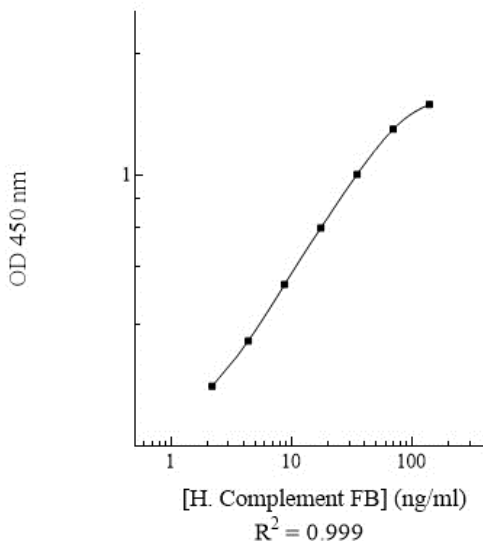
## 14. CALCULATIONS

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.

## 15. TYPICAL DATA

**TYPICAL STANDARD CURVE** – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

Human Complement Factor B Standard Curve



## 16. TYPICAL SAMPLE VALUES

### SENSITIVITY –

The minimum detectable dose of Human Factor B is typically ~0.71 ng/mL.

### RECOVERY –

Standard Added Value: 5 – 70 ng/mL

Recovery %: 88 – 110.

Average Recovery %: 98

### REFERENCE VALUES –

On average, normal Human Factor B plasma and serum levels is approximately 296 µg/mL.

### LINEARITY OF DILUTION –

Plasma Dilution	Average % Expected Value
1:8,000	102
1:16,000	100
1:32,000	97

Serum Dilution	Average % Expected Value
1:8,000	92
1:16,000	103
1:32,000	105

### PRECISION –

	Intra-Assay	Inter-Assay
% CV	5.3	10.2

## 17. ASSAY SPECIFICITY

Species	% Cross Reactivity
Canine	None
Equine	None
Mouse	None
Monkey	<5%
Bovine	None
Rat	None
Swine	None
Rabbit	None
Proteins	% Cross Reactivity
Factor I	None
Factor D	None
Factor P	None
Factor H	None

No significant cross-reactivity observed with complement factor D, factor H, factor I, factor P, C1, C2, C3, C4, C5, C6, C7, C8, and C9.

10% FBS in culture media will not affect this assay.

## 18. TROUBLESHOOTING

Problem	Cause	Solution
Poor standard curve	Improper standard dilution	Confirm dilutions made correctly
	Standard improperly reconstituted (if applicable)	Briefly spin vial before opening; thoroughly resuspend powder (if applicable)
	Standard degraded	Store sample as recommended
	Curve doesn't fit scale	Try plotting using different scale
Low signal	Incubation time too short	Try overnight incubation at 4°C
	Target present below detection limits of assay	Decrease dilution factor; concentrate samples
	Precipitate can form in wells upon substrate addition when concentration of target is too high	Increase dilution factor of sample
	Using incompatible sample type (e.g. serum vs. cell extract)	Detection may be reduced or absent in untested sample types
	Sample prepared incorrectly	Ensure proper sample preparation/dilution
Large CV	Bubbles in wells	Ensure no bubbles present prior to reading plate
	All wells not washed equally/thoroughly	Check that all ports of plate washer are unobstructed wash wells as recommended
	Incomplete reagent mixing	Ensure all reagents/master mixes are mixed thoroughly
	Inconsistent pipetting	Use calibrated pipettes and ensure accurate pipetting

## RESOURCES

Problem	Cause	Solution
	Inconsistent sample preparation or storage	Ensure consistent sample preparation and optimal sample storage conditions (eg. minimize freeze/thaws cycles)
Large CV	Improperly sealed microplate	Check the microplate pouch for proper sealing and has no punctures
High background/ Low sensitivity	Wells are insufficiently washed	Wash wells as per protocol recommendations
	Contaminated wash buffer	Make fresh wash buffer
	Waiting too long to read plate after adding STOP solution	Read plate immediately after adding STOP solution
	Improper storage of ELISA kit	Store all reagents as recommended. Please note all reagents may not have identical storage requirements.



## Technical Support

Copyright © 2023 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)