



ELISA PRODUCT INFORMATION & MANUAL

Fibrinogen *NBP2-60462*

Enzyme-linked Immunosorbent Assay for quantitative detection of Canine Fibrinogen. For research use only.
Not for diagnostic or therapeutic procedures.

Assay Summary

Step 1. Add 25 μ l of Standard or Sample and 25 μ l of Biotinylated Protein per well.
Incubate 2 hours.

Step 2. Wash, then add 50 μ l of SP Conjugate per well.
Incubate 30 minutes.

Step 3. Wash, then add 50 μ l of Chromogen Substrate per well.
Incubate 10 minutes.

Step 4. Add 50 μ l of Stop Solution per well.
Read at 450 nm immediately.

Assay Template

[illegible]

Canine Fibrinogen (FBG) ELISA Kit

Catalog No. NBP2-60462

Introduction

Fibrinogen (FBG) is a homodimer (340 kDa) that is made up of two sets of alpha, beta, and gamma polypeptide chains. FBG is synthesized in the parenchymal cell of the hepatocyte and in the megakaryocyte (1). FBG plays a major role in coagulation. Upon cleavage by thrombin in the initial stages of coagulation activation, FBG self-assembles to yield a fibrin clot matrix that subsequently is crosslinked by factor XIIIa to form an insoluble network. FBG also binds to the platelet glycoprotein IIb/IIIa receptor to form bridges between platelets, thus facilitating aggregation (2).

Principle of the Assay

The Canine Fibrinogen ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for detection of canine FBG in **plasma samples**. This assay employs a quantitative **competitive enzyme immunoassay** technique that measures canine FBG in less than 3 hours. A polyclonal antibody specific for canine FBG has been pre-coated onto a 96-well microplate with removable strips. Canine FBG in standards and samples is competed with a biotinylated canine FBG sandwiched by the immobilized antibody and streptavidin-peroxidase conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

Caution and Warning

- This product is for **Research Use Only** and is Not For Use In Diagnostic Procedures.
- Prepare all reagents (working diluent buffer, wash buffer, standard, biotinylated protein, 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 insert. However, the user should determine the optimal dilution factor.
- Spin down the SP conjugate vial before opening and using contents.
- The Stop Solution is an acidic solution.
- The kit should not be used beyond the expiration date.

Reagents

- **Canine FBG Microplate:** A 96-well polystyrene microplate (12 strips of 8 wells) coated with a polyclonal antibody against canine FBG.
- **Sealing Tapes:** Each kit contains 3 precut, pressure sensitive sealing tapes that can be cut to fit the format of the individual assay.
- **Canine FBG Standard:** Canine FBG in a buffered protein base (20 µg, lyophilized).
- **Biotinylated Canine FBG:** 2 vials, lyophilized.
- **EIA Diluent Concentrate (10x):** A 10-fold concentrated buffered protein base (30 ml).
- **Wash Buffer Concentrate (20x):** A 20-fold concentrated buffered surfactant (30 ml).
- **Streptavidin-Peroxidase Conjugate (SP Conjugate):** A 100-fold concentrate (80 µl).
- **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 Condition

- Upon arrival, immediately store components of the kit at recommended temperatures up to the expiration date.
- Store SP Conjugate at -20°C.
- Store Microplate, Diluent Concentrate (10x), Wash Buffer, Stop Solution, and Chromogen Substrate at 2-8°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°C.
- Store Standard and Biotinylated Protein at 2-8°C before reconstituting with Diluent and at -20°C after reconstituting with Diluent.

Other Supplies Required

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

Sample Collection, Preparation, and Storage

- 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 and use supernatants. Dilute samples 1:1000 into EIA 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 can also be used as an anticoagulant).

Refer to Sample Dilution Guidelines below for further instruction.

Guidelines for Dilutions of 1:100 or Greater <i>(for reference only; please follow the insert for specific dilution suggested)</i>	
1:100	1:10000
A) 4 µl sample : 396 µl buffer (100x) = 100 fold dilution <i>Assuming the needed volume is less than or equal to 400 µl.</i>	A) 4 µl sample : 396 µl buffer (100x) B) 4 µl of A : 396 µl buffer (100x) = 10000 fold dilution <i>Assuming the needed volume is less than or equal to 400 µl.</i>
1:1000	1:100000
A) 4 µl sample : 396 µl buffer (100x) B) 24 µl of A : 216 µl buffer (10x) = 1000 fold dilution <i>Assuming the needed volume is less than or equal to 240 µl.</i>	A) 4 µl sample : 396 µl buffer (100x) B) 4 µl of A : 396 µl buffer (100x) C) 24 µl of B : 216 µl buffer (10x) = 100000 fold dilution <i>Assuming the needed volume is less than or equal to 240 µl.</i>

Reagent Preparation

- Freshly dilute all reagents and bring all reagents to room temperature before use.
- EIA Diluent Concentrate (10x):** If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the EIA Diluent Concentrate 1:10 with reagent grade water. Store for up to 30 days at 2-8°C.
- Standard Curve:** Reconstitute the 20 µg of Canine FBG Standard with 1 ml of EIA Diluent to generate a 20 µg/ml standard stock 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 stock solution (20 µg/ml) 1:2 with EIA Diluent to produce 10, 5, 2.5, 1.25, and 0.625 µg/ml solutions. EIA Diluent serves as the zero standard (0 µg/ml). Any remaining solution should be frozen at -20°C and used within 5 days.

Standard Point	Dilution	[Canine FBG] (µg/ml)
P1	1 part Standard (20 µg/ml)	20.00
P2	1 part P1 + 1 part EIA Diluent	10.00
P3	1 part P2 + 1 part EIA Diluent	5.000
P4	1 part P3 + 1 part EIA Diluent	2.500
P5	1 part P4 + 1 part EIA Diluent	1.250
P6	1 part P5 + 1 part EIA Diluent	0.625
P7	EIA Diluent	0.000

- **Biotinylated Canine FBG (1x):** Reconstitute Biotinylated Canine FBG with 2 ml EIA Diluent to produce a working solution. Allow the biotin to sit for 10 minutes with gentle agitation prior to use. Any remaining solution should be frozen at -20°C and used within 5 days.
- **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 EIA Diluent. Any remaining solution should be frozen at -20°C.

Assay Procedure

- 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).
- Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccants inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
- Add 25 µl of Canine FBG Standard or sample per well and immediately add 25 µl of Biotinylated Canine FBG to each well (on top of the standard or sample) and tap plate to mix gently. Cover wells with a sealing tape and incubate for 2 hours. Start the timer after the last addition.
- 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, decanting the contents; hit 4-5 times on absorbent material to completely remove the liquid.
- Add 50 µl of Streptavidin-Peroxidase Conjugate to each well and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
- Wash the microplate as described above.

- Add 50 µl of Chromogen Substrate per well and incubate for 10 minutes or till the optimal color density develops. Gently tap plate to ensure thorough mixing and break the bubbles in the well with pipette tip.
- Add 50 µl of Stop Solution to each well. The color will change from blue to yellow.
- 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 low concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

Data Analysis

- Calculate the mean value of the duplicate or 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 (OD) 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.

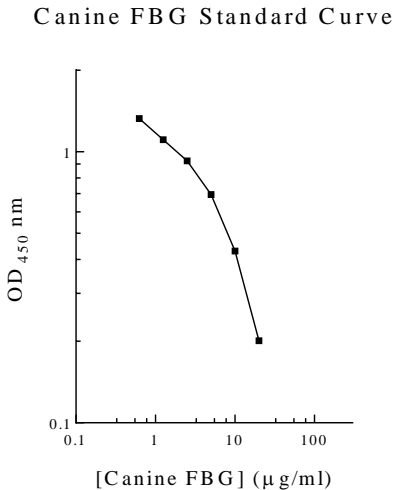
Typical Data

- The typical data is provided for reference only. Individual laboratory means may vary from the values listed. Variations between laboratories may be caused by technique differences.

Standard Point	µg/ml	Average OD
P1	20.00	0.201
P2	10.00	0.430
P3	5.000	0.695
P4	2.500	0.926
P5	1.250	1.108
P6	0.625	1.324
P7	0.000	1.601
Sample: Normal, Sodium Citrate Plasma (1000x)		1.031

Standard Curve

- 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 canine FBG as calculated by 2SD from the mean of a zero standard was established to be 0.4 µg/ml.
- Intra-assay precision was determined by testing replicates of three plasma samples in one assay.
- Inter-assay precision was determined by testing three plasma samples in twenty assays.

	Intra-Assay Precision			Inter-Assay Precision		
Sample	1	2	3	1	2	3
n	20	20	20	20	20	20
CV (%)	4.6%	5.0%	4.4%	8.6%	8.5%	9.0%
Average CV (%)	4.7%			8.7%		

Recovery

Standard Added Value	1 – 10 µg/ml
Recovery %	92 – 110%
Average Recovery %	96%

Linearity

- Plasma samples were serially-diluted to test for linearity.

Average Percentage of Expected Value (%)	
Sample Dilution	Plasma
1:500	104%
1:1000	96%
1:2000	101%

Cross-Reactivity

Species	Cross Reactivity (%)
Canine	100%
Bovine	None
Monkey	None
Mouse	<5%
Rat	<5%
Swine	10%
Rabbit	None
Human	None

Troubleshooting

Issue	Causes	Course of Action
Low Precision	Use of expired components	<ul style="list-style-type: none">Check the expiration date listed before use.Do not interchange components from different lots.
	Improper wash step	<ul style="list-style-type: none">Check that the correct wash buffer is being used.Check that all wells are dry after aspiration.Check that the microplate washer is dispensing properly.If washing by pipette, check for proper pipetting technique.
	Splashing of reagents while loading wells	<ul style="list-style-type: none">Pipette properly in a controlled and careful manner.
	Inconsistent volumes loaded into wells	<ul style="list-style-type: none">Pipette properly in a controlled and careful manner.Check pipette calibration.Check pipette for proper performance.
	Insufficient mixing of reagent dilutions	<ul style="list-style-type: none">Thoroughly agitate the lyophilized components after reconstitution.Thoroughly mix dilutions.

	Improperly sealed microplate	<ul style="list-style-type: none"> • Check the microplate pouch for proper sealing. • Check that the microplate pouch has no punctures. • Check that three desiccants are inside the microplate pouch prior to sealing.
Unexpectedly Low or High Signal Intensity	Microplate was left unattended between steps	<ul style="list-style-type: none"> • Each step of the procedure should be performed uninterrupted.
	Omission of step	<ul style="list-style-type: none"> • Consult the provided procedure for complete list of steps.
	Steps performed in incorrect order	<ul style="list-style-type: none"> • Consult the provided procedure for the correct order.
	Insufficient amount of reagents added to wells	<ul style="list-style-type: none"> • Check pipette calibration. • Check pipette for proper performance.
	Wash step was skipped	<ul style="list-style-type: none"> • Consult the provided procedure for all wash steps.
	Improper wash buffer	<ul style="list-style-type: none"> • Check that the correct wash buffer is being used.
	Improper reagent preparation	<ul style="list-style-type: none"> • Consult reagent preparation section for the correct dilutions of all reagents.
	Insufficient or prolonged incubation periods	<ul style="list-style-type: none"> • Consult the provided procedure for correct incubation time.
Deficient Standard Curve Fit	Non-optimal sample dilution	<ul style="list-style-type: none"> • Sandwich ELISA: If samples generate OD values higher than the highest standard point (P1), dilute samples further and repeat the assay. • Competitive ELISA: If samples generate OD values lower than the highest standard point (P1), dilute samples further and repeat the assay. • User should determine the optimal dilution factor for samples.
	Contamination of reagents	<ul style="list-style-type: none"> • A new tip must be used for each addition of different samples or reagents during the assay procedure.
	Contents of wells evaporate	<ul style="list-style-type: none"> • Verify that the sealing film is firmly in place before placing the assay in the incubator or at room temperature.
	Improper pipetting	<ul style="list-style-type: none"> • Pipette properly in a controlled and careful manner. • Check pipette calibration. • Check pipette for proper performance.
	Insufficient mixing of reagent dilutions	<ul style="list-style-type: none"> • Thoroughly agitate the lyophilized components after reconstitution. • Thoroughly mix dilutions.

Version 2.2R