



TF ELISA Kit

Catalog Number KA2047

96 assays

Version: 04

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	5
Precautions for Use	5
Assay Protocol	6
Reagent Preparation	6
Sample Preparation	6
Assay Procedure	7
Data Analysis.....	9
Calculation of Results.....	9
Resources.....	10
Plate Layout	10

Introduction

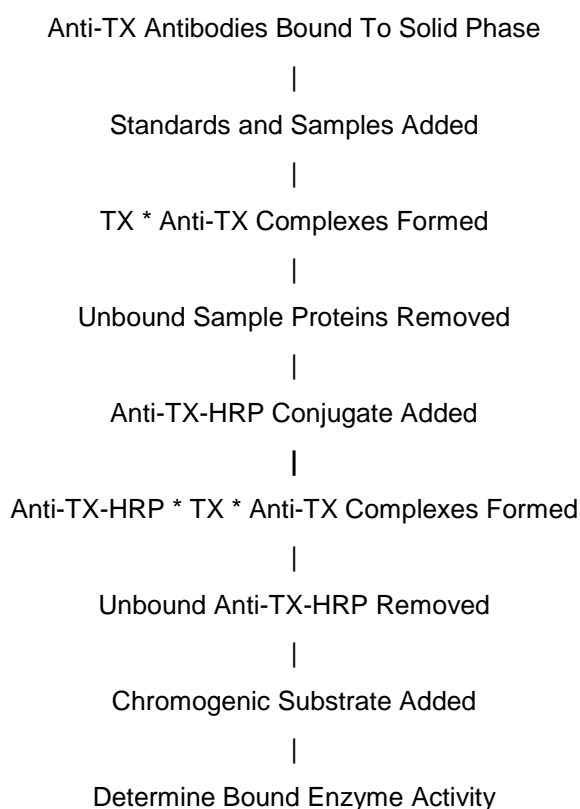
Intended Use

The TF (Bovine) ELISA Kit is a highly sensitive two-site enzyme-linked immunoassay (ELISA) for measuring transferrin (TX) in bovine biological samples. If the ELISA is to be used outside the intended use, the user may need to optimize for said use.

Principle of the Assay

The principle of the double antibody sandwich ELISA is represented in Figure 1. In this assay the transferrin (TX) present in samples reacts with the anti-TX antibodies, which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-TX antibodies conjugated with horseradish peroxidase (HRP) are added. These enzyme-labeled antibodies form complexes with the previously bound TX. Following another washing step, the enzyme bound to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of TX in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of TX in the test sample. The quantity of TX in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution

Figure 1.



General Information

Materials Supplied

List of component

Component	Description	Amount
ELISA Micro Plate, antibody coated	One plate of 12 removable 8 well strips, antibody coated. Ready to use as supplied.	96 (8x12) wells
Enzyme Conjugated Detection Antibody	One vial of 100X Horseradish Peroxidase Conjugated antibody in a stabilizing buffer. Dilute 1/100 immediately prior to use.	150 μ L
Calibrator	One vial of calibrator. Refer to Certificate of Analysis (CoA).	1 vial
Diluent Concentrate	One bottle of 20X diluent buffer. Dilute 1/20 to make 1X working solution.	50 mL
Wash Solution Concentrate	One bottle of 20X wash solution. Dilute 1/20 to make 1X working solution.	50 mL
Chromogen-Substrate Solution	One bottle of 3,3',5,5'- tetramethylbenzidine (TMB) and hydrogen peroxide in citric acid buffer at pH 3.3. Ready to use as supplied.	12 mL
STOP Solution	One bottle of 0.3 M sulfuric acid. <i>WARNING: Avoid Contact with Skin.</i> Ready to use as supplied.	12 mL

Storage Instruction

The expiration date for the kit and its components is stated on the box label. All components should be stable up to the expiration date if stored and used per this kit protocol insert.

Component	Storage	Stability
ELISA Micro Plate, antibody coated	4-8°C, in sealed foil bag with desiccant.	With proper storage the plate strips are stable until the expiration date.
Enzyme Conjugated Detection Antibody	4-8°C in the dark.	The working conjugate solution should be diluted immediately prior to use. The 100X conjugate is stable until the expiration date.
Calibrator	4-8°C for lyophilized calibrator. Aliquoted and frozen if reconstituted. Avoid multiple freeze-thaw cycles.	The working standard solutions should be prepared immediately prior to use.
Diluent Concentrate	4-8°C for both 1X working solution and 20X concentrate.	The 1X working solution is stable for at least one week from the date of preparation. The 20X concentrate is stable until the expiration date.

Wash Solution Concentrate	4-8°C for both 1X working solution and 20X wash concentrate.	The 1X working solution is stable for at least one week from the date of preparation. The 20X concentrate is stable until the expiration date.
Chromogen-Substrate Solution	4-8°C in the dark.	Protect from light. The Substrate Solution is stable until the expiration date.
STOP Solution	4-8°C.	The Stop Solution is stable until the expiration date.

Materials Required but Not Supplied

- ✓ Precision pipette (2 µL to 100 µL) for making and dispensing dilutions
- ✓ Test tubes
- ✓ Squirt bottler or Microtitre washer/aspirator
- ✓ Distilled or Deionized H₂O
- ✓ Microtitre Plate reader
- ✓ Assorted glassware for the preparation of reagents and buffer solutions
- ✓ Centrifuge for sample collection
- ✓ Anticoagulant for plasma collection
- ✓ Timer

Precautions for Use

Please read this protocol completely before using this product.

For Research Use Only, Not for Diagnostic Purposes. Not for Human and Animal Consumption. For In Vitro Laboratory Use Only.

- ✓ Limitation of the procedure
Reliable and reproducible results will be obtained when the assay procedure is carried out with a complete understanding of the information contained in the package insert instructions and with adherence to good laboratory practice. Factors that might affect the performance of the assay include instrument function, cleanliness of glassware, quality of distilled or deionized water, and accuracy of reagent and sample pipetting, washing technique, incubation time or temperature. Do not mix or substitute reagents with those from other lots or sources.

Assay Protocol

Reagent Preparation

Bring all reagents to room temperature (16°C to 25°C) before use.

- Diluent Concentrate

The Diluent solution supplied is a 20X Concentrate and must be diluted 1/20 with distilled or deionized water (1 part buffer concentrate, 19 parts dH₂O).

- Wash Solution Concentrate.

The Wash Solution supplied is a 20X concentrate and must be diluted 1/20 with distilled or deionized water (1 part buffer concentrate, 19 parts dH₂O). Crystal formation in the concentrate may occur when storage temperatures are low. Warming of the concentrate to 30-35°C before dilution can dissolve crystals.

- Enzyme-Antibody Conjugate

Calculate the required amount of working conjugate solution for each microtiter plate test strip by adding 10 µL Enzyme-Antibody Conjugate to 990 µL of 1X Diluent for each test strip to be used for testing. Dilute immediately before use and protect from light. Mix uniformly, but gently. Avoid foaming.

- Pre-coated ELISA Micro Plate

Ready to use as supplied. Unseal foil pouch and remove plate from pouch. Remove all strips and wells that will not be used in the assay and place back in pouch and re-seal along with desiccant.

- Bovine TX Calibrator

Prepare according to the lot specific Certificate of Analysis.

Sample Preparation

- Specimen Collection and Handling

All blood components and biological materials should be handled as potentially hazardous. Follow universal precautions when handling and disposing.

If blood samples are clotted, grossly hemolyzed, lipemic, or the integrity of the sample is of concern, make a note and interpret results with caution.

The sample collection and storage conditions listed below are intended as general guidelines. Sample stability has not been evaluated.

- Serum samples

Blood should be collected by venipuncture. The serum should be separated from the cells after clot formation by centrifugation. Remove serum and assay immediately or aliquot and store samples at -80°C (preferably) or -20°C. Avoid repeated freeze-thaw cycles.

- Plasma samples

Blood should be collected into a container with an anticoagulant and then centrifuged. Assay immediately or aliquot and store samples at -80°C (preferably) or -20°C. Avoid repeated freeze-thaw cycles.

- Urine samples

Collect mid-stream using sterile or clean urine collector. Centrifuge to remove cell debris. Assay immediately or aliquot and store samples at -80°C (preferably) or -20°C . Avoid repeated freeze-thaw cycles.

- Known interfering substances

Azide and thimerosal at concentrations higher than 0.1% inhibits the enzyme reaction.

- Dilution of Samples

The assay requires that each test sample be diluted before use. All samples should be assayed in duplicate each time the assay is performed. The recommended dilutions are only suggestions. Dilutions should be based on the expected concentration of the unknown sample such that the diluted sample falls within the dynamic range of the standard curve. If unsure of sample level, a serial dilution with one or two representative samples before running the entire plate is highly recommended.

- Serum samples

Recommended starting dilution is 1/40,000. To prepare a 1/40,000 dilution of a sample, transfer 5 μL of sample to 495 μL of 1X diluent. This gives you a 1/100 dilution. Next, dilute the 1/100 by transferring 2 μL into 798 μL of 1X diluent. This gives you a 1/40,000 dilution Mix thoroughly at each stage.

- Plasma samples

Recommended starting dilution is 1/40,000. To prepare a 1/40,000 dilution of a sample, transfer 5 μL of sample to 495 μL of 1X diluent. This gives you a 1/100 dilution. Next, dilute the 1/100 by transferring 2 μL into 798 μL of 1X diluent. This gives you a 1/40,000 dilution Mix thoroughly at each stage.

Assay Procedure

1. All samples and standards should be assayed in duplicates.
2. The Standards and the test sample(s) should be loaded into the ELISA wells as quickly as possible to avoid a shift in OD readings. Using a multichannel pipette would reduce this occurrence.

Pipette 100 μL of

Standard 0 (0.0 ng/mL) in duplicate

Standard 1 (18.75 ng/mL) in duplicate

Standard 2 (37.5 ng/mL) in duplicate

Standard 3 (75 ng/mL) in duplicate

Standard 4 (150 ng/mL) in duplicate

Standard 5 (300 ng/mL) in duplicate

Standard 6 (600 ng/mL) in duplicate

3. Pipette 100 μL of sample (in duplicate) into pre designated wells.
4. Incubate the Microtiter Plate at room temperature for thirty (30 ± 2) minutes. Keep plate covered and level during incubation.

5. Following incubation, aspirate the contents of the wells.
6. Completely fill each well with appropriately diluted Wash Solution and aspirate. Repeat three times, for a total of four washes. If washing manually: completely fill wells with wash buffer, invert the plate then pour/shake out the contents in a waste container. Follow this by sharply striking the wells on absorbent paper to remove residual buffer. Repeat 3 times for a total of four washes.
7. Pipette 100 μ L of appropriately diluted Enzyme-Antibody Conjugate to each well. Incubate at room temperature for twenty (20 ± 2) minutes. Keep plate covered in the dark and level during incubation.
8. Wash and blot the wells as described in Steps 5/6.
9. Pipette 100 μ L of TMB Substrate Solution into each well.
10. Incubate in the dark at room temperature for precisely ten (10) minutes.
11. After ten minutes, add 100 μ L of Stop Solution to each well.
12. Determine the absorbance (450 nm) of the contents of each well within 30 minutes. Calibrate the plate reader to manufacturer's specifications.

Data Analysis

Calculation of Results

1. Subtract the average background value (Average absorbance reading of Standard zero) from the test values for each sample.
2. Average the duplicate readings for each standard and use the results to construct a Standard Curve. Construct the standard curve by reducing the data using computer software capable of generating a four parameter logistic curve fit. A second order polynomial (quadratic) or other curve fits may also be used; however, they will be a less precise fit of the data.
3. Interpolate test sample values from standard curve. Correct for sera dilution factor to arrive at the TX concentration in original samples.

Resources

Plate Layout

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