



## HUMAN HEMOPEXIN ELISA Assay Kit

Immunoperoxidase Assay for Determination of Hemopexin in Human Samples

HPX39-K01

DIRECTIONS FOR USE

For Research Use Only, NOT for Diagnostic Purposes

***Please Read this Package Insert Completely Before Using This Product***

### INTENDED USE

The Human Hemopexin ELISA Assay Kit is a highly sensitive two-site enzyme linked immunoassay (ELISA) for measuring Hemopexin in serum, plasma of Human.

### INTRODUCTION

Hemopexin (HX) is the plasma protein with the highest binding affinity to heme among known proteins. It is mainly expressed in liver and belongs to acute phase reactants, the synthesis of which is induced after inflammation. This ELISA kit can be used to measure Hemopexin in serum, tissue extracts and other biological fluids.

### PRINCIPLE OF THE ASSAY

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

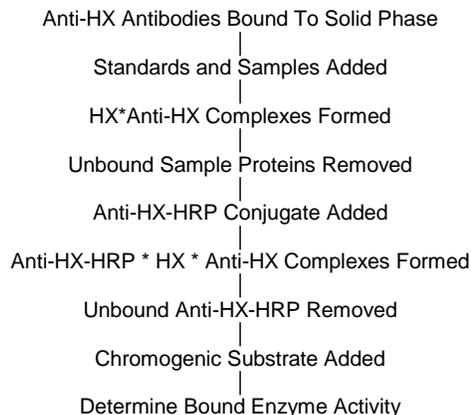


Figure 1.

### REAGENTS (Quantities sufficient for 96 determinations)

- 1. DILUENT CONCENTRATE (Running Buffer)**  
One bottle containing 50 ml of a 5X concentrated diluent running buffer.
- 2. WASH SOLUTION CONCENTRATE**  
One bottle containing 50 ml of a 20X concentrated wash solution.
- 3. ENZYME-ANTIBODY CONJUGATE 100X**  
One vial containing 150  $\mu$ L of affinity purified anti-Human Hemopexin antibody conjugated with horseradish peroxidase in a stabilizing buffer.
- 4. CHROMOGEN-SUBSTRATE SOLUTION**  
One vial containing 12 mL of 3,3',5,5'-tetramethylbenzidine (TMB) and hydrogen peroxide in citric acid buffer at pH 3.3.
- 5. STOP SOLUTION**  
One vial containing 12 ml 0.3 M sulfuric acid.

**WARNING: Avoid contact with skin.**

- 6. ANTI-HUMAN HEMOPEXIN ELISA MICRO PLATE**  
Twelve removable eight (8) well micro well strips in well holder frame. Each well is coated with affinity purified anti-Human HX.

## 7. HUMAN HEMOPEXIN CALIBRATOR

One vial containing a lyophilized Human Hemopexin calibrator.

### FOR RESEARCH USE ONLY

## REAGENT PREPARATION

### 1. DILUENT CONCENTRATE

The Diluent Solution supplied is a 5X Concentrate and must be diluted 1/5 with distilled or deionized water (1 part buffer concentrate, 4 parts dH<sub>2</sub>O).

### 2. 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 dH<sub>2</sub>O). Crystal formation in the concentrate is not uncommon when storage temperatures are low. Warming of the concentrate to 30-35°C before dilution can dissolve crystals.

### 3. ENZYME-ANTIBODY CONJUGATE

Calculate the required amount of working conjugate solution for each microtitre 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. Mix uniformly, but gently. Avoid foaming.

### 4. CHROMOGEN-SUBSTRATE SOLUTION

Ready to use as supplied.

### 5. STOP SOLUTION

Ready to use as supplied.

### 6. ANTI-HUMAN HEMOPEXIN ELISA MICRO PLATE

Ready to use as supplied. Unseal Microtiter 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.

### 7. HUMAN HEMOPEXIN CALIBRATOR

Add 1.0 ml of distilled or de-ionized water to the Human Hemopexin Calibrator and mix gently until dissolved. The calibrator is now at a concentration of 10.1 µg/ml (**the reconstituted calibrator should be aliquoted and frozen if future use is intended**). Human Hemopexin standards need to be prepared immediately prior to use (see the following chart). Mix well between each step. Avoid foaming.

Standard	ng/ml	Volume added to 1x Diluent	Volume of 1x Diluent
6	200	12 µl Human HX Calibrator	594 µl
5	100	300 µl standard 6	300 µl
4	50	300 µl standard 5	300 µl
3	25	300 µl standard 4	300 µl
2	12.5	300 µl standard 3	300 µl
1	6.25	300 µl standard 2	300 µl
0	0		600 µl

## STORAGE AND STABILITY

The expiration date for the package is stated on the box label.

### 1. DILUENT

The 5X Diluent Concentrate is stable until the expiration date. The 1X working solution is stable for at least one week from the date of preparation. Both solutions should be stored at 4-8°C.

### 2. WASH SOLUTION

The 20X Wash Solution Concentrate is stable until the expiration date. The 1X working solution is stable for at least one week from the date of preparation. Both solutions can be stored at room temperature (16-25°C) or at 4-8°C.

### 3. ENZYME-ANTIBODY CONJUGATE

Undiluted horseradish peroxidase anti-HX conjugate should be stored at 4-8°C and **diluted immediately prior to use**. The working conjugate solution is stable for up to 1 hour when stored in the dark.

### 4. CHROMOGEN-SUBSTRATE SOLUTION

The Substrate Solution should be stored at 4-8°C and is stable until the expiration date.

### 5. STOP SOLUTION

The Stop Solution should be stored at 4-8°C and is stable until the expiration date.

### 6. ANTI-HUMAN HEMOPEXIN ELISA MICRO PLATE

Anti-Human HX coated wells are stable until the expiration date, and should be stored at 4-8°C in sealed foil pouch with desiccant pack.

### 7. HUMAN HEMOPEXIN CALIBRATOR

**The lyophilized Human Hemopexin calibrator should be stored at 4°C or frozen until reconstituted. The reconstituted calibrator should be aliquoted out and stored frozen (Avoid multiple freeze-thaw cycles).** The working standard solutions should be prepared immediately prior to use and are stable for up to 8 hours

## INDICATIONS OF INSTABILITY

If the test is performing correctly, the results observed with the standard solutions should be within 20 % of the expected values.

## SPECIMEN COLLECTION AND HANDLING

Blood should be collected by venipuncture. The serum should be separated from the cells after clot formation by centrifugation. For plasma samples, blood should be collected into a container with an anticoagulant and then centrifuged. Care should be taken to minimize hemolysis, excessive hemolysis can impact your results. Assay immediately or aliquot and store samples at -20°C. Avoid repeated freeze-thaw cycles.

### 1. Precautions

For any sample that might contain pathogens, care must be taken to prevent contact with open wounds.

### 2. Additives and Preservatives

No additives or preservatives are necessary to maintain the integrity of the specimen. Avoid azide contamination.

### 3. Known interfering substances

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

## MATERIAL PROVIDED

See "REAGENTS"

## MATERIALS REQUIRED BUT NOT PROVIDED

- Precision pipette (2  $\mu$ L to 200  $\mu$ L) for making and dispensing dilutions
- Test tubes
- Microtitre washer/aspirator
- Distilled or Deionized H<sub>2</sub>O
- Microtitre Plate reader
- Assorted glassware for the preparation of reagents and buffer solutions
- Timer

## ASSAY PROTOCOL

### DILUTION OF SAMPLES

The assay for quantification of HX requires that each test sample be diluted before use. For a single step determination a dilution of serum/plasma at 1/40,000 is appropriate for most samples. For absolute quantification, samples that yield results outside the range of the standard curve, a lesser or greater dilution might be required. **If unsure of sample level, a serial dilution with one or two representative samples before running the entire plate is highly recommended.**

1. To prepare a 1/40,000 dilution of sample, transfer 5  $\mu$ L of sample to 495 $\mu$ L of 1X diluent. This gives you a 1/100 dilution. Next, dilute the 1/100 samples by transferring 2  $\mu$ L, to 798  $\mu$ L of 1X diluent. You now have a 1/40,000 dilution of your sample. Mix thoroughly at each stage.

### PROCEDURE

1. **Bring all reagents to room temperature before use.**

2. Pipette 100  $\mu$ L of  
Standard 0 (0.0 ng/ml) in duplicate  
Standard 1 (6.25 ng/ml) in duplicate  
Standard 2 (12.5 ng/ml) in duplicate  
Standard 3 (25 ng/ml) in duplicate  
Standard 4 (50 ng/ml) in duplicate  
Standard 5 (100 ng/ml) in duplicate  
Standard 6 (200 ng/ml) in duplicate

3. Pipette 100  $\mu$ L of sample (in duplicate) into pre designated wells.

4. Incubate the micro titer plate at room temperature for thirty (30  $\pm$  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  $\mu$ L of appropriately diluted Enzyme-Antibody Conjugate to each well. Incubate at room temperature for thirty (30  $\pm$  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  $\mu$ 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  $\mu$ L of Stop Solution to each well.
12. Determine the absorbance (450 nm) of the contents of each well. Calibrate the plate reader to manufacturer's specifications.

#### **STABILITY OF THE FINAL REACTION MIXTURE**

The absorbance of the final reaction mixture can be measured up to 2 hours after the addition of the Stop Solution. However, good laboratory practice dictates that the measurement be made as soon as possible.

#### **RESULTS**

1. Subtract the average background value from the test values for each sample.
2. Using the results observed for the standards construct a Standard Curve. The appropriate curve fit is that of a four-parameter logistics curve. A second order polynomial (quadratic) or other curve fits may also be used.

#### **LIMITED WARRANTY**

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or noninfringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.

***THANK YOU FOR USING EAGLE BIOSCIENCES!***

3. Interpolate test sample values from standard curve. Correct for sera dilution factor to arrive at the Hemopexin concentration in original samples.

#### **LIMITATION OF THE PROCEDURE**

1. 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.
2. Factors that might affect the performance of the assay include proper instrument function, cleanliness of glassware, quality of distilled or deionized water, and accuracy of reagent and sample pipettings, washing technique, incubation time or temperature.
3. Do not mix or substitute reagents with those from other lots or sources.

EAGLE BIOSCIENCES, INC.

20A NW BLVD, SUITE 112 NASHUA, NH 03063  
P: 617-419-2019 F: 617-419-1110

WWW.EAGLEBIO.COM — INFO@EAGLEBIO.COM



EAGLE  
BIOSCIENCES