



## **ELISA PRODUCT INFORMATION & MANUAL**

### **Human Erythropoietin ELISA Kit (Colorimetric)**

***NBP2-31082***

Enzyme-linked Immunosorbent Assay for quantitative  
detection. For research use only.

Not for diagnostic or therapeutic procedures.

[www.novusbio.com](http://www.novusbio.com) - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - [technical@novusbio.com](mailto:technical@novusbio.com)

Novus kits are guaranteed for 6 months from date of receipt

## TABLE OF CONTENTS

1	Intended Use	3
2	Summary	3
3	Principles of the Test	5
4	Reagents Provided	7
5	Storage Instructions – ELISA Kit	9
6	Specimen Collection and Storage Instructions	9
7	Materials Required But Not Provided	10
8	Precautions for Use	11
9	Preparation of Reagents	13
10	Test Protocol	17
11	Calculation of Results	22
12	Limitations	25
13	Performance Characteristics	26
14	Reagent Preparation Summary	33
15	Test Protocol Summary	34

## 1 Intended Use

The human Erythropoietin ELISA is an enzyme-linked immunosorbent assay for the quantitative detection of human Erythropoietin. **The human Erythropoietin ELISA is for research use only. Not for diagnostic or therapeutic procedures.**

## 2 Summary

Erythropoietin (EPO) is a hormone produced by the kidney that promotes the formation of red blood cells in the bone marrow. EPO is a glycoprotein with a molecular weight of 34,000.

The kidney cells that make EPO are specialized and sensitive to low oxygen levels in the blood. These cells release EPO when the oxygen level is low in the kidney. EPO then stimulates the bone marrow to produce more red cells and thereby increase the oxygen-carrying capacity of the blood.

EPO is the prime regulator of red blood cell production. Its major functions are to promote the differentiation and development of red blood cells and to initiate the production of hemoglobin, the molecule within red cells that transports oxygen.

The EPO gene has been found on human chromosome 7 (in band 7q21). EPO is produced not only in the kidney but also, to a lesser extent, in the liver. Different DNA sequences flanking the EPO gene act to control kidney versus liver production of EPO.

The measurement of EPO in the blood is useful in the study of bone marrow disorders and kidney disease. Elevated levels of EPO can be seen in polycythemia, a disorder in which there is an excess of red blood cells. Lower than normal levels of EPO are seen in chronic renal failure.

EPO plays an important role in the brain's response to neuronal injury. EPO is also involved in the wound healing process.

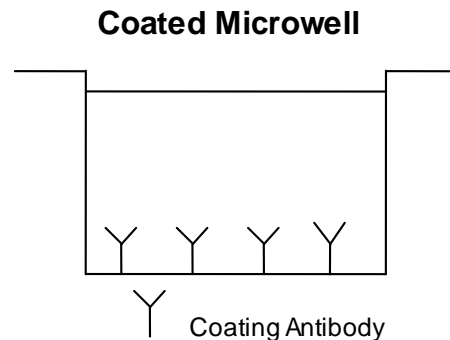
Measurement of serum immunoreactive erythropoietin has shown that overproduction of EPO can be an adaptive response to conditions producing tissue hypoxia, such as smoking chronic obstructive pulmonary disease, renal hypoxia or cyanotic heart disease. Elevated serum levels have further been found in patients suffering from various neoplastic diseases, such as renal carcinomas and benign renal tumors, liver carcinomas and hepatomas, cerebellar hemangioblastomas and other neoplastic disorders.

Serum levels of EPO lower than normal are found in various forms of anemias. Proinflammatory cytokines hereby act in lowering erythropoietin production. Mainly IL-1 and TNF alpha seem to be responsible for the defect in EPO production in severe systemic and renal inflammatory diseases. These include anemia of renal failure, anemias of chronic disorders like in AIDS or rheumatoid arthritis, anemias due to hypothyroidism, malnutrition, prematurity. In case EPO is not primarily involved in the cause of the anemia, elevated levels of EPO are found in serum of the patients suffering from diseases like aplastic anemias, iron deficiency, megaloblastic anemias, thalassemias and myelodysplastic syndromes.

### 3 Principles of the Test

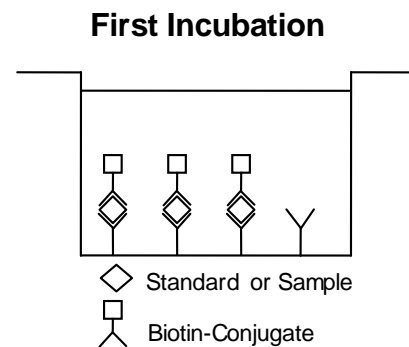
An anti-human Erythropoietin coating antibody is adsorbed onto microwells.

Figure 1



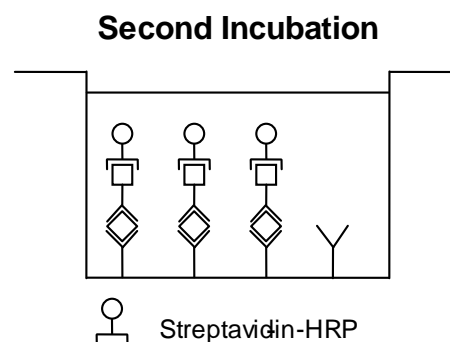
Human Erythropoietin present in the sample or standard binds to antibodies adsorbed to the microwells. A biotin-conjugated anti-human Erythropoietin antibody is added and binds to human Erythropoietin captured by the first antibody.

Figure 2



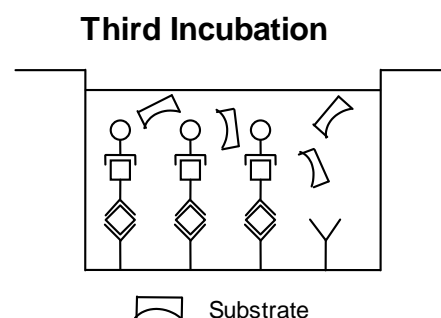
Following incubation unbound biotin-conjugated anti-human Erythropoietin antibody is removed during a wash step. Streptavidin-HRP is added and binds to the biotin-conjugated anti-human Erythropoietin antibody.

Figure 3



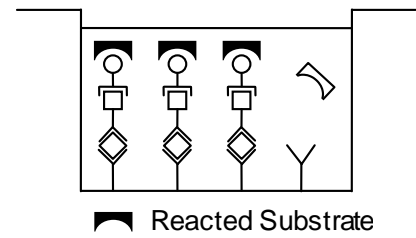
Following incubation unbound Streptavidin-HRP is removed during a wash step, and substrate solution reactive with HRP is added to the wells.

Figure 4



A coloured product is formed in proportion to the amount of human Erythropoietin present in the sample or standard. The reaction is terminated by addition of acid and absorbance is measured at 450 nm. A standard curve is prepared from 7 human Erythropoietin standard dilutions and human Erythropoietin sample concentration determined.

Figure 5



## 4 Reagents Provided

### 4.1 Reagents for human Erythropoietin ELISA NBP2-31082 (96 tests)

- 1 aluminium pouch with a **Microwell Plate coated** with monoclonal antibody to human Erythropoietin
- 1 vial (70 µl) **Biotin-Conjugate** monoclonal anti- Erythropoietin antibody
- 1 vial (150 µl) **Streptavidin-HRP**
- 2 vials human Erythropoietin **Standard** lyophilized, 200 mIU/ml upon reconstitution
- 1 vial **Control high**, lyophilized
- 1 vial **Control low**, lyophilized
- 1 vial (12 ml) **Sample Diluent**
- 1 vial (5 ml) **Assay Buffer Concentrate 20x**  
(PBS with 1% Tween 20, 10% BSA)
- 1 bottle (50 ml) **Wash Buffer Concentrate 20x**  
(PBS with 1% Tween 20)
- 1 vial (15 ml) **Substrate Solution** (tetramethyl-benzidine)
- 1 vial (15 ml) **Stop Solution** (1M Phosphoric acid)
- 4 **Adhesive Films**

## 4.2 Reagents for human Erythropoietin ELISA NBP2-31082 (10x96 tests)

- 10 aluminium pouches with a **Microwell Plate coated** with monoclonal antibody to human Erythropoietin
- 10 vials (70 µl) **Biotin-Conjugate** monoclonal anti-Erythropoietin antibody
- 10 vials (150 µl) **Streptavidin-HRP**
- 10 vials human Erythropoietin **Standard** lyophilized, 200 mIU/ml upon reconstitution
- 10 vials **Control high**, lyophilized
- 10 vials **Control low**, lyophilized
- 7 vials (12 ml) **Sample Diluent**
- 2 vials (5 ml) **Assay Buffer Concentrate 20x**  
(PBS with 1% Tween 20, 10% BSA)
- 6 bottles (50 ml) **Wash Buffer Concentrate 20x**  
(PBS with 1% Tween 20)
- 10 vials (15 ml) **Substrate Solution** (tetramethyl-benzidine)
- 1 vial (100 ml) **Stop Solution** (1M Phosphoric acid)
- 20 **Adhesive Films**



## 5 Storage Instructions – ELISA Kit

Store kit reagents between 2° and 8°C except controls. Store lyophilized controls at -20°C.

Immediately after use remaining reagents should be returned to cold storage (2° to 8°C), or to -20°C, respectively. Expiry of the kit and reagents is stated on labels.

Expiry of the kit components can only be guaranteed if the components are stored properly, and if, in case of repeated use of one component, this reagent is not contaminated by the first handling.

## 6 Specimen Collection and Storage Instructions

Cell culture supernatant, serum and plasma (EDTA, citrate, heparin), were tested with this assay. Other biological samples might be suitable for use in the assay. Remove serum or plasma from the clot or cells as soon as possible after clotting and separation.

Pay attention to a possible “**Hook Effect**” due to high sample concentrations (see chapter 11).

Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic specimens.

Samples should be aliquoted and must be stored frozen at -20°C to avoid loss of bioactive human Erythropoietin. If samples are to be run within 24 hours, they may be stored at 2° to 8°C (for sample stability refer to 13.5).

Avoid repeated freeze-thaw cycles. Prior to assay, the frozen sample should be brought to room temperature slowly and mixed gently.

## **7 Materials Required But Not Provided**

- 5 ml and 10 ml graduated pipettes
- 5  $\mu$ l to 1000  $\mu$ l adjustable single channel micropipettes with disposable tips
- 50  $\mu$ l to 300  $\mu$ l adjustable multichannel micropipette with disposable tips
- Multichannel micropipette reservoir
- Beakers, flasks, cylinders necessary for preparation of reagents
- Device for delivery of wash solution (multichannel wash bottle or automatic wash system)
- Microplate shaker
- Microwell strip reader capable of reading at 450 nm (620 nm as optional reference wave length)
- Glass-distilled or deionized water
- Statistical calculator with program to perform regression analysis

## 8 Precautions for Use

- All chemicals should be considered as potentially hazardous. We therefore recommend that this product is handled only by those persons who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. Wear suitable protective clothing such as laboratory overalls, safety glasses and gloves. Care should be taken to avoid contact with skin or eyes. In the case of contact with skin or eyes wash immediately with water. See material safety data sheet(s) and/or safety statement(s) for specific advice.
- Reagents are intended for research use only and are not for use in diagnostic or therapeutic procedures.
- Do not mix or substitute reagents with those from other lots or other sources.
- Do not use kit reagents beyond expiration date on label.
- Do not expose kit reagents to strong light during storage or incubation.
- Do not pipette by mouth.
- Do not eat or smoke in areas where kit reagents or samples are handled.
- Avoid contact of skin or mucous membranes with kit reagents or specimens.
- Rubber or disposable latex gloves should be worn while handling kit reagents or specimens.
- Avoid contact of substrate solution with oxidizing agents and metal.
- Avoid splashing or generation of aerosols.
- In order to avoid microbial contamination or cross-contamination of reagents or specimens which may invalidate the test use disposable pipette tips and/or pipettes.
- Use clean, dedicated reagent trays for dispensing the conjugate and substrate reagent.

- Exposure to acid inactivates the conjugate.
- Glass-distilled water or deionized water must be used for reagent preparation.
- Substrate solution must be at room temperature prior to use.
- Decontaminate and dispose specimens and all potentially contaminated materials as they could contain infectious agents. The preferred method of decontamination is autoclaving for a minimum of 1 hour at 121.5°C.
- Liquid wastes not containing acid and neutralized waste may be mixed with sodium hypochlorite in volumes such that the final mixture contains 1.0% sodium hypochlorite. Allow 30 minutes for effective decontamination. Liquid waste containing acid must be neutralized prior to the addition of sodium hypochlorite.

## 9 Preparation of Reagents

**Buffer Concentrates** should be brought to room temperature and should be diluted before starting the test procedure.

If crystals have formed in the **Buffer Concentrates**, warm them gently until they have completely dissolved.

### 9.1 Wash Buffer (1x)

Pour entire contents (50 ml) of the **Wash Buffer Concentrate (20x)** into a clean 1000 ml graduated cylinder. Bring to final volume of 1000 ml with glass-distilled or deionized water.

Mix gently to avoid foaming.

Transfer to a clean wash bottle and store at 2° to 25°C. Please note that Wash Buffer (1x) is stable for 30 days.

Wash Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Wash Buffer Concentrate (20x) (ml)	Distilled Water (ml)
1 - 6	25	475
1 - 12	50	950

### 9.2 Assay Buffer (1x)

Pour the entire contents (5 ml) of the **Assay Buffer Concentrate (20x)** into a clean 100 ml graduated cylinder. Bring to final volume of 100 ml with distilled water. Mix gently to avoid foaming.

Store at 2° to 8°C. Please note that the Assay Buffer (1x) is stable for 30 days.

Assay Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Assay Buffer Concentrate (20x) (ml)	Distilled Water (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

### 9.3 Biotin-Conjugate

**Please note that the Biotin-Conjugate should be used within 30 minutes after dilution.**

Make a 1:100 dilution of the concentrated **Biotin-Conjugate** solution with Assay Buffer (1x) in a clean plastic tube as needed according to the following table:

Number of Strips	Biotin-Conjugate (ml)	Assay Buffer (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

### 9.4 Streptavidin-HRP

**Please note that the Streptavidin-HRP should be used within 30 minutes after dilution.**

Make a 1:100 dilution of the concentrated **Streptavidin-HRP** solution with Assay Buffer (1x) in a clean plastic tube as needed according to the following table:

Number of Strips	Streptavidin-HRP (ml)	Assay Buffer (1x) (ml)
1 - 6	0.06	5.94
1 - 12	0.12	11.88

## 9.5 Human Erythropoietin Standard

Reconstitute **human Erythropoietin standard** by addition of distilled water. Reconstitution volume is stated on the label of the standard vial. Swirl or mix gently to insure complete and homogeneous solubilization (concentration of reconstituted standard = 200.0 mIU/ml). Allow the standard to reconstitute for 10-30 minutes. Mix well prior to making dilutions.

After usage remaining standard cannot be stored and has to be discarded.

**Standard dilutions** can be prepared directly on the microwell plate (see 10.c) or alternatively in tubes (see 9.5.1).

### 9.5.1 External Standard Dilution

Label 7 tubes, one for each standard point.

S1, S2, S3, S4, S5, S6, S7

Then prepare 1:2 serial dilutions for the standard curve as follows:  
Pipette 225 µl of Sample Diluent into each tube.

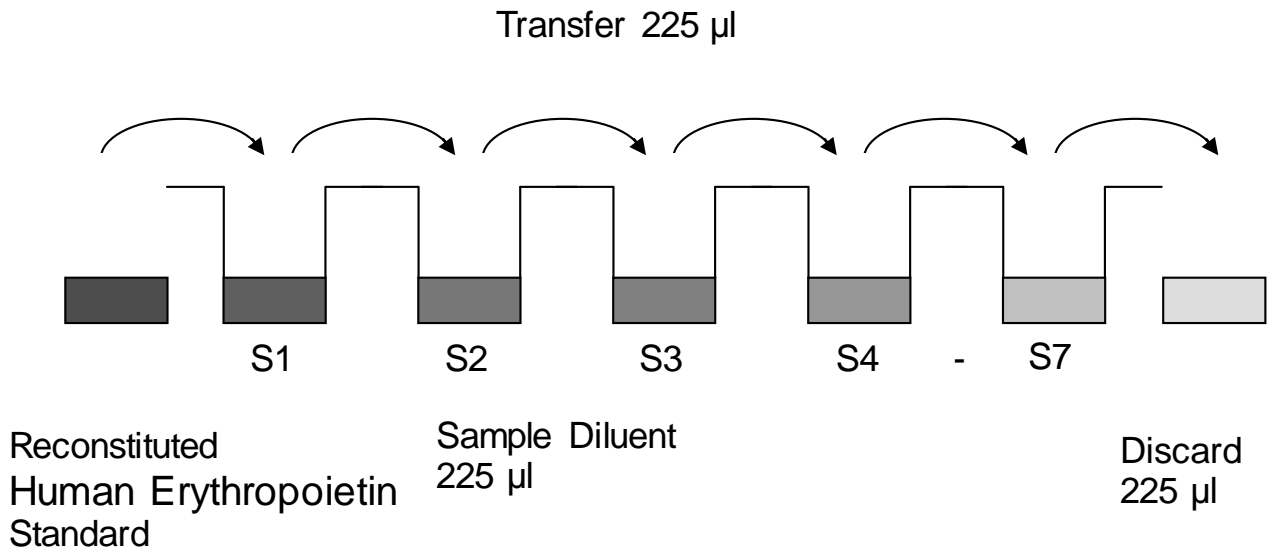
Pipette 225 µl of reconstituted standard (concentration of standard = 200.0 mIU/ml) into the first tube, labelled S1, and mix (concentration of standard 1 = 100.0 mIU/ml).

Pipette 225 µl of this dilution into the second tube, labelled S2, and mix thoroughly before the next transfer.

Repeat serial dilutions 5 more times thus creating the points of the standard curve (see Figure 6).

Sample Diluent serves as blank.

Figure 6



## 9.6 Controls

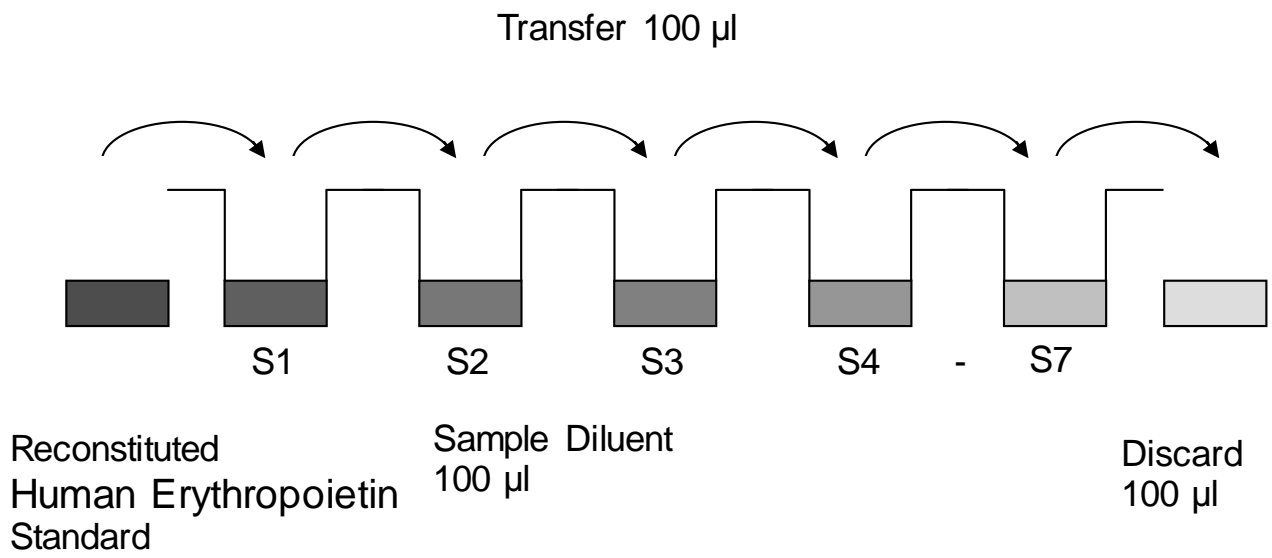
Reconstitute by adding 250  $\mu$ l distilled water to lyophilized controls (10-30 minutes). Swirl or mix gently to ensure complete and homogeneous solubilization. Further treat the controls like your samples in the assay. For control range please refer to certificate of analysis or vial label. Store reconstituted controls aliquoted at  $-20^{\circ}\text{C}$ . Avoid repeated freeze and thaw cycles



## 10 Test Protocol

- a. Determine the number of microwell strips required to test the desired number of samples plus appropriate number of wells needed for running blanks and standards. Each sample, standard, blank and optional control sample should be assayed in duplicate. Remove extra microwell strips from holder and store in foil bag with the desiccant provided at 2°-8°C sealed tightly.
- b. Wash the microwell strips twice with approximately 400 µl **Wash Buffer** per well with thorough aspiration of microwell contents between washes. Allow the Wash Buffer to sit in the wells for about **10 – 15 seconds** before aspiration. Take care not to scratch the surface of the microwells.  
After the last wash step, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess Wash Buffer. Use the microwell strips immediately after washing. Alternatively microwell strips can be placed upside down on a wet absorbent paper for not longer than 15 minutes. **Do not allow wells to dry.**
- c. **Standard dilution on the microwell plate** (Alternatively the standard dilution can be prepared in tubes - see 9.5.1):  
Add 100 µl of Sample Diluent in duplicate to all **standard wells**.  
Pipette 100 µl of prepared **standard** (see Preparation of Standard 9.5, concentration = 200.0 mIU/ml) in duplicate into well A1 and A2 (see Table 1). Mix the contents of wells A1 and A2 by repeated aspiration and ejection (concentration of standard 1, S1 = 100 mIU/ml), and transfer 100 µl to wells B1 and B2 (see Figure 7). Take care not to scratch the inner surface of the microwells.  
Continue this procedure 5 times, creating two rows of human Erythropoietin standard dilutions ranging from 100.0 to 1.6 mIU/ml. Discard 100 µl of the contents from the last microwells (G1, G2) used.

Figure 7



In case of an **external standard dilution** (see 9.5.1), pipette 100 µl of these standard dilutions (S1 - S7) in the standard wells according to Table 1.

Table 1

Table depicting an example of the arrangement of blanks, standards and samples in the microwell strips:

	<b>1</b>	<b>2</b>	<b>3</b>	<b>4</b>
<b>A</b>	Standard 1 (100.0 mIU/ml)	Standard 1 (100.0 mIU/ml)	Sample 1	Sample 1
<b>B</b>	Standard 2 (50.0 mIU/ml)	Standard 2 (50.0 mIU/ml)	Sample 2	Sample 2
<b>C</b>	Standard 3 (25.0 mIU/ml)	Standard 3 (25.0 mIU/ml)	Sample 3	Sample 3
<b>D</b>	Standard 4 (12.5 mIU/ml)	Standard 4 (12.5 mIU/ml)	Sample 4	Sample 4
<b>E</b>	Standard 5 (6.3 mIU/ml)	Standard 5 (6.3 mIU/ml)	Sample 5	Sample 5
<b>F</b>	Standard 6 (3.1 mIU/ml)	Standard 6 (3.1 mIU/ml)	Sample 6	Sample 6
<b>G</b>	Standard 7 (1.6 mIU/ml)	Standard 7 (1.6 mIU/ml)	Sample 7	Sample 7
<b>H</b>	Blank	Blank	Sample 8	Sample 8

- d. Add 100 µl of **Sample Diluent** in duplicate to the **blank wells**.
- e. Add 50 µl of **Sample Diluent** to the **sample wells**.
- f. Add 50 µl of each **sample** in duplicate to the **sample wells**.
- g. Prepare **Biotin-Conjugate** (see Preparation of Biotin-Conjugate 9.3).
- h. Add 50 µl of **Biotin-Conjugate** to all wells.
- i. Cover with an adhesive film and incubate at room temperature (18 to 25°C) for 1 hour, on a microplate shaker set at 400 rpm. (**Shaking is absolutely necessary for an optimal test performance**).
- j. Prepare **Streptavidin-HRP** (refer to Preparation of Streptavidin-HRP 9.4).
- k. Remove adhesive film and empty wells. **Wash** microwell strips 6 times according to point b. of the test protocol. Proceed immediately to the next step.
- l. Add 100 µl of diluted **Streptavidin-HRP** to all wells, including the blank wells.
- m. Cover with an adhesive film and incubate at room temperature (18° to 25°C) for 15 minutes, if available on a microplate shaker set at 400 rpm. (**Shaking is absolutely necessary for an optimal test performance**).
- n. Remove adhesive film and empty wells. **Wash** microwell strips 6 times according to point b. of the test protocol. Proceed immediately to the next step.
- o. Pipette 100 µl of **TMB Substrate Solution** to all wells.
- p. Incubate the microwell strips at room temperature (18° to 25°C) for about 10 min. Avoid direct exposure to intense light.

**The colour development on the plate should be monitored and the substrate reaction stopped (see next point of this protocol) before positive wells are no longer properly recordable. Determination of the ideal time period for colour development has to be done individually for each assay.**

It is recommended to add the stop solution when the highest standard has developed a dark blue colour. Alternatively the colour development can be monitored by the ELISA reader at 620 nm. The substrate reaction should be stopped as soon as Standard 1 has reached an OD of 0.9 – 0.95.

- q. Stop the enzyme reaction by quickly pipetting 100 µl of **Stop Solution** into each well. It is important that the Stop Solution is spread quickly and uniformly throughout the microwells to completely inactivate the enzyme. Results must be read immediately after the Stop Solution is added or within one hour if the microwell strips are stored at 2 - 8°C in the dark.
- r. Read absorbance of each microwell on a spectro-photometer using 450 nm as the primary wave length (optionally 620 nm as the reference wave length; 610 nm to 650 nm is acceptable). Blank the plate reader according to the manufacturer's instructions by using the blank wells. Determine the absorbance of both the samples and the standards.

**Shaking is absolutely necessary for an optimal test performance.**

## 11 Calculation of Results

- Calculate the average absorbance values for each set of duplicate standards and samples. Duplicates should be within 20 per cent of the mean value.
- Create a standard curve by plotting the mean absorbance for each standard concentration on the ordinate against the human Erythropoietin concentration on the abscissa. Draw a best fit curve through the points of the graph (a 5-parameter curve fit is recommended).
- To determine the concentration of circulating human Erythropoietin for each sample, first find the mean absorbance value on the ordinate and extend a horizontal line to the standard curve. At the point of intersection, extend a vertical line to the abscissa and read the corresponding human Erythropoietin concentration.
- **If instructions in this protocol have been followed, samples have been diluted 1:2 (50 µl sample + 50 µl Sample) and the concentration read from the standard curve must be multiplied by the dilution factor (x 2).**
- **Calculation of samples with a concentration exceeding standard 1 will result in incorrect, low human Erythropoietin levels (Hook Effect). Such samples require further external predilution according to expected human Erythropoietin values with Sample Diluent in order to precisely quantitate the actual human Erythropoietin level.**
- It is suggested that each testing facility establishes a control sample of known human Erythropoietin concentration and runs this additional control with each assay. If the values obtained are not within the expected range of the control, the assay results may be invalid.
- A representative standard curve is shown in Figure 8. This curve cannot be used to derive test results. Each laboratory must prepare a standard curve for each group of microwell strips assayed.

Figure 8

Representative standard curve for human Erythropoietin ELISA. Human Erythropoietin was diluted in serial 2-fold steps in Sample Diluent. Do not use this standard curve to derive test results. A standard curve must be run for each group of microwell strips assayed.

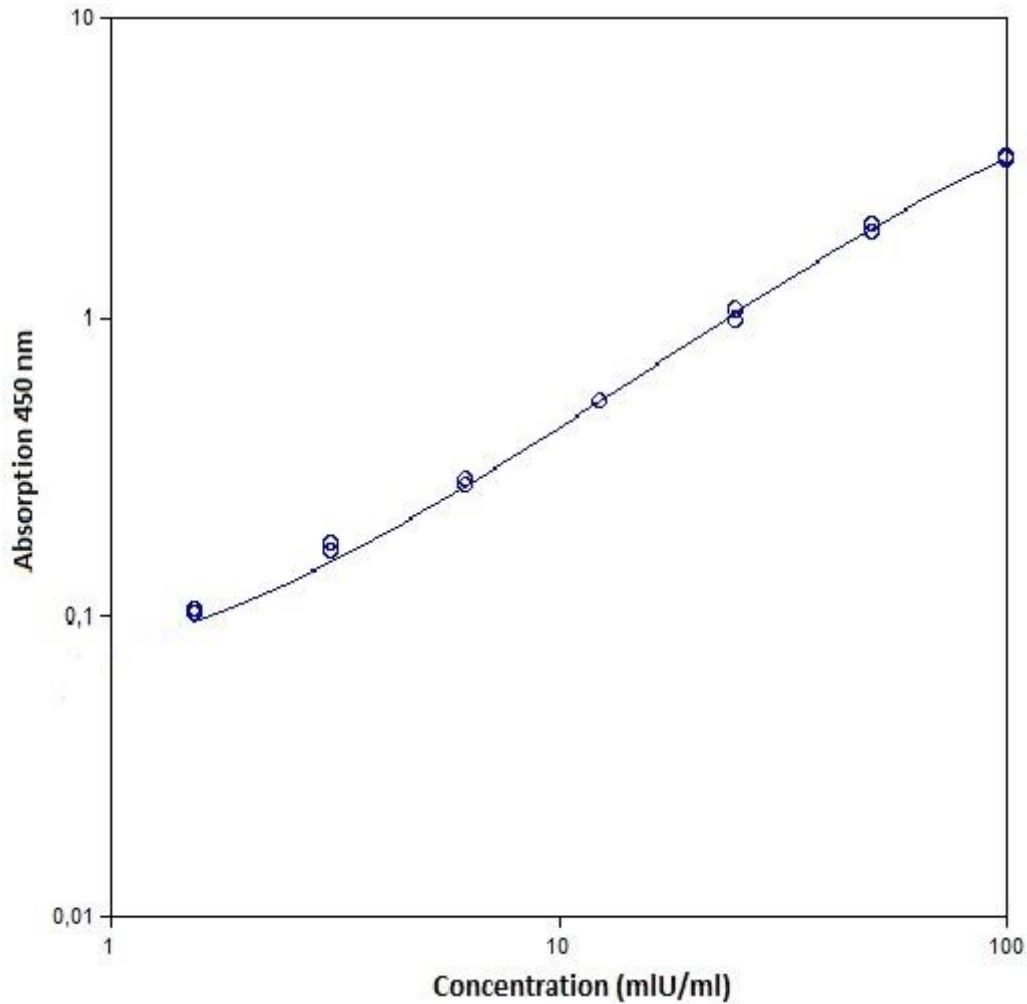


Table 2

Typical data using the human Erythropoietin ELISA

Measuring wavelength: 450 nm

Reference wavelength: 620 nm

Standard	Human Erythropoietin Concentration (mIU/ml)	O.D. at 450 nm	Mean O.D. at 450 nm	C.V. (%)
1	100.0	3.333 3.355	3.344	0.3%
2	50.0	1.881 2.014	1.948	3.4%
3	25.0	0.968 1.052	1.010	4.2%
4	12.5	0.514 0.515	0.515	0.1%
5	6.3	0.271 0.283	0.277	2.1%
6	3.1	0.162 0.174	0.168	3.5%
7	1.6	0.100 0.104	0.102	2.0%
Blank	0	0.028 0.029	0.028	2.0%

The OD values of the standard curve may vary according to the conditions of assay performance (e.g. operator, pipetting technique, washing technique or temperature effects). Furthermore shelf life of the kit may affect enzymatic activity and thus colour intensity. Values measured are still valid.



## 12 Limitations

- Since exact conditions may vary from assay to assay, a standard curve must be established for every run.
- Bacterial or fungal contamination of either screen samples or reagents or cross-contamination between reagents may cause erroneous results.
- Disposable pipette tips, flasks or glassware are preferred, reusable glassware must be washed and thoroughly rinsed of all detergents before use.
- Improper or insufficient washing at any stage of the procedure will result in either false positive or false negative results. Empty wells completely before dispensing fresh wash solution, fill with Wash Buffer as indicated for each wash cycle and do not allow wells to sit uncovered or dry for extended periods.
- The use of radioimmunotherapy has significantly increased the number of patients with human anti-mouse IgG antibodies (HAMA). HAMA may interfere with assays utilizing murine monoclonal antibodies leading to both false positive and false negative results. Serum samples containing antibodies to murine immunoglobulins can still be analysed in such assays when murine immunoglobulins (serum, ascitic fluid, or monoclonal antibodies of irrelevant specificity) are added to the sample.

## **13 Performance Characteristics**

### **13.1 Sensitivity**

The limit of detection of human Erythropoietin defined as the analyte concentration resulting in an absorbance significantly higher than that of the dilution medium (mean plus 2 standard deviations) was determined to be 0.17 mIU/ml (mean of 8 independent assays).

### **13.2 Reproducibility**

#### **13.2.1 Intra-assay**

Reproducibility within the assay was evaluated in 3 independent experiments. Each assay was carried out with replicates of 8 serum samples containing different concentrations of human Erythropoietin. 2 standard curves were run on each plate. Data below show the mean human Erythropoietin concentration and the coefficient of variation for each sample (see Table 3). The calculated overall intra-assay coefficient of variation was 3,4%.

Table 3

The mean human Erythropoietin concentration and the coefficient of variation for each sample

Sample	Experiment	Mean Human Erythropoietin Concentration (mIU/ml)	Coefficient of Variation (%)
1	1	25.5	2.5
	2	23.3	4.5
	3	24.6	3.0
2	1	25.1	3.4
	2	24.7	0.7
	3	25.0	3.3
3	1	58.3	2.2
	2	54.5	1.7
	3	55.3	2.8
4	1	65.4	4.0
	2	59.1	3.5
	3	60.1	3.7
5	1	8.6	3.3
	2	7.1	5.3
	3	8.3	7.2
6	1	11.0	2.1
	2	9.5	3.3
	3	10.2	5.0
7	1	16.6	4.8
	2	14.5	4.8
	3	15.0	4.6
8	1	7.5	2.1
	2	6.3	1.7
	3	7.7	2.8

### 13.2.2 Inter-assay

Assay to assay reproducibility within one laboratory was evaluated in 3 independent experiments. Each assay was carried out with replicates of 8 serum samples containing different concentrations of human Erythropoietin. 2 standard curves were run on each plate. Data below show the mean human Erythropoietin concentration and the coefficient of variation calculated on 18 determinations of each sample (see Table 4). The calculated overall inter-assay coefficient of variation was 6.2%.

Table 4

The mean human Erythropoietin concentration and the coefficient of variation of each sample

Sample	Mean Human Erythropoietin Concentration (mIU/ml)	Coefficient of Variation (%)
1	24	4.4
2	25	0.9
3	56	3.6
4	62	5.6
5	8	9.7
6	10	7.0
7	15	7.1
8	7	10.9

### 13.3 Spike Recovery

The spike recovery was evaluated by spiking 3 levels of human Erythropoietin into serum, plasma and cell culture supernatant samples. Recoveries were determined with 2 replicates each. The amount of endogenous human Erythropoietin in unspiked serum was subtracted from the spike values. For recovery data see Table 5.

Table 5

Sample matrix	Spike high		Spike medium		Spike low	
	Mean (%)	Range (%)	Mean (%)	Range (%)	Mean (%)	Range (%)
Serum	110	105 – 117	109	98 – 120	101	84 – 114
Plasma (EDTA)	128	122 – 133	106	104 – 107	99	85 – 113
Plasma (Heparin)	106		114		119	
Plasma (citrate)	97	97 – 97	101	99 – 104	109	107 – 112
Cell culture supernatant	136	128 – 144	135	127 - 144	139	128 – 150

### 13.4 Dilution Parallelism

Serum samples with different levels of human Erythropoietin were analysed at serial 2 fold dilutions with 4 replicates each.

The recovery ranged from 97.4% to 119.2% with an overall recovery of 105.2% (see Table 6).

Table 6

Sample matrix	Recovery of Exp. Val.		
	Dilution	Mean (%)	Range (%)
Serum	1:4	87	82 - 91
	1:8	90	89 - 92
	1:16	99	92 - 105
Plasma (EDTA)	1:4	85	82 - 89
	1:8	88	82 - 101
	1:16	92	71 - 122
Plasma (heparin)	1:4	89	85 - 92
	1:8	85	81 - 90
	1:16	83	78 - 91
Plasma (citrate)	1:4	98	95 - 102
	1:8	96	92 - 103
	1:16	99	94 - 103
Cell culture supernatant	1:4	79	75 - 83
	1:8	74	71 - 77
	1:16	71	66 - 75

## **13.5 Sample Stability**

### **13.5.1 Freeze-Thaw Stability**

Aliquots of serum samples (spiked or unspiked) were stored at -20°C and thawed 3 times, and the human Erythropoietin levels determined. There was no significant loss of human Erythropoietin immunoreactivity detected by freezing and thawing.

### **13.5.2 Storage Stability**

Aliquots of serum samples (spiked or unspiked) were stored at 2-8°C, room temperature (RT) and at 37°C, and the human Erythropoietin level determined after 24 h. There was no significant loss of human Erythropoietin immunoreactivity detected during storage under above conditions.

## **13.6 Specificity**

The assay detects both natural and recombinant human Erythropoietin. The interference of circulating factors of the immune system was evaluated by spiking these proteins at physiologically relevant concentrations into serum. There was no crossreactivity detected.

### 13.7 Expected Values

Panels of 40 serum as well as EDTA, citrate and heparin plasma samples from randomly selected apparently healthy donors (males and females) were tested for human Erythropoietin.

The levels measured may vary with the sample collection used. For detected human Erythropoietin levels see Table 7.

Table 7

Sample Matrix	Number of Samples Evaluated	Range (mIU/ml)	Mean (mIU/ml)	Standard Deviation (mIU/ml)
Serum	40	2.7 – 40.7	12.9	9.6
Plasma (EDTA)	40	1.6 – 57.2	8.5	10.3
Plasma (Citrate)	40	0.9 – 28.6	4.7	5.1
Plasma (Heparin)	40	2.1 – 41.9	13.7	9.0



## 14 Reagent Preparation Summary

### 14.1 Wash Buffer (1x)

Add **Wash Buffer Concentrate 20x** (50 ml) to 950 ml distilled water.

Number of Strips	Wash Buffer Concentrate (ml)	Distilled Water (ml)
1 - 6	25	475
1 - 12	50	950

### 14.2 Assay Buffer (1x)

Add **Assay Buffer Concentrate 20x** (5 ml) to 95 ml distilled water.

Number of Strips	Assay Buffer Concentrate (ml)	Distilled Water (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

### 14.3 Biotin-Conjugate

Make a 1:100 dilution of **Biotin-Conjugate** in Assay Buffer (1x):

Number of Strips	Biotin-Conjugate (ml)	Assay Buffer (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

### 15.4 Streptavidin-HRP

Make a 1:100 dilution of **Streptavidin-HRP** in Assay Buffer (1x):

Number of Strips	Streptavidin-HRP (ml)	Assay Buffer (1x) (ml)
1 - 6	0.06	5.94
1 - 12	0.12	11.88

### 14.5 Human Erythropoietin Standard

Reconstitute lyophilized **human Erythropoietin standard** with distilled water. (Reconstitution volume is stated on the label of the standard vial.)

### 14.6 Controls

Add 250 µl distilled water to lyophilized controls.

## 15 Test Protocol Summary

1. Determine the number of microwell strips required.
2. Wash microwell strips twice with Wash Buffer.
3. Standard dilution on the microwell plate: Add 100 µl Sample Diluent, in duplicate, to all standard. Pipette 100 µl prepared standard into the first wells and create standard dilutions by transferring 100 µl from well to well. Discard 100 µl from the last wells.  
Alternatively external standard dilution in tubes (see 9.5.1): Pipette 100 µl of these standard dilutions in the microwell strips.
4. Add 100 µl Sample Diluent in duplicate, to the blank wells.
5. Add 50 µl Sample Diluent to sample wells.
6. Add 50 µl sample in duplicate, to designated sample wells.
7. Prepare Biotin-Conjugate.
8. Add 50 µl Biotin-Conjugate to all wells.
9. Cover microwell strips and incubate 1 hour at room temperature (18° to 25°C).
10. Prepare Streptavidin-HRP.
11. Empty and wash microwell strips 6 times with Wash Buffer.
12. Add 100 µl diluted Streptavidin-HRP to all wells.
13. Cover microwell strips and incubate 15 minutes at room temperature (18° to 25°C).
14. Empty and wash microwell strips 6 times with Wash Buffer.
15. Add 100 µl of TMB Substrate Solution to all wells.
16. Incubate the microwell strips for about 10 minutes at room temperature (18° to 25°C).
17. Add 100 µl Stop Solution to all wells.
18. Blank microwell reader and measure colour intensity at 450 nm.

**Note: If instructions in this protocol have been followed, samples have been diluted 1:2 (50 µl sample + 50 µl Sample Diluent) and the concentration read from the standard curve must be multiplied by the dilution factor (x 2).**

