



Version 1.3.0

**EPO ELISA KIT**Immunoassay for Human  
Erythropoietin

Catalog #01630

1 kit

**PRODUCT DESCRIPTION:**

This kit is meant for the quantitative measurement of natural and recombinant human erythropoietin (EPO) in biological fluids, such as serum, plasma and tissue culture supernatants. The EPO ELISA utilizes two monospecific monoclonal antibodies raised against human urinary EPO. These antibodies bind two non-overlapping epitopes on the EPO polypeptide and show high-affinity binding to both natural and recombinant EPO.

*The ELISA is intended for measurement of human EPO only. It has not been validated for detection of EPO from other species.*

**COMPONENTS:**

The following is a list of EPO ELISA kit components:

COMPONENT	DESCRIPTION	SIZE	CATALOG #
<b>Buffer A</b>	Phosphate buffered saline containing protein, detergent and 0.1% (v/v) Proclin	3 mL	01600
<b>Buffer B</b>	Phosphate buffered saline containing protein, detergent and 0.1% (v/v) Proclin	50 mL	01601
<b>Biotinylated Anti-EPO Antibody</b>	Diluted in 7 mL Buffer B.	7 mL	01602
<b>Streptavidin-Peroxidase Conjugate</b>	Diluted in HRP stabilizing buffer.	12.5 mL	01603
<b>TMB Substrate Solution</b>	Tetramethylbenzidine and hydrogen peroxide solution	12.5 mL	01604
<b>Stop Solution</b>	0.5N sulphuric acid*	12 mL	01605
<b>10X Wash Buffer</b>	Concentrate of phosphate buffered saline containing detergent	100 mL	01606
<b>EPO Standard</b>	100 mU/mL	1 mL	01607
<b>EPO Standard</b>	50 mU/mL	1 mL	01608
<b>EPO Standard</b>	25 mU/mL	1 mL	01609
<b>EPO Standard</b>	12.5 mU/mL	1 mL	01610
<b>EPO Standard</b>	6.25 mU/mL	1 mL	01611
<b>EPO Standard</b>	3.12 mU/mL	1 mL	01612
<b>EPO Standard</b>	1.56 mU/mL	1 mL	01613
<b>EPO Standard</b>	0 mU/mL	1 mL	01614
<b>Microtiter Plate</b>	96 wells (12 strips of 8 wells) pre-coated with anti-EPO monoclonal antibody	1 plate	01615
<b>Adhesive Covers</b>	Plate covers	4 covers	01616
<b>Technical Manual</b>	Assay Procedure	1 manual	28379

**STABILITY AND STORAGE:**

Each component is stable for at least 1 year after the date of manufacture indicated on each vial when stored at 2-8°C.

**Do not freeze.**

**DIRECTIONS FOR USE:**

Please refer to the **Human Erythropoietin ELISA Assay Procedure Technical Manual** (Catalog #28379) for a detailed description of the assay principle, intended use and procedure. The manual is available on our website at [www.stemcell.com/technical/manuals.asp](http://www.stemcell.com/technical/manuals.asp).

**THIS REAGENT IS FOR RESEARCH ONLY.  
IT IS NOT TO BE ADMINISTERED TO HUMANS.**

*\*Warning: Sulphuric acid is corrosive. Handle with care!*

**StemCell Technologies**

In North America  
Tel: 1.604.877.0713  
Fax: 1.604.877.0704  
Toll Free Tel: 1.800.667.0322  
Toll Free Fax: 1.800.567.2899  
e-mail: [info@stemcell.com](mailto:info@stemcell.com)  
[www.stemcell.com](http://www.stemcell.com)

In the United Kingdom  
Tel: +44.(0).20.7691.3561  
Fax: +33.(0).4.76.18.99.63  
Toll Free within United Kingdom:  
Tel: 0800.731.27.14  
Fax: 0800.731.27.13  
e-mail: [info@stemcellgb.com](mailto:info@stemcellgb.com)

In Europe  
Tel: +33.(0).4.76.04.75.30  
Fax: +33.(0).4.76.18.99.63  
e-mail: [info@stemcellfrance.com](mailto:info@stemcellfrance.com)

Revised:  
August 2007