



**EAGLE**  
BIOSCIENCES

# **C-Reactive Protein HS ELISA**

Catalog Number: CRP31-K01

*For Research Use Only. Not for use in diagnostic procedures.*

*v. 1 (13 FEB 24)*

---

EAGLE BIOSCIENCES, INC.  
20A Northwest Blvd., Suite 112, Nashua, NH 03063  
Phone: 617-419-2019 Fax: 617-419-1110  
[WWW.EAGLEBIO.COM](http://WWW.EAGLEBIO.COM)



## INTENDED USE

The C-Reactive Protein HS ELISA is intended for the quantitative determination of C-reactive protein (CRP) concentration in serum .

*For further information about this kit, its application, or the procedures in this insert, please contact the Technical Service Team at Eagle Biosciences, Inc at [www.EagleBio.com](http://www.EagleBio.com) or at 866-411-8023.*

## ASSAY BACKGROUND

The C-Reactive Protein HS ELISA is based on the principle of a solid phase enzyme-linked immunosorbent assay. The assay system utilizes a unique monoclonal antibody directed against a distinct antigenic determinant on the on the CRP molecule. This mouse monoclonal anti-CRP antibody is used for solid phase immobilization (on the microtiter wells). A goat anti-CRP antibody is in the antibody-enzyme (horseradish peroxidase) conjugate solution. The test sample is allowed to react simultaneously with the two antibodies, resulting in the CRP molecules being sandwiched between the solid phase and enzyme-linked antibodies. After a 45-minute incubation at room temperature, the wells are washed with water to remove unbound labeled antibodies. A tetramethylbenzidine (TMB) reagent is added and incubated for 20 minutes, resulting in the development of blue color. The color development is stopped with the addition of 1 N HCl changing the color to yellow.

The concentration of CRP is directly proportional the color intensity of the test sample. Absorbance is measured spectrophotometrically at 450 nm.

## MATERIALS PROVIDED

- **Antibody-coated wells** ( 1 plate, 96 wells)  
Microtiter wells coated with mouse monoclonal anti-CRP
- **Reference Standard Set** (1.0 mL/vial)  
Contains 0 – 0.005 – 0.010 – 0.025 – 0.050 – and 1.00 mg/L CRP in serum based buffer with preservatives
- **HsCRP Sample Diluent** (50 mL/vial)  
Contains phosphate buffer-BSA solution with preservatives.
- **CRP Enzyme Conjugate Reagent** ( 12 mL/vial)  
Contains goat anti-CRP conjugated t horseradish peroxidase with preservatives
- **TMB Reagent** (11 mL/Bottle)  
Contains one-step TMB solution
- **Stop Solution** (11mL/bottle)  
Contains diluted hydrochloric acid (1N HCl)

## MATERIALS REQUIRED BUT NOT PROVIDED

- Precision pipettes 5  $\mu$ L, 10 $\mu$ L, 50 $\mu$ L, 100 $\mu$ L and 1.0 mL
- Disposable pipette tips
- Microtiter well reader capable of reading absorbance at 450nm
- Vortex mixer, or equivalent
- Absorbent paper



- Graph paper

### STORAGE CONDITONS

- Store the unopened kit at 2-8C upon receipt and when it is not in use, until the expiration shown on the kit label. Refer to the package label for expiration date.
- Keep microtiter plate in sealed bag with desiccant to minimize exposure to damp air.

### ASSAY PREPARATIONS

- All reagents should be allowed to reach room temperature ( 18°C – 25°C )before use.
- Serum should be diluted 100-fold prior to use. Prepare a series of small tubes (i.e 1.5mL microcentrifuge tubes) and mix 5µL of serum with 495 µL (0.495mL) Sample Diluent.

#### **DO NOT DILUTE THE STANDARDS**

- Samples with expected CRP concentration over 10mg/L may be quantitated by further dilution (10 fold) of the 100-fold diluted solution with sample diluent. (i.e., 10 µL of the 100-fold diluted sample to 90µL sample diluent)

### INSTRUMENTATION

A microtiter well reader with a bandwidth of 10nm or less and an optical density range of 0 to 3 OD or greater at 450nm wavelength is acceptable for absorbance measurement.

### ASSAY PROCEDURE

1. Serum and control serum should be diluted **100 fold** prior to use. See Reagent Preparation. **PLEASE DO NOT DILUTE THE STANDARDS**
2. Secure the desired number of coated wells in the holder
3. Dispense **10µL** of CRP Standards, **DILUTED** Specimens, and **DILUTED** controls into appropriate wells.
4. Dispense **100 µL** of CRP Enzyme Conjugate Reagent into each well.
5. Thoroughly mix for 30 seconds. It is very important to mix completely.
6. Incubate at room temperature (18°C-25°C) for 45 minutes
7. Remove the incubation mixture by flicking plate contents into a waste container. Rinse and flick the microtiter wells 5 times with deionized water. **DO NOT USE TAP WATER**
8. Strike the wells sharply onto absorbent paper or paper towels to remove all residual water droplets.
9. Dispense **100µL** TMB solution into each well. Gently mix for 10 seconds.
10. Incubate at room temperature for 20 minutes.
11. Stop the reaction by adding **100 µL** of Stop solution to each well.
12. Gently mix for 30 seconds. **It is important to make sure that all the blue color changes to yellow color completely.**
13. Read absorbance at 450nm with microtiter well reader **within 15 minutes.**



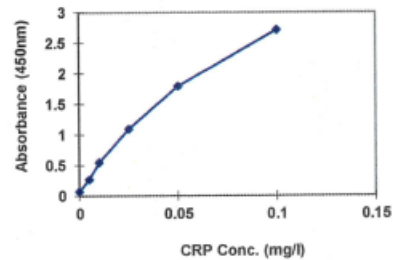
## CALCULATION OF RESULTS

1. Calculate the mean absorbance value ( $OD_{450}$ ) for each set of reference standards, controls and samples.
2. Construct a standard curve by plotting the mean absorbance obtained for each reference standard against its concentration in mg/L on graph paper, with absorbance on the vertical (y) axis and concentration on the horizontal (x) axis
3. Using the mean absorbance value for each sample, determine the corresponding concentration of CRP (mg/L) from the standard curve. Depending on the experience and/or the availability of computer capability, other methods of data reduction may be employed.
4. The obtained values of the samples and control sera should be multiplied by the dilution factor of 100 to obtain CRP results in mg/L
5. Samples with CRP concentration greater than 10 mg/L should be further diluted 10-fold after the initial 100-fold dilution (total dilution 1:1000), and the final CRP values should be multiplied by 1,000 to obtain CRP results in mg/L.
6. NOTE: Samples with CRP concentration less than 0.1 mg/L should be reported as "<0.1mg/L CRP".

## EXAMPLE STANDARD CURVE

Results of a typical standard run with absorbance readings at 450nm shown on the Y axis against CRP concentrations shown on the x axis. **NOTE:** This standard curve is for the purpose of illustration only, and should not be used to calculate unknowns. Each laboratory must provide its own data and standard curve in each experiment.

CRP (mg/L)	Absorbance (450 nm)
0	0.066
0.005	0.264
0.010	0.457
0.025	1.092
0.050	1.788
0.100	2.710





## **WARRANTY INFORMATION**

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 non-infringement 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, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.*