



EAGLE
BIOSCIENCES

Sheep IgG ELISA Assay Kit

Catalog Number:

IGS99-K01 (1 x 96 wells)

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

v. 1.0

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



Introduction

The Eagle Biosciences Sheep IgG ELISA Assay Kit is designed for the quantitative determination of sheep IgG in complex samples (serum or other biological samples). The Sheep IgG ELISA Assay Kit is for research use only and should not be used in diagnostic procedures.

Principle of the Assay

The determination of sheep IgG is carried out as direct sandwich ELISA. An antibody specific for sheep IgG has been pre-coated onto a microplate. Standards and samples are pipetted into the wells and any IgG present is bound. After washing away any unbound substances, an enzyme-linked antibody is added. Following a wash, a substrate solution is added to the wells and color develops in proportion to the amount of antibody conjugate. The absorption at 450 nm is proportional to the IgG concentration.

Precautions

- Store the Sheep IgG ELISA Assay Kit at 2-8 °C.
- The Sheep IgG ELISA Assay Kit is for research use only. Not for use in diagnostic procedures.
- Do not use the reagents beyond the expiration date marked on box label.
- Please read the instructions carefully before using the Sheep IgG ELISA Assay Kit.
- Do not mix reagents from different lots.
- Some components of this Sheep IgG ELISA Assay Kit contain Thimerosal, a mercury containing compound. Follow routine precautions for handling hazardous chemicals. Avoid contact with skin and mucous membranes when handling reagents, which contain preservatives (see materials provided). Wash thoroughly with water in case of contact and possibly look up a doctor.
- The stop solution contains 0.5 M sulphuric acid. Wash thoroughly with water in case of contact with skin. In case of contact with eyes rinse with much water and look up a doctor.
- Do not allow the wells to become dry once the assay has begun.

Other supplies required

- Deionized or distilled water
- Graduated cylinder
- Micropipettes, multipipette
- Microplate reader

Preparation of reagents and samples

- Bring all reagents of the Sheep IgG ELISA Assay Kit to room temperature before use. If crystals have formed, mix gently until the crystals have completely dissolved.

- The microplate strips are ready to use. Remove excess strips (breakable) from the frame, reseal in the bag with the desiccant and store at 2-8 °C

- Dilute the wash buffer with deionized or distilled water **1:10** (e. g. 40 ml + 360 ml water). The diluted solution is stable for 30 days at 2-8 °C.



- Use the Standard concentrate to produce a 1:2-dilution series with diluent (e. g. 250 μ l + 250 μ l diluent):

Standard	Preparation	Conc. (ng/ml)
S 7	standard conc. Undiluted	300
S 6	S 7 1:2 diluted	150
S 5	S 6 1:2 diluted	75
S 4	S 5 1:2 diluted	37.5
S 3	S 4 1:2 diluted	18.75
S 2	S 3 1:2 diluted	9.38
S 1	S 2 1:2 diluted	4.69

- Dilute the HRP conjugate with diluent **1:100** (e. g. 50 μ l + 4950 μ l diluent). The required amount of conjugate solution should be prepared freshly.

Assay procedure

It is recommended that all samples and standards be assayed in duplicate.

1. Prepare all reagents, standard curve and samples as directed in the previous section.
2. Pipette 100 μ l of samples, standards, or diluent (as negative control) into the wells.
3. Seal wells with adhesive strip and incubate for 1 hour at room temperature with shaking.
4. Aspirate fluid from wells and wash three times with 300 μ l wash buffer. After the last wash, invert the plate and tap on a clean paper towel.
5. Add 100 μ l of HRP conjugate to each well.
6. Seal wells with adhesive strip and incubate for 1 hour at room temperature with shaking.
7. Repeat the wash as in step 4.
8. Dispense 100 μ l of TMB substrate solution into each well.
9. Incubate for 30 minutes at room temperature in the dark.
10. Add 100 μ l of stop solution to each well.
11. Determine the absorbance within 30 minutes at 450 nm. A reference wavelength of 620 nm/690 nm is recommended.

Calculation of results

Create a standard curve for the Sheep IgG ELISA Assay Kit using computer software capable of generating a curve fit (e.g. x-axis: log, IgG concentration; y-axis: linear, absorbance; four parameter curve fit). As an alternative, draw a standard curve on semi-log paper (x-axis: log, IgG concentration; y-axis: linear, absorbance). The IgG concentrations can be calculated from the standard curve. The calculated concentrations must be multiplied by the sample dilution factor.



If the absorbance of some samples is outside the standard curve a subsequent determination with changed dilutions will provide a proper result.

Materials provided:

Number of determinations	1x96 determinations
Microplate strips, antibody coated	12 x 8
Wash buffer, 10fold conc. ◆	50 ml
Diluent, ready to use ◆	100 ml
Standard concentrate ◆	1 ml
Anti-IgG(sheep)-Ab., HRP conjugate, 100fold conc. ◆	0.2 ml
TMB substrate, ready to use	12 ml
Stop solution, ready to use (0.5 M sulphuric acid)	12 ml

◆: contains Thimerosal

Assay procedure summary:

A. Preparation

1. Bring all reagents to room temperature
2. Dilute wash buffer 1:10
3. Prepare the standard curve from a 1:2-dilution series of standard concentrate with diluent
4. Dilute samples with diluent
5. Dilute freshly HRP conjugate 1:100 with diluent

B. Performance

1. Pipette 100 μ l of samples, standards, controls into the wells
2. Incubate for 1 hour at room temperature with shaking
3. Wash three times with 300 μ l of wash buffer
4. Add 100 μ l of HRP conjugate to each well
5. Incubate for 1 hour at room temperature with shaking
6. Wash three times with 300 μ l of wash buffer
7. Dispense 100 μ l of TMB substrate solution
8. Incubate for 30 minutes at room temperature in the dark
9. Add 100 μ l of stop solution
10. Measure absorption at 450 nm



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 or at 866-411-8023.