

# Bovine IgG ELISA Kit ab205078

[2 References](#) [1 Image](#)

### Overview

**Product name** Bovine IgG ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

**Sample type** Cell culture supernatant, Milk, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 1.245 ng/ml

**Range** 3.125 ng/ml - 100 ng/ml

**Recovery** > 85 %

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	3.125ng/ml - 100ng/ml

**Assay time** 0h 50m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Cow

**Does not react with:** Mouse, Rat, Rabbit, Human

**Product overview** IgG in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IgG protein in bovine cell culture supernatant, milk, plasma and serum.

**Notes** In this assay the IgG present in samples reacts with the anti-IgG antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-IgG antibodies conjugated with horseradish peroxidase (HRP), are added. These

enzyme-labeled antibodies form complexes with the previously bound IgG. 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 IgG in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of IgG in the test sample. The quantity of IgG in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X HRP Conjugated Enzyme Antibody	1 x 150µl
20X Diluent Concentration	1 x 50ml
20X Wash Buffer Concentrate	1 x 50ml
Anti-Bovine IgG EIISA Microplate	1 unit
Bovine IgG Calibrator (Lyophilized)	1 vial
Chromogen Substrate Solution	1 x 12ml
Stop Solution	1 x 12ml

**Relevance** Immunoglobulin G (IgG) are antibody molecules. Each IgG is composed of four peptide chains - two heavy chains and two light chains. Each IgG has two antigen binding sites. Other Immunoglobulins may be described in terms of polymers with the IgG structure considered the monomer.

## Images



Representative standard curve using ab205078 Bovine IgG ELISA Kit.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors