

Horse IgG ELISA Kit ab190531

1 Image

Overview

Product name Horse 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 Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 2.217 ng/ml

Range 6.25 ng/ml - 400 ng/ml

Recovery

Sample specific recovery

| Sample type | Average % | Range |
|-------------|-----------|-------|
| Serum | > 85 | % - % |

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Horse

Product overview

Abcam's IgG Horse ELISA kit (ab190531) is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of IgG in horse serum and plasma.

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.

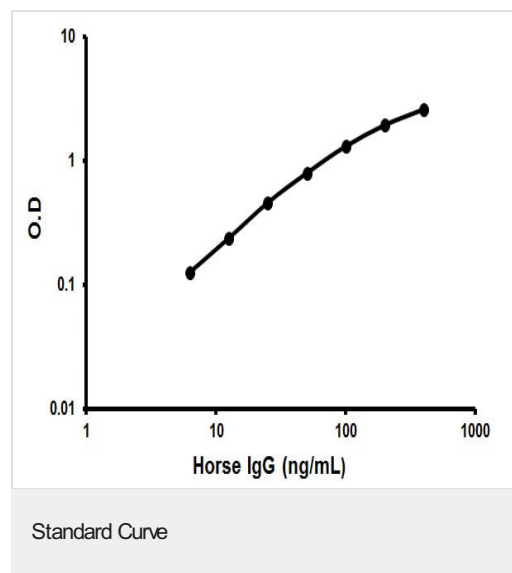
Platform Microplate

Properties

| Components | 1 x 96 tests |
|--------------------------------------|--------------|
| 20X Wash Buffer Concentrate | 1 x 50ml |
| 5X Diluent Concentrate | 1 x 50ml |
| Chromogen Substrate Solution | 1 x 12ml |
| IgG Horse Antibody coated microwells | 1 unit |
| IgG Horse Calibrator (Lyophilized) | 1 vial |
| IgG Horse HRP Conjugate | 1 vial |
| Stop Solution | 1 x 12ml |

Cellular localization Secreted

Images



Representative standard curve using ab190531 IgG Horse 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