

Product datasheet

Guinea Pig IgG ELISA Kit ab190525

1 Image

Overview

Product name Guinea Pig IgG ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Serum				< 10%

Sample type

Serum

Assay type

Sandwich

Sensitivity

= 1.069 ng/ml

Range

6.25 ng/ml - 400 ng/ml

Recovery

> 85 %

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	6.25ng/ml - 400ng/ml

Assay duration

Multiple steps standard assay

Species reactivity

Reacts with: Guinea pig

Product overview

Abcam's IgG Guinea pig ELISA Kit (ab190525) is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of IgG in serum.

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.

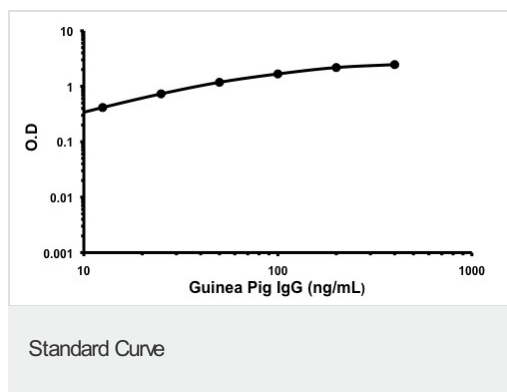
Platform Microplate

Properties

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
IgG Guinea Pig Antibody Coated Microplate	1 x 96 tests
IgG Guinea Pig Calibrator (lyophilized)	1 vial
IgG Guinea Pig HRP Conjugate	1 vial
Stop Solution	1 x 12ml
Chromogen Substrate Solution	1 x 12ml

Cellular localization Secreted

Images



Representative standard curve using ab190525 IgG Guinea pig 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