

2 References

Overview

Product name	Chicken IgY 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	1.5305 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:** Chicken

Product overview Abcam's IgY Chicken ELISA kit is a highly sensitive two-site enzyme linked immunoassay (ELISA) for the quantitative measurement of IgY in chicken serum or plasma samples.

In this assay the IgY present in samples reacts with the anti-IgY antibodies which have been adsorbed to the surface of polystyrene microtiter wells. After the removal of unbound proteins by washing, anti- IgY antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound IgY. 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 IgY in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of IgY in the test sample. The quantity of IgY in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Get results in 90 minutes with Chicken IgY ELISA Kit ([**ab189577**](#)) from our SimpleStep ELISA[®] range.

Platform

Microplate

Properties

Storage instructions

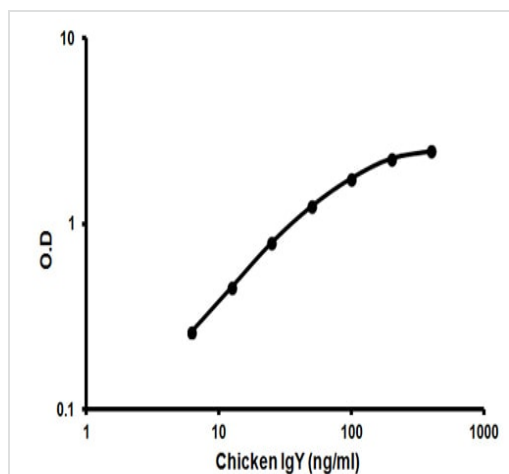
Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X HRP-conjugated anti-chicken IgY antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chicken IgY Calibrator (Lyophilized)	1 vial
Chicken IgY ELISA Microplate	1 unit
Chromogen Substrate Solution	1 x 12ml
Stop Solution	1 x 12ml

Relevance

Secreted immunoglobulins have two separable functions: one is to bind specifically to pathogenic molecules, the other is to recruit cells and molecules to destroy the opsonized pathogen. IgY has a tetrameric structure consisting of two identical light chains ($M_r = 24$ kDa) and two identical heavy chains ($M_r = 73$ kDa).

Images



Representative standard curve using ab157693 IgY Chicken ELISA Kit.

Standard Curve

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