

Mouse IgG1 ELISA Kit ab133045

★★★★☆ [1 Abreviews](#) [8 References](#) [1 Image](#)

Overview

Product name Mouse IgG1 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Low	19	49.9ng/ml		1.2%
Medium	19	99.6ng/ml		2.7%
High	19	160.4ng/ml		3.6%

Inter-assay

Sample	n	Mean	SD	CV%
Low	16	48.5ng/ml		3.1%
Medium	16	98.3ng/ml		4.4%
High	16	148.7ng/ml		8.1%

Sample type Cell culture supernatant, Serum

Assay type Sandwich (quantitative)

Sensitivity 0.116 ng/ml

Range 7.81 ng/ml - 250 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	102.8	% - %
Tissue Culture Media	105.3	% - %

Assay time 0h 90m

Assay duration

Multiple steps standard assay

Species reactivity**Reacts with:** Mouse**Product overview**

Abcam's Mouse IgG₁ *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Mouse IgG₁ in Culture supernatants and Serum.

IgG₁ specific antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and along with an HRP-conjugated IgG₁ detection antibody and the microplate is then incubated at room temperature. After the removal of unbound proteins by washing, TMB is used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a colored product that changes after adding acidic stop solution. The density of coloration is directly proportional to the IgG₁ amount of sample captured in plate.

Notes

IgG is divided into four subclasses; IgG1, IgG2, IgG3, and IgG4. IgG₁ is the most abundant immunoglobulin found in the blood. It is a glycoprotein which consists of two identical heavy chains (50 kDa each) and two identical light chains (25 kDa each), to give a combined mass of approximately 150 kDa. The chains are held in place by covalent disulfide bonds. Each light chain contains two immunoglobulin (Ig) domains, while the heavy chains contain four Ig domains each. In the middle of each heavy chain is a relative varying portion called the "hinge region" which is unique to each IgG. This region allows for molecular flexibility and sets IgG₁ apart from its IgG counterparts. IgG1 properties and functions include neutralization, opsonization, activation of the complement system, diffusion into extravascular sites and crossing the placenta.

Cross Reactivity

Compound	Cross Reactivity
Rat IgG1	0.9%
Mouse IgG2b	0.21
Human IgG1	<0.01%
Mouse IgG2a	<0.01%
Mouse IgG3	<0.01%
Mouse IgM	<0.01%

Recommended dilution of serum is 1:20,000

Platform

Microplate

Properties**Storage instructions**

Please refer to protocols.

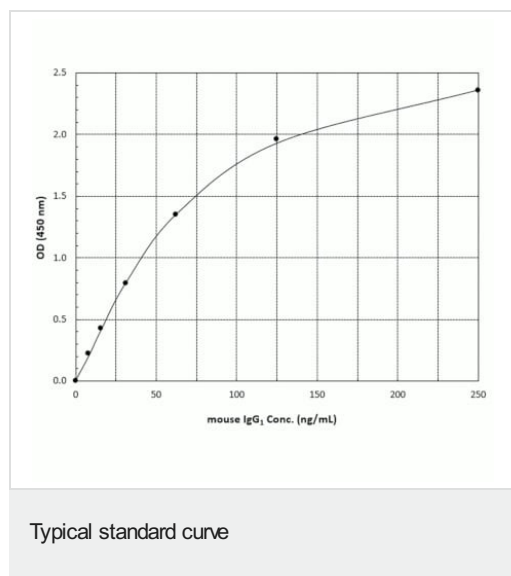
Components	Identifier	1 x 96 tests
20X Wash Buffer Concentrate		1 x 100ml

Components	Identifier	1 x 96 tests
Assay Buffer 13 Concentrate		1 x 50ml
Goat anti-mouse IgG Microplate (12 x 8 wells)		1 unit
Mouse IgG ₁ Conjugate Concentrate		1 x 0.07ml
Mouse IgG ₁ Standard		1 x 0.25ml
Mouse IgG1 Conjugate Diluent	Blue	1 x 6ml
Plate Sealer		2 units
Stop Solution 2		1 x 11ml
TMB Substrate		1 x 12ml

Relevance

There are four IgG subclasses (IgG1, 2, 3 and 4) in humans, named in order of their abundance in serum (IgG1 being the most abundant).

Images



Representative Standard Curve using ab133045.

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