

Mouse IgG2c ELISA Kit ab157720

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

Overview

Product name Mouse IgG2c 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 0.38 ng/ml

Range 3.13 ng/ml - 200 ng/ml

Recovery > 85 %

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	% - %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's IgG2c Mouse ELISA kit is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of IgG2c in mouse samples.

In this assay the IgG2c present in samples reacts with the anti-IgG2c antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-IgG2c antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound IgG2c. 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 IgG2c in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of IgG2c in the test sample. The quantity of IgG2c in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Platform

Microplate

Properties

Storage instructions

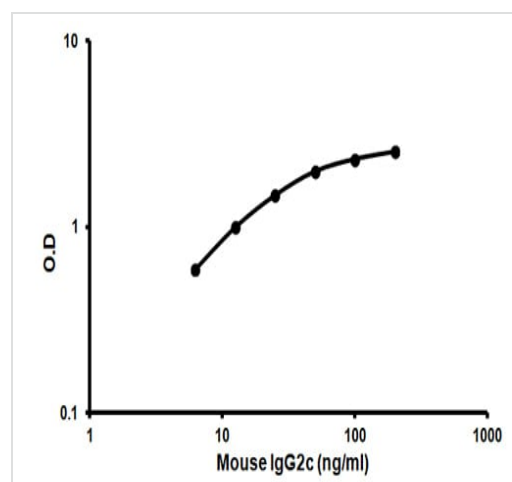
Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X HRP-conjugated anti-mouse IgG2c antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Mouse IgG2c Calibrator (lyophilized)	1 vial
Mouse IgG2c ELISA Microplate	1 unit
Stop Solution	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). IgG2 is the only IgG subclass which passes through the placenta at a level generally lower than that found in the mother. A deficiency of IgG2 indicates a poor antibody response to bacterial polysaccharides and can lead to increased susceptibility to infections caused by encapsulated bacteria.

Images



Representative standard curve using ab157720 IgG2c Mouse 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