

# Human IgG3 ELISA Kit ab137981

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

### Overview

**Product name** Human IgG3 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall				6.5%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				9.5%

**Sample type** Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Cell Lysate, Tissue Lysate, Cerebral Spinal Fluid

**Assay type** Sandwich (quantitative)

**Sensitivity** > 0.52 ng/ml

**Range** 0.391 ng/ml - 25 ng/ml

**Recovery** 102 %

**Assay time** 4h 0m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's IgG3 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IgG3 in plasma, serum, milk, urine, saliva, CSF, cell culture, cell lysate, and tissue samples.

An IgG3 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an IgG3 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of IgG3 captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

**Platform** Microplate

## Properties

---

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human IgG3 Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
IgG3 Microplate (12 x 8 well strips)	1 unit
IgG3 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

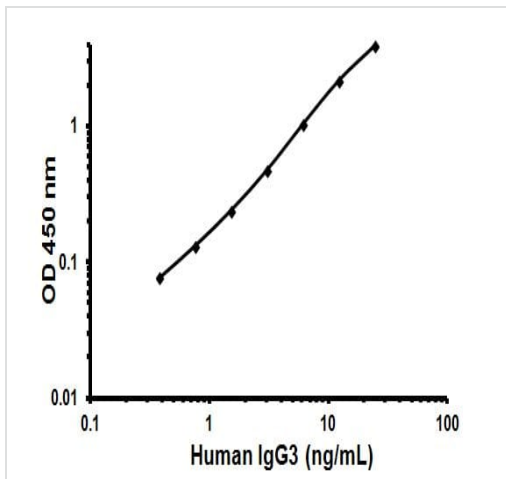
## Relevance

IgG3, like IgG1, activates complement and has an affinity for Fc receptors; thus IgG1 and IgG3 are the more effective of the IgG subclasses. An IgG3 deficiency is associated with Wiscott-Aldrich disease, systemic lupus erythematosus, juvenile diabetes mellitus, and periodontal infections. IgG anti-viral antibodies are restricted to IgG subclasses 1 and 3, with IgG3 antibodies being the first to appear in response to infection.

## Images

---

Representative Standard Curve using ab137981



Typical 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