



ELISA PRODUCT INFORMATION & MANUAL

IgG ELISA Development Kit ***NBP3-11757***

Enzyme-linked Immunosorbent Assay for quantitative
detection. For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

IgG ELISA DEVELOPMENT KIT

ELISA Development Kit for quantitative determination of native human IgG in solution, e.g. serum/plasma samples or cell supernatants.

The kit includes		NBP3-11757 for 6 plates
Capture mAb: MT145 (0.5 mg/ml)		300 µl
Detection mAb: MT78, biotinylated (0.5 mg/ml)		50 µl
Streptavidin-HRP		80 µl
Purified human IgG ELISA standard		1 vial
Standard reconstitution buffer A5		1 ml

To ensure total recovery of the stated quantity, vials have been overfilled.

Shipping and storage

Shipped at ambient temperature. All reagents should be stored at 4-8 °C upon receipt, except the standard which should be stored at -20 °C. Antibodies are supplied in sterile-filtered PBS with sodium azide (0.02%). Streptavidin-HRP is supplied in PBS with 0.002% Kathon CG. The expiry date indicates how long unopened products, stored according to instructions, are recommended for use.

General and Preparations

Specificity

The kit contains a matched pair of monoclonal antibodies (mAbs) specific for the Fc part of human IgG. The mAbs cross-react with IgG from non-human primates (NHP).

Standard range

0.1-10 ng/ml

Calibration

The ELISA standard has been calibrated against an international standard from the National Institute of Biological Standards and Control (NIBSC), Potters Bar, Hertfordshire EN6 3QG, UK. One µg of supplied standard equals 11 mU NIBSC-standard. Please note that the calibration is batch specific.

Analysis of serum and plasma samples

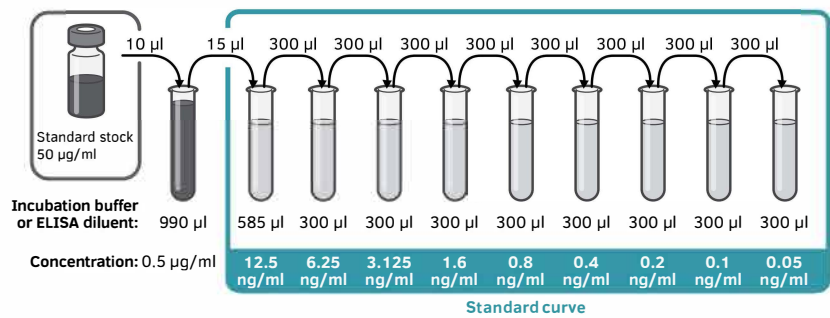
Analysis of serum/plasma requires the use of ELISA diluent. The ELISA diluent blocks heterophilic antibodies, commonly found in serum/plasma, from cross-linking the assay antibodies, thereby preventing false positive read-outs. The ELISA diluent should be used for dilution of standard, samples, and detection antibody.

Reconstitution of ELISA standard

Reconstitute the ELISA standard to a stock solution of 50 µg/ml by adding 0.5 ml of the standard reconstitution buffer. Allow the standard to dissolve for 5 minutes and mix thoroughly. The standard should be kept in aliquots at -20 °C. Avoid repeated freeze-thaw cycles.

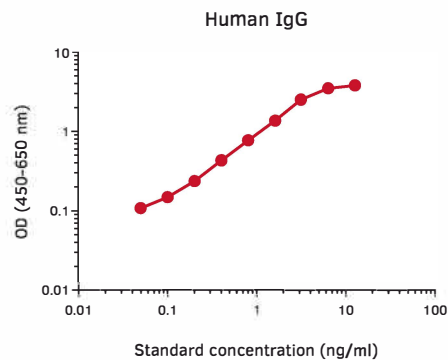
Preparation of standard curve

Prepare within 30 minutes of use. Volumes are sufficient for duplicates.



Protocol

- Day 1**
1. Add 100 μl /well of capture mAbs MT145 diluted to 2 $\mu\text{g}/\text{ml}$ in PBS, pH 7.4. Use high protein binding ELISA plates. Incubate overnight at 4–8 $^{\circ}\text{C}$.
- Day 2**
2. Empty the plate and add 200 μl /well of PBS with 0.05% Tween 20 and 0.1% BSA (incubation buffer) to block the plate. Incubate for 1 hour at room temperature.
 3. Wash the plate 5 times with PBS containing 0.05% Tween 20 (300 μl /well).
 4. Add 100 μl /well of samples or standards diluted in incubation buffer or ELISA diluent. Include assay background control, i.e. wells without standard. Incubate for 2 hours at room temperature.
 5. Wash as above.
 6. Add 100 μl /well of detection mAb MT78-biotin diluted to 0.25 $\mu\text{g}/\text{ml}$ in incubation buffer or ELISA diluent. Incubate for 1 hour at room temperature.
 7. Wash as above.
 8. Add 100 μl /well of Streptavidin-HRP diluted 1:1000 in incubation buffer. Incubate for 1 hour at room temperature. Please note that sodium azide used in buffers will inhibit HRP activity.
 9. Wash as above.
 10. Add 100 μl /well of TMB substrate and incubate for 15 minutes.
 11. Add 100 μl /well of 0.2 M H_2SO_4 to stop the reaction.
 12. Measure the optical density in an ELISA reader at 450 nm within 15 min. Preferably use a reader capable of subtracting a reference wavelength of between 570 and 650 nm. Representative standard curve shown below.



Quality management system complies with
the standards
ISO 9001:2015 & ISO 13485:2016.

The products are for research use only.

