

## ab285253 – Cytomegalovirus (CMV) IgG ELISA Kit

For detecting the presence of Cytomegalovirus IgG in human serum or plasma  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab285253>

### Storage and Stability

On receipt entire assay kit should be stored at 4°C, protected from light. Upon opening, use kit within 12 months. Keep microwells sealed in a dry bag with desiccants.

### Materials Supplied

Item	Quantity	Storage Condition
CMV Antigen pre-coated Microplate	8 wells × 12 strips	4°C
Sample Diluent	22 mL	4°C
Calibrator	1 mL	4°C
Positive Control	1 mL	4°C
Negative Control	1 mL	4°C
Enzyme Conjugate	12 mL	4°C
TMB Substrate	12 mL	4°C
Stop Solution	12 mL	4°C
20X Wash Concentrate	25 mL	4°C

### Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Distilled or deionized water
- Adjustable Precision pipettes
- Disposable pipette tips
- Microplate reader capable of measuring absorbance at 450 nm.
- Absorbent paper

### Reagent Preparation

- Before using the kit, spin the tubes prior to opening.

**Wash Buffer:** Prepare 1X Wash buffer by adding the contents of the bottle (25 mL, 20X) to 475 mL of distilled or deionized water. Store at room temperature after dilution.

### Sample Preparation

- Collect blood specimens and separate the serum.  
**Δ Note:** Specimens may be refrigerated at 2-8° C for up to seven days or frozen for up to six months.
- Avoid repeated freeze-thaw cycles.
- Lipemic or haemolyzed samples may cause erroneous results.

### Assay Protocol

- Bring all reagents and samples to room temperature and mix gently 30 minutes prior to the assay.

1. Place the desired number of coated strips into the holder.
2. Negative control, positive control, and calibrator are ready to use. Prepare 1:21 dilution of test samples by adding 10 µL of the sample to 200 µL of sample diluent. Mix well.

3. Dispense 100 µL of diluted sera, calibrator and controls into the appropriate wells. For the reagent blank, dispense 100µL sample diluent in 1A well position. Tap the holder to remove air bubbles from the liquid and mix well. Incubate at room temperature for 20 minutes.
4. Remove liquid from all wells. Wash wells three times with 300 µL of 1X wash buffer. Blot on absorbance paper or paper towel to remove residual wash buffer.
5. Dispense 100 µL of enzyme conjugate to each well and incubate for 20 minutes at room temperature.
6. Remove enzyme conjugate from all wells. Wash wells three times with 300 µL of 1X wash buffer. Blot on absorbance paper or paper towel to remove residual wash buffer.
7. Dispense 100 µL of TMB substrate and incubate for 10 minutes at room temperature.
8. Add 100 µL of stop solution.
9. Read O.D. at 450 nm using ELISA reader within 15 min. A dual wavelength is recommended with reference filter of 600-650 nm.

### Calculation

- Check Calibrator Factor (CF) value on the calibrator bottle. This value might vary depending on the lot number of the kit.
- Calculate the cut-off value: Calibrator OD x Calibrator Factor (CF).
- Calculate the Antibody (Ab) Index: Dividing the O.D. value of each sample by cut-off value.

**Δ Note:** Background subtraction for each reading is optional for calculating the sample Cortisol concentration, and will not change the final results.

### Quality Control:

- The test run may be considered valid provided the following criteria are met:
  1. The O.D. of the Calibrator should be greater than 0.25
  2. The Ab index for Negative control should be less than 0.9
  3. The Ab Index for Positive control should be greater than 1.2

### Final Result Interpretation:

Ab Index	Result
< 0.9	No detectable antibody
0.9 – 1.1	Borderline Positive
> 1.1	Detectable antibody

### **Technical Support**

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)