



Product datasheet

Human Anti-Cytomegalovirus IgG ELISA Kit (CMV) ab108724

★★★★★ [2 Abreviews](#) [14 References](#)

概述

产品名称	人Anti-Cytomegalovirus IgG ELISA试剂盒(CMV)			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Pos.Serum	12		1.9%
	批次间			
	样品	n	Mean	SD
	Pos.Serum	4		6.2%
样品类型	Serum, Hep Plasma, Cit plasma			
检测类型	Indirect			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	Abcam's anti-Cytomegalovirus (CMV) IgG Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the qualitative measurement of IgG class antibodies against Cytomegalovirus in Human serum and plasma.			

A 96-well plate has been precoated with Cytomegalovirus antigens to bind cognate antibodies. Controls or test samples are added to the wells and incubated. Following washing, a horseradish peroxidase (HRP) labelled anti-Human IgG conjugate is added to the wells, which binds to the immobilized Cytomegalovirus-specific antibodies. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The density of yellow coloration is directly proportional to the amount of Cytomegalovirus IgG sample captured in plate.

平台 Microplate

性能

存放说明

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

组件	标识符	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Cover foil		1 unit
Cytomegalovirus antigens Coated Microplate (12 x 8 wells)		1 unit
Cytomegalovirus anti-IgG HRP Conjugate	Black cap	1 x 20ml
Cytomegalovirus IgG Cut-off Control	Green cap	1 x 3ml
Cytomegalovirus IgG Negative Control	Blue cap	1 x 2ml
Cytomegalovirus IgG Positive Control	Red cap	1 x 2ml
IgG Sample Diluent	White cap	1 x 100ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors