

Product datasheet

Human Anti-Herpes simplex virus Type 2 IgM ELISA Kit (HSV2) ab108740

[1 Image](#)

Overview

Product name Human Anti-Herpes simplex virus Type 2 IgM ELISA Kit (HSV2)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Pos.Serum	24			6.9%
Pos.Serum	20			7.7%

Inter-assay

Sample	n	Mean	SD	CV%
Pos.Serum	14			5.6%
Pos.Serum	14			4.2%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Indirect

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's anti-Herpes simplex virus Type 2 (HSV2) IgM Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Herpes simplex virus Type 2 in Human serum and plasma.

A 96-well plate has been precoated with Herpes simplex virus Type 2 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 IgM conjugate is added to the wells, which binds to the immobilized Herpes simplex virus Type 2-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 Herpes simplex virus Type 2 IgM sample captured in plate.

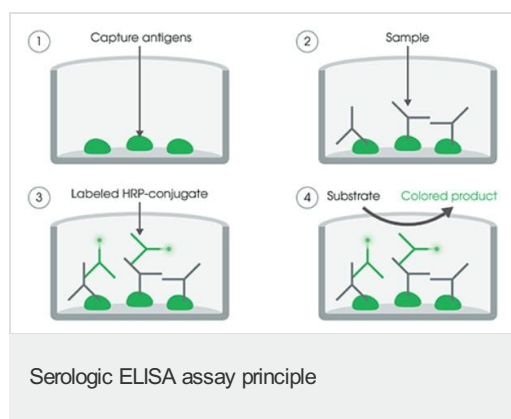
Platform

Microplate

Properties**Storage instructions**

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Cover foil		1 unit
Herpes simplex virus Type 2 (IgM) Coated Microplate (12 x 8 wells)		1 unit
Herpes simplex virus Type 2 anti-IgM HRP Conjugate	Red , Black cap	1 x 20ml
Herpes simplex virus Type 2 IgM Cut-off Control	Yellow , Green cap	1 x 3ml
Herpes simplex virus Type 2 IgM Negative Control	Yellow , Blue cap	1 x 2ml
Herpes simplex virus Type 2 IgM Positive Control	Yellow , Red cap	1 x 2ml
IgM Sample Diluent	Green , White cap	1 x 100ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

Images

Specific antigens are coated on the 96-well plate, controls or test samples are added to the well and incubated. The wells are washed to remove any unbound Human anti-antigen antibodies (Ig). A horseradish peroxidase (HRP) labelled anti-Human Ig conjugate is added to the wells. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The intensity of yellow coloration is directly proportional to the amount of Human anti-antigen Ig captured on the plate.

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