

Human Anti-Toxoplasma gondii IgM ELISA Kit ab108778

1 References

Overview

Product name Human Anti-Toxoplasma gondii IgM ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum 1	24			2.9%
Serum 2	24			1.5%

Inter-assay

Sample	n	Mean	SD	CV%
Serum 2	12			6.5%
Serum 3	12			4.5%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Sandwich (qualitative)

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview ab108778 Toxoplasma gondii IgM Human ELISA Kit is intended for the qualitative determination of IgM class antibodies against Toxoplasma gondii in human serum or plasma (citrate).

Toxoplasma gondii is a small intracellular parasite, whose live cycle has a sexual and an asexual phase. Sexual development is restricted to the intestinal cells of (probably exclusively) cats; the oocysts formed are excreted and due to their resistant cell walls they may be infectious under advantageous circumstances for at least 1 year. Animals and man are intermediate hosts for the asexual proliferation of T. gondii: the ingested parasites will proliferate explosively within the host cells lysing them eventually. They disseminate throughout the body via circulation and lymphatic system and may infect any cell type. In muscle and brain cells cysts are formed which are spheroidal and about 5-100 µm in diameter. Cysts are virtually immortal in the intermediate host.

Toxoplasma gondii is the most common parasite in humans, but its abundance (7-80%) is highly dependent on the geographic area, the socio-economic status and the nutritional customs. Infection only rarely causes toxoplasmosis and usually clinical symptoms are absent, but may produce severe problems in immunosuppressed persons and fetus.

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
Anti-Human (IgM) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Cover foil		1 unit
IgM Sample Diluent	Colored yellow	1 x 100ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml
Toxoplasma gondii HRP Conjugate		1 x 15ml
Toxoplasma gondii IgM Cut-off Control	Colored yellow; Green cap	1 x 3ml
Toxoplasma gondii IgM Negative Control	Colored yellow; Blue cap	1 x 2ml
Toxoplasma gondii IgM Positive Control	Colored yellow; Red cap	1 x 2ml

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