

# Human Anti-Toxoplasma gondii IgM ELISA Kit ab108778

**概述**

<b>产品名称</b>	人Anti-Toxoplasma gondii IgM ELISA试剂盒																														
<b>检测方法</b>	Colorimetric																														
<b>精确度</b>	<p style="text-align: right;">批次内</p> <table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th style="width: 20%;">样品</th> <th style="width: 10%;">n</th> <th style="width: 20%;">Mean</th> <th style="width: 10%;">SD</th> <th style="width: 30%;">CV%</th> </tr> </thead> <tbody> <tr> <td>Serum 1</td> <td>24</td> <td></td> <td></td> <td>2.9%</td> </tr> <tr> <td>Serum 2</td> <td>24</td> <td></td> <td></td> <td>1.5%</td> </tr> </tbody> </table> <p style="text-align: right;">批次间</p> <table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th style="width: 20%;">样品</th> <th style="width: 10%;">n</th> <th style="width: 20%;">Mean</th> <th style="width: 10%;">SD</th> <th style="width: 30%;">CV%</th> </tr> </thead> <tbody> <tr> <td>Serum 2</td> <td>12</td> <td></td> <td></td> <td>6.5%</td> </tr> <tr> <td>Serum 3</td> <td>12</td> <td></td> <td></td> <td>4.5%</td> </tr> </tbody> </table>	样品	n	Mean	SD	CV%	Serum 1	24			2.9%	Serum 2	24			1.5%	样品	n	Mean	SD	CV%	Serum 2	12			6.5%	Serum 3	12			4.5%
样品	n	Mean	SD	CV%																											
Serum 1	24			2.9%																											
Serum 2	24			1.5%																											
样品	n	Mean	SD	CV%																											
Serum 2	12			6.5%																											
Serum 3	12			4.5%																											
<b>样品类型</b>	Serum, Plasma																														
<b>检测类型</b>	Sandwich (qualitative)																														
<b>实验步骤</b>	Multiple steps standard assay																														
<b>种属反应性</b>	<b>与反应:</b> Human																														
<b>产品概述</b>	<p>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).</p> <p>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.</p> <p>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.</p>																														

经测试应用

适用于: Sandwich ELISA

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
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		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

## 应用

Our [Abpromise guarantee](#) covers the use of **ab108778** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### 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 <http://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