

## Product datasheet

# anti-Omalizumab ELISA Kit ab237664

### Overview

---

<b>Product name</b>	anti-Omalizumab ELISA Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Serum, Plasma
<b>Assay type</b>	Sandwich (qualitative)
<b>Assay duration</b>	Multiple steps standard assay
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	anti-Omalizumab ELISA Kit (ab237664) is designed to quantify/measure the antibody against Omalizumab with high specificity and sensitivity in biological matrices.
<b>Notes</b>	This product is manufactured by BioVision, an Abcam company and was previously called E4395 BioSim™ anti-Omalizumab (Human) ELISA Kit. E4395-100 is the same size as the 96 test size of ab237664.
<b>Platform</b>	Microplate

### Properties

---

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
Assay Buffer	1 x 12ml
Micro ELISA Plate	1 unit
Negative Control	1 x 0.5ml
Peroxidase Conjugate	1 x 12ml
Plate sealers	2 units
Positive Control	1 x 0.25ml
Stop Solution	1 x 12ml
TMB substrate	1 x 12ml

Components	1 x 96 tests
Wash buffer (20X)	1 x 50ml

### Relevance

Omalizumab is a recombinant DNA-derived humanized IgG1k monoclonal antibody that selectively binds to human immunoglobulin E (IgE). Xolair inhibits the binding of IgE to the high-affinity IgE receptor (FcεRI) on the surface of mast cells and basophils. Reduction in surface-bound IgE on FcεRI-bearing cells limits the degree of release of mediators of the allergic response. Xolair is used to treat severe, persistent asthma.

**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