

Datasheet: MCA2475KZZ

Description:	MOUSE OVALBUMIN SPECIFIC IgE ELISA ASSAY KIT
Name:	MOUSE OVALBUMIN SPECIFIC IgE ELISA ASSAY KIT
Format:	Kit
Product Type:	Kits
Quantity:	96 TESTS

Product Details

Applications

This product has been reported to work in the following applications. This information is derived from testing within our laboratories, peer-reviewed publications or personal communications from the originators. Please refer to references indicated for further information. For general protocol recommendations, please visit www.bio-rad-antibodies.com/protocols.

	Yes	No	Not Determined	Suggested Dilution
ELISA	■			

Where this product has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the product for use in their own system using appropriate negative/positive controls.

Target Species

Mouse

Product Information

Mouse Ovalbumin specific IgE ELISA Assay Kit accurately assays Ovalbumin specific IgE levels in a two-step direct ELISA using a rat anti-mouse IgE monoclonal antibody and HRP-conjugated Ovalbumin.

Reagents In The Kit

Assay Buffer:	1 bottle (30 ml)
Detection Conjugate:	1 bottle (15 ml) HRP-labeled Ovalbumin, ready-to-use
Antibody-coated assay plate:	1 plate (96 wells) coated with rat anti-mouse IgE monoclonal antibody
Microplate for predispensing samples:	1 plate (96 wells)
Wash Buffer Concentrate:	1 bottle (90 ml)
HRP Substrate:	1 bottle (15 ml), ready-to-use
Stop Solution:	1 bottle (15 ml), ready-to-use
Standards:	6 vials

Instructions For Use

Instructions for use can be found at www.bio-rad-antibodies.com/uploads/IFU/MCA2475KZZ.pdf

References

1. Hajek, A.R. *et al.* (2008) 12/15-Lipoxygenase deficiency protects mice from allergic airways inflammation and increases secretory IgA levels. [J Allergy Clin Immunol. 122 \(3\): 633-9.e3.](#)
2. Aranzamendi, C. *et al.* (2012) Protection against allergic airway inflammation during the chronic and acute phases of *Trichinella spiralis* infection. [Clin Exp Allergy. 43: 103-15.](#)
3. Naura, A.S. *et al.* (2013) Minocycline blocks asthma-associated inflammation in part by interfering with the T Cell receptor-NF- κ B-GATA-3-IL-4 axis without a prominent effect on PARP. [J Biol Chem. 288: 1458-68.](#)
4. Lee, M.Y. *et al.* (2013) Pinellia ternata Breitenbach attenuates ovalbumin-induced allergic airway inflammation and mucus secretion in a murine model of asthma. [Immunopharmacol Immunotoxicol. 35 \(3\): 410-8.](#)
5. He, C. *et al.* (2013) Measles virus-derived peptide/food antigen adducts facilitate the establishment of antigen specific oral tolerance. [J Physiol Pharmacol. 64 \(1\): 95-102.](#)
6. Toomer, O.T. *et al.* (2014) Maternal and postnatal dietary probiotic supplementation enhances splenic regulatory T helper cell population and reduces ovalbumin allergen-induced hypersensitivity responses in mice. [Immunobiology. 219 \(5\): 367-76.](#)
7. Ghonim, M.A. *et al.* (2015) DNA-dependent protein kinase inhibition blocks asthma in mice and modulates human endothelial and CD4⁺ T-cell function without causing severe combined immunodeficiency. [J Allergy Clin Immunol. 135 \(2\): 425-40.](#)
8. Chung, S.H. *et al.* (2015) The C-C Chemokine receptor 6 (CCR6) is crucial for Th2-driven allergic conjunctivitis. [Clin Immunol. pii: S1521-6616\(15\)30021-8.](#)
9. Kwon, J.Y. *et al.* (2016) TRPV1 Antagonist Suppresses Allergic Conjunctivitis in a Murine Model. [Ocul Immunol Inflamm. Oct 11 \[Epub ahead of print\]](#)
10. Lew, D.B. *et al.* (2017) Beneficial Effects of Prebiotic *Saccharomyces cerevisiae* Mannan on Allergic Asthma Mouse Models. [J Immunol Res. 2017: 3432701.](#)

Storage

Store in a cool place(+4°C).
Avoid light and freezing.

Shelf Life

6 months from the date of despatch.

Health And Safety Information

Material Safety Datasheet documentation #10294 #10310 #10311 #10312 #10313 available at:
Antibody (10294): <https://www.bio-rad-antibodies.com/uploads/MSDS/10294.pdf>
Ovalbumin:HRP (10310): <https://www.bio-rad-antibodies.com/uploads/MSDS/10310.pdf>
Buffer/Wash Buffer Concentrate (10311): <https://www.bio-rad-antibodies.com/uploads/MSDS/10311.pdf>
TMB Substrate (10312): <https://www.bio-rad-antibodies.com/uploads/MSDS/10312.pdf>
Stop Solution (10313): <https://www.bio-rad-antibodies.com/uploads/MSDS/10313.pdf>

Regulatory

For research purposes only

North & South America

Tel: +1 800 265 7376

Fax: +1 919 878 3751

Email: antibody_sales_us@bio-rad.com

Worldwide

Tel: +44 (0)1865 852 700

Fax: +44 (0)1865 852 739

Email: antibody_sales_uk@bio-rad.com

Europe

Tel: +49 (0) 89 8090 95 21

Fax: +49 (0) 89 8090 95 50

Email: antibody_sales_de@bio-rad.com

'M277975:151208'

Printed on 14 May 2018