

anti-Ipilimumab ELISA Kit ab237667

Overview

Product name	anti-Ipilimumab ELISA Kit
Detection method	Colorimetric
Sample type	Serum, Plasma
Assay type	Sandwich (qualitative)
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	anti-Ipilimumab ELISA Kit (ab237667) is designed to quantify/measure the antibody against Ipilimumab with high specificity and sensitivity in biological matrices.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called E4398 BioSim™ anti-Ipilimumab (Yervoy®) (Human) ELISA Kit. E4398-100 is the same size as the 96 test size of ab237667.
Platform	Microplate

Properties

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

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

Relevance

Ipilimumab is a fully human IgG1 kappa antibody that binds to CTLA-4 (cytotoxic T lymphocyte associated antigen 4), a molecule on T cells that is indicated for unresectable or metastatic melanoma. The absence or presence of CTLA-4 can augment or suppress the immune system's T-cell response in fighting disease. Ipilimumab is designed to block the activity of CTLA-4, thereby sustaining an active immune response in its attack on cancer cells. The proposed mechanism of action is indirect, and may be through T-cell - mediated anti-tumor immune responses.

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