

Product datasheet

anti-Etanercept ELISA Kit ab237661

Overview

Product name	anti-Etanercept 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-Etanercept ELISA Kit (ab237661) is designed to quantify/measure the antibody against Etanercept with high specificity and sensitivity in biological matrices.
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 1ml
Peroxidase Conjugate	1 x 12ml
Plate sealers	2 units
Positive Control	1 x 0.3ml
Stop Solution	1 x 12ml
TMB substrate	1 x 12ml
Wash buffer (20X)	1 x 50ml

Relevance Etanercept is a therapeutic fusion protein specific for Tumor Necrosis Factor-Alpha (TNF- α) and

is used to treat rheumatic arthritis, intestinal disorders, dermatological diseases and cancer. Etanercept specifically binds to TNF alpha and blocks its interaction with cell surface TNF receptors and reduces the inflammation and subsequently improves the patient's health. Drug level quantification can be important to adapt patient prescription or to switch to an alternative TNF inhibitor drug.

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