

Product datasheet

anti-Golimumab ELISA Kit ab237662

Overview

Product name	anti-Golimumab 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-Golimumab ELISA Kit (ab237662) is designed to quantify/measure the antibody against Golimumab with high specificity and sensitivity in biological matrices.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called E4393 BioSim™ anti-Golimumab (Simponi®) (Human) ELISA Kit. E4393-100 is the same size as the 96 test size of ab237662.
Platform	Microplate

Properties

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

Components	1 x 96 tests
Micro ELISA Plate	1 unit
Positive Control	1 x 0.25ml
Negative Control	1 x 0.5ml
Assay Buffer	1 x 12ml
Peroxidase Conjugate	1 x 12ml
TMB substrate	1 x 12ml
Stop Solution	1 x 12ml
Wash buffer (20X)	1 x 50ml

Components	1 x 96 tests
Plate sealers	1 unit

Relevance

Golimumab (Simponi®) is a human immunoglobulin G1? monoclonal antibody which is specific for pro-inflammatory cytokine, tumor necrosis factor- α (TNF α). In 2009, it was approved by FDA for the treatment of rheumatoid arthritis, psoriatic arthritis and ankylosing spondylitis in adult patients. Elevated levels of TNF are found in the synovial fluid of rheumatoid arthritis, including juvenile idiopathic arthritis, psoriatic arthritis, and ankylosing spondylitis patients and play an important role in both the pathologic inflammation and the joint destruction that are hallmarks of these diseases. Increased levels of TNF are also found in psoriasis (Ps) plaques. Golimumab binds to both the soluble and transmembrane bioactive forms of human TNF and prevent TNF from binding to its receptors and finally inhibits biological activity of TNF.

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