

2 References

Overview

Product name	Adalimumab ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 15%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 15%
Sample type	Serum, Plasma				
Assay type	Quantitative				
Sensitivity	10 ng/ml				
Range	30 ng/ml - 1000 ng/ml				
Recovery	Sample specific recovery				
	Sample type	Average %		Range	
	Serum			85% - 115%	
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Note: Adalimumab ELISA kit ab237641 is intended for Research Use Only, not for use in diagnostic procedures.				

This kit is a sandwich based ELISA kit. The color developed is proportional to the amount of Adalimumab in the sample or standard. Results of samples can be determined directly using the standard curve.

Notes

This product is manufactured by BioVision, an Abcam company and was previously called E4372 BioSim™ Adalimumab (Human) ELISA Kit. E4372-100 is the same size as the 96 test size of

ab237641.

Platform

Microplate

Properties

Storage instructions

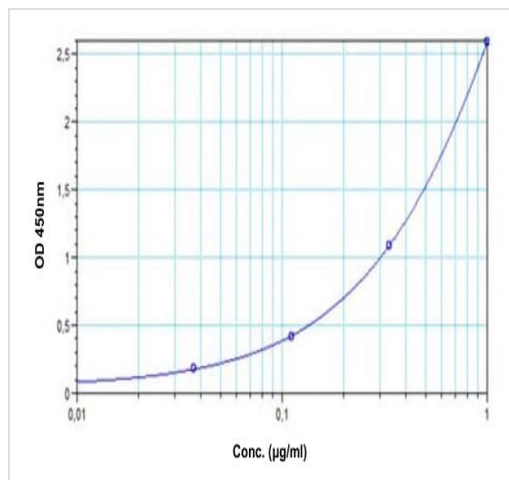
Store at +4°C. Please refer to protocols.

Components	96 tests
Adalimumab Standard S1	1 x 1ml
Adalimumab Standard S2	1 x 1ml
Adalimumab Standard S3	1 x 1ml
Adalimumab Standard S4	1 x 1ml
Adalimumab Standard S5	1 x 1ml
Adalimumab Standard S6	1 x 1ml
Adalimumab Standard S7	1 x 1ml
Assay Buffer	1 x 50ml
HRP-conjugate Probe	1 x 12ml
Micro ELISA Plate	1 unit
Plate sealers	2 units
Stop Solution	1 x 12ml
TMB substrate	1 x 12ml
Wash buffer (20X)	1 x 50ml

Relevance

Adalimumab is a recombinant human IgG1 monoclonal antibody specific for Tumor Necrosis Factor-Alpha (TNF- α) and is used to treat rheumatic arthritis, intestinal disorders, dermatological diseases and cancer. Adalimumab specifically binds to TNF α and blocks its interaction with p55 and p75 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.

Images



Typical Standard Curve: This standard curve is for demonstration only. A standard curve must be run with each assay.

Example of Adalimumab standard curve

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