

### Anti-Cetuximab ELISA Kit (qualitative) ab282902

#### Overview

**Product name** Anti-Cetuximab ELISA Kit (qualitative)

**Detection method** Colorimetric

**Precision** Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 30%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 30%

**Sample type** Serum, Plasma

**Assay type** Sandwich (qualitative)

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Note: Anti-Cetuximab ELISA kit ab282902 is intended for Research Use Only, not for use in diagnostic procedures.

Anti-Cetuximab ELISA kit (qualitative) is a sandwich ELISA assay for the qualitative detection of antibodies to Cetuximab in human serum and plasma samples. Controls and samples (serum and plasma) are incubated in the microtiter plate coated with the drug Cetuximab. After incubation, the wells are washed and the enzymatic activity is detected by the addition of chromogen-substrate. The enzyme-substrate reaction is terminated with an acidic stop solution. The color developed is proportional to the amount of antibodies specific for Cetuximab present in the samples and controls. The qualitative test results can be evaluated using cut-off value.

Cross Reactivity: Except for Cetuximab, there is no cross reaction with other therapeutic antibodies and native serum immunoglobins.

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called E4873 BioSim™ anti-Cetuximab II (Human) ELISA Kit. E4873-100 is the same size as the 1 x 96 tests size of ab282902.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

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

Components	1 x 96 tests
Anti-Cetuximab Negative Control	1 x 1ml
Anti-Cetuximab Positive Control	1 x 300µl
Assay Buffer	1 x 12ml
Cetuximab coated microtitre plate	1 unit
HRP-conjugate Probe	1 x 12ml
Plate sealers	2 x 1ml
Stop Solution	1 x 12ml
TMB substrate	1 x 12ml
Wash Buffer (20X)	1 x 50ml

**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