

### Bovine IL-10 ELISA Kit ab277386

[1 Image](#)

#### Overview

**Product name** Bovine IL-10 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

**Sample type** Cell culture supernatant, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 420 pg/ml

**Range** 420 pg/ml - 7000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	130.5	120% - 139%
Plasma	110.3	71% - 139%
Cell culture media	109.5	79% - 140%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Cow

**Product overview** Bovine IL-10 ELISA Kit (ab277386) is an *in-vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Bovine IL-10 serum (bovine IL-10 concentration is low in normal serum/plasma, and may not be detectable in this assay), plasma and cell culture supernatants.

This assay employs an antibody specific for Bovine IL-10 coated on a 96-well plate. Standards

and samples are pipetted into the wells and Bovine IL-10 present in a sample is bound to the wells by the immobilized antibody. The wells are washed, and biotinylated anti-Bovine IL-10 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Bovine IL-10 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Biotinylated Anti-Bovine IL-10 Antibody	2 vials
Anti-Bovine IL-10 coated Microplate	1 unit
Bovine IL-10 standard protein (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

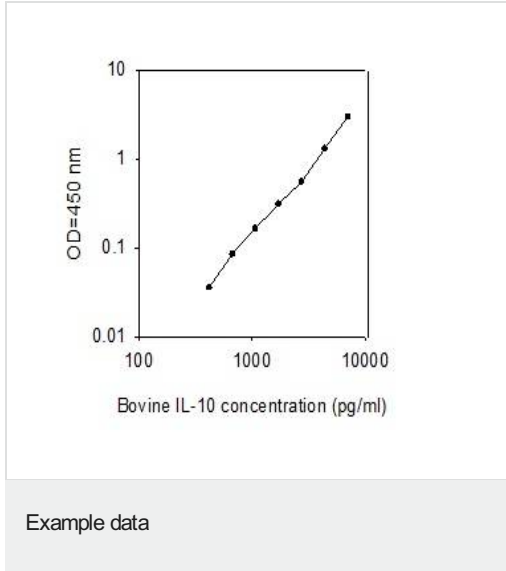
**Function** Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

**Tissue specificity** Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types.

**Sequence similarities** Belongs to the IL-10 family.

**Cellular localization** Secreted.

## Images



These standard curves are for demonstration only. A standard curve must be run with each assay.

**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