

Product datasheet

Equine IL-10 ELISA Kit ab155466

1 Image

Overview

Product name Equine 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 < 90 pg/ml

Range 0.1 ng/ml - 25 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	104.9	96% - 112%
Plasma	94.01	86% - 102%
Cell culture media	97.3	90% - 105%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Horse

Product overview

Abcam's IL-10 Equine ELISA Kit is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of equine IL-10 in serum, plasma and cell culture supernatants.

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

and samples are pipetted into the wells and IL-10 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-equine IL-10 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted into 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 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 Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Detection Antibody IL-10 : biotinylated anti-equine IL-10	2 vials
IL-10 Microplate (12 strips x 8 wells)	1 unit
IL-10 Standard equine(recombinant) (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

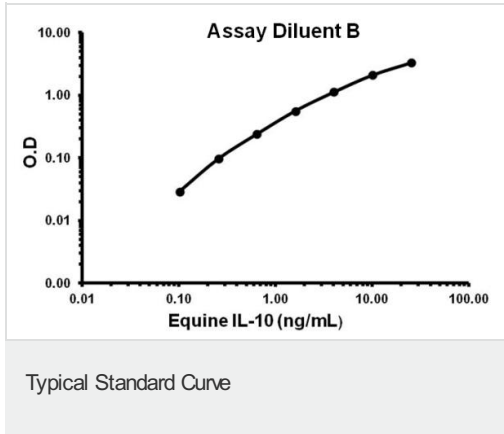
Function Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

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



Representative standard curve using ab155466 - Assay Diluent B.

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