

Product datasheet

Ovine IL-12 p70 ELISA Kit ab273255

1 Image

Overview

Product name Ovine IL-12 p70 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, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity 2.5 pg/ml

Range 2.05 pg/ml - 200 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
-------------	-----------	-------

Serum 72.35 68% - 79%

Plasma 74.38 68% - 79%

Cell culture media 94.78 80% - 114%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Sheep

Product overview

Ovine IL-12 p70 ELISA Kit (ab273255) is an *in-vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Ovine IL-12 p70 in serum, plasma (Collect plasma using EDTA as an anticoagulant. **Heparin and citrate are not recommended**) and cell culture supernatants.

This assay employs an antibody specific for Ovine IL-12 p70 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-12 p70 present in a sample is bound to the wells by the immobilized antibody. The wells are washed, and biotinylated anti-Ovine IL-12 p70 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 IL-12 p70 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Tested applications

Suitable for: Sandwich ELISA

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
150X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Anti-Ovine IL-12 p70 coated Microplate (12 x 8 wells)	1 unit
Assay Diluent C	1 x 30ml
Biotinylated Anti-Ovine IL-12 p70 Antibody	2 vials
Ovine IL-12 p70 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Relevance

IL-12 p70 is a disulfide bonded heterodimer of the IL-12 40 kDa and 35 kDa subunits.

Cellular localization

Secreted

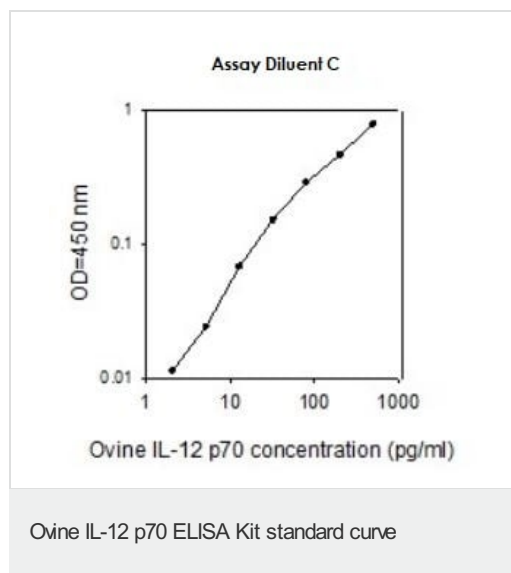
Applications

Our [Abpromise guarantee](#) covers the use of **ab273255** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



This standard curve is 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