

Porcine IL-17F ELISA Kit ab273224

1 Image

Overview

Product name Porcine IL-17F 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, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 6 pg/ml

Range 4.1 pg/ml - 1000 pg/ml

Recovery Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	76.04	73% - 84%
Serum	74.84	63% - 71%
Plasma	82.75	75% - 91%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rat, Pig

Product overview Porcine IL-17F ELISA Kit (ab273224) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of Porcine IL-17F in serum, plasma (Collect plasma using EDTA and citrate as an anti-coagulant. **Heparin is not recommended as anti-coagulant for use in this assay**) and cell culture supernatants.

This assay employs an antibody specific for Porcine IL-17F coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-17F present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Porcine IL-17F 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-17F 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
150X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Anti-Porcine IL-17F coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Porcine IL-17F Antibody	2 vials
Porcine IL-17F Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function Ligand for IL17RA and IL17RC (PubMed:17911633). The heterodimer formed by IL17A and IL17F is a ligand for the heterodimeric complex formed by IL17RA and IL17RC (PubMed:18684971). Involved in stimulating the production of other cytokines such as IL6, IL8 and CSF2, and in regulation of cartilage matrix turnover (PubMed:11591732, PubMed:11591768, PubMed:11574464). Also involved in stimulating the proliferation of peripheral blood mononuclear cells and T-cells and in inhibition of angiogenesis (PubMed:11591732). Plays a role in the induction of neutrophilia in the lungs and in the exacerbation of antigen-induced pulmonary allergic inflammation.

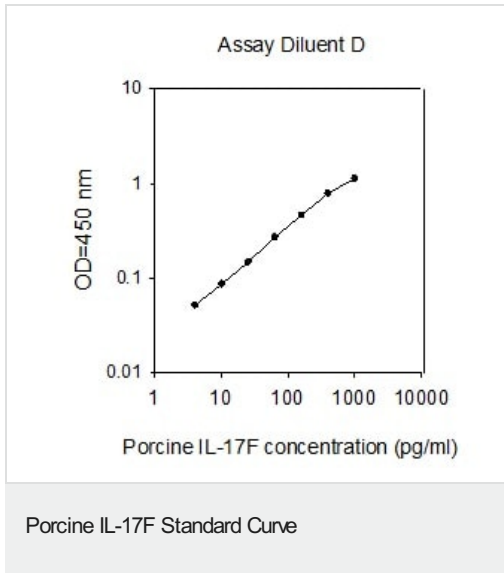
Tissue specificity Expressed in activated, but not resting, CD4+ T-cells and activated monocytes.

Involvement in disease Candidiasis, familial, 6

Sequence similarities Belongs to the IL-17 family.

Cellular localization Secreted.

Images



Typical standard curve – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.

Porcine IL-17F ELISA kit (ab273224) 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