

Feline IL-4 ELISA Kit ab273226

1 Image

Overview

Product name	Feline IL-4 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	0.3 ng/ml				
Range	0.33 ng/ml - 80 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		126.7		109% - 143%
	Plasma		115.2		107% - 124%
	Cell culture media		86.89		77% - 91%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Cat				
Product overview	Feline IL-4 ELISA Kit (ab273226) is an <i>in-vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Feline IL-4 in serum, plasma and cell culture supernatants.				

This assay employs an antibody specific for Feline IL-4 coated on a 96-well plate. Standards and samples are pipetted into the wells and Feline IL-4 present in a sample is bound to the wells by

the immobilized antibody. The wells are washed, and biotinylated anti-Feline IL-4 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 Feline IL-4 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
20X Wash Buffer	1 x 25ml
500X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Anti-Feline IL-4 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Feline IL-4 Antibody	2 vials
Feline IL-4 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

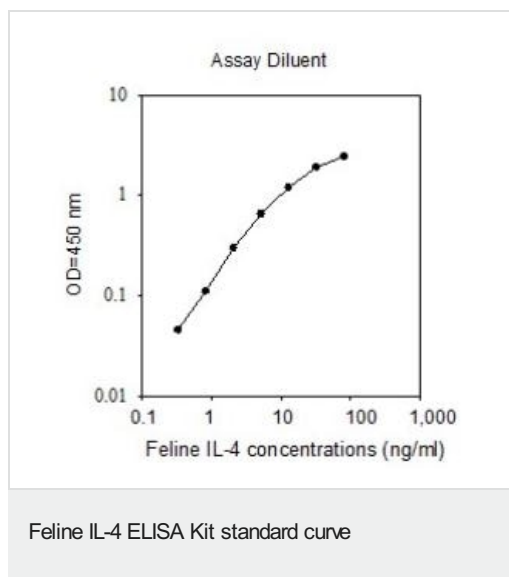
Function Participates in at least several B-cell activation processes as well as of other cell types. It is a costimulator of DNA-synthesis. It induces the expression of class II MHC molecules on resting B-cells. It enhances both secretion and cell surface expression of IgE and IgG1. It also regulates the expression of the low affinity Fc receptor for IgE (CD23) on both lymphocytes and monocytes.

Involvement in disease Genetic variations in IL4 may be a cause of susceptibility to ischemic stroke (ISCHSTR) [MIM:601367]; also known as cerebrovascular accident or cerebral infarction. A stroke is an acute neurologic event leading to death of neural tissue of the brain and resulting in loss of motor, sensory and/or cognitive function. Ischemic strokes, resulting from vascular occlusion, is considered to be a highly complex disease consisting of a group of heterogeneous disorders with multiple genetic and environmental risk factors.

Sequence similarities Belongs to the IL-4/IL-13 family.

Cellular localization Secreted.

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