

Canine IL8 ELISA Kit (Interleukin-8) ab155465

1 References 1 Image

Overview

Product name Canine IL8 ELISA Kit (Interleukin-8)

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 < 6 pg/ml

Range 6.1 pg/ml - 1500 pg/ml

Recovery	Sample specific recovery		
	Sample type	Average %	Range
	Serum	110.2	102% - 118%
	Plasma	103.2	84% - 122%
	Cell culture media	128.2	122% - 132%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Dog

Product overview Abcam's Interleukin-8 (IL-8) Canine ELISA Kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of dog IL-8 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for dog IL-8 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-8 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-dog IL-8 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-8 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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Detection Antibody IL-8 (biotinylated anti-Canine IL-8)	2 vials
IL-8 Microplate (12 strips x 8 wells) coated with anti canine IL-8	1 unit
IL-8 Standard canine (recombinant) (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

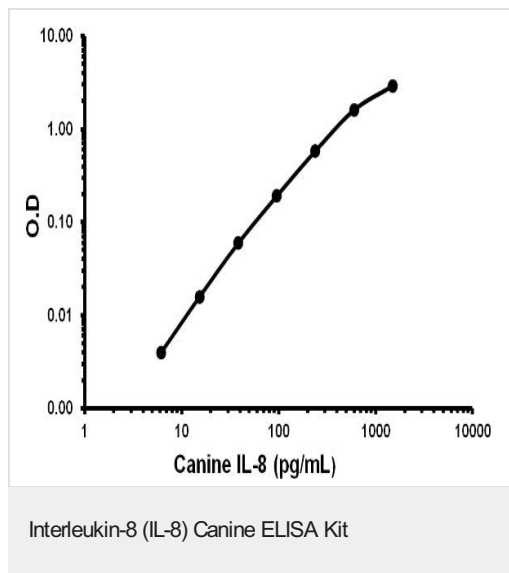
Function IL-8 is a chemotactic factor that attracts neutrophils, basophils, and T-cells, but not monocytes. It is also involved in neutrophil activation. It is released from several cell types in response to an inflammatory stimulus. IL-8(6-77) has a 5-10-fold higher activity on neutrophil activation, IL-8(5-77) has increased activity on neutrophil activation and IL-8(7-77) has a higher affinity to receptors CXCR1 and CXCR2 as compared to IL-8(1-77), respectively.

Sequence similarities Belongs to the intercrine alpha (chemokine CxC) family.

Post-translational modifications Several N-terminal processed forms are produced by proteolytic cleavage after secretion from at least peripheral blood monocytes, leukocytes and endothelial cells. In general, IL-8(1-77) is referred to as interleukin-8. IL-8(6-77) is the most prominent form.

Cellular localization Secreted.

Images



Representative standard curve using ab155465

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