

## Product datasheet

# Canine IL-6 ELISA Kit ab193686

[1 References](#) [2 Images](#)

### Overview

**Product name** Canine IL-6 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** 0.1 ng/ml

**Range** 0.1 ng/ml - 25 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	79.27	68% - 90%
Plasma	78.06	69% - 91%
Cell culture media	108	99% - 118%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Dog

**Product overview** Abcam's IL-6 Canine ELISA Kit (ab193686) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of canine IL-6 in serum, plasma and cell culture supernatant.

This assay employs an antibody specific for canine IL-6 coated on a 96-well plate. Standards and

samples are pipetted into the wells and the immobilized antibody captures IL-6 present in the samples. The wells are washed and biotinylated anti-canine IL-6 antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of IL-6 bound in each well. Addition of the Stop Solution will change 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 Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated canine IL-6 detection antibody (lyophilized)	2 vials
Canine IL-6 Standards (lyophilized)	2 vials
Pre-coated canine IL-6 microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Cytokine with a wide variety of biological functions. It is a potent inducer of the acute phase response. Plays an essential role in the final differentiation of B-cells into Ig-secreting cells. Involved in lymphocyte and monocyte differentiation. It induces myeloma and plasmacytoma growth and induces nerve cells differentiation. Acts on B-cells, T-cells, hepatocytes, hematopoietic progenitor cells and cells of the CNS. Also acts as a myokine. It is discharged into the bloodstream after muscle contraction and acts to increase the breakdown of fats and to improve insulin resistance.

**Involvement in disease** Genetic variations in IL6 are associated with susceptibility to rheumatoid arthritis systemic juvenile (RASJ) [MIM:604302]. An inflammatory articular disorder with systemic-onset beginning before the age of 16. It represents a subgroup of juvenile arthritis associated with severe extraarticular features and occasionally fatal complications. During active phases of the disorder, patients display a typical daily spiking fever, an evanescent macular rash, lymphadenopathy, hepatosplenomegaly, serositis, myalgia and arthritis.  
Note=A IL6 promoter polymorphism is associated with a lifetime risk of development of Kaposi sarcoma in HIV-infected men.

**Sequence similarities** Belongs to the IL-6 superfamily.

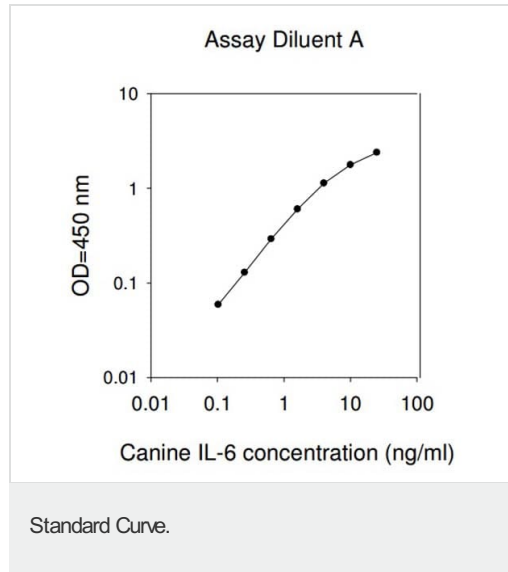
**Post-translational** N- and O-glycosylated.

## modifications

### Cellular localization

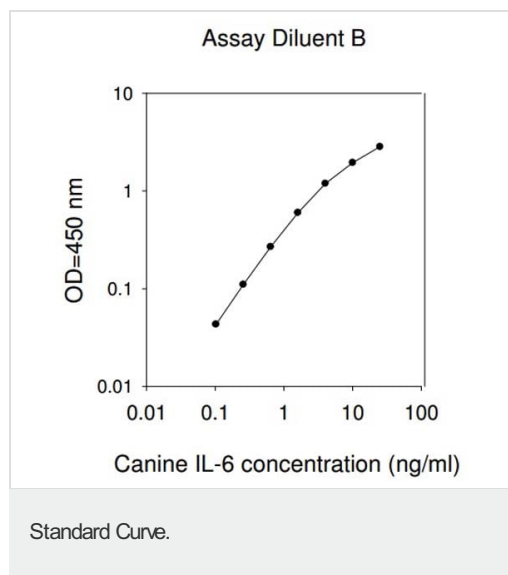
Secreted.

## Images



Standard Curve in Assay Diluent A using ab193686 IL-6 Canine ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Standard Curve in Assay Diluent B using ab193686 IL-6 Canine ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

**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