

Bovine IL-6 ELISA Kit ab205080

1 References 1 Image

Overview

Product name	Bovine 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				< 10%
Sample type	Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	0.03 ng/ml				
Range	15.63 pg/ml - 1000 pg/ml				
Recovery	> 85 %				
	Sample specific recovery				
	Sample type	Average %		Range	
	Serum	> 85		0.063ng/ml - 4ng/ml	
Assay time	2h 10m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Cow				
Product overview	IL-6 in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-6 protein in bovine plasma and serum.				
Notes	In this assay the Interleukin 6 present in samples reacts with the anti-Interleukin 6 antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, the Detection antibody, biotin conjugated anti-IL-6 is added and complexes are formed. Following a wash step, the horseradish peroxidase (HRP) conjugated Streptavidin is				

added and complexes are formed. After another washing step, the complexes are assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of IL-6 in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of IL-6 in the test sample. The quantity of IL-6 in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X Detection Antibody Conjugate	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentration	1 x 50ml
Anti-Bovine Interleukin 6 ELISA Microplate	1 unit
Bovine Interleukin 6 Calibrator (Lyophilized)	1 vial
Chromogen Substrate Solution	1 x 12ml
HRP-STREPTAVIDIN 100X	1 x 150µl
Stop Solution	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.

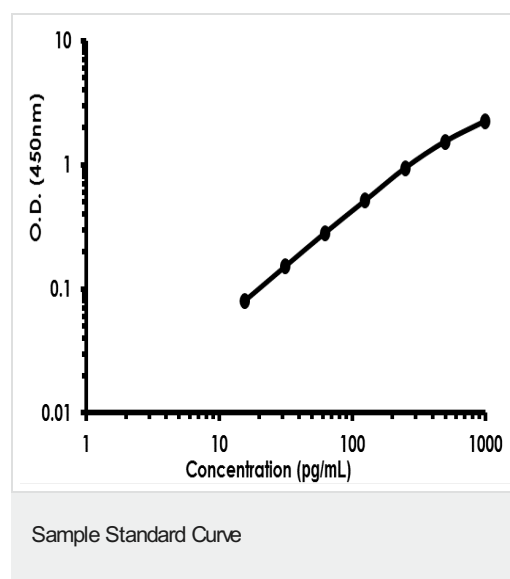
Involvement in disease Genetic variations in IL6 are associated with susceptibility to rheumatoid arthritis systemic juvenile (RASJ) [MIM:604302]. An inflammatory articular disorder with systemic-onset beginning before the age of 16. It represents a subgroup of juvenile arthritis associated with severe 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 modifications N- and O-glycosylated.

Cellular localization Secreted.

Images



Representative standard curve using ab205080 Bovine IL-6 ELISA Kit.

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