

2 References

Overview

Product name	Canine C Reactive Protein 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, Other biological fluids
--------------------	--

Assay type	Sandwich (quantitative)
-------------------	-------------------------

Sensitivity 0.6467 ng/ml

Range 3.125 ng/ml - 200 ng/ml

Recovery	Sample specific recovery
----------	--------------------------

Sample type	Average %	Range
Serum	> 85	% - %

Assay duration	Multiple steps standard assay
-----------------------	-------------------------------

Species reactivity Reacts with: Dog

Product overview Abcam's C Reactive Protein Dog ELISA kit is a highly sensitive two-site enzyme linked immunoassay (ELISA) for the quantitative measurement of C Reactive Protein in dog serum and plasma samples.

In this assay the C Reactive Protein present in samples reacts with the anti-C Reactive Protein antibodies which have been adsorbed to the surface of polystyrene microtiter wells. After the removal of unbound proteins by washing, anti-C Reactive Protein antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound C Reactive Protein. Following another washing step, the enzyme bound to the immunosorbent is 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 C Reactive Protein in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of C Reactive Protein in the test sample. The quantity of C Reactive Protein in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Platform Microplate

Properties

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

Components	1 x 96 tests
100X HRP-conjugated anti-dog C Reactive Protein antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Dog C Reactive Protein Calibrator (lyophilized)	1 vial
Dog C Reactive Protein ELISA Microplate	1 unit
Stop Solution	1 x 12ml

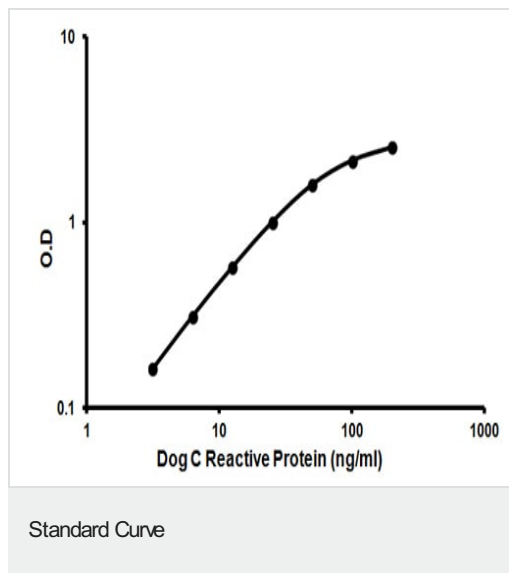
Function Displays several functions associated with host defense: it promotes agglutination, bacterial capsular swelling, phagocytosis and complement fixation through its calcium-dependent binding to phosphorylcholine. Can interact with DNA and histones and may scavenge nuclear material released from damaged circulating cells.

Tissue specificity Found in plasma.

Sequence similarities Belongs to the pentaxin family.
Contains 1 pentaxin domain.

Cellular localization Secreted.

Images



Representative standard curve using ab157698 C Reactive Protein
Dog 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