



# **Canine (Dog) IgA ELISA**

Catalog Number:

**IGA49-K01 (1 x 96 wells)**

*For Research Use Only. Not for use in diagnostic procedures.*

*v. 1.0*

---

EAGLE BIOSCIENCES, INC.  
20A Northwest Blvd., Suite 112, Nashua, NH 03063  
Phone: 617-419-2019 Fax: 617-419-1110  
[WWW.EAGLEBIO.COM](http://WWW.EAGLEBIO.COM)



## Introduction

The Eagle Biosciences Canine (Dog) IgA ELISA Assay Kit is designed for the quantitative determination of canine (dog) IgA in complex samples (serum or other biological samples). The Dog IgA ELISA Assay Kit is for research use only and should not be used in diagnostic procedures.

## Principle of the Assay

The determination of dog IgA is carried out as direct sandwich ELISA. An antibody specific for dog IgA has been pre-coated onto a microplate. Standards and samples are pipetted into the wells and any IgA present is bound. After washing away any unbound substances, an enzyme-linked antibody is added. Following a wash, a substrate solution is added to the wells and color develops in proportion to the amount of antibody conjugate. The absorption at 450 nm is proportional to the IgA concentration.

## Assay Procedure

Incubation of samples/standards	100 µl/well	1 hr with shaking or 2 hrs without shaking
Wash		
Incubation of detection antibody	100 µl/well	1 hr with shaking or 2 hrs without shaking
Wash		
Substrate incubation	100 µl/well	10 min, room temperature
Add stop solution	100 µl/well	
Read at 450 nm		



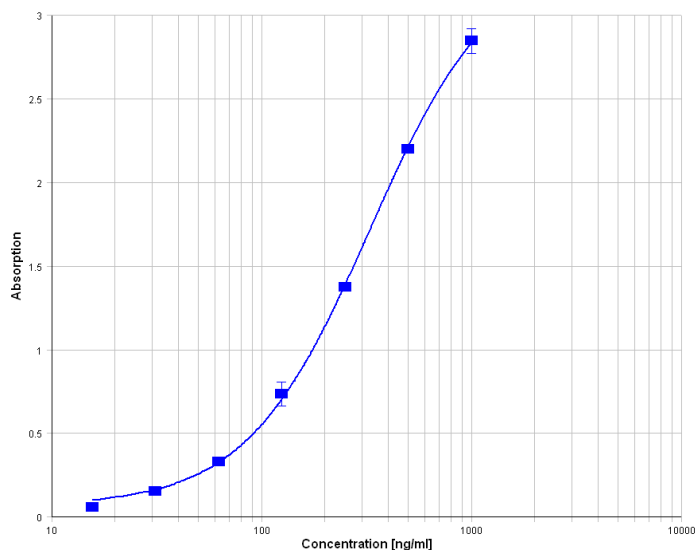
## Performance Characteristics

Standard curve:  
7 standards between 15.6 ng/ml and 1000 ng/ml

Sensitivity:  
156 ng/ml (sample dilution 1:10)

Sample materials:  
Serum, Biological Fluids, Protein solutions

## Typical Standard Curve



## Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

*For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.*