

Human IgD ELISA Kit ab157708

1 Image

Overview

Product name	Human IgD ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			CV%
				< 10%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			CV%
				< 10%
Sample type	Serum, Plasma, Other biological fluids			
Assay type	Sandwich (quantitative)			
Sensitivity	4.4545 ng/ml			
Range	6.25 ng/ml - 400 ng/ml			
Recovery	> 85 %			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	<p>Abcam's IgD Human ELISA kit is an <i>in vitro</i> enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of IgD in Human samples.</p> <p>In this assay the IgD present in samples reacts with the anti-IgD antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-IgD antibodies conjugated with horseradish peroxidase (HRP) are added. These enzyme-labeled antibodies form complexes with the previously bound IgD. 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 IgD in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of IgD in the test sample. The quantity of IgD in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.</p>			

Properties

Storage instructions

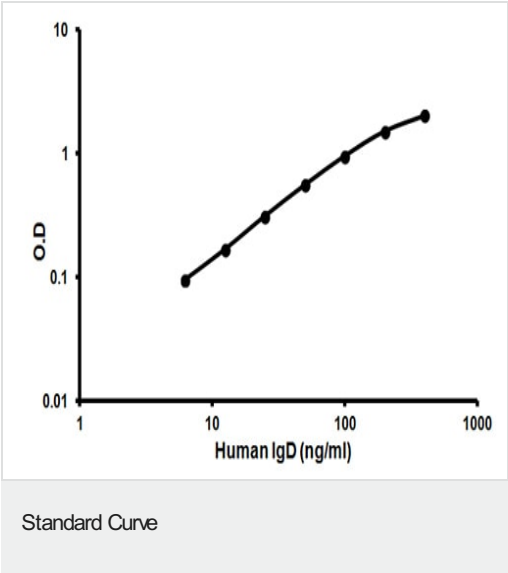
Store at +4°C. Please refer to protocols.

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

Relevance

Little is known about the normal function of IgD, and few clinical signs or symptoms are associated with its absence. Individuals with low or absent levels of IgD do not appear unusually predisposed to infections. In mice, IgD may substitute for some functions of IgM when IgM is absent. IgD is expressed on peripheral B cells and is used as a pan B cell marker.

Images



Representative standard curve using ab157708 IgD Human 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