

Human CD200 ELISA Kit ab155429

1 Image

Overview

Product name	Human CD200 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	Serum, Plasma, Cell culture media				
Assay type	Sandwich (quantitative)				
Sensitivity	< 20 pg/ml				
Range	0.025 ng/ml - 6 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		= 86.14		77% - 95%
	Plasma		= 78.35		67% - 86%
	Cell culture media		= 81.85		74% - 89%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's CD200 Human ELISA (Enzyme-Linked Immunosorbent Assay) Kit is an <i>in vitro</i> enzyme-linked immunosorbent assay designed for the quantitative measurement of Human in serum, plasma and cell culture supernatants.				

the immobilized antibody. The wells are washed and biotinylated anti-Human CD200 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated Streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of CD200 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Notes** Optimisation may be required with urine samples.

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

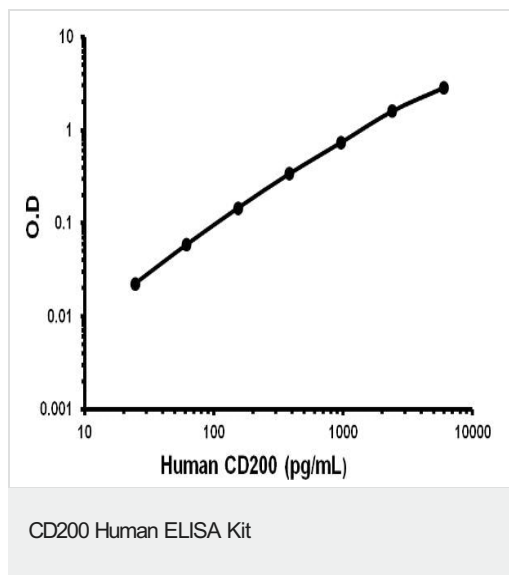
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
240X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Human CD200	2 vials
CD200 Microplate (12 x 8 wells)	1 unit
Recombinant Human CD200 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Costimulates T-cell proliferation. May regulate myeloid cell activity in a variety of tissues.

**Sequence similarities** Contains 1 Ig-like C2-type (immunoglobulin-like) domain.  
Contains 1 Ig-like V-type (immunoglobulin-like) domain.

**Cellular localization** Cell membrane.

## Images



Representative standard curve using ab155429

**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