

## Product datasheet

# Human CD163 ELISA Kit (M130) $\alpha$ b155428

[2 Images](#)

### Overview

**Product name** Human CD163 ELISA Kit (M130)

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
overall				< 12%

**Sample type** Cell culture supernatant, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** < 30 pg/ml

**Range** 0.032 ng/ml - 8 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	= 104.8	97% - 113%
Plasma	= 102.1	94% - 110%
Cell culture media	= 99.57	93% - 106%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Does not react with:** Mouse

**Product overview**

Abcam's CD163 (M130) Human ELISA (Enzyme-Linked Immunosorbent Assay) Kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of Human CD163 in serum, plasma (EDTA plasma is not suitable for use in this assay due to its chelating

property) and cell culture supernatants.

This assay employs an antibody specific for Human CD163 (M130) coated on a 96-well plate. Standards and samples are pipetted into the wells and CD163 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CD163 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 CD163 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.

**Tested applications** **Suitable for:** Sandwich ELISA

**Platform** Microplate

## Properties

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

Components	1 x 96 tests
200x HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CD163	2 vials
CD163 Microplate (12 x 8 wells)	1 unit
Recombinant Human CD163 Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Cellular localization** Cell Membrane

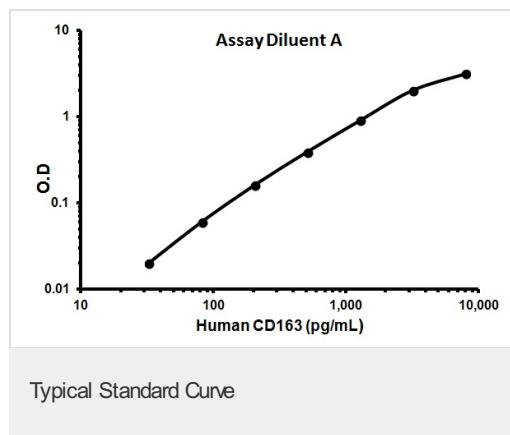
## Applications

Our [Abpromise guarantee](#) covers the use of **ab155428** in the following tested applications.

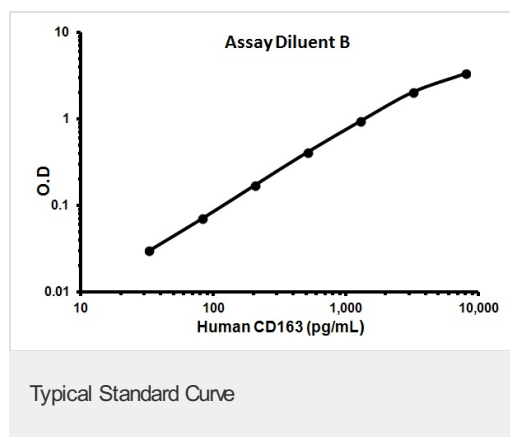
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

## Images



Representative standard curve using ab155428 - Assay Diluent A.



Representative standard curve using ab155428 - Assay Diluent B

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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 <http://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