

Human CD14 ELISA Kit ab45523

1 References 2 Images

Overview

Product name Human CD14 ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Sample A	8	210.5ng/ml	2.6	1.23%
	Sample B	8	23.4ng/ml	1.14	4.6%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Sample A	8	210.5ng/ml	1.55	0.73%
	Sample B	8	24ng/ml	2	8.3%

Sample type Cell culture supernatant, Milk, Serum, Plasma, Cell Lysate

Assay type Sandwich (quantitative)

Sensitivity < 6 ng/ml

Range 6.89 ng/ml - 220 ng/ml

Assay time 3h 45m

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview Abcam's CD14 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of CD14 in Human serum, plasma, buffered solutions or cell culture media.

A monoclonal antibody specific for CD14 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known CD14 concentrations and unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for CD14 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB

substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of CD14 present in the samples.

This kit will recognize both endogenous and recombinant Human CD14.

Get higher sensitivity in only 90 minutes with Human CD14 ELISA Kit ([ab208983](#)) from our SimpleStep ELISA® range.

Platform Microplate

Properties

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

Components	Identifier	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Biotinylated anti-CD14		2 x 0.3ml	1 x 0.3ml
CD14 Microplate (12 x 8 well strips)		2 units	1 unit
CD14 Standard (lyophilized)		4 vials	2 vials
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5ml	2 x 5ml

Function Cooperates with MD-2 and TLR4 to mediate the innate immune response to bacterial lipopolysaccharide (LPS). Acts via MyD88, TIRAP and TRAF6, leading to NF-kappa-B activation, cytokine secretion and the inflammatory response. Up-regulates cell surface molecules, including adhesion molecules.

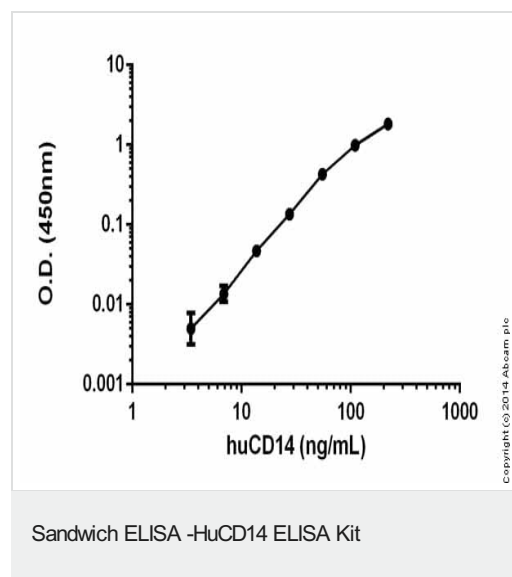
Tissue specificity Expressed strongly on the surface of monocytes and weakly on the surface of granulocytes; also expressed by most tissue macrophages.

Sequence similarities Contains 11 LRR (leucine-rich) repeats.

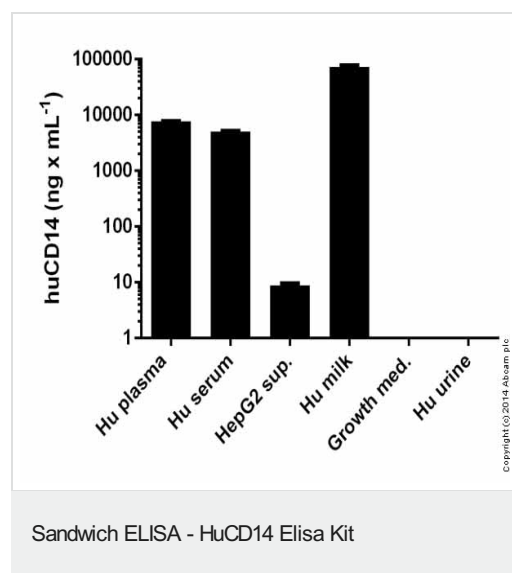
Post-translational modifications N- and O- glycosylated. O-glycosylated with a core 1 or possibly core 8 glycan.

Cellular localization Cell membrane.

Images



Colorimetric standard curve: mean of duplicates (+/-SD) with background readings subtracted



hu CD14 measured in biologicals showing quantity (ng) per mL of tested sample

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