

Human TLR2 ELISA Kit (CD282) ab131556

[1 References](#) [3 Images](#)

Overview

Product name Human TLR2 ELISA Kit (CD282)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	369pg/ml	18.81	= 4.3%
2	16	2018pg/ml	133.18	= 6.6%
3	16	5303pg/ml	212.12	= 4%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	392pg/ml	22.34	= 5.7%
2	24	1927pg/ml	150.3	= 7.8%
3	24	5643pg/ml	242.64	= 4.3%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

Assay type Sandwich (qualitative)

Sensitivity < 10 pg/ml

Range 156 pg/ml - 10000 pg/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Human TLR2 (CD282) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human soluble TLR2 in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

A TLR2 (CD282) specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection

polyclonal antibody from goat, specific for TLR2 is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human TLR2 amount of sample captured in plate.

Platform Microplate

Properties

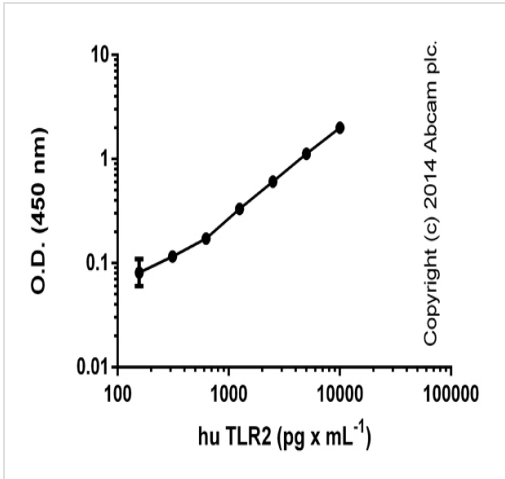
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human TLR2 coated antibody Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human TLR2 antibody		1 x 100µl
Lyophilized recombinant Human TLR2 standard		2 x 10ng
Plate Seal		1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

Relevance Cooperates with LY96 to mediate the innate immune response to bacterial lipoproteins and other microbial cell wall components. Cooperates with TLR1 to mediate the innate immune response to bacterial lipoproteins or lipopeptides. Acts via MYD88 and TRAF6, leading to NF-kappa-B activation, cytokine secretion and the inflammatory response. May also promote apoptosis in response to lipoproteins. Recognizes mycoplasmal macrophage-activating lipopeptide-2kD (MALP-2), soluble tuberculosis factor (STF), phenol-soluble modulin (PSM) and B.burgdorferi outer surface protein A lipoprotein (OspA-L) cooperatively with TLR6.

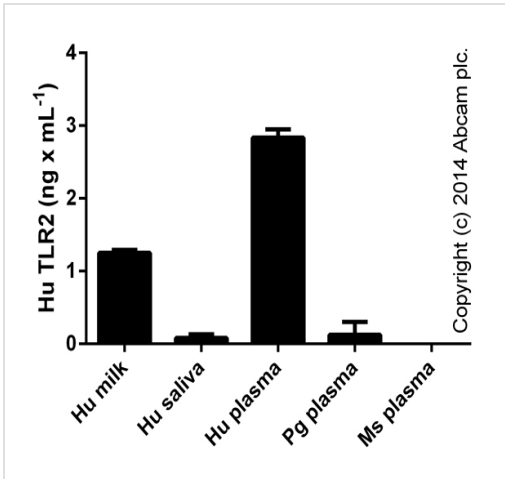
Cellular localization Membrane; Single-pass type I membrane protein.

Images



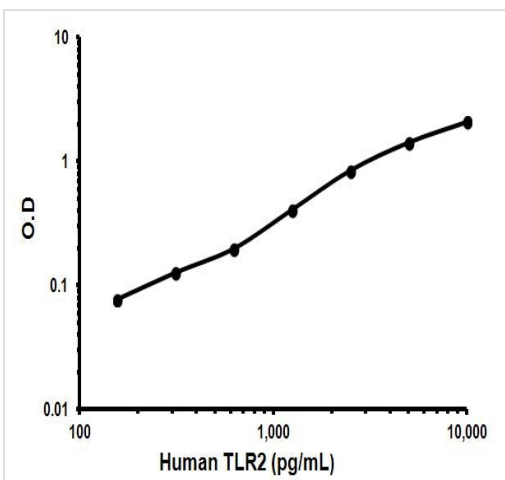
Standard curve with background signal subtracted (duplicates; +/- SD).

ELISA: TLR2 (CD282) Human ELISA Kit (ab131556)



TLR2 measured in undiluted biological fluids (duplicates +/- SD).

ELISA: TLR2 (CD282) Human ELISA Kit (ab131556)



Representative standard curve using ab131556

Typical standard curve

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