

Product datasheet

Human TLR3 ELISA Kit (CD283) ab131557

[1 Image](#)

Overview

Product name	Human TLR3 ELISA Kit (CD283)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, 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	<p>Abcam's Human TLR3 (CD283) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human TLR3 (CD283) in cell culture supernatants, serum and plasma (heparin, EDTA).</p> <p>A TLR3 (CD283) 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 TLR3 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 TLR3 amount of sample captured in plate.</p>
Tested applications	Suitable for: Sandwich ELISA
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 TLR3 antibody Microplate (12 x 8 wells)		1 unit

Components	Identifier	1 x 96 tests
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 130µl
Biotinylated anti-Human TLR3 antibody		1 x 130µl
Lyophilized recombinant Human TLR3 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

Relevance

Key component of innate and adaptive immunity. TLRs (Toll-like receptors) control host immune response against pathogens through recognition of molecular patterns specific to microorganisms. TLR3 is a nucleotide-sensing TLR which is activated by double-stranded RNA, a sign of viral infection. Acts via MYD88 and TRAF6, leading to NF-kappa-B activation, cytokine secretion and the inflammatory response.

Cellular localization

Endoplasmic reticulum membrane; Single-pass type I membrane protein. Endosome membrane.

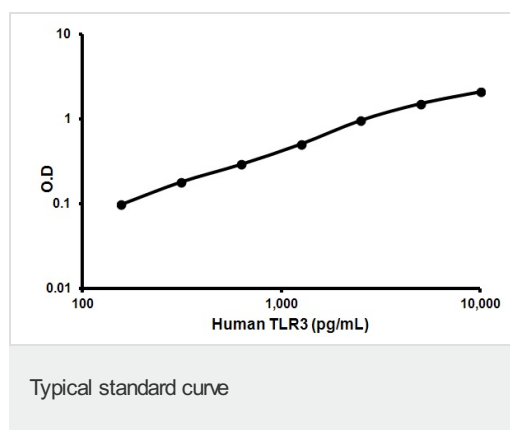
Applications

Our [Abpromise guarantee](#) covers the use of **ab131557** 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 ab131557

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