

Human CD105 ELISA Kit (Endoglin) ab100507

3 References 1 Image

Overview

Product name	Human CD105 ELISA Kit (Endoglin)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	12.35 pg/ml - 3000 pg/ml
Recovery	> 96 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	96.92	87% - 110%
Serum	120.1	111% - 130%
Plasma	116.7	94% - 132%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's CD105 (Endoglin) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human CD105 in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human CD105 coated on a 96-well plate. Standards and samples are pipetted into the wells and CD105 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CD105 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 CD105 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
Notes	Optimisation may be required with urine samples.
Platform	Microplate

Properties

Storage instructions

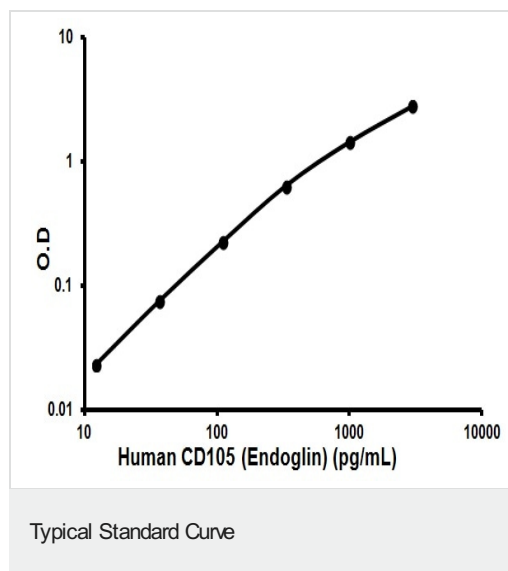
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated anti-Human CD105	2 vials
CD105 Microplate (12 x 8 well strips)	1 unit
Recombinant Human CD105 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance

CD105 or Endoglin is a Type I transmembrane protein, which is highly expressed on human vascular endothelial cells. It exists on an O- and N-glycosylated homodimer. Up regulation of endoglin expression has been demonstrated in tumor vasculature and proliferating cells, suggesting that it is a proliferation associated endothelial marker. CD105 binds to TGF beta 1 and 3 with high affinity but not to TGF beta 2.

Images



Representative Standard Curve using ab100507.

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