

Human Siglec 9 ELISA Kit (CD329) ab113339

[2 Images](#)

Overview

Product name	Human Siglec 9 ELISA Kit (CD329)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 130 pg/ml
Range	0.123 ng/ml - 30 ng/ml
Recovery	85 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	95.43	88% - 102%
Serum	82.7	72% - 91%
Plasma	78.16	70% - 91%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Siglec 9 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Siglec 9 in serum, plasma, cell culture supernatants.

This assay employs an antibody specific for Human Siglec 9 coated on a 96-well plate. Standards and samples are pipetted into the wells and Siglec 9 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Siglec 9 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 Siglec 9 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.

Platform Microplate

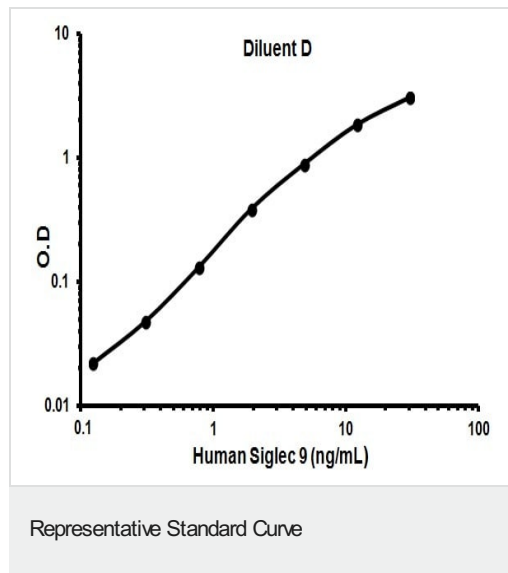
Properties

Storage instructions

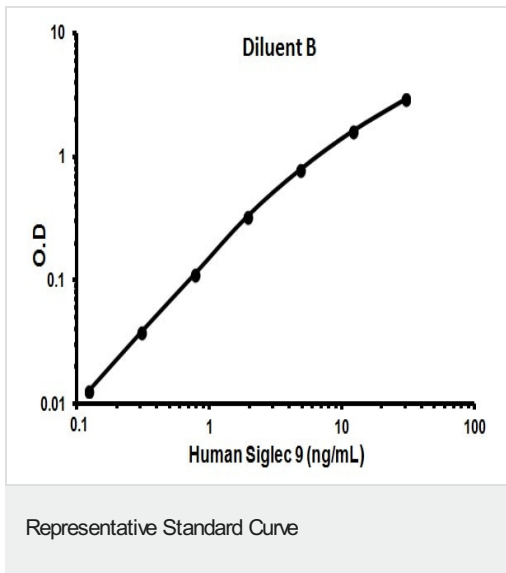
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent D	1 x 15ml
Biotinylated anti-Human Siglec 9	2 vials
Recombinant Human Siglec 9 Standard (lyophilized)	2 vials
Siglec 9 Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Images



Representative Standard Curve using ab113339



Representative Standard Curve using ab113339

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