

Human Siglec 9 ELISA Kit (CD329) ab113339

2 图像

概述

产品名称	人Siglec 9 ELISA试剂盒(CD329)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 130 pg/ml
范围	0.123 ng/ml - 30 ng/ml
回收率	85 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	95.43	88% - 102%
Serum	82.7	72% - 91%
Plasma	78.16	70% - 91%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 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.

说明 Optimization may be required with urine samples.

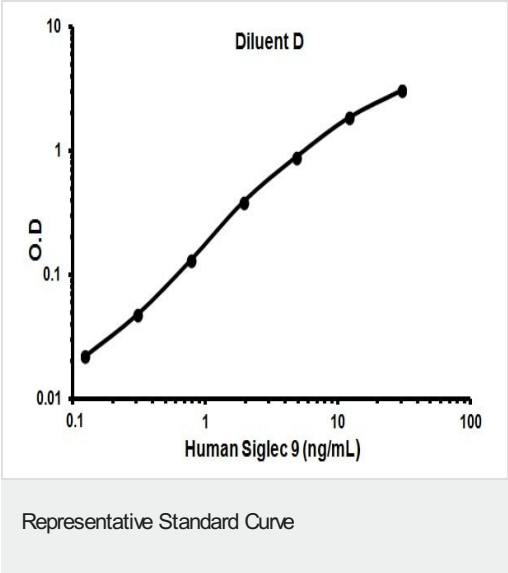
平台 Microplate

性能

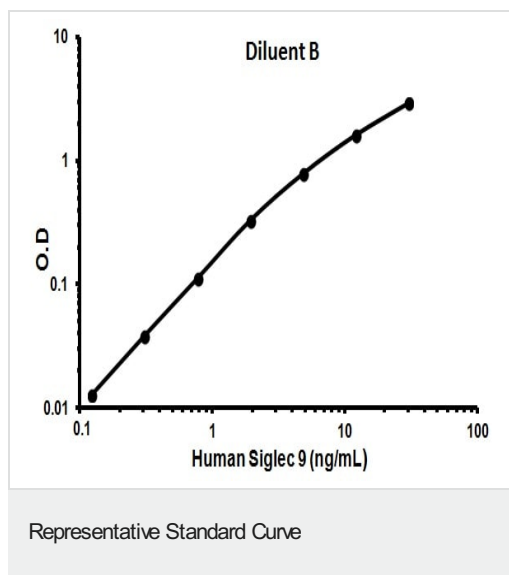
存放说明Store at -20°C. Please refer to protocols.

组件	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

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors