

## Product datasheet

# Human SIGLEC5 ELISA Kit ab100641

[2 Images](#)

### Overview

<b>Product name</b>	Human SIGLEC5 ELISA Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 8 pg/ml
<b>Range</b>	6.86 pg/ml - 5000 pg/ml
<b>Recovery</b>	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	101.3	91% - 108%
Serum	106.3	93% - 114%
Plasma	102.4	91% - 111%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's SIGLEC5 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human SIGLEC5 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human SIGLEC5 coated on a 96-well plate. Standards and samples are pipetted into the wells and SIGLEC5 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human SIGLEC5 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 SIGLEC5 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.

**Tested applications** **Suitable for:** Sandwich ELISA

**Platform** Microplate

## Properties

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

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human SIGLEC5	2 vials
Recombinant Human SIGLEC5 Standard	2 vials
SIGLEC5 Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 25ml
TMB One-Step Substrate Reagent	1 x 12ml

## Relevance

Putative adhesion molecule that mediates sialic-acid dependent binding to cells. Binds equally to alpha2,3-linked and alpha2,6-linked sialic acid. The sialic acid recognition site may be masked by cis interactions with sialic acids on the same cell surface. This molecule is expressed by monocytic/myeloid lineage cells. Found at high levels in peripheral blood leukocytes, spleen, bone marrow and at lower levels in lymph node, lung, appendix, placenta, pancreas and thymus. Expressed by monocytes and neutrophils but absent from leukemic cell lines representing early stages of myelomonocytic differentiation.

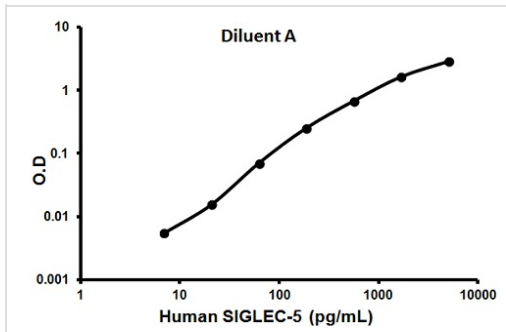
## Applications

Our [Abpromise guarantee](#) covers the use of **ab100641** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

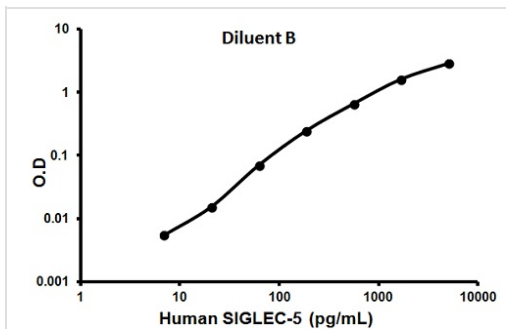
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

## Images



Representative standard curve using  
ab100641

Typical standard curve



Representative standard curve using  
ab100641

Typical standard curve

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### 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 <http://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