

Human SDF1 alpha + SDF1 beta ELISA Kit ab100637

★★★★★ [2 Abreviews](#) [14 References](#) [2 Images](#)

Overview

Product name	Human SDF1 alpha + SDF1 beta ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	= 42 pg/ml				
Range	42 pg/ml - 6000 pg/ml				
Recovery	> 80 %				
	Sample specific recovery				
	Sample type		Average %		Range
	Cell culture supernatant		94.63		86% - 104%
	Plasma		89.34		85% - 104%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Human SDF1 alpha ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab100637) is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human SDF1 alpha in plasma and cell culture supernatants. (Human SDF1 alpha concentration is pretty low in normal plasma, it may not be detected in this assay).				

This assay employs an antibody specific for Human SDF1 alpha coated on a 96-well plate.

Standards and samples are pipetted into the wells and SDF1 alpha present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human SDF1 alpha 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 SDF1 alpha bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect the endogenous Human SDF-1 alpha in normal serum with ab100637, only in serum spiked with Human SDF-1 alpha.

Notes

Optimization may be required with urine samples.

Platform

Microplate

Properties

Storage instructions

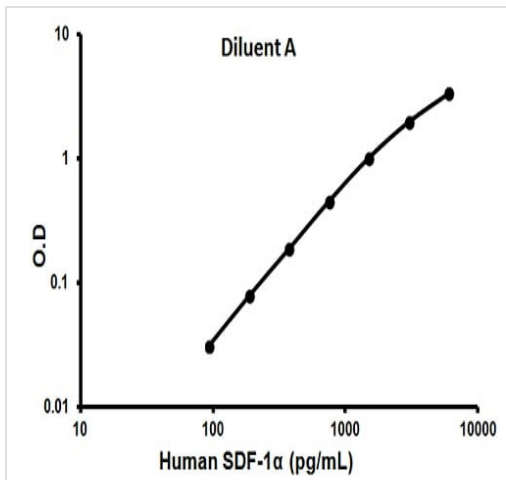
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human SDF1 alpha	2 vials
Recombinant Human SDF1 alpha Standard (lyophilized)	2 vials
SDF1 alpha Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Cellular localization

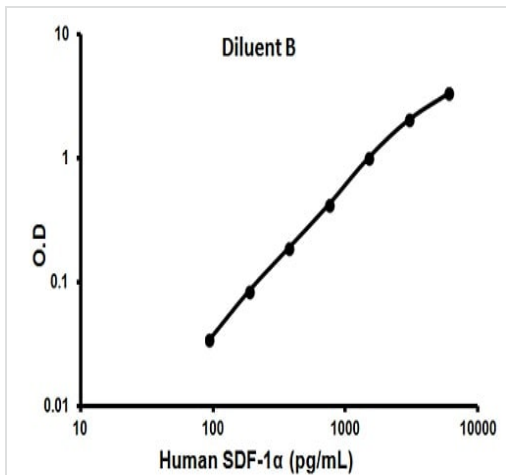
SDF1 beta: Secreted.

Images



Representative standard curve using ab100637

Typical standard curve



Representative standard curve using ab100637

Typical standard curve

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