

Human SDF1 alpha + SDF1 beta ELISA Kit ab100637

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

概述

产品名称 人SDF1 alpha + SDF1 beta ELISA试剂盒

检测方法 Colorimetric

精确度	批次内				
	样品	n	Mean	SD	CV%
	Overall				< 10%

	批次间				
	样品	n	Mean	SD	CV%
	Overall				< 12%

样品类型 Cell culture supernatant, Plasma

检测类型 Sandwich (quantitative)

灵敏度 = 42 pg/ml

范围 42 pg/ml - 6000 pg/ml

回收率 > 80 %

	特定样本回收率		
	样品类型	平均%	范围
	Cell culture supernatant	94.63	86% - 104%
	Plasma	89.34	85% - 104%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Human SDF1 alpha ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab100637) is an *in vitro* 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.

说明

Optimization may be required with urine samples.

平台

Microplate

性能

存放说明

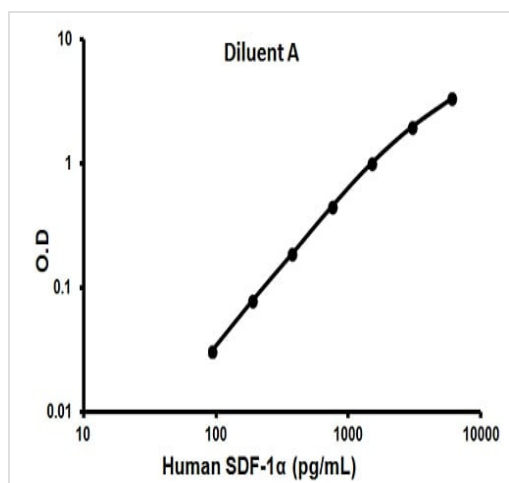
Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
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

细胞定位

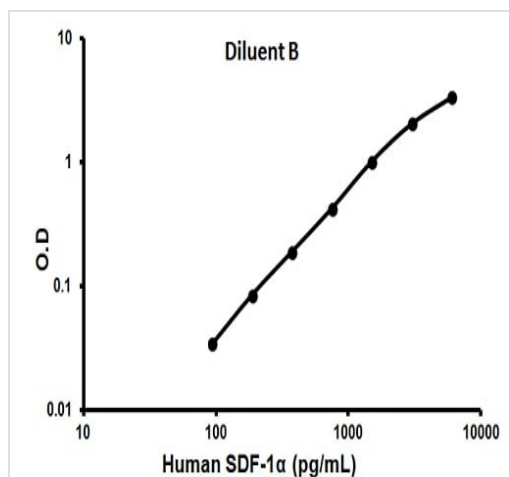
SDF1 beta: Secreted.

图片



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.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