



Product datasheet

Human SCF ELISA Kit ab176109

SimpleStep ELISA

1 References 5 图像

概述

产品名称	人SCF ELISA试剂盒																				
检测方法	Colorimetric																				
精确度	<div>批次内</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Pooled Serum</td><td>24</td><td></td><td></td><td>6%</td></tr></table> <div>批次间</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Pooled Serum</td><td>6</td><td></td><td></td><td>2%</td></tr></table>	样品	n	Mean	SD	CV%	Pooled Serum	24			6%	样品	n	Mean	SD	CV%	Pooled Serum	6			2%
样品	n	Mean	SD	CV%																	
Pooled Serum	24			6%																	
样品	n	Mean	SD	CV%																	
Pooled Serum	6			2%																	
样品类型	Serum, Plasma, Cell culture media, Hep Plasma, Cit plasma																				
检测类型	Sandwich (quantitative)																				
灵敏度	0.51 pg/ml																				
范围	15.63 pg/ml - 1000 pg/ml																				
回收率	<div>特定样本回收率</div> <table><tr><th>样品类型</th><th>平均%</th><th>范围</th></tr><tr><td>Serum</td><td>102</td><td>92% - 111%</td></tr><tr><td>Cell culture media</td><td>106</td><td>100% - 113%</td></tr><tr><td>Hep Plasma</td><td>105</td><td>91% - 115%</td></tr><tr><td>Cit plasma</td><td>89</td><td>79% - 103%</td></tr></table>	样品类型	平均%	范围	Serum	102	92% - 111%	Cell culture media	106	100% - 113%	Hep Plasma	105	91% - 115%	Cit plasma	89	79% - 103%					
样品类型	平均%	范围																			
Serum	102	92% - 111%																			
Cell culture media	106	100% - 113%																			
Hep Plasma	105	91% - 115%																			
Cit plasma	89	79% - 103%																			
检测时间	1h 30m																				
实验步骤	One step assay																				
种属反应性	与反应: Human																				

产品概述

Human Stem Cell Factor (SCF) ELISA kit (ab176109) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of SCF protein in human plasma, serum and cell culture supernatant samples. It uses our proprietary SimpleStep ELISA® technology. Quantitate human SCF with 0.51 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

SCF is the ligand for the receptor-type protein-tyrosine kinase KIT. SCF is expressed in fibroblasts, liver cells, Sertoli cells, endothelial cells, neurons, macrophages, oocytes, Schwann cells and numerous carcinoma cell lines. SCF can be found as a cell membrane protein or as a soluble protein after secretion due to proteolytic processing in the extracellular domain. Both forms have growth factor activities; however the membrane form is also important for cell-cell adhesion and interaction.

Soluble SCF is a heavily N- and O-glycosylated protein that exists as a non-covalently linked homodimer or as a heterotetramer with its receptor. Binding to the receptor, results in KIT activation by autophosphorylation. This leads to multiple signaling cascades (via AKT, GRB2, RAS, MAPK amongst others) that are important in the regulation of cell survival, proliferation, hematopoiesis, stem cell maintenance, gametogenesis, mast cell development, migration, melanogenesis and neurite-inductive activity of dorsal root ganglia neurons.

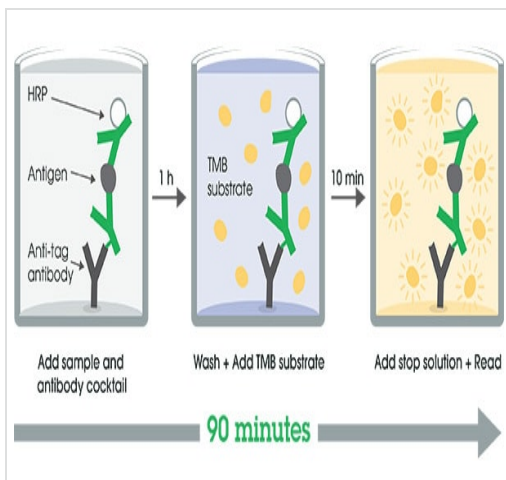
Genetic defects in SCF are the cause of familial progressive hyperpigmentation (melanosis universalis hereditaria). SCF has also been involved in other conditions such as mastocytosis, mast cell leukemia, myeloid leukemia, aplastic anemia and gastrointestinal stromal tumor.

平台 Microplate

性能

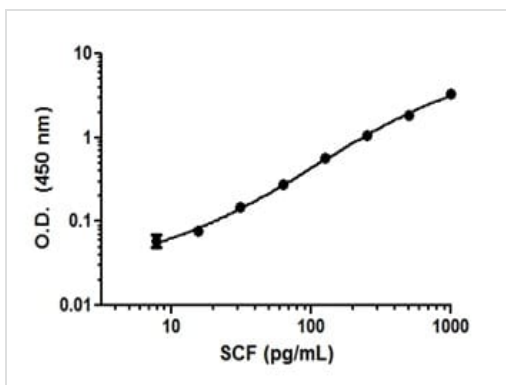
存放说明 Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Human SCF Capture Antibody	1 x 600µl
10X Human SCF Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml



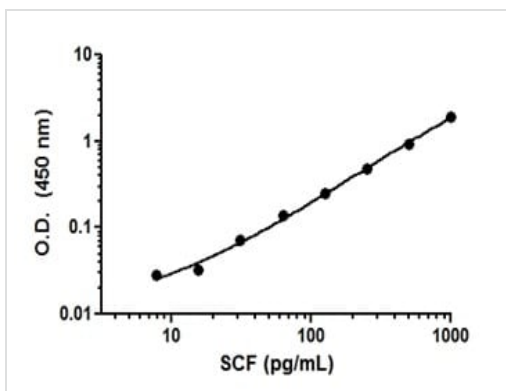
Other - Human SCF ELISA Kit (ab176109)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



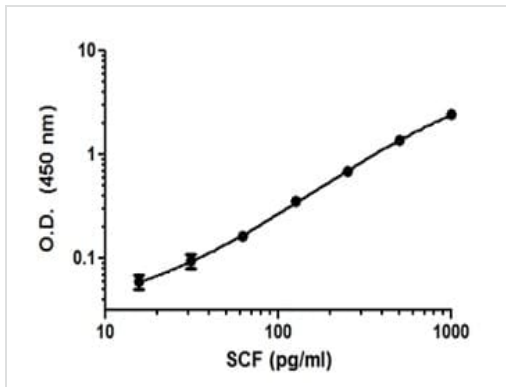
Example of SCF standard curve in Sample Diluent NS.

Background-subtracted data values (mean \pm SD) are graphed.



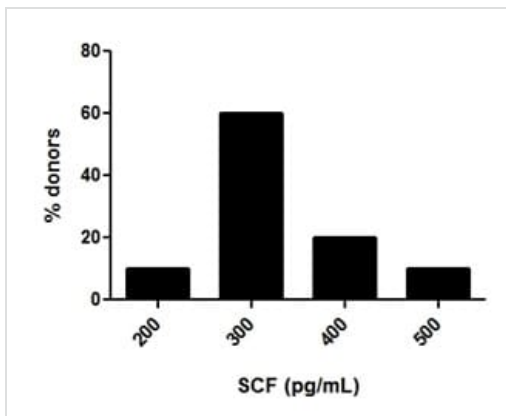
Example of SCF standard curve in Sample Diluent 50BP.

Background-subtracted data values (mean \pm SD) are graphed.



Background-subtracted data values (mean \pm SD) are graphed.

Example of SCF standard curve in 100% cell culture media.



The levels of SCF in serum samples were tested from ten individual healthy donors. Levels were interpolated from the standard curve in Sample diluent 50BP and corrected for sample dilution. The levels of SCF are shown for the percentage of individuals within each 100 pg/mL bin center of the distribution. The mean level of SCF was 330 pg/mL with a range of 185 to 498 pg/mL and a standard deviation of 82 pg/mL.

Frequency histogram of SCF levels in serum of individual normal healthy donors.

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