

Human IL-23 ELISA Kit ab64708

3 References 1 图像

概述

产品名称 人IL-23 ELISA试剂盒

检测方法 Colorimetric

精确度	批次内				
	样品	n	Mean	SD	CV%
	A	6	4637pg/ml	72	1.6%
	B	6	3908pg/ml	36	1%
	C	6	3016pg/ml	72	2.4%

	批次间				
	样品	n	Mean	SD	CV%
	A	9	4439pg/ml	276	6.2%
	B	9	3807pg/ml	353	9.2%
	C	9	2570pg/ml	388	15%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 20 pg/ml

范围 156.25 pg/ml - 5000 pg/ml

回收率 89 %

特定样本回收率		
样品类型	平均%	范围
Serum	118	103% - 134%

检测时间 2h 45m

实验步骤 Multiple steps standard assay

种属反应性

产品概述

平台

性能

存放说明

		与反应: Human	
		Abcam's IL-23 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-23 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.	
		A monoclonal antibody specific for IL-23 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-23 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-23 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-23 present in the samples.	
		This kit will recognize both endogenous and recombinant Human IL-23.	
		Microplate	
		Store at +4°C. Please refer to protocols.	
组件	标识符	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Biotinylated anti-IL-23		1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-23 Microplate(12 x 8 well strips)		1 unit	2 units
IL-23 Standard (Lyophilized)	Yellow	2 vials	4 vials
Plastic Plate Covers		2 units	4 units
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

功能

Associates with IL12B to form the IL-23 interleukin, an heterodimeric cytokine which functions in innate and adaptive immunity. IL-23 may constitute with IL-17 an acute response to infection in peripheral tissues. IL-23 binds to an heterodimeric receptor complex composed of IL12RB1 and IL23R, activates the Jak-Stat signaling cascade, stimulates memory rather than naive T-cells and promotes production of proinflammatory cytokines. IL-23 induces autoimmune inflammation and

thus may be responsible for autoimmune inflammatory diseases and may be important for tumorigenesis.

组织特异性

Secreted by activated dendritic and phagocytic cells and keratinocytes. Also expressed by dermal Langerhans cells (at protein level).

序列相似性

Belongs to the IL-6 superfamily.

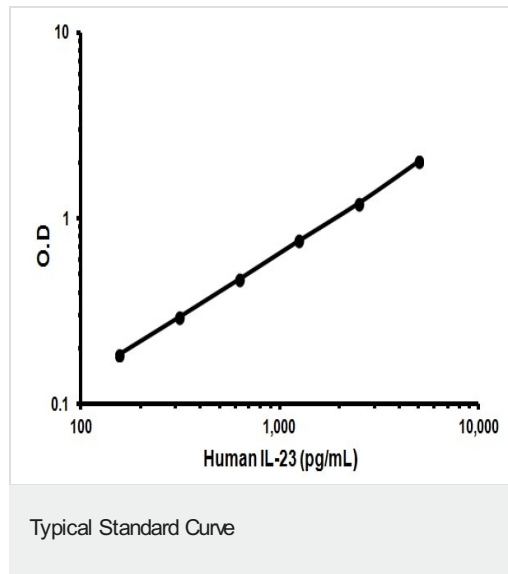
发展阶段

Expressed by newborns dendritic cells.

细胞定位

Secreted. Secreted upon association with IL12B.

图片



Representative Standard Curve using ab64708

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