

Human IL-23 ELISA Kit ab64708

3 References 画像数 1

カルタヘナ法

製品の概要

製品名	Human IL-23 ELISA Kit
検出方法	Colorimetric
再現性	Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
A	6	4637pg/ml	72	1.6%
B	6	3908pg/ml	36	1%
C	6	3016pg/ml	72	2.4%

Inter-Assay (日差再現性)				
サンプル	N	平均値	SD	CV%
A	9	4439pg/ml	276	6.2%
B	9	3807pg/ml	353	9.2%
C	9	2570pg/ml	388	15%

サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 20 pg/ml
検出範囲	156.25 pg/ml - 5000 pg/ml
添加回収試験	89 %

特定サンプルでの回収試験		
サンプルの種類	平均 %	測定範囲
Serum	118	103% - 134%

全工程の試験時間	2h 45m
----------	--------

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>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.</p> <p>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.</p> <p>This kit will recognize both endogenous and recombinant Human IL-23.</p>
試験プラットフォーム	Microplate

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	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

機能 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.

組織特異性

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

配列類似性

Belongs to the IL-6 superfamily.

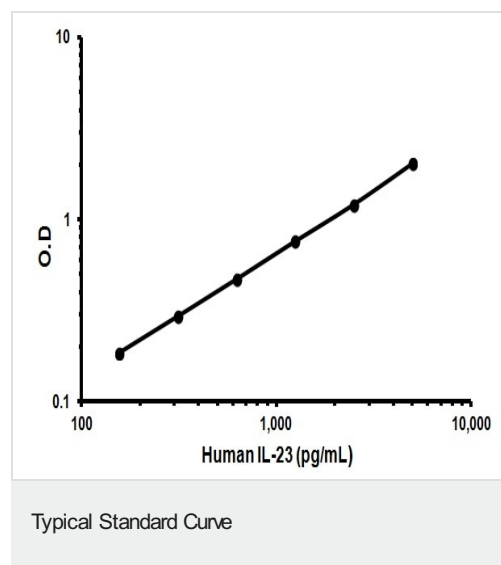
発生段階

Expressed by newborns dendritic cells.

細胞内局在

Secreted. Secreted upon association with IL12B.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors