

Human CCL16 ELISA Kit ab243673

リコンビナント SimpleStep ELISA

画像数 5

製品の概要

製品名	Human CCL16 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%
	sample	8			1.9%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	sample	3			4.7%
サンプルの種類	Cell culture supernatant, Serum, Cit plasma				
アッセイタイプ	Sandwich (quantitative)				
検出感度	1 pg/ml				
検出範囲	2.34 pg/ml - 150 pg/ml				
添加回収試験	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	
	Cell culture supernatant	110		103% - 121%	
	Serum	110		81% - 128%	
	Cit plasma	97		80% - 117%	
全工程の試験時間	1h 30m				
ステップ	One step assay				
種交差性	交差種: Human 非交差種: Cow				
製品の概要	Human CCL16 ELISA Kit (ab243673) is a single-wash 90 min sandwich ELISA designed for the				

quantitative measurement of CCL16 protein in cell culture supernatant, cit plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human CCL16 with 1 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.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

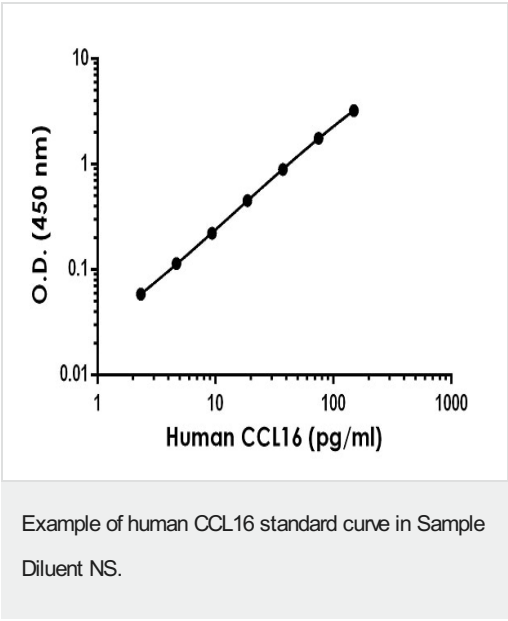
内容	1 x 96 tests
10X Human CCL16 Capture Antibody	1 x 600µl
10X Human CCL16 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 5BI	1 x 6ml
Human CCL16 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

機能

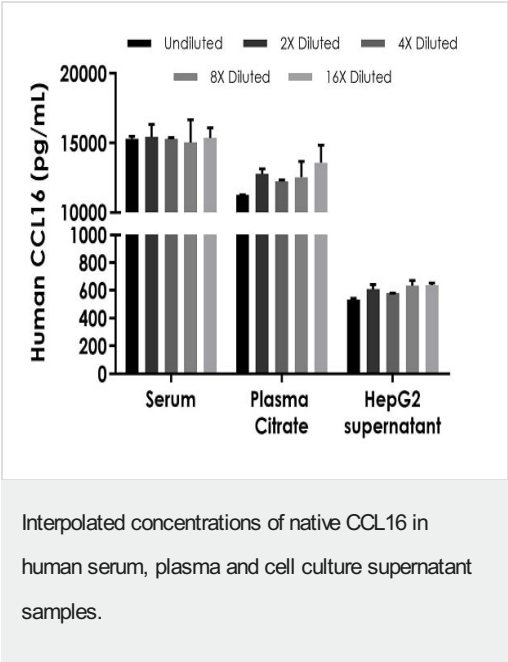
Shows chemotactic activity for lymphocytes and monocytes but not neutrophils. Also shows potent myelosuppressive activity, suppresses proliferation of myeloid progenitor cells. Recombinant SCYA16 shows chemotactic activity for monocytes and THP-1 monocytes, but not for resting lymphocytes and neutrophils. Induces a calcium flux in THP-1 cells that were desensitized by prior

	expression to RANTES.
組織特異性	Mainly expressed in liver, also found in spleen and thymus. Highly expressed in LPS- and IFN-gamma-activated monocytes, weakly in some lymphocytes, including natural killer cells, gamma-delta T-cells, and some T-cell clones.
配列類似性	Belongs to the intercrine beta (chemokine CC) family.
細胞内局在	Secreted.

画像



The CCL16 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



The concentrations of CCL16 were measured in duplicates, interpolated from the CCL16 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1%, plasma (citrate) 1.5%, and HepG2 supernatant 30%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CCL16 concentration was determined to be 15284.9 pg/mL in serum, 12472.9 pg/mL in plasma (citrate) and 599.3 pg/mL in HepG2 supernatant.

Powered by
recombinant antibodies



Research with
confidence
Consistent and
reproducible results



Long-term and
scalable supply
Recombinant
technology



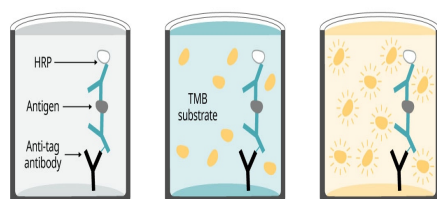
Success from the
first experiment
Confirmed
specificity



Ethical standards
compliant
Animal-free
production

Sandwich ELISA - Human CCL16 ELISA Kit
(ab243673)

To learn more about the advantages of recombinant antibodies see [here](#).



90 minutes

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.

Sandwich ELISA - Human CCL16 ELISA Kit
(ab243673)

**Get more done with
SimpleStep ELISA**



Easy to use
Single-wash 90-minute
protocol



Flexible
Matched antibody pairs
available



Precision antibodies
High sensitivity, specificity
and reproducibility



Scalable
Now in 10-pack and
384-well formats

Sandwich ELISA - Human CCL16 ELISA Kit
(ab243673)

To learn more about the advantages of SimpleStep ELISA® kits
see [here](#).

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