

Monkey IL-6 ELISA Kit ab242233

リコンビナント SimpleStep ELISA

2 References 画像数 7

製品の概要

製品名	Monkey IL-6 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
	cell culture	8		CV%
				5.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	cell culture	3		CV%
				5%
サンプルの種類	Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	3.2 pg/ml			
検出範囲	12.5 pg/ml - 800 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	90	86% - 94%
Serum	76	75% - 77%
Cell culture media	101	95% - 109%
Hep Plasma	103	94% - 114%
EDTA Plasma	85	78% - 92%
Cit plasma	78	76% - 81%

全工程の試験時間	1h 30m
ステップ	One step assay
種交差性	交差種: Monkey, Cynomolgus monkey, Rhesus monkey
製品の概要	<p>Monkey IL-6 ELISA Kit (ab242233) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IL-6 protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Monkey IL-6 with 3.2 pg/ml sensitivity.</p>

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

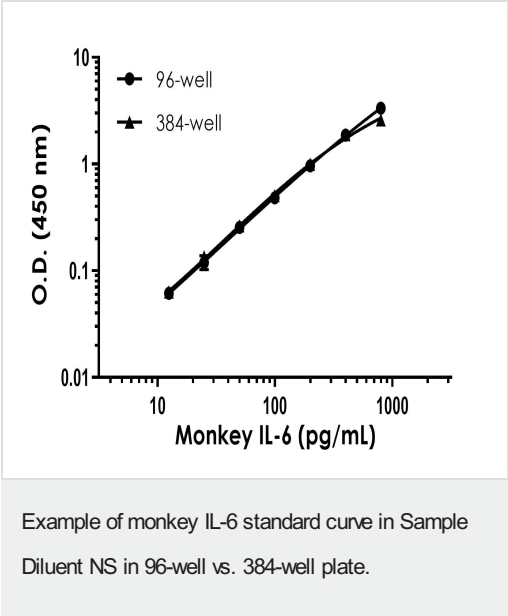
保存方法	Store at +4°C. Please refer to protocols.
------	---

内容	1 x 96 tests	10 x 96 tests	1 x 384 tests
10X Monkey IL-6 Capture Antibody	1 x 600µl	1 x 6000µl	1 x 600µl
10X Monkey IL-6 Detector Antibody	1 x 600µl	1 x 6000µl	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 200ml	1 x 20ml
384 well CaptSure™ microplates	0 x 0 unit	0 x 0 unit	1 unit
Antibody Diluent 5BI	1 x 6ml	10 x 6ml	1 x 6ml
Monkey IL-6 Lyophilized recombinant Protein	2 vials	2 x 10 vials	2 vials
Plate Seals	1 unit	1 x 10 units	1 unit
Sample Diluent NS (ab193972)	1 x 50ml	2 x 250ml	2 x 20ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	1 x 10 units	0 x 0 unit
Stop Solution	1 x 12ml	1 x 120ml	2 x 12ml

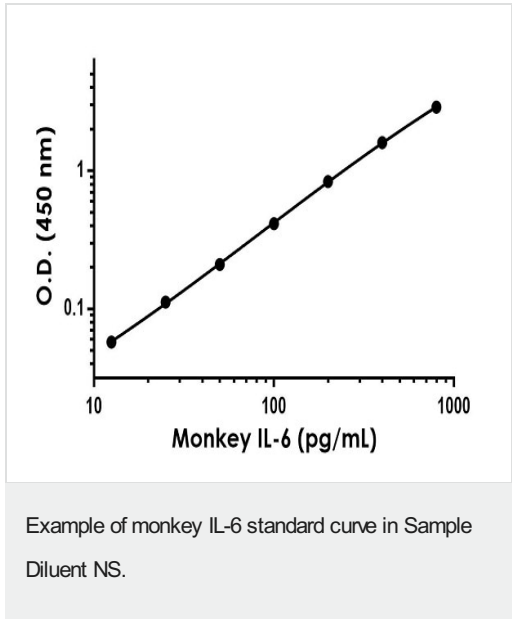
内容	1 x 96 tests	10 x 96 tests	1 x 384 tests
TMB Development Solution	1 x 12ml	1 x 120ml	2 x 12ml

機能	Cytokine with a wide variety of biological functions. It is a potent inducer of the acute phase response. Plays an essential role in the final differentiation of B-cells into Ig-secreting cells Involved in lymphocyte and monocyte differentiation. It induces myeloma and plasmacytoma growth and induces nerve cells differentiation Acts on B-cells, T-cells, hepatocytes, hematopoietic progenitor cells and cells of the CNS. Also acts as a myokine. It is discharged into the bloodstream after muscle contraction and acts to increase the breakdown of fats and to improve insulin resistance.
関連疾患	Genetic variations in IL6 are associated with susceptibility to rheumatoid arthritis systemic juvenile (RASJ) [MIM:604302]. An inflammatory articular disorder with systemic-onset beginning before the age of 16. It represents a subgroup of juvenile arthritis associated with severe extraarticular features and occasionally fatal complications. During active phases of the disorder, patients display a typical daily spiking fever, an evanescent macular rash, lymphadenopathy, hepatosplenomegaly, serositis, myalgia and arthritis. Note=A IL6 promoter polymorphism is associated with a lifetime risk of development of Kaposi sarcoma in HIV-infected men.
配列類似性	Belongs to the IL-6 superfamily.
翻訳後修飾	N- and O-glycosylated.
細胞内局在	Secreted.

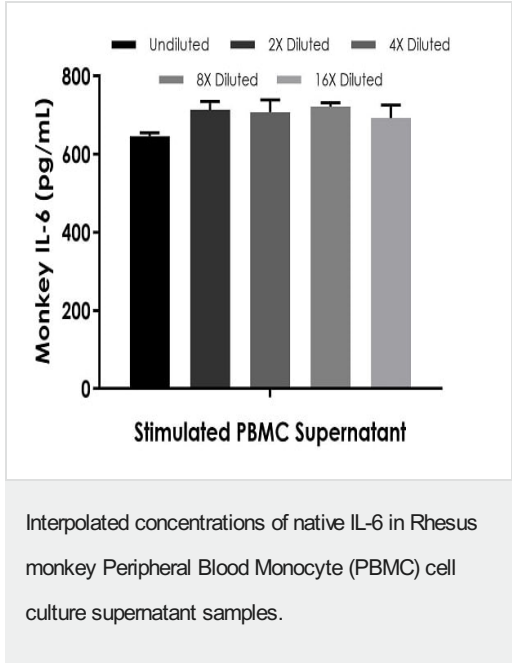
画像



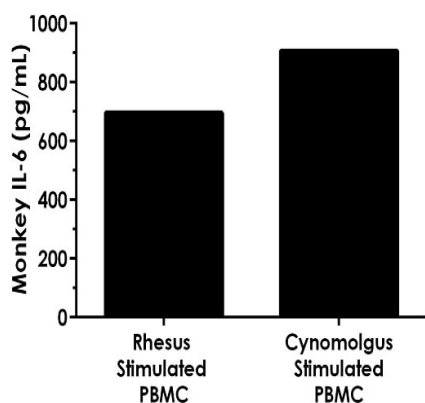
Example of monkey IL-6 standard curve in 96-well vs. 384-well plate. Background-subtracted data values (mean +/- SD) are graphed.



The IL-6 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.

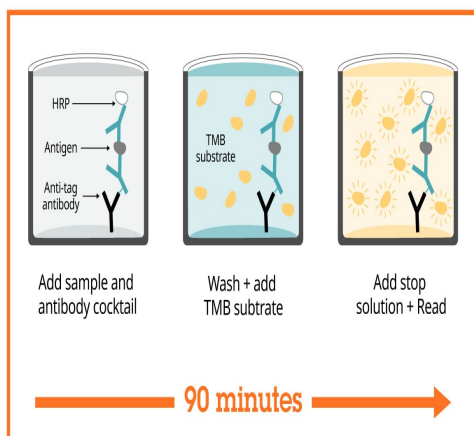


PBMCs were cultured in the presence (stimulated) or absence of 50 ng/mL PMA + 1 µg/mL Ionomycin for 24 hours. The concentrations of IL-6 were measured in duplicates, interpolated from the IL-6 standard curves and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean IL-6 concentration was determined to be 696.3 pg/mL in neat stimulated PBMC cell culture supernatant, and undetectable in unstimulated PBMC cell culture supernatant, and naive cell culture media controls (not shown).



Comparison of native IL-6 in rhesus and cynomolgus stimulated peripheral blood mononuclear cell (PBMC) supernatant samples.

PBMCs were cultured in the presence or absence of 50 ng/mL PMA and 1 µg/mL Ionomycin for 24 hours. The concentrations of IL-6 were measured in duplicates, interpolated from the IL-6 standard curves and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean IL-6 concentration was determined to be 696.3 pg/mL in neat stimulated rhesus PBMC supernatant and 907.3 pg/mL in neat stimulated cynomolgus PBMC supernatant. Values for unstimulated supernatants were undetectable in both rhesus and cynomolgus samples (not shown).



Sandwich ELISA - Monkey IL-6 ELISA Kit
(ab242233)

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.

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 - Monkey IL-6 ELISA Kit
(ab242233)

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

**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 - Monkey IL-6 ELISA Kit
(ab242233)

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