

## Human FAS ELISA Kit ab183360

リコンビナント SimpleStep ELISA

画像数 8

### 製品の概要

製品名 Human FAS 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%
Media	8			3.8%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Media	3			9.5%

サンプルの種類 Cell culture supernatant, Serum, Cell culture media, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 26 pg/ml

検出範囲 46.88 pg/ml - 3000 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	97	91% - 108%
Serum	97	88% - 107%
Cell culture media	100	94% - 105%
Hep Plasma	105	98% - 111%
EDTA Plasma	92	81% - 102%
Cit plasma	98	90% - 105%

**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

**製品の概要** **As of October 15, 2019, Human FAS ELISA kit has been re-developed. We have identified new recombinant monoclonal antibodies to provide improved performance and consistency.**

Human FAS ELISA kit (ab183360) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of FAS protein in human cell culture supernatant, serum, and plasma samples. It uses our proprietary SimpleStep ELISA® technology. Quantitate human FAS with 2 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.

**特記事項**

FAS is a plasma membrane receptor for FASLG ligand (TNFSF6/FASLG). Upon ligand binding, the FASLG Receptor is activated and the adapter molecule FADD recruits caspase-8 to the activated receptor. The resulting death-inducing signaling complex (DISC) performs caspase-8 proteolytic activation which initiates the subsequent cascade of caspases (aspartate-specific cysteine proteases) activation mediating apoptosis. FASLG-mediated apoptosis may have a role in the induction of peripheral tolerance, in the antigen-stimulated suicide of mature T-cells, or both. The secreted isoforms 2 to 6 of the FASLG Receptor block apoptosis as they sequester the Fas ligand.

This kit uses a recombinant protein fragment corresponding to the extracellular portion (amino acid 17-172) of the full length Human Fas Receptor protein as standard sample.

**試験プラットフォーム**

Microplate

**製品の特性**

**保存方法**

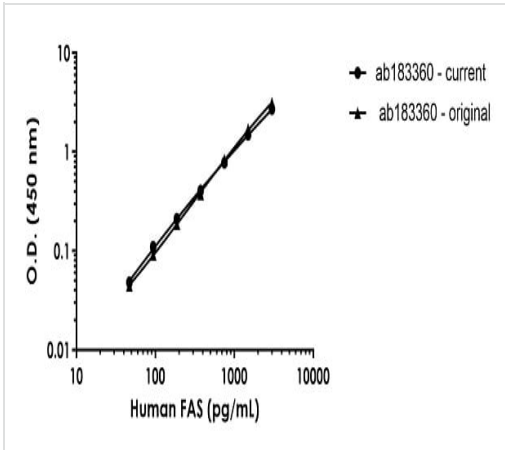
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Human FAS Capture Antibody	1 x 600µl
10X Human FAS Detector Antibody	1 x 600µl

内容	1 x 96 tests
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml
Human FAS 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

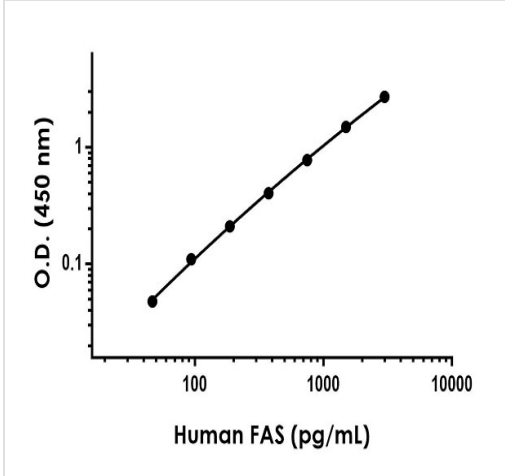
<b>機能</b>	Receptor for TNFSF6/FASLG. The adapter molecule FADD recruits caspase-8 to the activated receptor. The resulting death-inducing signaling complex (DISC) performs caspase-8 proteolytic activation which initiates the subsequent cascade of caspases (aspartate-specific cysteine proteases) mediating apoptosis. FAS-mediated apoptosis may have a role in the induction of peripheral tolerance, in the antigen-stimulated suicide of mature T-cells, or both. The secreted isoforms 2 to 6 block apoptosis (in vitro).
<b>組織特異性</b>	Isoform 1 and isoform 6 are expressed at equal levels in resting peripheral blood mononuclear cells. After activation there is an increase in isoform 1 and decrease in the levels of isoform 6.
<b>関連疾患</b>	Defects in FAS are the cause of autoimmune lymphoproliferative syndrome type 1A (ALPS1A) [MIM:601859]; also known as Canale-Smith syndrome (CSS). ALPS is a childhood syndrome involving hemolytic anemia and thrombocytopenia with massive lymphadenopathy and splenomegaly.
<b>配列類似性</b>	Contains 1 death domain. Contains 3 TNFR-Cys repeats.
<b>ドメイン</b>	Contains a death domain involved in the binding of FADD, and maybe to other cytosolic adapter proteins.
<b>細胞内局在</b>	Secreted and Cell membrane.

## 画像



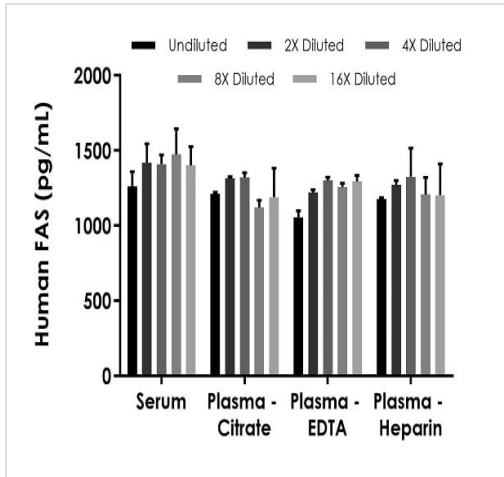
Human FAS standard curve comparison

Standard Curve comparison between the original Human FAS SimpleStep ELISA kit and current Human FAS SimpleStep ELISA kit.



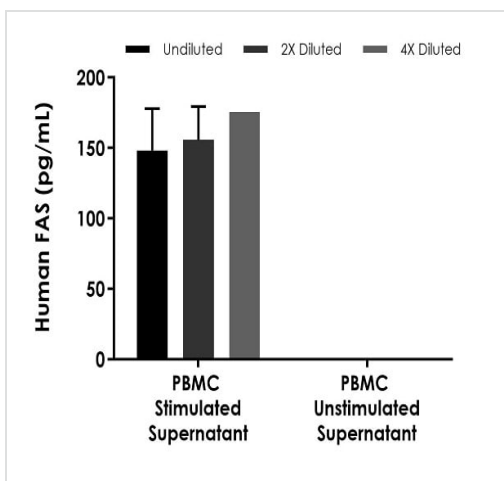
Example of human FAS standard curve in Sample Diluent NS.

The FAS 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.



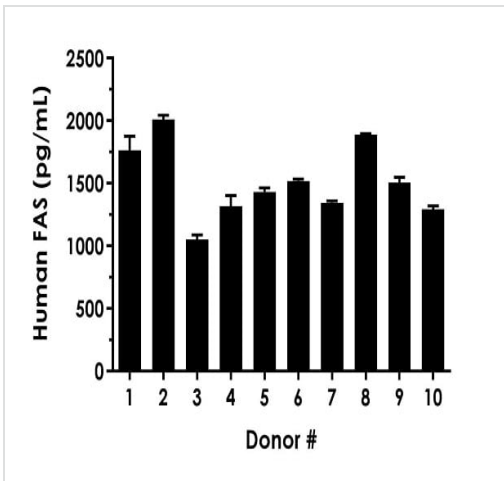
Interpolated concentrations of native FAS in human serum and plasma samples.

The concentrations of FAS were measured in duplicates, interpolated from the FAS standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate, EDTA, heparin) 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean FAS concentration was determined to be 1,393 pg/mL in neat serum, 1,231 pg/mL in neat plasma (citrate) and 1,226 pg/mL in neat plasma (EDTA) and 1,237pg/mL in neat plasma (heparin).



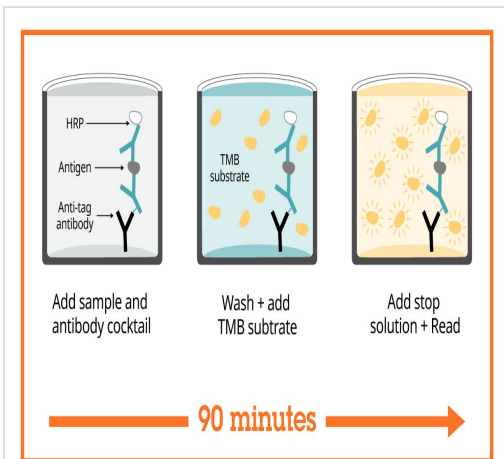
Interpolated concentrations of native FAS in human PBMC cell culture supernatant derived from cells seeded at 107 cells/mL and cultured in the presence or absence of 1.5% PHA for 36 hours.

The concentrations of FAS were measured in duplicate and interpolated from the FAS standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean FAS concentration was determined to be 157 pg/mL in neat stimulated PBMC cell culture supernatant and undetectable in unstimulated PBMC cell culture supernatant.



Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean FAS concentration was determined to be 1,513 pg/mL with a range of 1,050 – 2,009 pg/mL.

Serum from ten individual healthy human male donors was measured in duplicate.



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 FAS ELISA Kit  
(ab183360)

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 FAS ELISA Kit  
(ab183360)

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 - Human FAS ELISA Kit  
(ab183360)

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

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

**Our Promise 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