

## Product datasheet

# Human Fas ELISA Kit (Apo-1) ab100513

2 Images

### Overview

<b>Product name</b>	Human Fas ELISA Kit (Apo-1)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 5 pg/ml
<b>Range</b>	2.74 pg/ml - 2000 pg/ml
<b>Recovery</b>	> 85 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	91.24	82% - 105%
Serum	84.46	78% - 102%
Plasma	86.37	80% - 104%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's Fas Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Fas in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human Fas coated on a 96-well plate. Standards and samples are pipetted into the wells and Fas present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Fas antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Fas bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human FASLG Receptor ELISA Kit (CD95) ([ab183360](#)) from our SimpleStep ELISA® range.

**Notes** Optimization may be required with urine samples

**Tested applications** **Suitable for:** Sandwich ELISA

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
640X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Fas (lyophilized)	2 vials
Fas Microplate (12 strips x 8 wells)	1 unit
Recombinant Human Fas Standard	2 vials
Stop Solution	1 x 8ml
TMB Substrate Reagent	1 x 12ml

**Function** 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).

**Tissue specificity** 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.

**Involvement in disease** 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.

**Sequence similarities** Contains 1 death domain.  
Contains 3 TNFR-Cys repeats.

**Domain** Contains a death domain involved in the binding of FADD, and maybe to other cytosolic adapter proteins.

**Cellular localization** Secreted and Cell membrane.

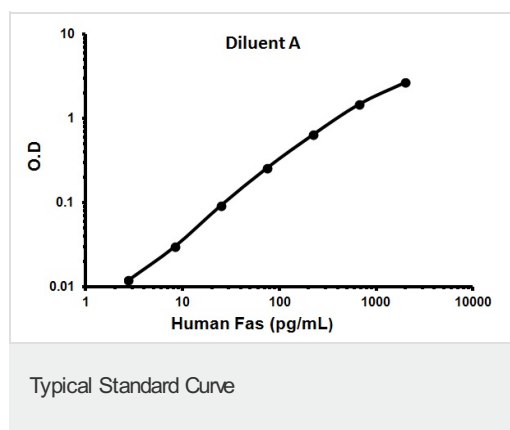
## Applications

Our [Abpromise guarantee](#) covers the use of **ab100513** in the following tested applications.

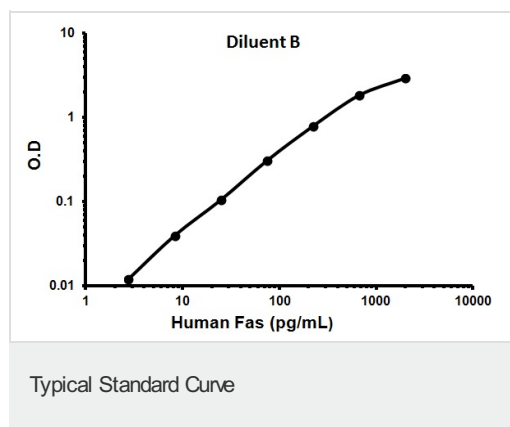
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

## Images



Representative Standard Curve using ab100513.



Representative Standard Curve using ab100513.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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 <http://www.abcam.com/abpromise> or contact our technical team.

## Terms and conditions

---

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