

# Human FAS Ligand ELISA Kit (CD95L) ab45892

[1 References](#)   [1 Image](#)

Overview

Product name	Human FAS Ligand ELISA Kit (CD95L)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Sample A	40	588.4pg/ml	22.6	3.8%
	Sample B	35	302.2pg/ml	14.6	4.5%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	< 12 pg/ml				
Range	62.5 pg/ml - 2000 pg/ml				
Assay time	3h 45m				
Assay duration	Multiple steps standard assay				
Species reactivity	<b>Reacts with:</b> Human				
Product overview	<p>Abcam's Human FAS Ligand <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of FAS Ligand in Human serum, plasma, buffered solutions and cell culture media.</p> <p>A monoclonal antibody specific for FAS Ligand has been coated onto the wells of the microtiter strips provided. Samples, including standards of known FAS Ligand 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 FAS Ligand 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 FAS Ligand present in the samples.</p> <p>This kit will recognize both endogenous and recombinant Human FAS Ligand.</p>				
Platform	Microplate				

**Storage instructions**

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

Components	Identifier	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 7.5ml	1 x 13ml
Biotinylated anti-FAS Ligand	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
Control	Silver	2 vials	4 vials
FAS Ligand Microplate (12 x 8 well strips)		1 unit	2 units
FAS Ligand Standard (Lyophilized)	Yellow	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

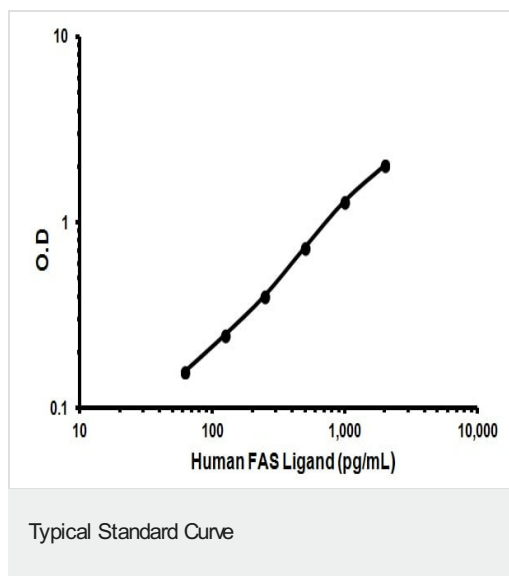
**Relevance**

Fas / APO1 (CD95) is an important member of the tumour necrosis factor (TNF) superfamily involved in membrane-mediated apoptosis. Ligation of Fas by Fas Ligand or an anti Fas cross linking antibody, triggers activation of the caspase cascade. Functional impairment of the Fas / Fas Ligand system is associated with the development and progression of malignancies. Fas gene mutations have been suggested to have a role in testicular germ cell tumours. Tumour cells frequently exhibit de novo expression of Fas Ligand, which plays a significant role in local tissue destruction, metastatic spread, and immune escape of the tumor cells. The apoptosis of lymphocytes, which occurs in autoimmune diseases, is usually induced by the Fas/Fas Ligand system.

**Cellular localization**

Cell Membrane (single-pass type II membrane protein) and Secreted

**Images**



Representative standard curve using ab45892.

**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.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