

Product datasheet

Human Fas Ligand ELISA Kit (APTL) ab100515

★★★★☆ [2 Abreviews](#) [2 References](#) [2 Images](#)

Overview

Product name	Human Fas Ligand ELISA Kit (APTL)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 2 pg/ml
Range	1.37 pg/ml - 1000 pg/ml
Recovery	75 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	76.99	70% - 90%
Serum	76.36	70% - 87%
Plasma	73.42	67% - 86%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

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

This assay employs an antibody specific for Human Fas Ligand coated on a 96-well plate. Standards and samples are pipetted into the wells and Fas Ligand present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human Fas Ligand 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 Ligand bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples

Platform Microplate

Properties

Storage instructions

Store at -20°C. Please refer to protocols.

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

Function

Cytokine that binds to TNFRSF6/FAS, a receptor that transduces the apoptotic signal into cells. May be involved in cytotoxic T-cell mediated apoptosis and in T-cell development. TNFRSF6/FAS-mediated apoptosis may have a role in the induction of peripheral tolerance, in the antigen-stimulated suicide of mature T-cells, or both. Binding to the decoy receptor TNFRSF6B/DcR3 modulates its effects.

Involvement in disease

Defects in FASLG are the cause of autoimmune lymphoproliferative syndrome type 1B (ALPS1B) [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

Belongs to the tumor necrosis factor family.

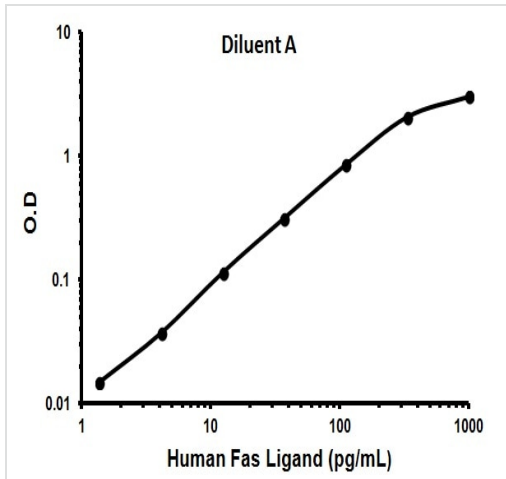
Post-translational modifications

N-glycosylated.
The soluble form derives from the membrane form by proteolytic processing.

Cellular localization

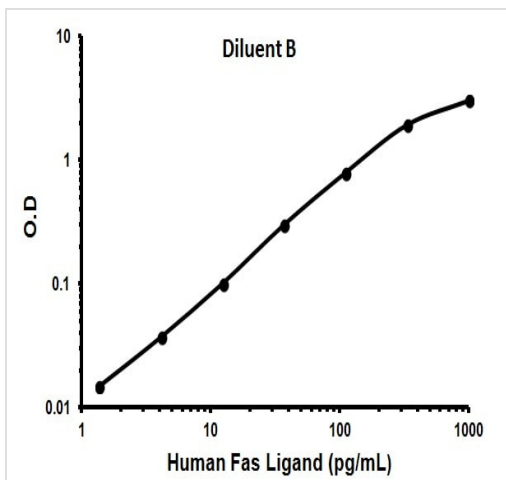
Cell membrane. Secreted. May be released into the extracellular fluid, probably by cleavage from the cell surface.

Images



Representative Standard Curve using ab100515.

Typical Standard Curve



Representative Standard Curve using ab100515.

Typical Standard Curve

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.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