

Human Fas Ligand ELISA Set (without plates) ab83703

[1 Image](#)

Overview

Product name	Human Fas Ligand ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 44 pg/ml
Range	62.5 pg/ml - 2000 pg/ml
Assay time	3h 45m
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's Fas Ligand (Fas-L) Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Fas-L in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The Fas L ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>This ELISA Set is intended to measure Fas L in cell -culture supernatants, as well as in serum samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>SPECIFICITY: Recognizes both natural and recombinant soluble human Fas-Ligand RANGE: 62.5pg/ml - 2000pg/ml SENSITIVITY: <12pg/ml</p>
Tested applications	Suitable for: Sandwich ELISA
Platform	Reagents

Properties

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

Components	Identifier	5 x 96 tests	10 x 96 tests
Detection biotinylated anti-Fas Ligand antibody		1 x 0.55ml	2 x 0.55ml
Fas Ligand Capture Antibody		1 x 0.5ml	2 x 0.5ml
FAS Ligand Standard (Lyophilized)	Yellow	5 vials	10 vials
Ready-to-use TMB		2 vials	4 vials
Streptavidin-HRP		1 vial	2 vials

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 tumors. Tumor 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. The full length of FasL contains 281 amino acids, consisting of the transmembrane domain, extracellular domain, and the cytoplasmic region. A metalloproteinase cleaves the membrane-bound FasL to produce soluble Fas (sFasL), which exists as a homotrimer. The sFasL shows lower cytotoxicity than membrane bound FasL.

Cellular localization

Cell membrane: single pass type II membrane protein. Secreted protein: may be released into the extracellular fluid, probably by cleavage from the cell surface.

Applications

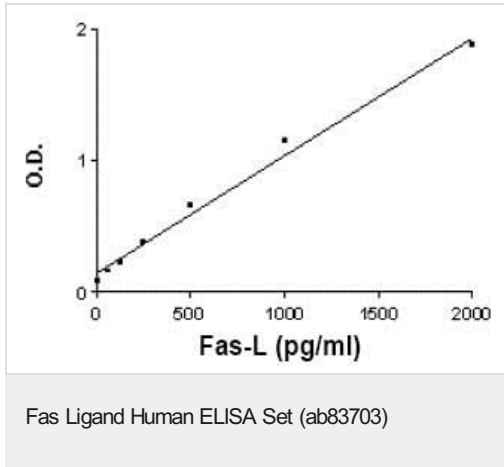
The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab83703 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



Fas Ligand standard curve ranging from 62.5 to 2000 pg/mL

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