

Human Fas Ligand ELISA Set (without plates) ab83703

画像数 1

製品の概要

製品名	Human Fas Ligand ELISA Set (without plates)
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 44 pg/ml
検出範囲	62.5 pg/ml - 2000 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<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>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	5 x 96 tests	10 x 96 tests
Detection biotinylated anti-Fas Ligand antibody		1 x 0.55ml	2 x 0.55ml

内容	ラベル	5 x 96 tests	10 x 96 tests
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

関連性

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.

細胞内局在

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

アプリケーション

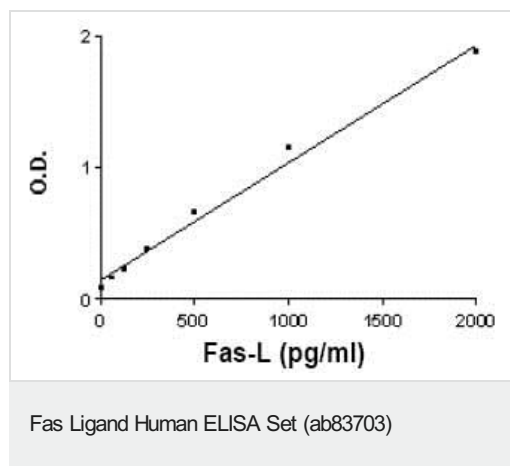
The Abpromise guarantee

Abpromise保証は、次のテスト済みアプリケーションにおけるab83703の使用に適用されます

アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご確認ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



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.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