

Human DR4 ELISA Kit (CD261) ab46510

画像数 1

製品の概要

製品名	Human DR4 ELISA Kit (CD261)				
検出方法	Colorimetric				
再現性	Intra-Assay (同時再現性)				
	サンプル	N	平均値	SD	CV%
	A	8	475pg/ml	13	2.76%
	B	8	54pg/ml	4	7.57%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	A	12	498pg/ml	40	8.11%
	B	12	53pg/ml	7	13.87%
サンプルの種類	Cell culture supernatant, Serum, Cell Lysate				
アッセイタイプ	Sandwich (quantitative)				
検出感度	< 8 pg/ml				
検出範囲	31.25 pg/ml - 1000 pg/ml				
添加回収試験	特定サンプルでの回収試験				
	サンプルの種類	平均 %	測定範囲		
	Serum	93.4	250pg/ml - 1000pg/ml		
全工程の試験時間	2h 45m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's DR4 (CD261) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of DR4 in serum, plasma, buffered solutions and cell culture media.				

A monoclonal antibody specific for DR4 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known DR4 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 DR4 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 DR4 present in the samples.

This kit will recognize both endogenous and recombinant Human DR4.

特記事項

The standard is affinity purified.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

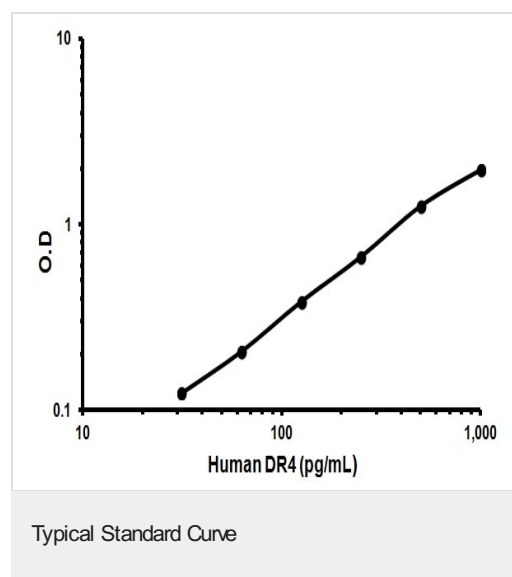
内容	ラベル	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Biotinylated anti-DR4	Red	2 x 400µl	1 x 400µl
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
DR4 Microplate (12 x 8 well strips)		2 units	1 unit
DR4 Standard (lyophilized)	Yellow	4 vials	2 vials
HRP Diluent	Red	1 x 23ml	1 x 12ml
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

関連性

Human Trail R1, CD261 or DR4 (Death Receptor 4) is a TNF receptor, which is a receptor for Trail (APO2 ligand). DR4 is involved in apoptosis. In the Trail receptor family, DR4, but also Trail R2 (DR5) transduce an apoptosis signal, whereas Trail R3 (DcR1) and Trail R4 (DcR2) antagonize TRAIL-induced apoptosis. Binding of trimeric Trail to DR4 induces apoptosis by oligomerization (likely to be trimerization) of the receptor. DR4 cDNA encodes a 468 amino acid residue precursor protein containing extracellular cysteine-rich domains, a transmembrane domain and a cytoplasmic death domain. DR4 shares 55% of its amino acid sequence with Trail R2. DR4 is widely expressed in normal tissue.

細胞内局在

Cell Membrane; single-pass type I membrane protein



Representative Standard Curve using ab46510

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