

Human DR4 ELISA Kit (CD261) ab46510

1 图像

概述

产品名称 人DR4 ELISA试剂盒(CD261)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
A	8	475pg/ml	13	2.76%
B	8	54pg/ml	4	7.57%

批次间

样品	n	Mean	SD	CV%
A	12	498pg/ml	40	8.11%
B	12	53pg/ml	7	13.87%

样品类型 Cell culture supernatant, Serum, Cell Lysate

检测类型 Sandwich (quantitative)

灵敏度 < 8 pg/ml

范围 31.25 pg/ml - 1000 pg/ml

回收率 特定样本回收率

样品类型	平均%	范围
Serum	93.4	250pg/ml - 1000pg/ml

检测时间 2h 45m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's DR4 (CD261) Human *in vitro* 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.

说明 The standard is affinity purified.

平台 Microplate

性能

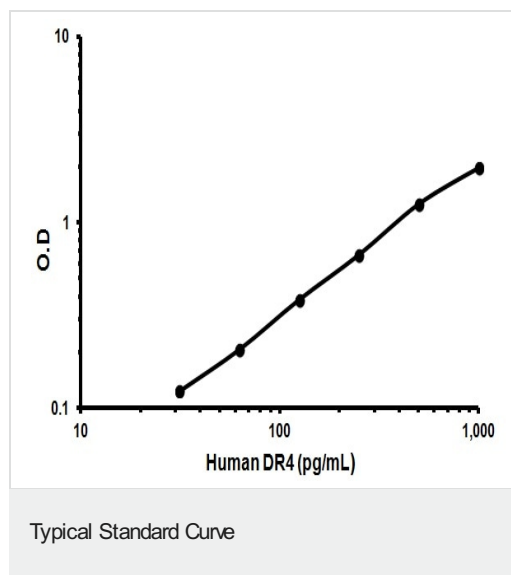
存放说明 Store at +4°C. Please refer to protocols.

组件	标识符	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

相关性 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.

细胞定位 Cell Membrane; single-pass type I membrane protein

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors