



Product datasheet

Human DcR1 ELISA Kit (CD263) αb46016

1 References 1 Image

Overview

Product name Human DcR1 ELISA Kit (CD263)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Sample A	8	7171pg/ml	192	2.7%
Sample B	8	3584pg/ml	244	6.8%

Inter-assay

Sample	n	Mean	SD	CV%
Sample A	9	7311pg/ml	515	7%
Sample B	9	2654pg/ml	356	9.7%

Sample type Serum, Cell Lysate

Assay type Sandwich (quantitative)

Sensitivity < 147 pg/ml

Range 312.5 pg/ml - 10000 pg/ml

Assay time 3h 45m

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview Abcam’s DcR1 (CD263) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of DcR1 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

A monoclonal antibody specific for DcR1 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known DcR1 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 DcR1 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 DcR1 present in the samples.

This kit will recognize both endogenous and recombinant Human DcR1.

#### Platform

Microplate

#### Properties

#### Storage instructions

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

Components	Identifier	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Biotinylated anti-DcR1	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
DcR1 Microplate (12 x 8 well strips)		1 unit	2 units
DcR1 Standard (Lyophilized)	Yellow	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

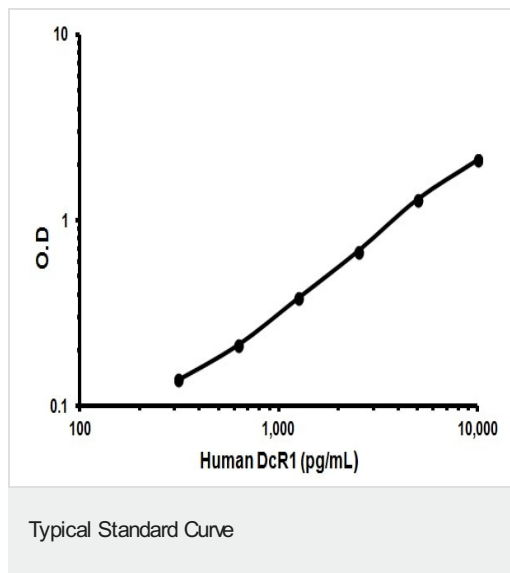
#### Relevance

DcR1 is a member of the TNF-receptor superfamily. It contains an extracellular TRAIL-binding domain and a transmembrane domain, but no cytoplasmic death domain. It is not capable of inducing apoptosis, and is thought to function as an antagonistic receptor that protects cells from TRAIL-induced apoptosis. This gene was found to be a p53-regulated DNA damage-inducible gene. The expression of this gene was detected in many normal tissues but not in most cancer cell lines, which may explain the specific sensitivity of cancer cells to the apoptosis-inducing activity of TRAIL.

#### Cellular localization

Cell membrane; lipid-anchor; GPI-anchor

#### Images



Representative Standard Curve using ab46016

**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