

Human IL-2 Receptor ELISA Kit ab46036

[2 References](#) [1 Image](#)

Overview

Product name Human IL-2 Receptor ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				4.5%

Inter-assay

Sample	n	Mean	SD	CV%
1	6	1594pg/ml	184	11.6%
2	6	1528pg/ml	135	8.8%
3	6	1705pg/ml	204	12%
4	6	1124pg/ml	171	15.2%
5	6	2305pg/ml	165	7.2%
6	6	1824pg/ml	126	69%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity = 32.5 pg/ml

Range 68.75 pg/ml - 2200 pg/ml

Recovery 87 %

Sample specific recovery

Sample type	Average %	Range
Serum	99	86% - 122%

Assay time 3h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's IL-2 Receptor Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-2 Receptor in supernatants, buffered solutions or serum and plasma samples. This assay will recognise both natural and recombinant Human IL-2 Receptor.

A monoclonal antibody specific for IL-2 Receptor has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-2 Receptor 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 IL-2 Receptor 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 IL-2 Receptor present in the samples.

This kit will recognize both endogenous and recombinant Human IL-2 Receptor.

Platform Microplate

Properties

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

Components	Identifier	2 x 96 tests	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml	1 x 25ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml	1 x 13ml
Biotinylated anti-IL-2 Receptor	Red	2 x 0.4ml	1 x 0.4ml	2 x 0.4ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 23ml	1 x 12ml	1 x 23ml
IL-2 Receptor Control	Silver	4 vials	2 vials	4 vials
IL-2 Receptor Microplate (12 x 8 well strips)		2 units	1 unit	2 units
Plastic Plate Covers		4 units	2 units	4 units
Stop Reagent	Black	2 x 11ml	1 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl	4 x 5µl

Function Receptor for interleukin-2.

Involvement in disease Genetic variations in IL2RA are associated with susceptibility to diabetes mellitus insulin-

dependent type 10 (IDDM10) [MIM:601942]. A multifactorial disorder of glucose homeostasis that is characterized by susceptibility to ketoacidosis in the absence of insulin therapy. Clinical features are polydipsia, polyphagia and polyuria which result from hyperglycemia-induced osmotic diuresis and secondary thirst. These derangements result in long-term complications that affect the eyes, kidneys, nerves, and blood vessels.

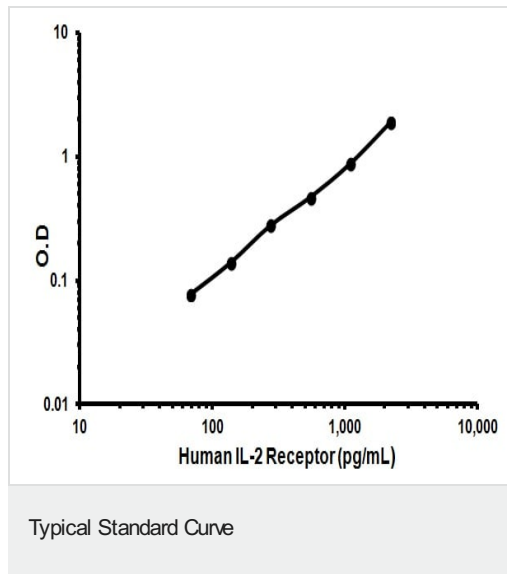
Sequence similarities

Contains 2 Sushi (CCP/SCR) domains.

Cellular localization

Membrane.

Images



Representative Standard Curve using ab46036

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