

Rat IL-4 ELISA Kit ab46073

1 Image

Overview

Product name	Rat IL-4 ELISA Kit																																	
Detection method	Colorimetric																																	
Precision	<div>Intra-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>A</td><td>18</td><td>379.3pg/ml</td><td>13.5</td><td>2.5%</td></tr><tr><td>B</td><td>19</td><td>532.5pg/ml</td><td>5.99</td><td>1.6%</td></tr></table> <div>Inter-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>A</td><td>6</td><td>128.6pg/ml</td><td>7.6</td><td>5.9%</td></tr><tr><td>B</td><td>6</td><td>532.2pg/ml</td><td>20.6</td><td>3.8%</td></tr></table>				Sample	n	Mean	SD	CV%	A	18	379.3pg/ml	13.5	2.5%	B	19	532.5pg/ml	5.99	1.6%	Sample	n	Mean	SD	CV%	A	6	128.6pg/ml	7.6	5.9%	B	6	532.2pg/ml	20.6	3.8%
Sample	n	Mean	SD	CV%																														
A	18	379.3pg/ml	13.5	2.5%																														
B	19	532.5pg/ml	5.99	1.6%																														
Sample	n	Mean	SD	CV%																														
A	6	128.6pg/ml	7.6	5.9%																														
B	6	532.2pg/ml	20.6	3.8%																														
Sample type	Cell culture supernatant, Serum																																	
Assay type	Sandwich (quantitative)																																	
Sensitivity	< 15 pg/ml																																	
Range	31.25 pg/ml - 1000 pg/ml																																	
Assay time	4h 45m																																	
Assay duration	Multiple steps standard assay																																	
Species reactivity	Reacts with: Rat																																	
Product overview	Abcam's IL-4 Rat <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-4 in cell culture supernatants, buffered solutions and plasma.																																	

A monoclonal antibody specific for IL-4 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-4 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-4 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-4 present in the samples.

This kit will recognize both endogenous and recombinant rat IL-4.

Platform Microplate

Properties

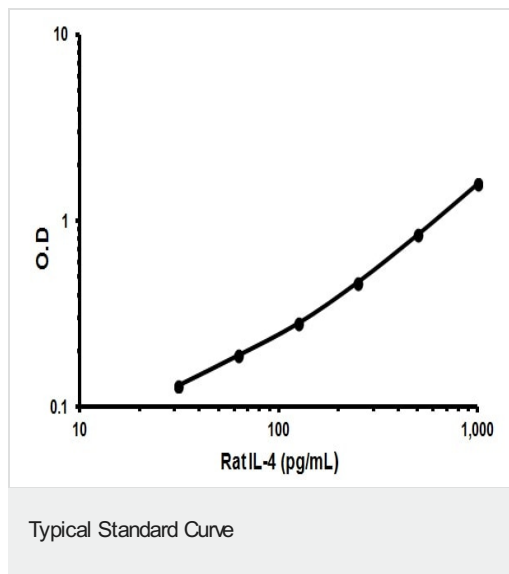
Storage instructions Store at +4°C. Please refer to protocols.

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 7.5ml	1 x 13ml
Biotinylated anti-IL-4	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-4 Microplate (12 x 8 well strips)		1 unit	2 units
Rat IL-4 Standard (Lyophilized)	Yellow	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

Relevance IL-4 is a pleiotropic cytokine produced by activated T cells. It is a ligand for interleukin 4 receptor. The interleukin 4 receptor also binds to IL13, which may contribute to many overlapping functions of this cytokine and IL13. STAT6, a signal transducer and activator of transcription, has been shown to play a central role in mediating the immune regulatory signal of IL-4. IL-4, IL3, IL5, IL13, and CSF2 form a cytokine gene cluster on chromosome 5q, with this gene particularly close to IL13. IL-4, IL13 and IL5 are found to be regulated coordinately by several long-range regulatory elements in an over 120 kilobase range on the chromosome. Two alternatively spliced transcript variants of IL-4 encoding distinct isoforms have been reported.

Cellular localization Secreted

Images



Representative Standard Curve using ab46073

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