

Mouse IL-2 ELISA Kit ab46096

3 References 1 Image

Overview

Product name	Mouse IL-2 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>10</td><td>170pg/ml</td><td>13</td><td>7.8%</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>B</td><td>8</td><td>137pg/ml</td><td>8.9</td><td>6.4%</td></tr></table>				Sample	n	Mean	SD	CV%	A	10	170pg/ml	13	7.8%	Sample	n	Mean	SD	CV%	B	8	137pg/ml	8.9	6.4%
Sample	n	Mean	SD	CV%																				
A	10	170pg/ml	13	7.8%																				
Sample	n	Mean	SD	CV%																				
B	8	137pg/ml	8.9	6.4%																				
Sample type	Cell culture supernatant, Serum																							
Assay type	Sandwich (quantitative)																							
Sensitivity	< 3.4 pg/ml																							
Range	15.6 pg/ml - 500 pg/ml																							
Assay time	3h 45m																							
Assay duration	Multiple steps standard assay																							
Species reactivity	Reacts with: Mouse																							
Product overview	<p>Abcam's IL-2 mouse <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-2 in mouse serum, buffered solutions and cell culture medium.</p> <p>A monoclonal antibody specific for IL-2 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-2 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 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 present in the samples.</p> <p>This kit will recognize both endogenous and recombinant mouse IL-2.</p>																							

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

Function

Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

Involvement in disease

Note=A chromosomal aberration involving IL2 is found in a form of T-cell acute lymphoblastic leukemia (T-ALL). Translocation t(4;16)(q26;p13) with involves TNFRSF17.

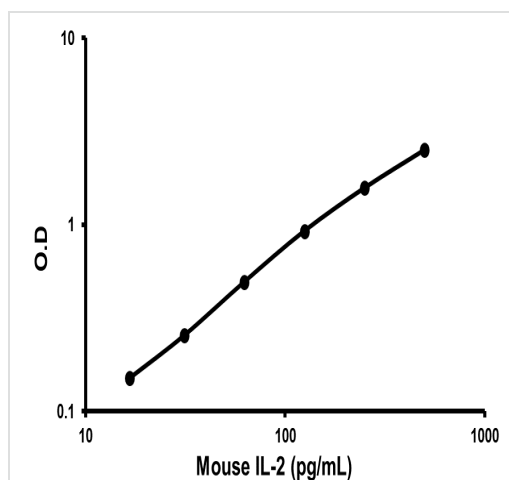
Sequence similarities

Belongs to the IL-2 family.

Cellular localization

Secreted.

Images



Representative Standard Curve using ab46096.

Typical Standard Curve

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