

IL-6 (mouse), ELISA kit

Catalog #: ADI-900-045

96 Well Kit

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

Description	2
Introduction	2
Precautions	3
Materials Supplied	4
Storage	5
Other Materials Needed	5
Sample Handling	5
Procedural Notes	6
Reagent Preparation	7
Assay Procedure	8
Calculation of Results	9
Typical Results	10
Typical Standard Curve	10
Performance Characteristics	11
Sample Dilution Recommendations	12
References	13
Contact Information	14

DESCRIPTION

The IL-6 (mouse), ELISA kit is a complete kit for the quantitative determination of mouse IL-6 in biological fluids. Please read the complete kit insert before performing this assay. The kit uses a monoclonal antibody to mouse IL-6 immobilized on a microtiter plate to bind the mouse IL-6 in the standards or sample. A recombinant mouse IL-6 Standard is provided in the kit. After a short incubation the excess sample or standard is washed out and a biotinylated monoclonal antibody to mouse IL-6 is added. This antibody binds to the mouse IL-6 captured on the plate. After a short incubation the excess antibody is washed out and Streptavidin conjugated to Horseradish peroxidase is added, which binds to the biotinylated mouse IL-6 antibody. Excess conjugate is washed out and substrate is added. After a short incubation, the enzyme reaction is stopped and the color generated is read at 450nm. The measured optical density is directly proportional to the concentration of mouse IL-6 in either standards or samples. For further explanation of the principles and practices of immunoassays please see the excellent books by Chard¹ or Tijssen².

INTRODUCTION

Interleukin-6 (IL-6) is a cytokine critical to the regulation of immune and hematopoietic systems. It has been called interferon β_2 , hybridoma/plasmacytoma growth factor, B-cell stimulatory factor 2, 26kDa inducible protein, hepatocyte stimulating factor, hematopoietic colony stimulating factor, monocyte granulocyte inducer type 2 and cytotoxic T-cell differentiation factor^{3-5,10}. A 211-212 amino acid protein, IL-6 is expressed by cell types such as T-cells, mast cells, monocytes, macrophages, fibroblasts, endothelial cells, keratinocytes, and many tumor cell lines. It appears to take part in acute phase reactions and response to injury and inflammation^{3-5,9}. IL-6 also stimulates the differentiation of B-cells for antibody production, promotes expansion of activated T-cells, expands hematopoietic cell production, and induces the expression of acute phase proteins⁵⁻¹⁰. Rodent IL-6 is 42-58% homologous with human IL-6 at the amino acid level. Rodent IL-6 is not N-glycosylated.

SAFETY WARNINGS & PRECAUTIONS



Handle
with care



Avoid
freeze /
thaw cycles

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

- Stop Solution is a 1 normal (1N) hydrochloric acid solution. This solution is caustic; care should be taken in use.
- The activity of the Horseradish peroxidase conjugate is affected by nucleophiles such as azide, cyanide and hydroxylamine.
- We test this kit's performance with a variety of samples, however it is possible that high levels of interfering substances may cause variation in assay results.
- The mouse IL-6 Standard provided, Catalog No. 80-0656, should be handled with care because of the known and unknown effects of mouse IL-6.
- The mouse IL-6 Standard should be stored at or below -20°C. Do not repeatedly freeze-thaw.

MATERIALS SUPPLIED

- 1. mouse IL-6 Microtiter Plate, One Plate of 96 Wells, Catalog No. 80-0590**
A plate using break-apart strips coated with monoclonal antibody specific to mouse IL-6.
- 2. mouse IL-6 ELISA Antibody, 5ml, Catalog No. 80-0441**
A yellow solution of biotinylated rat monoclonal antibody to mouse IL-6.
- 3. Assay Buffer 13, 55ml, Catalog No. 80-1500**
Tris buffered saline containing proteins and detergents.
- 4. mouse IL-6 ELISA Conjugate 5ml, Catalog No. 80-1626**
A blue solution of Streptavidin conjugated to Horseradish peroxidase.
- 5. Wash Buffer Concentrate, 100ml, Catalog No. 80-1287**
Tris buffered saline containing detergents.
- 6. mouse IL-6 Standard, each, Catalog No. 80-0656**
Two vials containing 2,000pg each of lyophilized recombinant mouse IL-6. Avoid repeated freeze/thaw cycles.
- 7. TMB Substrate, 5ml, Catalog No. 80-0615**
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide. Ready to use. Protect from prolonged exposure to light.
- 8. Stop Solution 2, 10ml, Catalog No. 80-0377**
A 1N solution of hydrochloric acid in water. Keep tightly capped. Caution: Caustic.
- 9. mouse IL-6 Assay Layout Sheet, 1 each, Catalog No. 30-0077**
- 10. Plate Sealer, 3 each, Catalog No. 30-0012**



Reagents
require
separate
storage
conditions.

STORAGE

All components of this kit, except the Standard, are stable at 4°C until the kit's expiration date. The Standard must be stored at or below -20°C.

OTHER MATERIALS NEEDED

1. Deionized or distilled water.
2. Precision pipets for volumes between 100µl and 1,000µl.
3. Disposable test tubes for dilution of samples and standards.
4. Repeater pipets for dispensing 100µl.
5. Disposable beakers for diluting buffer concentrates.
6. Graduated cylinders.
7. A 37°C Incubator.
8. Adsorbent paper for blotting.
9. Microplate reader capable of reading at 450nm, preferably with correction between 570nm and 590nm.
10. Graph paper for plotting the standard curve.

SAMPLE HANDLING

The IL-6 (mouse), ELISA kit is compatible with mouse IL-6 samples in a wide range of matrices. Samples diluted sufficiently into the proper diluent can be read directly from a standard curve. Please refer to the Sample Recovery recommendations on page 12 for details of suggested dilutions. Culture fluids and serum are suitable for use in the assay. Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic specimens. Samples in the majority of tissue culture media, including those containing fetal bovine serum, can also be read in the assay, provided the standards have been diluted into the tissue culture media instead of Assay Buffer 13. There will be a small change in binding associated with running the standards and samples in media. Users should only use standard curves generated in media or buffer to calculate concentrations of mouse IL-6 in the appropriate matrix.

Samples must be stored frozen to avoid loss of bioactive mouse IL-6. If samples are to be run within 24 hours, they may be stored at 4°C. Otherwise, samples must be stored frozen at -70°C to avoid loss of bioactive mouse IL-6. Excessive freeze/thaw cycles should be avoided. Prior to assay, frozen sera should be brought to room temperature slowly and gently mixed by hand. Do not thaw samples in a 37°C incubator. Do not vortex or sharply agitate samples.

PROCEDURAL NOTES

1. Do not mix components from different kit lots or use reagents beyond the kit expiration date.
2. Allow all reagents to warm to room temperature for at least 30 minutes before opening.
3. Standards can be made up in either glass or plastic tubes.
4. Pre-rinse the pipet tip with reagent, use fresh pipet tips for each sample, standard and reagent.
5. Pipet standards and samples to the bottom of the wells.
6. Add the reagents to the side of the well to avoid contamination.
7. This kit uses break-apart microtiter strips, which allow the user to measure as many samples as desired. Unused wells must be kept desiccated at 4°C in the sealed bag provided. The wells should be used in the frame provided.
8. Prior to addition of substrate, ensure that there is no residual wash buffer in the wells. Any remaining wash buffer may cause variation in assay results.
9. It is important that the matrix for the standards and samples be as similar as possible. Mouse IL-6 samples diluted with Assay Buffer 13 should be run with a standard curve diluted in the same buffer. Serum and plasma samples should be evaluated against a standard curve run in Assay Buffer 13, while tissue culture samples should be read against a standard curve diluted in the same complete but non-conditioned media. See Reagent Preparation, step #2.

REAGENT PREPARATION

1. Wash Buffer

Prepare the Wash Buffer by diluting 50ml of the supplied concentrate with 950ml of deionized water. This can be stored at room temperature until the kit expiration, or for 3 months, whichever is earlier.

2. mouse IL-6 Standards

Allow the 2,000pg mouse IL-6 standard vial to warm to room temperature. Reconstitute one vial of mouse IL-6 Standard with 2mls standard diluent (Assay Buffer 13 or Tissue Culture Media). Label seven 12 x 75mm glass tubes #2 through #8. Label vial of reconstituted mouse IL-6 Std #1. Pipet 500 μ l of standard diluent into tubes #2 through #8. Remove 500 μ l from vial #1 and add to tube #2. Vortex thoroughly. Remove 500 μ l from tube #2 and add to tube #3 and vortex. Continue this for tubes #4 through #8. The concentration of mouse IL-6 in tubes #1 through #8 will be 1,000, 500, 250, 125, 62.5, 31.25, 15.63 and 7.81pg/ml respectively. See mouse IL-6 Assay Layout Sheet for dilution details. **STORE STANDARD AT -20°C**, avoid repeated freeze-thaws.

ASSAY PROCEDURE

Bring all reagents to room temperature for at least 30 minutes prior to opening. Plates will require shaking on an orbital rotor at 500rpm. All standards, controls and samples should be run in duplicate.

1. Refer to the Assay Layout Sheet to determine the number of wells to be used and put any remaining wells with the desiccant back into the pouch and seal the ziploc. Store unused wells at 4°C.
2. Pipet 50µl of standard diluent (Assay Buffer 13 or tissue culture media) into the S0 (0 pg/ml standard) wells.
3. Pipet 50µl of Standards #1 through #8 into the appropriate wells.
4. Pipet 50µl of the Samples into the appropriate wells.
5. Tap the plate gently to mix the contents.
6. Seal the plate and incubate at room temperature on a plate shaker for 1 hour at ~500rpm*.
7. Empty the contents of the wells and wash by adding 400µl of wash solution to every well. Repeat the wash 3 more times for a total of 4 washes. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
8. Pipet 50µl of yellow Antibody into each well, except the Blank.
9. Seal the plate and incubate at room temperature on a plate shaker for 1 hour at ~500rpm*.
10. Empty the contents of the wells and wash by adding 400µl of wash solution to every well. Repeat the wash 3 more times for a total of 4 washes. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
11. Add 50µl of blue Conjugate to each well, except the Blank.
12. Seal the plate and incubate at room temperature on a plate shaker for 30 minutes at ~500rpm*.
13. Empty the contents of the wells and wash by adding 400µl of wash solution to every well. Repeat the wash 3 more times for a total of 4 washes. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
14. Pipet 50µl of Substrate Solution into each well.

15. Incubate for 15 minutes at room temperature on a plate shaker for 15 minutes at ~500rpm*.
16. Pipet 50µl Stop Solution 2 to each well. This stops the reaction and the plates should be read immediately.
17. Blank the plate reader against the Blank wells, read the optical density at 450nm, preferably with correction between 570 and 590nm. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the Blank wells from all the readings.

* The plate shaker speed was based on a BellCo Mini Orbital Shaker (mod no. 7744-08096). The actual speed of the plate shaker should be such that the liquid in the plate wells mixes thoroughly, but does not splash out of the well.

CALCULATION OF RESULTS

Several options are available for the calculation of the concentration of mouse IL-6 in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic curve fitting program. If data reduction software is not readily available, the concentration of mouse IL-6 can be calculated as follows:

1. Calculate the average net Optical Density (OD) bound for each standard and sample by subtracting the average Blank OD from the average OD for each standard and sample.

Average Net OD = Average OD - Average Blank OD

2. Using linear graph paper, plot the Average Net OD for each standard versus mouse IL-6 concentration in each standard. Approximate a straight line through the points. The concentration of mouse IL-6 in the unknowns can be determined by interpolation.

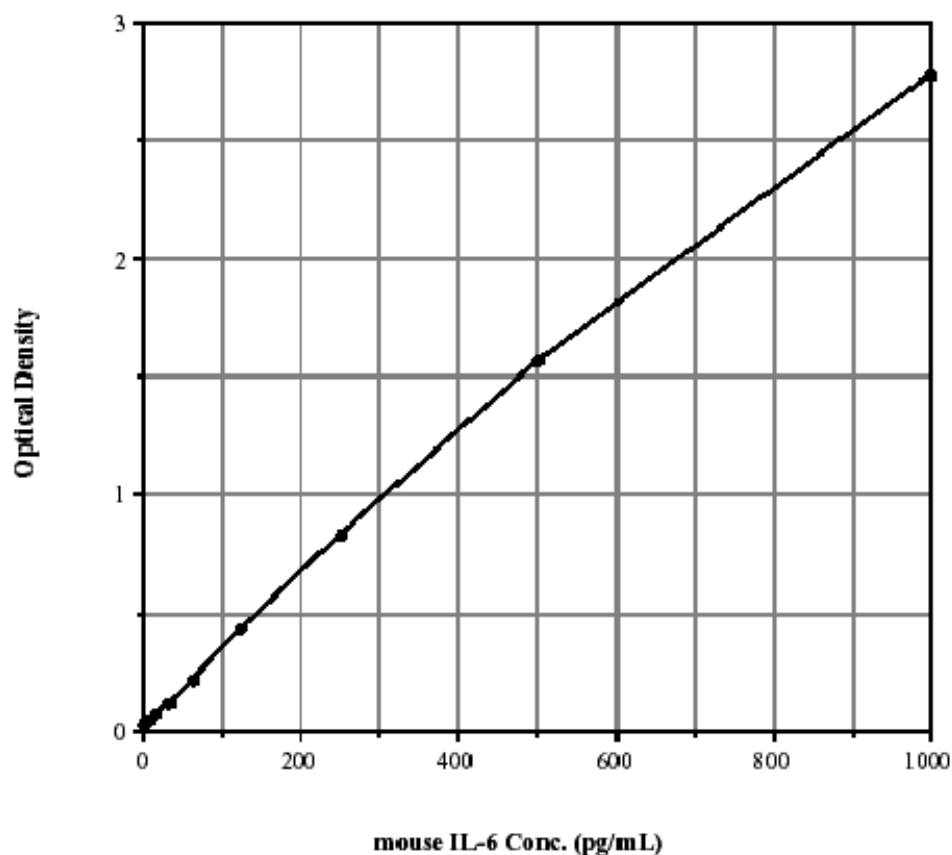
TYPICAL RESULTS

The results shown below are for illustration only and should not be used to calculate results from another assay.

Sample	Average OD	Net OD	mouse IL-6 (pg/mL)
Blank	0.041		
S0	0.067	0.0	
S1	2.815	2.774	1,000
S2	1.605	1.564	500
S3	0.868	0.827	250
S4	0.473	0.432	125
S5	0.262	0.221	62.5
S6	0.157	0.116	31.25
S7	0.109	0.068	15.63
S8	0.088	0.047	7.81
Unknown #1	2.382	2.341	821.1
Unknown #2	1.346	1.305	412.1

TYPICAL STANDARD CURVES

A typical standard curve is shown below. This curve must not be used to calculate mouse IL-6 concentrations; each user must run a standard curve for each assay.



PERFORMANCE CHARACTERISTICS

The following parameters for this kit were determined using the guidelines listed in the National Committee for Clinical Laboratory Standards (NCCLS) Evaluation Protocols¹¹

Sensitivity

Sensitivity was calculated by determining the average optical density bound for sixteen (16) wells run with 0 pg/ml Standard, and comparing to the average optical density for sixteen (16) wells run with Standard #8. The detection limit was determined as the concentration of mouse IL-6 measured at two (2) standard deviations from the 0 pg/ml Standard along the standard curve.

Mean OD for S0 = 0.029 ± 0.002 (7.4%)

Mean OD for Standard #8 = 0.060 ± 0.005 (8.0%)

Delta Optical Density = $(7.81 - 0 \text{ pg/ml}) = 0.060 - 0.029 = 0.031$

2 SD's of 0 pg/ml Standard = $2 \times 0.002 = 0.004$

Sensitivity = $\frac{0.004}{0.031} \times 7.81 \text{ pg/ml} = 1.01 \text{ pg/ml}$

Linearity

A sample containing 821.07 pg/ml mouse IL-6 was serially diluted 7 times 1:2 in the Assay Buffer 13 supplied in the kit and measured in the assay. The data was plotted graphically as actual mouse IL-6 concentration versus measured mouse IL-6 concentration.

The line obtained had a slope of 1.0216 with a correlation coefficient of 0.9999.

Precision

Intra-assay precision was determined by taking samples containing low, medium and high concentrations of mouse IL-6 and running these samples multiple times (n=16) in the same assay. Inter-assay precision was determined by measuring three samples with low, medium and high concentrations of mouse IL-6 in multiple assays (n=6). The precision numbers listed below represent the percent coefficient of variation for the concentrations of mouse IL-6 determined in these assays as calculated by a 4 parameter logistic curve fitting program.

	mouse IL-6 (pg/ml)	Intra-assay %CV	Inter-assay %CV
Low	207.7	8.9	
Medium	340.5	11.2	
High	505.9	10.0	
Low	249.0		10.3
Medium	414.6		8.2
High	651.4		7.1

Cross Reactivities

The IL-6 (mouse), ELISA kit is specific for bioactive mouse IL-6. It is unaffected by the presence of the following recombinant molecules: mouse IL-1 α , mouse IL-1 β , mouse IL-2, mouse IL-3, mouse IL-4, mouse IL-5, mouse IL-7, mouse IL-10, mouse IFN- γ , mouse GM-CSF, mouse TNF- α and human IL-6.

SAMPLE RECOVERIES

Please refer to pages 5-7 for Sample Handling recommendations and Standard preparation. Mouse IL-6 concentrations were measured in mouse serum and tissue culture media. Mouse IL-6 was spiked into the undiluted samples of these matrices which were then diluted with the appropriate diluent and assayed in the kit. The following results were obtained:

Sample	% Recovery*	Recommended Dilution*
mouse Serum	91.6	$\geq 1:8$
Tissue Culture Media	93.0	None

* See Sample Handling instructions on page 5 for details.

REFERENCES

1. T. Chard, in "An Intro. to Radioimmunoassay & Related Tech.", (1990), 4th Ed., Elsevier, Amsterdam.
2. P. Tijssen, in "Practice & Theory of Enz. Immunoassays", (1985), Elsevier, Amsterdam.
3. P. Carmeliet, Interleukin-6: a putative mediator of anterior pituitary hormone secretion during the stress of infection. *Trends in Endocrinology and Metabolism* 237 May/June (1990).
4. T. Hirano et al., Biological and clinical aspects of interleukin-6. *Immunol. Today* 11:443 (1990).
5. Kishimoto, The biology of Interleukin-6. *Blood* 74:1 (1989).
6. T. Kishimoto and T. Hirano, Molecular regulation of β lymphocyte response. *Annu. Rev. Immunol.* 6:485 (1988).
7. Y. Takai et al., β cell stimulatory factor-2 is involved in the differentiation of cytotoxic T lymphocytes. *J.Immunol.* 140:508 (1988).
8. G. Tosato et al., Monocyte-derived human B-cell growth factor identified as interferon-B2 (BSF-2, IL-6). *Science* 239:502 (1988).
9. J. Van Snick, Interleukin-6: an overview. *Annu. Rev. Immunol.* 8:253 (1990).
10. Wong, G. and Clark, S. Multiple actions of Interleukin-6 within a cytokine network. *Immunol. Today* 9:137 (1988).
11. National Committee for Clinical Laboratory Standards Evaluation Protocols, SC1, 1989, NCCLS, Villanova, PA, 19085



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com