



Human IL-6 Conferma™ ELISA

96-Well Plate Assay

**Cat. # EZIL6-98K
(EZIL6-98K5PK,
EZIL6-98K10PK)**

HUMAN IL-6 CONFERMA™ ELISA KIT
96-Well Plate

Cat. # EZIL6-98K
(EZIL6-98K5PK, EZIL6-98K10PK)

<u>TABLE OF CONTENTS</u>	<u>PAGE</u>
Intended Use	2
Principles Of Assay	2
Reagents Supplied	3
Storage and Stability	3
Reagent Precautions	4
Materials Required But Not Provided	6
Sample Collection And Storage	7
Reagent Preparation	8
Assay Procedure	9
Microtiter Plate Arrangement	12
Graph of Typical Reference Curve	13
Assay Characteristics	13
Quality Controls	17
Troubleshooting Guide	18
Ordering Information	18

Millipore Sigma Corporation • 3050 Spruce Street • St. Louis, Missouri 63103 USA
Toll Free U.S: 1-800 645-5476 • Fax: 1-800-645-5439
MilliporeSigma.com

The M logo is a trademark of Merck KGaA, Darmstadt, Germany.
© 2013 EMD Millipore Corporation, Billerica, MA 01821 USA.

EZIL6-98K

INTENDED USE

This Human IL-6 Conferma™ ELISA kit is used for the non-radioactive quantification of Human IL-6 in serum and plasma samples. One kit is sufficient to measure 36 unknown samples in duplicate. ***This kit is for Research Use Only. Not for Use in Diagnostic Procedures.***

PRINCIPLES OF ASSAY

This assay is a Sandwich ELISA which uses in-house developed critical reagents, including the Monoclonal Antibodies (mAb) and calibration material, to detect endogenous IL-6 in biological fluids such as Human serum or plasma.

The Sandwich ELISA first binds IL-6 using a specific capture Mouse anti-human IL-6 monoclonal antibody bound to the wells of a 96 well microtitre plate. Following the addition of the sample, the assay is incubated for two hours, during which time endogenous or recombinant antigen (depending on the well) is bound by the mAb. The unbound material is washed off post-incubation, and a biotinylated mouse anti-human IL-6 monoclonal antibody is added to complete the "Sandwich". After an incubation period, the unbound material is washed off. The next step is a final incubation step during which a streptavidin-horseradish peroxidase conjugate binds to the immobilized biotinylated antibodies, and following a final wash horseradish peroxidase substrate, 3,3',5,5'-tetramethylbenzidine is added. The enzyme activity is measured spectrophotometrically by the increased absorbance at 450-590 nm after acidification of formed products by the addition of Stop Solution. The increase in absorbance is directly proportional to the amount of captured Human IL-6. Quantitation of the analyte is derived by interpolation from a reference curve comprised of standard points of known concentrations of recombinant human IL-6.

REAGENTS SUPPLIED

Each kit is sufficient to run one 96-well plate and contains the following reagents:

Note: Store all reagents at 2-8°C

Reagents Supplied	Catalog Number	Volume	Quantity
Human IL-6 ELISA plate with 2 sealers	EP98	-----	1 plate 2 sealers
Human IL-6 Standard	E8098-K	lyophilized	1 vial
Human IL-6 Quality Controls 1, 2 and 3	E6098-1-K E6098-2-K E6098-3-K	lyophilized	1 vial each
Serum Matrix	EMTX-98	lyophilized	1 vial
Assay Buffer	EAB098	10 mL	1 vial
10X Wash Buffer	EWB-HRP98	50 mL	2 bottles
Human IL-6 Detection Antibody	E1098	12 mL	1 bottle
Enzyme Solution (100X)	EHRP-98	150 µL	1 bottle
Enzyme Solution Diluent	ED-098	12 mL	1 bottle
Substrate Solution	ESS-TMB98	12 mL	1 bottle
Stop Solution	ET-TMB98	12 mL	1 bottle

STORAGE AND STABILITY

The recommended storage for kit components is 2-8°C.

All components are shipped and stored at 2-8°C. Reconstituted standards and controls can be frozen for future use, but repeated freeze/thaw cycles should be avoided.


10X Wash Buffer does not contain a preservative. After dilution, the 1X Wash Buffer may be filter sterilized (Stericup® filter, Millipore Sigma- Cat# SCGPU11RE) for storage of up to 1 month at 2 - 8°C. If not filter sterilized, all remaining 1X wash buffer should not be used after one week.







Refer to expiration dates on all reagents before use. Do not mix reagents from different kits unless they have the same lot numbers.

REAGENT PRECAUTIONS

Sodium Azide has been added to some reagents as a preservative. Although the concentration is low, Sodium Azide may react with lead and copper plumbing to form highly explosive metal azides. Dispose of unused contents and waste in accordance with international, federal, state, and local regulations.

FULL HAZARD LABEL:

Ingredient, Cat #		Full Label	
Human IL-6 Standard	E8098-K		<p>Danger: Harmful if swallowed or if inhaled. Toxic in contact with skin. Causes serious eye damage. May cause damage to the brain through prolonged or repeated exposure. Do not breathe dust/ fume/ gas/ mist/ vapors/ spray. Wash skin thoroughly after handling. Do not eat, drink, or smoke when using this product. Use only outdoors or in a well-ventilated area. Wear protective gloves/ eye protection/ face protection. IF SWALLOWED: Call a POISON CENTER/ doctor if you feel unwell. IF ON SKIN: Wash with plenty of soap and water. IF INHALED: Remove person to fresh air and keep comfortable for breathing. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present, and easy to do. Continue rinsing. Immediately call a POISON CENTER/ doctor. Specific measures (see supplemental first aid instructions on this label). Rinse mouth. Remove/ Take off immediately all contaminated clothing. Wash contaminated clothing before reuse. Store locked up. Dispose of contents/ container to an approved waste disposal plant.</p>

Human IL-6 Quality Control 1, 2 & 3	E6098-1-K E6098-2-K E6098-3-K	  	<p>Danger: Harmful if swallowed or if inhaled. Toxic in contact with skin. Causes serious eye damage. May cause damage to the brain through prolonged or repeated exposure. Do not breathe dust/ fume/ gas/ mist/ vapors/ spray. Wash skin thoroughly after handling. Do not eat, drink, or smoke when using this product. Use only outdoors or in a well-ventilated area. Wear protective gloves/ eye protection/ face protection. IF SWALLOWED: Call a POISON CENTER/ doctor if you feel unwell. IF ON SKIN: Wash with plenty of soap and water. IF INHALED: Remove person to fresh air and keep comfortable for breathing. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present, and easy to do. Continue rinsing. Immediately call a POISON CENTER/ doctor. Specific measures (see supplemental first aid instructions on this label). Rinse mouth. Remove/ Take off immediately all contaminated clothing. Wash contaminated clothing before reuse. Store locked up. Dispose of contents/ container to an approved waste disposal plant.</p>
Human IL-6 detection antibody	E1098		<p>Warning: Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present, and easy to do. Continue rinsing.</p>
Serum Matrix	EMTX-98	No label	Harmful to aquatic life with long lasting effects. Avoid release to the environment.
Assay Buffer	EAB098		<p>Warning: Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present, and easy to do. Continue rinsing.</p>
Stop Solution	ET-TMB98		<p>Warning: May be corrosive to metals.</p>

MATERIALS REQUIRED BUT NOT PROVIDED

1. Multi-channel Pipettes and pipette tips: 5-50 μL and 50-300 μL
2. Pipettes and pipette tips: 10 μL -20 μL or 20 μL -100 μL
3. Reagent Reservoirs
4. Polypropylene Microfuge Tubes
5. Vortex Mixer
6. De-ionized water
7. Microtiter Plate Reader capable of reading absorbency at 450 nm
8. Orbital Microtiter Plate Shaker
9. Absorbent Paper or Cloth

SAMPLE COLLECTION AND STORAGE

A. Preparation of Serum Samples:

- Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000xg. Remove serum and assay immediately or aliquot and store samples at $\leq -20^{\circ}\text{C}$.
- Avoid multiple >2 freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing, and centrifuge prior to use in the assay to remove particulates.
- Serum samples should be used neat.

B. Preparation of Plasma Samples:

- Plasma collection using EDTA as an anticoagulant is recommended. Centrifuge for 10 minutes at 1000xg within 30 minutes of blood collection. Remove plasma and assay immediately or aliquot and store samples at $\leq -20^{\circ}\text{C}$.
- Avoid multiple >2 freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing, and centrifuge prior to use in the assay to remove particulates.
- Plasma samples should be used neat.

C. Preparation of Tissue Culture Supernatant:

- Centrifuge the sample to remove debris and assay immediately or aliquot and store samples at $\leq -20^{\circ}\text{C}$.
- Avoid multiple (>2) freeze/thaw cycles.
- Tissue culture supernatant may require dilution with an appropriate control medium prior to assay. Tissue/cell extracts should be done in neutral buffers containing reagents and conditions that do not interfere with assay performance. Excess concentrations of detergent, salt, denaturants, high or low pH, etc. will negatively affect the assay. Organic solvents should be avoided. The tissue/cell extract samples should be free of particles such as cells or tissue debris.

NOTE:

- A maximum of 50 μL per well of neat serum or plasma can be used. Tissue culture or other media may also be used.
- All samples must be stored in polypropylene tubes. **DO NOT STORE SAMPLES IN GLASS.**
- Avoid debris, lipids, and cells when using samples with gross hemolysis or lipemia.
- Care must be taken when using heparin as an anticoagulant since an excess of heparin will provide falsely high values. Use no more than 10 IU heparin per mL of blood collected.

REAGENT PREPARATION

A. Human IL-6 Standard Preparation

Use care in opening the lyophilized Standard vial. Refer to the Standard reconstitution instructions provided on the Certificate of analysis to hydrate the stock standard vial to 1X concentration

- For dilution series, Label 7 polypropylene microfuge tubes as Std 7, Std 6, Std 5, Std 4, Std 3, Std 2, and Std 1. Add 200 μ L of Assay Buffer to each of the 7 tubes. Prepare serial dilutions by adding 200 μ L of the reconstituted standard to the Std 7 tube, mix well and transfer 200 μ L of the Std 7 to the Std 6 tube, mix well and transfer 200 μ L of the Std 6 to the Std 5 tube, mix well and transfer 200 μ L of the Std 5 to the Std 4 tube, mix well and transfer 200 μ L of the Std 4 to the Std 3 tube, mix well and transfer 200 μ L of the Std 3 to the Std 2 tube, mix well and transfer 200 μ L of the Std 2 to the Std 1 tube and mix well. The 0 pg/mL standard (Background) will be Assay Buffer.

Note: Change tip for every dilution. Wet tip with the standard before dispensing. Unused portions of the reconstituted standard should be stored in small aliquots at $\leq -20^{\circ}\text{C}$. Avoid multiple freeze/thaw cycles.

Tube #	Volume of Deionized Water to Add	Volume of Assay Buffer to Add	Standard Stock Concentration
Reconstituted standard	Refer to COA	Refer to COA	150 pg/mL

Tube #	Volume of Assay Buffer to Add	Volume of Standard to Add	Standard Concentration pg/mL
Standard 7	200 μ L	200 μ L of reconstituted standard	75
Standard 6	200 μ L	200 μ L of Standard 7	37.5
Standard 5	200 μ L	200 μ L of Standard 6	18.75
Standard 4	200 μ L	200 μ L of Standard 5	9.38
Standard 3	200 μ L	200 μ L of Standard 4	4.69
Standard 2	200 μ L	200 μ L of Standard 3	2.34
Standard 1	200 μ L	200 μ L of Standard 2	1.17

B. Human IL-6 Quality Control 1, 2 and 3 Preparation

Use care in opening the lyophilized Quality Control vials. Reconstitute each Human IL-6 Quality Control 1, 2, and 3 as per the instructions provided in the Certificate of Analysis. Once hydrated, controls can be stored in small aliquots at $\leq -20^{\circ}\text{C}$. Avoid further freeze/thaw cycles.

REAGENT PREPARATION (continued)

C. Preparation of Wash Buffer

Bring the 10X Wash Buffer to room temperature and mix to bring all salts into solution. Dilute 100 mL of 10X Wash Buffer (two bottles) with 900 mL of de-ionized water.

NOTE: 10X Wash Buffer does not contain a preservative. For storage of up to 1 month at 2 - 8°C, the 1X Wash Buffer may need to be filter sterilized (Stericup® filter, Millipore Sigma- Cat# SCGPU11RE)

D. Preparation of Serum Matrix

Add 1.5 mL distilled or de-ionized water to the bottle containing lyophilized Serum Matrix. Mix well. Allow at least 15 minutes for complete reconstitution. Leftover reconstituted Serum Matrix should be stored at $\leq -20^{\circ}\text{C}$ for up to one month.

E. Preparation of Enzyme Solution

Add 120 μL of 100X enzyme solution to the bottle containing 12 mL of enzyme solution diluent. Mix well. Store unused portion at 2-8°C for up to one month.

Human IL-6 Conferma™ ELISA ASSAY PROCEDURE

Warm all reagents to room temperature before setting up the assay.

1. Remove the required number of strips from the Microtiter Assay Plate. Unused strips should be resealed in the foil pouch and stored at 2-8°C. Assemble the strips in an empty plate holder. Add 300 μL diluted Wash Buffer to each well of the plate. Decant Wash Buffer and remove the residual volume by inverting the plate and tapping it smartly onto absorbent towels several times. **Do not let wells dry before proceeding to the next step.** If an automated machine is used for the assay, follow the manufacturer's instructions for all washing steps described in this protocol.
2. Add 50 μL of appropriate Matrix Solution to Blank, Standards, and Quality Control wells (refer to Microtiter Plate Arrangement section for suggested sample order placement). When assaying serum or plasma, use EMTX-98. When assaying tissue culture or other supernatants, use proper control culture medium as the matrix solution.
3. Add 50 μL Assay Buffer to each of the Blank and Sample wells.
4. Add 50 μL Standards or Controls to the appropriate wells.
5. Add 50 μL of neat sample to the appropriate wells.
6. Cover the plate with a plate sealer and incubate at room temperature for 2 hours on an orbital microtiter plate shaker set to rotate at moderate speed, about 400 to 500 rpm.

Human IL-6 Conferma™ ELISA ASSAY PROCEDURE (continued)

7. Remove plate sealer and decant reagents from the plate. Tap as before to remove residual volume in well. Wash wells 3 times with diluted Wash Buffer, 300 µL per well per wash. Decant and tap after each wash to remove residual buffer.
8. Add 100 µL Detection Antibody to each well. Re-cover plate with sealer and incubate at room temperature for 1 hour on an orbital microtiter plate shaker set to rotate at moderate speed, approximately 400 rpm.
9. Remove plate sealer and decant reagents from the plate. Tap as before to remove residual volume in well. Wash wells 3 times with diluted Wash Buffer, 300 µL per well per wash. Decant and tap after each wash to remove residual buffer.
10. Add 100 µL of 1X Enzyme Solution to each well. Cover the plate with sealer and incubate with moderate shaking at room temperature for 30 minutes on the microtiter plate shaker.
11. Remove sealer, decant reagents from the plate, and tap the plate to remove the residual volume. Wash wells 5 times with diluted Wash Buffer, 300 µL per well per wash. Decant and tap after each wash to remove residual buffer.
12. Add 100 µL of Substrate Solution to each well, cover plate with sealer, and shake on the plate shaker for approximately 15 minutes. Blue color should be formed in wells of the IL-6 standards with intensity proportional to increasing concentrations of IL-6.

Note: Please be aware that the color may develop more quickly or more slowly than the recommended incubation time, depending on the localized room temperature. Please visually monitor the color development to optimize the incubation time.

13. Remove sealer and add 100 µL Stop Solution [**CAUTION: CORROSIVE SOLUTION**] and gently shake plate by hand to ensure complete mixing of the solution in all wells. The blue color should turn to yellow after acidification. Wipe the bottom of the microtiter plate to remove any residue prior to reading on a plate reader. Read absorbance at 450 nm and 590 nm in a plate reader within 5 minutes and ensure that there are no air bubbles in any well. Record the difference of absorbance units. The absorbance of the highest IL-6 standard should be approximately 2.0 - 3.0, or not to exceed the capability of the plate reader used.

Note: When sample volumes assayed differ from 50 µL, an appropriate mathematical adjustment must be made to accommodate for the dilution factor (e.g., if 25 µL of sample is used, then calculated data must be multiplied by 2). When the sample volume assayed is less than 50 µL, compensate for the volume deficit with the matrix solution.

Assay Procedure for Human IL-6 Conferma™ ELISA Kit (Cat. # EZIL6-98K)

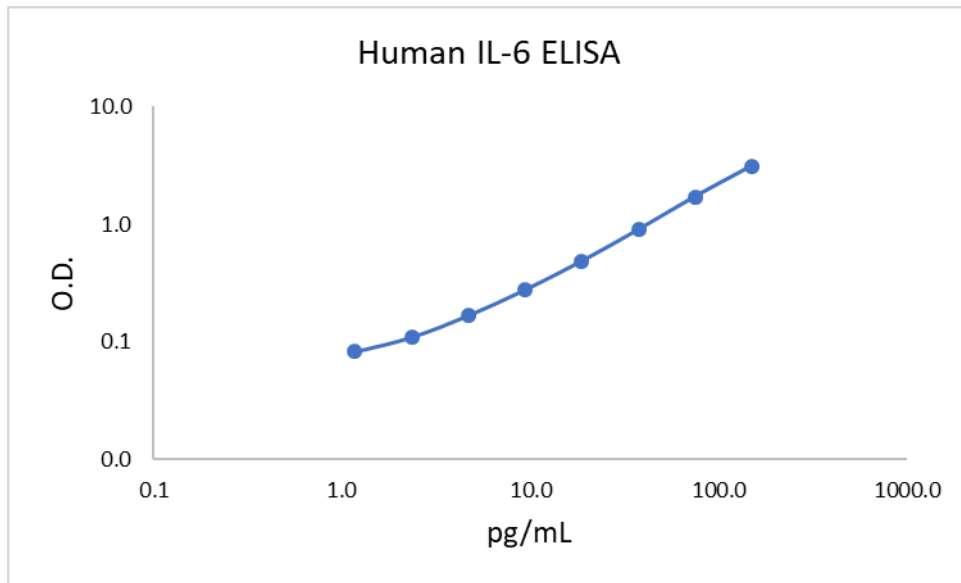
	Step 1	Step 2	Step 3	Step 4-5	Step 6-7	Step 8	Step 9	Step 10	Step 11	Step 12-13			
Well #	Wash plate 1X with 300 µL 1X Wash Buffer. Remove residual buffer by tapping smartly on absorbent towels.	Matrix Solution	Assay Buffer	Standards/ QCs/ Samples	Seal, Agitate, Incubate 2 hour at Room Temperature on a plate shaker. Wash 3X with 300 µL Wash Buffer.	Detection Antibody	Seal, Agitate, Incubate 1 hour at Room Temperature on a plate shaker. Wash 3X with 300 µL Wash Buffer.	Enzyme Solution	Seal, Agitate, Incubate 30 minutes at Room Temperature on a plate shaker. Wash 5X with 300 µL Wash Buffer.	Substrate	Seal, Agitate, Incubate 15 minutes at Room Temperature on a plate shaker.	Stop	Read Absorbance at 450 nm and 590 nm.
A1, B1		50 µL	50 µL	--		100 µL		100 µL		100 µL		100 µL	
C1, D1		50 µL	--	50 µL of Std1		↓		↓		↓		↓	
E1, F1		50 µL	--	50 µL of Std 2									
G1, H1		50 µL	--	50 µL of Std 3									
A2, B2		50 µL	--	50 µL of Std 4									
C2, D2		50 µL	--	50 µL of Std 5									
E2, F2		50 µL	--	50 µL of Std 6									
G2, H2		50 µL	--	50 µL of Std 7									
A3, B3		50 µL	--	50 µL of Reconstituted standard									
C3, D3		50 µL		50 µL of QC 1									
E3, F3		50 µL		50 µL of QC 2									
G3, H3		50 µL		50 µL of QC 3									
A4, B4 Etc.			50 µL	50 µL of sample									

MICROTITER PLATE ARRANGEMENT

Human IL-6 Conferma™ ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	Std 4	Reconstituted Standard	Sample [#]								
B	Blank	Std 4	Reconstituted Standard	Sample [#]								
C	Std 1	Std 5	QC1									
D	Std 1	Std 5	QC1									
E	Std 2	Std 6	QC2									
F	Std 2	Std 6	QC2									
G	Std 3	Std 7	QC3									
H	Std 3	Std 7	QC3									

GRAPH OF TYPICAL REFERENCE CURVE



Typical Standard Curve, not to be used to calculate data.

ASSAY CHARACTERISTICS

A. Sensitivity

The lower limit of quantitation (LLOQ) of IL-6 assay is 1.17 pg/mL using Belysa™ Immunoassay Analysis software from Millipore Sigma. LLOQ is calculated by back interpolation of the standard point that provides $CV \leq 20\%$ and recovery $\pm 20\%$ of the expected.

B. Specificity

The antibody pair used in this assay is specific to Human IL-6 and does not cross-react to the following molecules/hormones tested:

Mouse IL-6, Human IL-1a, IL-2, IL-3, IL-4, IL-5, IL-7, IL-8, IL-9, IL-10, IL-12, IL-13, IL-15, $TNF\alpha$, $TNF\beta$, GMCSF, GCSF

C. Precision

Mean Intra-assay precision is calculated from the results of twenty replicates each of the three different concentrations of human IL-6 in a single assay. The mean inter-assay precision is generated from the results of eight separate assays with duplicate samples in each assay for the three different concentrations of IL-6.

ASSAY CHARACTERISTICS (continued)

Intra-Assay Variation

	Mean IL-6 Levels pg/mL	Intra-Assay %CV
1	3.6	4.1
2	10.6	2.1
3	32.1	1.6

Inter-Assay Variation

	Mean IL-6 Levels pg/mL	Inter-Assay %CV
1	3.1	12.9
2	9.6	7.4
3	28.3	6.5

D. Spike Recovery of IL-6 in Blood Samples

Varying amounts of Human IL-6 were added to 10 individual human serum and plasma samples, and the resulting IL-6 content of each sample was assayed by Human IL-6 Conferma™ ELISA.

The recovery = [(observed- Basal / (spiked IL-6 concentration))] x 100%.

Sample	Spiked Concentration of IL-6 pg/mL	Concentration observed in the assay (pg/mL)	Recovery%
Serum 1	0	0	
	4.7	4.4	94
	9.4	9.0	96
	18.8	17.7	95
Serum 2	0	0.0	
	4.7	5.0	107
	9.4	9.9	105
	18.8	18.9	101
Serum 3	0	3.5	
	4.7	8.2	99
	9.4	12.6	97
	18.8	22.1	99

Sample	Spiked Concentration of IL-6 pg/mL	Concentration observed in the assay (pg/mL)	Recovery%
Serum 4	0	0.4	
	4.7	4.8	93
	9.4	9.2	94
	18.8	18.0	94
Serum 5	0	0.3	
	4.7	4.6	91
	9.4	8.9	91
	18.8	17.5	91
Average			93

Sample	Spiked Concentration of IL-6 pg/mL	Concentration observed in the assay (pg/mL)	Recovery%
Plasma 1	0	0.63	
	4.7	5.6	105
	9.4	10.0	100
	18.8	21.2	109
Plasma 2	0	0.0	
	4.7	4.5	96
	9.4	9.3	99
	18.8	19.7	105
Plasma 3	0	0.7	
	4.7	5.6	104
	9.4	10.1	99
	18.8	20.5	105
Plasma 4	0	0.4	
	4.7	5.3	104
	9.4	9.9	101
	18.8	19.2	100
Plasma 5	0	0.4	
	4.7	5.2	104
	9.4	10.1	104
	18.8	20.8	109
Average			103

ASSAY CHARACTERISTICS (continued)

E. Linearity of Dilution

10 spiked individual human serum and plasma samples were assayed for linearity studies. Sample volumes of 50 µL, 25 µL, 12.5 µL, and 6.25 µL represents dilution factors of neat, 2, 4, and 8, respectively. Required amounts of serum matrix were added to compensate for the lost volumes below 50 µL.

$$\text{Dilution linearity} = (\text{observed/expected}) \times 100\%$$

Observed = mean calculated dilution corrected concentration at each dilution

Expected = mean calculated concentration of the neat sample

Sample	Sample volume (µL)	Mean (pg/mL)	Dilution Corrected (pg/mL)	Linearity%
Serum 1	50	16.1	16.1	
	25	8.5	17.0	106
	12.5	4.4	17.8	111
	6.25	2.3	18.6	116
Serum 2	50	17.7	17.7	
	25	9.5	19.0	108
	12.5	4.9	19.6	111
	6.25	2.5	20.2	114
Serum 3	50	20.3	20.3	
	25	10.7	21.5	106
	12.5	5.7	22.7	112
	6.25	2.8	22.6	112
Serum 4	50	18.0	18.0	
	25	9.5	18.9	105
	12.5	4.9	19.7	110
	6.25	2.5	20.3	113
Serum 5	50	16.7	16.7	
	25	9.0	18.0	108
	12.5	4.7	18.8	112
	6.25	2.4	19.2	115
Average				110

Sample	Sample volume (µL)	Mean (pg/mL)	Dilution Corrected (pg/mL)	Linearity%
Plasma 1	50	20.0	20.0	
	25	10.0	19.9	100
	12.5	4.9	19.6	98
	6.25	2.2	17.7	88
Plasma 2	50	19.6	19.6	
	25	9.5	19.0	97
	12.5	4.5	18.2	93
	6.25	2.3	18.3	93
Plasma 3	50	19.8	19.8	
	25	9.8	19.7	100
	12.5	5.0	19.9	101
	6.25	2.3	18.7	95
Plasma 4	50	21.0	21.0	
	25	10.3	20.5	98
	12.5	5.2	20.7	98
	6.25	2.5	19.9	94
Plasma 5	50	19.7	19.7	
	25	9.8	19.6	100
	12.5	5.0	20.0	102
	6.25	2.6	20.6	105
Average				97

NOTE: More data related to assay characteristics can be found in Human IL-6 Conferma™ ELISA verification report.

QUALITY CONTROLS

The ranges for Quality Control 1, 2, and 3 are provided on the card insert or can be located at the MILLIPORE SIGMA website www.milliporesigma.com.

TROUBLESHOOTING GUIDE

1. To obtain reliable and reproducible results, the operator should carefully read this manual and fully understand all aspects of each assay step before attempting to run the assay.
2. Throughout the assay, the operator should adhere strictly to the procedures with good laboratory practice.
3. Have all necessary reagents and equipment ready on hand before starting. Once the assay has been started, all steps should be completed with precise timing and without interruption.
4. Avoid cross-contamination of any reagents or samples to be used in the assay.
5. Make sure all reagents and samples are added to the bottom of each well.
6. Careful and complete mixing of solutions in the well is critical. Poor assay precision will result from incomplete mixing or cross well contamination due to inappropriate mixing.
7. Remove any air bubbles formed in the well after acidification of the substrate solution because bubbles interfere with spectrophotometric readings.
8. High signal in the background or blank wells could be due to 1.) cross well contamination by standard solution or sample or 2.) inadequate washing of wells with Wash Buffer or 3.) overexposure to light after the substrate has been added.

ORDERING INFORMATION

To place an order or to obtain additional information about our immunoassay products, please contact your Customer Service or Technical Support Specialist. Contact information for each region can be found on our website:

emdmillipore.com/contact

Conditions of Sale

For Research Use Only. Not for Use in Diagnostic Procedures.

Safety Data Sheets (SDS)

Safety Data Sheets for EMD Millipore products may be ordered by fax or phone or through our website at emdmillipore.com/msds.