

Mouse Adiponectin
96-Well Plate Assay
Cat. # EZMADP-60K

# MOUSE ADIPONECTIN ELISA KIT 96-Well Plate

# (Cat. # EZMADP-60K)

TAE	BLE OF CONTENTS	<u>PAGE</u>
I.	Intended Use	2
II.	Principles Of Procedure	2
III.	Reagents Supplied	2
IV.	Storage and Stability	4
٧.	Reagent Precautions	4
VI.	Materials Required But Not Provided	6
VII.	Sample Collection And Storage	6
VIII.	Sample Preparation	6
IX.	Standard and Quality Controls Preparation	7
Χ.	Assay Procedure	8
XI.	Microtiter Plate Arrangement	11
XII.	Calculations	12
XIII.	Interpretation	12
XIV.	Standard Curve	13
XV.	Assay Characteristics	14
XVI.	Correlation Graph	17
XVII.	Quality Controls	17
(VIII.	Troubleshooting Guide	18
XIX.	Replacement Reagents	18
XX.	Ordering Information	19

# MOUSE ADIPONECTIN ELISA KIT 96-Well Plate (Cat. # EZMADP-60K)

#### I. INTENDED USE

This Mouse Adiponectin (ACRP30) ELISA kit is used for the non-radioactive quantification of Mouse Adiponectin in serum, plasma, and adipocyte extracts or cell culture media samples. This kit has 100% cross reactivity to Mouse Adiponectin. There is no binding to Mouse Adiponectin globular domain, or to Rat Adiponectin. One kit is sufficient to measure 38 unknown samples in duplicate. *For Research Use Only. Not for Use in Diagnostic Procedures.* 

#### II. PRINCIPLES OF PROCEDURE

This assay is a Sandwich ELISA based, sequentially, on: 1) concurrent capture of Mouse Adiponectin molecules from samples to the wells of a microtiter plate coated by a pre-titered amount of anti- mouse adiponectin monoclonal antibodies, and binding of a second biotinylated anti- mouse polyclonal antibody to the captured molecules, 2) wash away of unbound materials from samples, 3) conjugation of horseradish peroxidase to the immobilized biotinylated antibodies, 4) wash away of free enzyme conjugates, and 5) quantification of immobilized antibody-enzyme conjugates by monitoring horseradish peroxidase activities in the presence of the substrate 3,3',5,5'-tetramethylbenzidine. The enzyme activity is measured spectrophotometrically by the increased absorbency at 450 nm, corrected from the absorbency at 590nm, after acidification of formed products. Since the increase in absorbency is directly proportional to the amount of captured Mouse Adiponectin in the unknown sample, the latter can be derived by interpolation from a reference curve generated in the same assay with reference standards of known concentrations of Mouse Adiponectin.

#### III. REAGENTS SUPPLIED

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

# A. Mouse Adiponectin ELISA Plate

Coated with Rat Monoclonal anti-Mouse Adiponectin Antibodies

Quantity: 1 plate

Preparation: Ready to Use

Note: Unused strips should be resealed in the foil pouch with the desiccant provided and

stored at 2-8°C.

#### **B.** Adhesive Plate Sealer

Quantity: 1 sheet

Preparation: Ready to Use

#### III. REAGENTS SUPPLIED (continued)

#### C. 10X HRP Wash Buffer Concentrate

10X concentrate of 50mM Tris Buffered Saline containing Tween-20.

Quantity: Two bottles containing 50 mL each

Preparation: Dilute 1:10 with distilled or deionized water.

# D. Mouse Adiponectin Standard

Mouse Adiponectin, lyophilized. Quantity: 1mL upon hydration

Preparation: Contents Lyophilized. Reconstitute with 1 mL distilled or deionized water. The actual concentration of Mouse Adiponectin present in the vial will be lot dependent. Please refer to the analysis sheet for exact concentration present in a specific lot.

#### E. Quality Controls 1 and 2

Purified Recombinant Mouse Adiponectin in Assay Buffer, lyophilized.

Quantity: 1mL/vial upon hydration

Preparation: Contents Lyophilized. Reconstitute with 1 mL distilled or deionized water.

#### F. 10X Assay Buffer

Quantity: 50 mL

Preparation: Dilute 1:10 with distilled or deionized water to make 1X assay buffer (0.05M Phosphosaline containing 0.025M EDTA, 0.08% Sodium Azide, 1% BSA).

Note: Use 1X Assay Buffer to dilute samples (Section VIII, SAMPLE PREPARATION) and Standard Curve (SECTION IX, STANDARD AND QUALITY CONTROL PREPARATION) and in Assay Procedure (SECTION X, ASSAY PROCEDURE).

### G. Mouse Adiponectin Detection Antibody

Pre-titered Biotinylated Goat anti-Mouse Adiponectin Polyclonal Antibody

Quantity: 3.0 mL

Preparation: Ready to Use

### H. Enzyme Solution

Pre-titered Streptavidin-Horseradish Peroxidase Conjugate in Buffer

Quantity: 12 mL

Preparation: Ready to Use

# I. Substrate (Light sensitive, avoid unnecessary exposure to light)

3, 3', 5, 5'-tetramethylbenzidine in buffer

Quantity: 12 mL

Preparation: Ready to Use.

#### J. Stop Solution (Caution: Corrosive Solution)

0.3 M HCI

Quantity: 12 mL

Preparation: Ready to Use

#### IV. STORAGE AND STABILITY

Recommended storage for kit components is 2-8°C.

All components are shipped and stored at 2-8°C. Reconstituted standards and quality controls can be frozen for future use, but repeated freeze thaws should be avoided. Refer to expiration dates on all reagents prior to use. Do not mix reagents from different kits unless they have the same lot numbers

#### V. REAGENT PRECAUTIONS

#### A. Sodium Azide

Sodium Azide or Proclin has been added to some reagents as a preservative. Although the concentrations are low, Sodium Azide and Proclin 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

#### **B.** Hydrochloric Acid

Hydrochloric Acid is corrosive and can cause eye and skin burns. It is harmful if swallowed and can cause respiratory and digestive tract burns. Avoid contact with skin and eyes. Do not swallow or ingest.

Note: See Full Labels of Hazardous components on next page.

# Full Hazardous Labeling for components in this kit:

Ingredient, Cat #		Full Label			
Mouse Adiponectin Quality Controls 1 & 2	E6300-K	<b>1 1 1 1 1 1 1 1 1 1</b>	Danger. Toxic if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment. IF exposed or concerned: immediately call a POISON CENTER or doctor/physician.		
Mouse Adiponectin Standards	E8060-K		Danger. Toxic if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment. IF exposed or concerned: immediately call a POISON CENTER or doctor/physician.		
Stop Solution	ET-TMB		Warning. May be corrosive to metals.		
10X HRP Wash Buffer Concentrate	EWB-HRP	<u>(!)</u>	Warning. May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.		

#### VI. MATERIALS REQUIRED BUT NOT PROVIDED

- 1. Pipettes and Pipette Tips: 10μL 20 μL or 20μL 100 μL
- 2. Multi-Channel Pipettes and Pipette Tips: 5 ~ 50 μL and 50 ~ 300 μL
- 3. Buffer and Reagent Reservoirs
- 4. Vortex Mixer
- 5. Deionized Water
- 6. Microtiter Plate Reader capable of reading absorbency at 450 nm
- 7. Orbital Microtiter Plate Shaker
- 8. Absorbent Paper or Cloth

#### VII. SAMPLE COLLECTION AND STORAGE

1. To prepare serum samples, whole blood is directly drawn into a centrifuge tube that contains no anti-coagulant. Let blood clot at room temperature for 30 min.

Promptly centrifuge the clotted blood at 2,000 to 3,000 x g for 15 minutes at  $4 \pm 2^{\circ}$ C.

Transfer and store serum samples in separate tubes. Date and identify each sample.

Use freshly prepared serum or aliquot and store samples at  $\leq$  -20°C for later use. For long-term storage, keep at -70°C. Avoid freeze/thaw cycles.

- 2. To prepare plasma samples, whole blood should be collected into centrifuge tubes containing enough K<sub>3</sub>EDTA to achieve a final concentration of 1.735 mg/mL and centrifuged immediately after collection. Observe the same precautions in the preparation of serum samples.
- 3. If heparin is to be used as an anticoagulant, the effect on the assay outcome at the dose of heparin used should be pre-determined.
- 4. Avoid using samples with gross hemolysis or lipemia.

#### VIII. SAMPLE PREPARATION

- 1. Allow all the reagents to come to room temperature.
- 2. Dilute serum or plasma samples 1:1000 in 1X Assay Buffer (See Section III, F). Cellular extract and culture media dilutions will vary.
  - a. Make Dilution A by adding 10 µL sample to 990 µL Assay Buffer and vortex.
  - b. Make Dilution B by adding 100  $\mu$ L of Dilution A to 900  $\mu$ L Assay Buffer and vortexing. Use Dilution B (1:1000) for the assay procedure.

#### IX. STANDARD AND QUALITY CONTROLS PREPARATION

#### A. Mouse Adiponectin Standard Preparation

- 1. Use care in opening the lyophilized Standard vial. Using a pipette, reconstitute the Mouse Adiponectin Standard with 1.0 mL distilled or deionized water into the vial to give a concentration prescribed in the analysis sheet. Invert and mix gently, let sit for 5 minutes then mix well.
- 2. Label seven tubes 1, 2, 3, 4, 5, 6, and 7. Add 0.5 mL Assay Buffer to each of the seven tubes. Prepare serial dilutions by adding 0.5 mL of the reconstituted standard to tube 1, mix well and transfer 0.5 mL of tube 1 to tube 2, mix well and transfer 0.5 mL of tube 2 to tube 3, mix well and transfer 0.5 mL of tube 3 to tube 4, mix well and transfer 0.5 mL of tube 4 to tube 5, mix well and transfer 0.5 mL of tube 5 to tube 6, mix well and transfer 0.5 mL of tube 6 to tube 7 and mix well.

**Note:** Do not use a Repeater pipette. Change tip for every dilution. Wet tip with Standard before dispensing. Unused portions of reconstituted standard should be stored at  $\leq$  -20°C. Avoid multiple freeze/thaw cycles.

Volume of	Volume of Standard to	Standard
Deionized Water	Add	Concentration ng/mL
to Add		-
1 mL	0	X
		(Refer to analysis
		sheet for exact
		concentration)

	Volume of Assay	Volume of Standard to	Standard
Tube #	Buffer to Add	Add	Concentration ng/mL
1	0.5 mL	0.5 mL of reconstituted standard	X/2
2	0.5 mL	0.5 mL of tube 1	X/4
3	0.5 mL	0.5 mL of tube 2	X/8
4	0.5 mL	0.5 mL of tube 3	X/16
5	0.5 mL	0.5 mL of tube 4	X/32
6	0.5 mL	0.5 mL of tube 5	X/64
7	0.5 mL	0.5 mL of tube 6	X/128

#### B. Mouse Adiponectin Quality Control 1 and 2 Preparation

 Use care in opening the lyophilized Quality Control vials. Using a pipette, reconstitute each of the Mouse Adiponectin Quality Control 1 and Quality Control 2 with 1 mL distilled or deionized water into the vials. Invert and mix gently, let sit for 5 minutes then mix well.

#### X. ASSAY PROCEDURE

Pre-warm all reagents to room temperature prior to setting up the assay.

Note: Since the standard concentration may vary from lot to lot, please make sure to change the standard concentration in the template while reading absorbance in the spectrophotometer.

- 1. Dilute the 10X Wash Buffer concentrate 10 fold by mixing the entire content of each bottle of Wash Buffer with 450 mL deionized water (dilute both bottles with 900 mL deionized water).
- 2. 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 and wash each well 3 times with 300 μL of diluted Wash Buffer per wash. Decant Wash Buffer and remove the residual amount from all wells 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, use a gentle wash program for all washing steps described in this protocol.
- 3. Add 60 µL Assay Buffer to all wells.
- 4. Add in duplicate 20 μL Assay Buffer to the blank wells.
- 5. Add in duplicate 20  $\mu$ L Mouse Adiponectin Standards in the order of ascending concentration to the appropriate wells. Add in duplicate 20  $\mu$ L QC1 and 20  $\mu$ L QC2 to the appropriate wells. Add sequentially 20  $\mu$ L of the unknown samples in duplicate to the remaining wells
- 6. Add 20 µL Detection Antibody to all wells. For best result all additions should be completed within one hour. Cover the plate with plate sealer and incubate at room temperature for 2 hours on an orbital microtiter plate shaker set to rotate at moderate speed, approximately 400 to 500 rpm.
- 7. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in the wells.
- 8. Wash wells 5 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap firmly after each wash to remove residual buffer.
- 9. Add 100 µL Enzyme Solution to each well. Cover plate with sealer and incubate with moderate shaking at room temperature for 30 minutes on the microtiter plate shaker.
- Remove sealer, decant solutions from the plate, and tap plate to remove the residual fluid.

#### X. ASSAY PROCEDURE (continued)

- 11. Wash wells 5 times with diluted Wash Buffer, 300 µl per well per wash. Decant and tap firmly 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** 5 to 20 minutes. Blue color should be formed in wells of Adiponectin standards with intensity proportional to increasing concentrations of Adiponectin.

**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. One can monitor color development using 370 nm filter, if available on the spectrophotometer. When the absorbance is between 1.2 and 1.8 at 370 nm, the stop solution can be added to terminate the color development.

13. Remove sealer and add 100 µl Stop Solution **[CAUTION: CORROSIVE SOLUTION]** and shake plate by hand to ensure complete mixing of 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 plate reader. Read absorbance at 450 nm and 590nm 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 highest Adiponectin standard should be approximately 2.2-2.8 or not to exceed the capability of the plate reader used.

# Assay Procedure for Mouse Adiponectin ELISA kit (Cat. # EZMADP-60K)

	Step 1	Step 2	Step 3-4	Step 5	Step 6	Step 6-8	Step 9	Step 10-11	Step 12	Step 12	Step 13	Step 13
Well #			Assay Buffer	Standards/ Controls/ Samples	Detectio n Ab		Enzyme Solution		Substrate		Stop Solution	
A1, B1	ater.	<u>s</u>	80 µL		20 µL		100 μL		100 µL		100 μL	
C1, D1	zed Wa	nt towe	60 µL	20 μL of Tube 7		ure.		iture .		20 minutes at Room Temperature.		
E1, F1	Deioni	er. sorber	60 µL	20 μL of Tube 6		at Room Temperature. Wash Buffer		mpera		edwej		
G1, H1	0mL [	μL Wash Buffer. smartly on absor	60 µL	20 μL of Tube 5		ate, Incubate 2 hours at Room Terr Wash 5X with 300 µL Wash Buffer		om Te 3uffer		L moo!		nm.
A2, B2	vith 45	L Was nartly	60 µL	20 μL of Tube 4		at Rool Nash I		at Ro Nash I		es at R		nd 590
C2, D2	uffer v	300 μ Jing sr	60 µL	20 μL of Tube 3		hours a		inutes 00 µL \		minute		Read Absorbance at 450 nm and 590 nm.
E2, F2	/ash B	Wash plate 3X with 300 idual buffer by tapping s	60 µL	20 μL of Tube 2		ate 2 h with 30		e 30 m with 30				at 450
G2, H2	10X W	olate 3 uffer k	60 µL	20 μL of Tube 1		Incubash 5X v		cubate sh 5X v		ubate		bance
A3, B3	ttle of	Nash g	60 µL	20 μL of QC I		gitate, Was		ate, In Was		te, Inc		Absor
C3, D3	ach bo	ا ر	60 µL	20 μL of QC II		Seal, Agitate, Incubate Wash 5X with		Seal, Agitate, Incubate 30 minutes at Room Temperature Wash 5X with 300 µL Wash Buffer		Seal, Agitate, Incubate 5		Read
E3, F3	Dilute each bottle of 10X Wash Buffer with 450mL Deionized Water.	Wash plate 3X with 300 μL Wash Buffer. Remove residual buffer by tapping smartly on absorbent towels	60 µL	20 μL of Sample		Ø		Sea		Seal,		
G3, H3	Ö	<u> </u>	60 µL	20 μL of Sample								
A4, B4 ↓			60 µL	20 μL of Sample								

# XI. MICROTITER PLATE ARRANGEMENT

Mouse Adiponectin ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
А	Blank	Tube 4	QC 1	Sample								
В	Blank	Tube 4	QC 1	Sample								
С	Tube 7	Tube 3	QC 2	Etc.								
D	Tube 7	Tube 3	QC 2									
Е	Tube 6	Tube 2	Sample									
F	Tube 6	Tube 2	Sample									
G	Tube 5	Tube 1	Sample									
Н	Tube 5	Tube 1	Sample									

#### XII. CALCULATIONS

The dose-response curve of this assay fits best to a sigmoidal 4- or 5-parameter logistic equation. The results of unknown samples can be calculated with any computer program having a 4- or 5-parameter logistic function. Final results should be multiplied by a 1000 dilution factor.

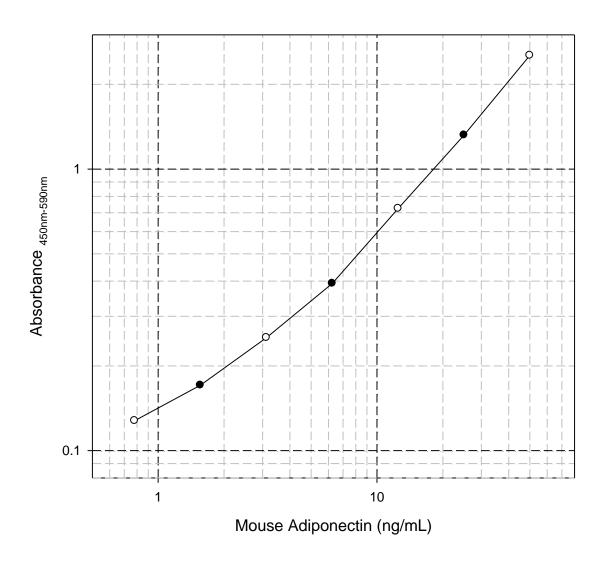
**Note:** When sample volumes assayed differ from 20  $\mu$ L, an appropriate mathematical adjustment must be made to accommodate for the dilution factor (e.g., if 10  $\mu$ L of sample is used, then calculated data must be multiplied by 2). When sample volume assayed is less than 20  $\mu$ L, compensate the volume deficit with assay buffer.

#### XIII. INTERPRETATION - Acceptance Criteria

- The assay will be considered accepted when all Quality Control values fall within the calculated Quality Control Range. If any QC's fall outside the control range, review results with a supervisor.
- 2. If the difference between duplicate results of a sample is >15% CV, repeat the sample.
- 3. The limit of sensitivity of this assay is 0.2 ng/mL mouse Adiponectin (20  $\mu$ L sample size).
- 4. The approximate dynamic range of this assay is 1 50 ng/mL mouse Adiponectin (20  $\mu$ l sample size). Any result greater than 50 ng/mL in a 20  $\mu$ L sample should be diluted using 1X assay buffer, and the assay repeated until the results fall within range.

# **XIV.STANDARD CURVE**

# **Mouse Adiponectin ELISA**



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

#### XV. ASSAY CHARACTERISTICS

#### A. Sensitivity

The lowest level of Adiponectin that can be detected by this assay is 0.2 ng/mL when using a  $20 \mu L$  sample size.

## B. Specificity

The antibody pair used in this assay is specific to mouse adiponectin and does not significantly cross-react with human adiponectin, globular domain of mouse adiponectin, and other cytokine or hormone molecules tested, as shown in the following table.

Analyte	Max. Conc.	Cross Reactivity
Human Adiponectin	200ng/mL	< 2%
Mouse gAcrp	1000ng/mL	n.d.
Mouse Endocrines:		
Insulin	10nM	n.d.
Amylin	10nM	n.d.
Leptin	10nM	n.d.
Glucagon	10nM	n.d.
GLP-1	10nM	n.d.
Mouse Cytokines:	40.14	
IL-1ß	10nM	n.d.
IL-2	10nM	n.d.
IL-4	10nM	n.d.
IL-5	10nM	n.d.
IL-6	10nM	n.d.
IL-9	10nM	n.d.
IL-10	10nM	n.d.
IL-12	10nM	n.d.
IL-13	10nM	n.d.
IFN-γ TNF-α	10nM 10nM	n.d.
GMCSF	10nM	n.d. n.d.
MIP-1α	10nM	n.d.
KC	10nM	n.d.
RANTES	10nM	n.d.
MCP-1	10nM	n.d.
IVIOI - I	TOTTIVI	n.u.

#### n.d. = not detectable

No cross reactivity to rat serum or plasma samples.

#### XV. ASSAY CHARACTERISTICS (continued)

#### C. Precision

Within and Between Assay Variation

Sample No.	Mean Adiponectin	Within% CV	Between% CV
	Levels (ng/mL)		
1	9.18	3.8	6.8
2	13.74	5.4	1.4
3	6.89	8.2	4.9
4	7.23	5.6	10.8

The assay variations of EMD Millipore Mouse Adiponectin ELISA kits were studied on four mouse serum samples with varying concentrations of endogenous Adiponectin. The mean within variation was calculated from results of six duplicate determinations in each assay of the indicated samples. The mean between variations of each sample was calculated from results of four separate assays with duplicate samples in each assay.

#### D. Recovery

Spike & Recovery of Mouse Adiponectin in Serum

Sample	Adiponectin	Expected	Observed	% of
No.	Added	(ng/mL)	(ng/mL)	Recovery
	(ng/mL)		, , ,	•
1	0	7.1	7.1	100
	5	12.1	12.6	104
	10	17.1	17	99
	20	27.1	26.7	99
2	0	7.2	7.2	100
	5	12.2	13.7	112
	10	17.5	17.8	103
	20	27.2	28.9	106
3	0	10.49	10.49	100
	5	15.49	15.76	102
	10	20.49	22.17	108
	20	30.49	31.07	102

Varying amounts of mouse Adiponectin were added to three mouse serum samples and the Adiponectin content was determined in three separate assays. The % of recovery = observed Adiponectin concentrations/expected Adiponectin concentrations x 100%.

# XV. ASSAY CHARACTERISTICS (continued)

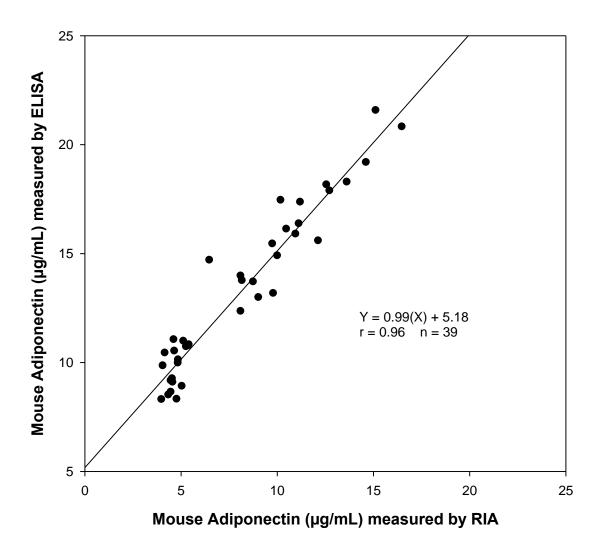
#### E. Linearity

#### Effect of Serum Dilution

Sample	Volume	Expected	Observed	% Of
No.	Sampled	(ng/mL)	(ng/mL)	Expected
	(µL)			
1	20	14.3	14.3	100
	10	7.15	7.3	102
	5	3.6	4.1	114
	3.33	2.4	2.9	121
2	20	14.99	14.99	100
	10	7.49	7.23	96
	5	3.75	4.37	116
	3.33	2.5	2.81	112
3	20	14.43	14.43	100
	10	7.2	7.17	99
	5	3.61	4.16	115
	3.33	2.41	2.58	107

Three mouse serum samples with the indicated sample volumes were assayed in three separate experiments. Required amounts of assay buffer were added to compensate for lost volumes below 20  $\mu$ L. The resulting dilution factors of 1.0, 2.0, 4.0, and 6.0 representing 20  $\mu$ L, 10  $\mu$ L, 5  $\mu$ L, and 3.3  $\mu$ L sample volumes assayed, respectively, were applied in the calculation of observed Adiponectin concentrations. % expected = observed/expected x 100%.

# Adiponectin Correlation RIA vs. ELISA in Mouse Serum Samples



Serum samples obtained from 39 mouse subjects were assayed for Adiponectin content using both EMD Millipore Mouse Adiponectin RIA Kit (Catalogue #MADP-60HK) and Mouse Adiponectin ELISA Kit (Catalogue EZMADP-60K). Correlation of the two kits is derived by linear regression analysis of paired results from each sample.

#### **XVII. QUALITY CONTROLS**

The ranges for Quality Control 1 and 2 are provided on the card insert or can be located at the EMD Millipore website <a href="mailto:emdmillipore.com">emdmillipore.com</a> using the catalog number as the keyword.

#### XVIII. 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 bubble formed in the well after acidification of substrate solution because bubbles interfere with spectrophotometric readings.
- 8. Do not let the absorbency reading of the highest standard reach 3.0 units or higher after acidification.
- High absorbance in background or blank wells could be due to 1) cross well contamination by standard solution or sample or 2) inadequate washing of wells with HRP Wash Buffer or 3) overexposure to light after substrate has been added.

#### XIX. REPLACEMENT REAGENTS

Reagents	Cat. #
Mouse Adiponectin ELISA Plate 10X HRP Wash Buffer Concentrate (50 mL) Mouse Adiponectin Standards Quality Controls 1 and 2 Assay Buffer (10X) Mouse Adiponectin Detection Antibody Enzyme Solution Substrate Stop Solution	EP60 EWB-HRP E8060-K E6300-K EAB-10XP E1060 EHRP-3 ESS-TMB ET-TMB
10-pack of Mouse Adiponectin (ACRP30) ELISA kits	EZMADP-60BK

#### XX. 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.