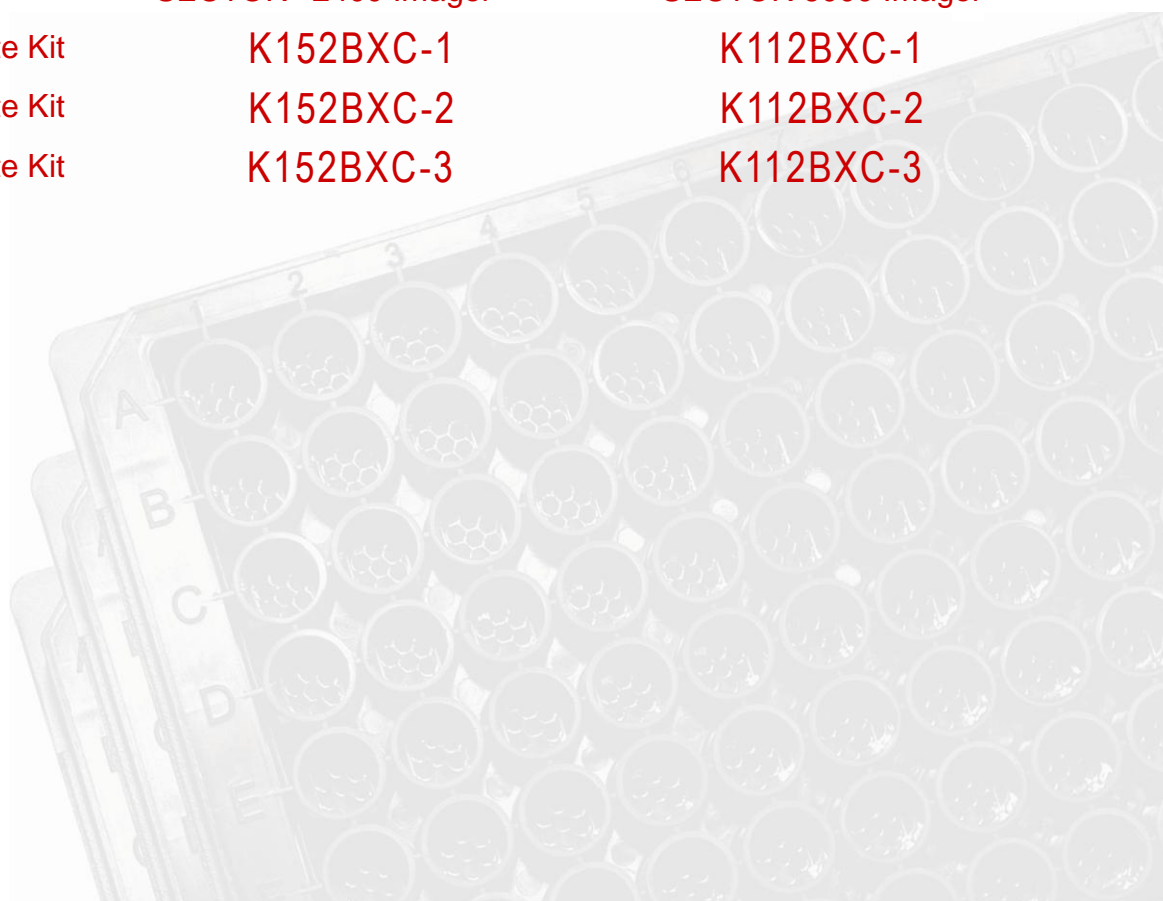


# Meso Scale Discovery<sup>®</sup>

## MULTI-ARRAY<sup>®</sup> Assay System

### Mouse Adiponectin Assay Kit

	SECTOR <sup>®</sup> 2400 Imager	SECTOR 6000 Imager
1-Plate Kit	K152BXC-1	K112BXC-1
5-Plate Kit	K152BXC-2	K112BXC-2
20-Plate Kit	K152BXC-3	K112BXC-3



Meso Scale Discovery Meso Scale Discovery Meso Scale Discovery Meso Scale Discovery Meso Scale Discovery Meso Scale Discovery Meso Scale Di



# MSD Metabolic Assays

## Mouse Adiponectin Assay Kit

*This package insert must be read in its entirety before using this product.*

**FOR RESEARCH USE ONLY.**

**NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

### Meso Scale Discovery

A division of Meso Scale Diagnostics, LLC.

9238 Gaither Road

Gaithersburg, MD 20877 USA

[www.mesoscale.com](http://www.mesoscale.com)

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, SECTOR, SECTOR HTS and SECTOR PR are trademarks and/or service marks of Meso Scale Diagnostics, LLC.  
© 2010 Meso Scale Diagnostics, LLC. All rights reserved.

# Table of Contents

table of contents

I. Introduction .....	4
II. Principle of the Assay .....	4
III. Reagents Supplied.....	5
IV. Required Material and Equipment – not supplied .....	5
V. Safety.....	5
VI. Reagent Preparation.....	6
VII. Assay Protocol.....	7
VIII. Analysis of Results.....	7
IX. Typical Standard Curve .....	8
X. Sensitivity.....	8
XI. Endogenous levels.....	9
XII. Spike Recovery.....	9
XIII. Linearity .....	10
XIV. Assay Components.....	10
XV. References .....	11
Summary Protocol .....	13
Plate Diagrams .....	15

## Ordering Information

ordering information

### MSD Customer Service

Phone: 1-301-947-2085  
Fax: 1-301-990-2776  
Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)

### MSD Scientific Support

Phone: 1-301-947-2025  
Fax: 1-240-632-2219 attn: Scientific Support  
Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)

# Introduction

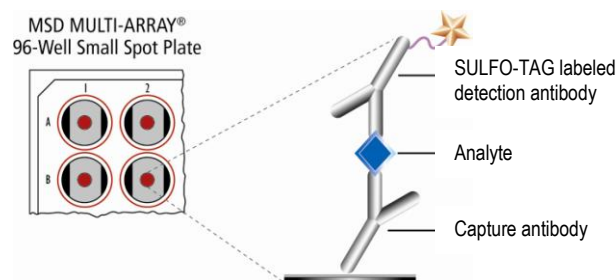
introduction

**Adiponectin** is a 30 kD protein secreted exclusively by adipocytes and is thought to play a role in lipid and glucose metabolism. Adiponectin enhances insulin action by activating glucose uptake and fatty acid oxidation. In addition, this adipokine has potent anti-inflammatory and anti-atherosclerotic properties. Adiponectin circulates at high levels (between 2-20  $\mu\text{g/mL}$ ) in plasma as trimers, hexamers and high molecular weight multimers. In contrast to most adipokines, plasma adiponectin is often negatively correlated with body mass index (BMI) in humans and rodents. However, expression in various adipose tissues and disease models may vary.

## Principle of the Assay

principle of the assay

MSD<sup>®</sup> metabolic assays provide rapid and convenient methods for measuring the levels of protein targets within single small-volume samples. The assays are available in both singleplex and multiplex formats. In a singleplex assay, an antibody for a specific protein target is coated on one electrode (or “spot”) per well. In a multiplex assay, an array of capture antibodies against different targets is patterned on distinct spots in the same well. Our Mouse Adiponectin Assay detects adiponectin in a sandwich immunoassay (Figure 1). MSD provides a plate that has been pre-coated with adiponectin antibody. The user adds the sample and a solution containing the labeled detection antibody—anti-adiponectin labeled with an electrochemiluminescent compound, MSD SULFO-TAG<sup>™</sup> label—over the course of one or more incubation periods. Adiponectin in the sample binds to capture antibody immobilized on the working electrode surface; recruitment of the labeled detection antibody by bound analyte completes the sandwich. The user adds an MSD read buffer that provides the appropriate chemical environment for electrochemiluminescence and loads the plate into an MSD SECTOR instrument for analysis. Inside the SECTOR instrument, a voltage applied to the plate electrodes causes the labels bound to the electrode surface to emit light. The instrument measures intensity of emitted light to afford a quantitative measure of adiponectin present in the sample.



**Figure 1.** Sandwich immunoassay on MSD platform



# Reagents Supplied

reagents supplied

Product Description	Storage	Quantity per Kit		
		K152BXC-1 K112BXC-1	K152BXC-2 K112BXC-2	K152BXC-3 K112BXC-3
MULTI-ARRAY 96-well Mouse Adiponectin Plate(s) L452BXB-1 (K152BXC) L412BXB-1 (K112BXC)	2-8°C	1 plate	5 plates	20 plates
SULFO-TAG Anti-mAdiponectin Antibody <sup>1</sup> (100X)	2-8°C	1 vial (40 µL)	1 vial (200 µL)	4 vials (200 µL ea)
Mouse Adiponectin Calibrator 50 µg/mL	≤-70°C	1 vial (15 µL)	5 vials (15 µL ea)	20 vials (15 µL ea)
Blocker A Kit R93AA-2 (250 mL)	RT	1 bottle (250 mL)	1 bottle (250 mL)	4 bottles (250 mL ea)
Diluent 100 R50AA-2 (200 mL) R50AA-3 (1 L)	2-8°C	1 bottle (200 mL)	2 bottles (200 mL ea)	2 bottles (200 mL & 1 L)
Read Buffer T (4X) R92TC-3 (50 mL) R92TC-2 (200 mL)	RT	1 bottle (50 mL)	1 bottle (50 mL)	1 bottle (200 mL)



## Required Materials and Equipment - not supplied

required materials and equipment - not supplied

- Deionized water for diluting concentrated buffers
- 50 mL tubes for reagent preparation
- 15 mL tubes for reagent preparation
- Microcentrifuge tubes for preparing serial dilutions
- Phosphate buffered saline plus 0.05% Tween-20 (PBS-T) for plate washing
- Appropriate liquid handling equipment for desired throughput, capable of dispensing 25 to 150 µL into a 96-well microtiter plate
- Plate washing equipment: automated plate washer or multichannel pipette
- Adhesive plate seals
- Microtiter plate shaker



## Safety

safety

Safe laboratory practices and personal protective equipment such as gloves, safety glasses, and lab coats should be used at all times during the handling of all kit components. All hazardous samples should be handled and disposed of properly, in accordance with local, state, and federal guidelines.

<sup>1</sup> Some SULFO-TAG labeled detection antibodies may be light-sensitive, so they should be stored in the dark.

# VI Reagent Preparation

## reagent preparation

Bring all reagents to room temperature and thaw the Calibrator stock on ice.

### Prepare Blocker A Solution

Follow instructions included with the Blocker A Kit.

### Prepare Calibrator and Control Solutions

Calibrator for the Mouse Adiponectin Assay is supplied at 50 µg/mL. For the assay, an 8-point standard curve is recommended with 5-fold serial dilution steps and a zero Calibrator. The table below shows the concentrations of the 8-point standard curve:

Standard	Adiponectin conc. (ng/mL)	Dilution Factor
Stock Cal. Vial	50000	
STD-01	1000	50
STD-02	200	5
STD-03	40	5
STD-04	8.0	5
STD-05	1.6	5
STD-06	0.32	5
STD-07	0.064	5
STD-08	0	n/a

To prepare this 8-point standard curve:

- 1) Prepare the highest Calibrator by transferring 10 µL of the Calibrator stock vial to 490 µL of Diluent 100.
- 2) Prepare the next Calibrator by transferring 20 µL of the diluted Calibrator to 80 µL of Diluent 100. Repeat 5-fold serial dilutions 5 additional times to generate 7 Calibrators.
- 3) The recommended 8<sup>th</sup> Standard is Diluent 100 (i.e. zero Calibrator).
- 4) Diluted Calibrators should be kept on ice prior to addition to the plate.

**Note:** The standard curve can be modified as necessary to meet specific assay requirements.

### Preparation of Serum and Plasma Samples

The assay format requires 10 µL of pre-diluted serum/plasma samples per well. An adequate volume of each sample should be prepared depending upon desired number of replicates. The serum/plasma samples should be pre-diluted with Diluent 100 one thousand fold (some samples may require larger dilutions). A two-step dilution is recommended; an initial 10-fold dilution followed by a second 100-fold dilution.

- 10 µL of sample + 90 µL of Diluent 100
- 10 µL of the above dilution + 990 µL of Diluent 100

### Prepare Detection Antibody Solution

The Detection Antibody is provided at 100X stock solution. The final concentration of the working Detection Antibody Solution should be at 1X. For each plate used, dilute a 50 µL aliquot of the stock Detection Antibody into 4.95 mL of Diluent 100.

## Prepare Read Buffer

The Read Buffer should be diluted 4-fold in deionized water to make a final concentration of 1X Read Buffer T. Add 5 mL of 4X Read Buffer T to 15 mL of deionized water for each plate.

## Prepare MSD Plate

This plate has been pre-coated with antibody for the analyte shown in Figure 1. The plate can be used as delivered; no additional preparation (e.g., pre-wetting) is required. The plate has also been exposed to a proprietary stabilizing treatment to ensure the integrity and stability of the immobilized antibodies.

# VII Assay Protocol

assay protocol

1. **Addition of Blocker A Solution:** Dispense 150  $\mu\text{L}$  of Blocker A Solution into each well. Seal the plate with an adhesive plate seal and incubate for 1 hour with vigorous shaking (300–1000 rpm) at room temperature.
2. **Wash and Addition of the Detection Antibody Solution followed by Sample or Calibrator:** Wash the plate 3 times with PBS-T. Dispense 40  $\mu\text{L}$  of 1X Detection Antibody Solution into each well of the MSD plate. Immediately add 10  $\mu\text{L}$  of sample or Calibrator into the appropriate wells of the MSD plate. Seal the plate with an adhesive plate seal and incubate for 2 hours with vigorous shaking (300–1000 rpm) at room temperature.
3. **Wash and Read:** Wash the plate 3 times with PBS-T. Add 150  $\mu\text{L}$  of 1X Read Buffer T to each well of the MSD plate. Analyze the plate on the SECTOR Imager. Plates may be read immediately after the addition of Read Buffer.

## Notes

*Shaking a 96-well MSD MULTI-ARRAY plate typically accelerates capture at the working electrode.*

*Bubbles in the fluid will interfere with reliable reading of MULTI-ARRAY plate. Use reverse pipetting techniques to insure bubbles are not created when dispensing the Read Buffer.*

# VIII Analysis of Results

analysis of results

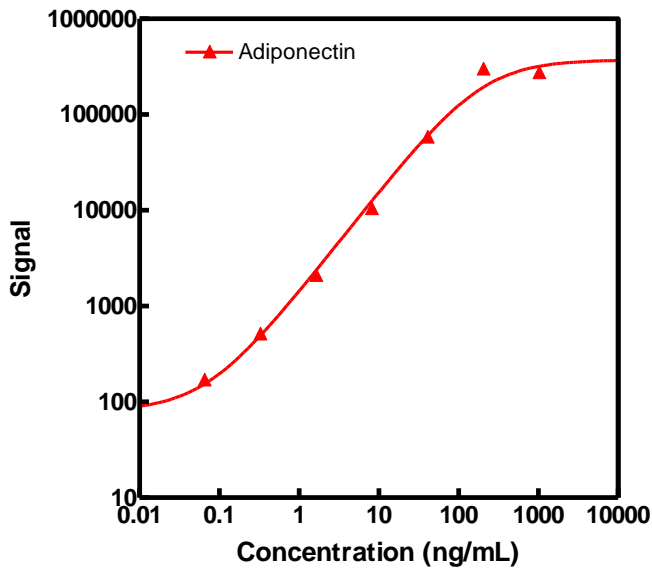
The Calibrators should be run in duplicate to generate a standard curve. The standard curve is modeled using least squares fitting algorithms so that signals from samples with known levels of the analyte of interest can be used to calculate the concentration of analyte in the sample. The assays have a wide dynamic range (3–4 logs) which allows accurate quantification in many samples without the need for dilution. The MSD DISCOVERY WORKBENCH<sup>®</sup> analysis software utilizes a 4-parameter logistic model (or sigmoidal dose-response) and includes a  $1/Y^2$  weighting function. The weighting functionality is important because it provides a better fit of data over a wide dynamic range, particularly at the low end of the standard curve.

# IX Typical Standard Curve

typical standard curve

The MSD Mouse Adiponectin Assay is designed for use with mouse serum and plasma samples.

The following standard curve is an example of the dynamic range of the assay. The actual signals may vary. A standard curve should be run for each set of samples and on each plate for the best quantification of unknown samples.



Adiponectin		
Conc. (ng/mL)	Average Signal	%CV
0	84	4.3
0.064	175	2.0
0.32	529	1.0
1.6	2165	4.3
8.0	10735	3.1
40	60188	2.9
200	308547	2.0
1000	283295	3.1

# X Sensitivity

sensitivity

The lower limit of detection (LLOD) is the calculated concentration of the signal that is 2.5 standard deviations over the zero Calibrator. The value below represents the average LLOD over multiple kit lots.

Adiponectin	
LLOD (ng/mL)	0.04

# XI

## Endogenous Levels

endogenous levels

Endogenous levels of mouse adiponectin in normal and obese mouse serum samples. Samples were diluted 1000 fold prior to measurement. The data below represents final concentrations which have been corrected for dilutions.

Sample	Serum (µg/mL)	Obese Serum (ob/ob) (µg/mL)
1	6.4	14.9
2	10.2	13.2
3	9.3	14.7
4	7.6	14
5	8.4	13.8
6	7.2	
7	6.5	
8	6.2	

# XII

## Spike Recovery

spike recovery

Diluted serum, EDTA plasma, and heparin plasma were spiked with the Calibrators at multiple values throughout the range of the assay. Measured analyte represents average spike recovery in 3 pooled mouse samples.

% Recovery = measured / expected x 100

	Spike Conc. (pg/mL)	% Recovery
Spiked Serum	5	96
	50	95
	500	104
Spiked EDTA Plasma	5	110
	50	112
	500	106
Spiked Heparin Plasma	5	113
	50	105
	500	109

# XIII Linearity

linearity

Linearity was measured by spiking Calibrator levels in 3 pooled mouse samples followed by subsequent dilution.

Percent recovery is calculated as the measured concentration divided by the concentration of the previous dilution (expected).

$$\% \text{ Recovery} = \text{measured} \times \text{dilution factor} / \text{expected} \times 100$$

	Fold Dilution	% Recovery
Serum	2	105
	4	101
	8	94
EDTA Plasma	2	111
	4	122
	8	117
Heparin Plasma	2	113
	4	96
	8	105

# XIV Assay Components

assay components

Calibrator	
Analyte	Mouse adiponectin
Source	Purified, recombinant mouse adiponectin with a FLAG-Tag expressed in HEK 293

Capture Antibody	
Analyte	Mouse adiponectin
Source	Rat monoclonal
Isoforms Recognized	Detects full-length mouse adiponectin (amino acids 18 - 247), does not cross-react with recombinant human adiponectin or the C-terminal globular domain of mouse adiponectin (amino acids 111 - 247)
Species cross-reactivity	Mouse

Detection Antibody	
Analyte	Mouse adiponectin
Source	Goat polyclonal
Isoforms Recognized	n/a
Species cross-reactivity	Mouse

1. Lihn AS, Pedersen SB, Richelsen B. Adiponectin: action, regulation and association to insulin sensitivity. *Obesity Reviews*. 2005 Feb;6(1):13-21
2. Tsao TS, Murry HE, Hug C et al.. Oligomerization state-dependent activation of NF-kappa B signaling pathway by adipocyte complement-related protein of 30 kDa (Acrp30). 2002 *J. Biol. Chem.* Aug;277(33): 29359-62
3. Combs TP, Pajvani UB, Berg AH et al. A transgenic mouse with a deletion in the collagenous domain of adiponectin displays elevated circulating adiponectin and improved insulin sensitivity. *Endocrinology*. 2004 Jan;145(1)::367-83
4. Fruebis J, Tsao TS, Jauorschi S et al. Proteolytic cleavage product of 30-kDa adipocyte complement-related protein increases fatty acid oxidation in muscle and causes weight loss in mice. 2001 *PNAS* Feb;98(4): 2005-10
5. Fasshauer M, Paschke R, Stumvoll M. Adiponectin, obesity, and cardiovascular disease. 2004 *Biochimie* Nov;86(11):779-86



## Summary Protocol

### MSD 96-well MULTI-ARRAY Mouse Adiponectin Assay Kit

MSD provides this summary protocol for your convenience.  
Please read the entire detailed protocol prior to performing the Mouse Adiponectin Assay.

#### Step 1 : Sample and Reagent Preparation

Bring all reagents to room temperature and thaw the Calibrator stock on ice.

Prepare Blocker A Solution.

Prepare serum or plasma samples.

Prepare an 8-point standard curve using supplied Calibrator:

- The Calibrator should be diluted in Diluent 100.
- Dilute the stock Calibrator 1:50 in Diluent 100 then perform a series of 5-fold dilution steps and a no Calibrator blank.
- Diluted Calibrators should be kept on ice until use.

**Note:** *The standard curve can be modified as necessary to meet specific assay requirements.*

Prepare Detection Antibody Solution by diluting the 100X Anti-mAdiponectin Antibody to 1X in 5.0 mL of Diluent 100 per plate.

Prepare 20 mL of 1X Read Buffer T by diluting 4X Read Buffer T with deionized water.

#### Step 2 : Add Blocker A Solution

Dispense 150  $\mu$ L/well Blocker A Solution.

Incubate at room temperature with vigorous shaking (300–1000 rpm) for 1 hour.

#### Step 3 : Wash and Add Detection Antibody Solution Followed by Sample or Calibrator

Wash plate 3 times with PBS-T.

Dispense 40  $\mu$ L/well 1X Detection Antibody Solution.

Immediately, dispense 10  $\mu$ L/well Calibrator or Sample.

- Samples should be diluted 1000-fold as described in the Reagent Preparation section.

Incubate at room temperature with vigorous shaking (300–1000 rpm) for 2 hours.

#### Step 4 : Wash and Read Plate

Wash plate 3 times with PBS-T.

Dispense 150  $\mu$ L/well 1X Read Buffer T.

Analyze plate on SECTOR instrument.



