

Mouse Adiponectin ELISA

Enzyme Immunoassay for quantitative Determination of
Mouse Adiponectin
English

For Research Use Only.
Not for use in diagnostic procedures.

distributed in the US/Canada by:
EAGLE BIOSCIENCES, INC.


20A NW Blvd, Suite 112 Nashua, NH 03063
Phone: 617-419-2019 FAX: 617-419-1110
www.EagleBio.com • info@eaglebio.com



REF **E091-M**



Gesellschaft für Forschung und Herstellung von Diagnostika GmbH

 : Aspenhastr. 25 • D-72770 Reutlingen / Germany
Telefon: + 49 - (0) 7121 51484-0 • Fax: + 49 - (0) 7121 51484-10
E-Mail: contact@mediagnost.de • <http://www.mediagnost.de>

Symbols/ Symbole /Symboles/ Simboli/ Símbolos/ Símbolos/ symbolen/ Symboler/ Symboler/Symbole/ Szimbólumok/ Symboly/ Symboly/ Символи/ Sümboldid/ Σύμβολα/ Simboluri/ Simboli/ Symbolit

according to DIN EN 980 and EDMA recommendations Standard News 6 2001



Expiry date/ Verfallsdatum/ Date de péremption/ Data di scadenza/ Fecha de caducidad/ Data de validade/ Uiterste gebruiksdatum/ Udløbsdato/ Bäst före-datum/ Termin ważności/ Lejárati idő/ Čas expirácie/ Doba expirace/ Срок на годност/ Αεγυμίσκωρπάεβ/ Ημερομηνία λήξης/ Data de expirare/ Rok uporabe/ Viimeinen käyttöpäivä



Consider instructions for use/ Bitte Gebrauchsanweisung beachten/ Consultez la notice d'utilisation/ Consultare le istruzioni per l'uso/ Consulte las instrucciones de uso/ Respeitar as instruções de utilização./ A.u.b de gebruiksaanwijzing volgen/ Se brugsanvisningen/ Läs anvisningarna före användning/ Proszę przeczytać instrukcję obsługi/ Vegye figyelembe a használati utasításban foglaltakat/ Postupujte podľa pokynov na použitie/ Dodržujte návod k použití/ Моля, спазвайте инструкцията за употреба/ Palun järgige kasutusjuhendit./ Λάβετε υπόψη σας τις οδηγίες χρήσης/ Vá rugám sã respectați instrucțiunile de utilizare/ Upoštevejate navodila za uporabo! Lue käyttöohje huolellisesti!



Lot-Batch Number/ Charge-Chargennummer/ Lot-Code du lot/ Lotto-Numero di lotto/ Lote-Código de lote/ Lote-Código do Lote/Lot-Partijnummer/ Lot-Batchkode/ Partisatskod/ Numer serii/ Tétel-sarzs szám/ Číslo šarže/ Číslo šarže/ Партиден номер/Partii – partii number/ Παρτίδα-αριθμός παρτίδας/ Lot-număr lot/ Številka serije/ erä



Manufactured by/ Hergestellt von/Fabriqué par/ Prodotto da/Fabricado por/ Fabricado por/ Vervaardigd doo/Fabrikation af /Tillverkad av/ Wyprodukowane przez/ Gyártotta / Vyrobené/ Vyrobeno v/ Производител/ Tootja/Κατασκευάζεται από/ Produs de/Proizvajalec/ Valmistaja



Catalogue Number/ Bestellnummer/ Numéro de référence/Numero di riferimento/Número de referencia/ Número de Referência/ Referentienummer/Referencenummer /Beställningsnummer/ Numer katalogowy/ Rendelési szám/Katalógovné číslo/ Objednací číslo/Καταλογος αριθμός/Tellimisnumber/ Αρ. παραγγελίας/Număr comandă/ Številka naročila/ viite tai tilausnumero



Store at between/ Lagerung bei zwischen/ Conserver à entre/ Conservare a tra/ Conservar a temp. Entre/ Armazenar entre/ Bewaar bij tussen/ Opbevares mellem/ Förvaras vid/ Przechowywać w/ Tárolási tartomány/ Skladujte v rozsahu / Skladujte v rozmezi/ Температурно ограничение/ Säilittä temperaturauidel/ Φύλαξη σε θερμοκρασία/ Depozitare între/ Skladiščenje med/ Säilytys x-y Celsiusasteen lämpötilassa



Contains sufficient for x tests/ Inhalt ausreichend für x Tests/ Contient suffisant pour x tests/ Contenido suficiente per x test/ Contenido suficiente para x pruebas/ Conteúdo suficiente para x testes/ Bevat voldoende voor x bepalingen/ Indeholder tilstrækkeligt til x prøver/Innehållet räcker till x analyser/ Zawartość na x testów/ Tartalma x teszt elvégzésére elegendő/ Obsahuje materiál pre x testov / Obsah dostačuje pro x testů/ Съдържание достатъчно за x тестове/ Sisust jätkub x katse jaoks/ Το περιεχόμενο επαρκεί για x δοκιμές/ Conținut suficient pentru x teste/Vsebina zadostuje za x preizkusov/ Sisältö riittää x testille



Keep away from sunlight/ Nicht dem Sonnenlicht aussetzen/ Conserver à l'abri de la lumière/ Conservare al riparo della luce solare/ No exponer a la luz solar/ Proteger da luz solar/ Niet aan zonlicht blootstellen/ Må ikke udsættes for sollys/ Utsatt inte för solljus/ Nie wystawiać na słońce/ Napfénytől távol tartandó/ Nevystavovat slnečnému svetlu/ Nevystavovat slunečnému světlu/ Да се предпазва от слънчева светлина/ Kaitsta otsese päikesekiirguse eest/ Κρατήστε το μακριά από την ηλιακή ακτινοβολία/ Τηνετή departe de lumina soarelui/ Ne izpostavljajte sončni svetlobi/ suojaa auringonvalolta



Incubation time/ Inkubationszeit/ Temps d'incubation/ Tempo d'incubazione/ Tiempo de incubación/ Tempo de incubação/ incubatietijd/Inkubationstid/ inkubationstid/ Czas inkubacji/ Inkubációs idő/ Inkubačná lehota/ Inkubační doba/ Инкубационен период/ Inkubatsiooniaeg/ Χρόνος επώασης/ Timp de incubare/Inkubacijska doba/ inkubaatioaika



incubate at / Inkubation bei/ Incuber à/Incubare a/incubar a/Incubar a/incubatietemperatuur/Inkubation ved/inkubation vid/Inkubacja przy/Inkubáció hőmérséklete/Inkubácia pri/Inkubace při/Инкубира се при/Inkubatsioon temperatuuril/Επώαση στους/Incubare la/Inkubacija pri/ inkubaatiolämpötila



Shaking/ Schütteln/ Mélanger/ Agitare/ Agitar/ Agitação/ Schudden/ Ryster/ Skaka/ Wstrząsanie/ Rázás/ Pretrepat/ Protřepat/ Разклатяване/ Raputada/ Ανακινήστε/ Vibrare/ Stresite/ Sekoita



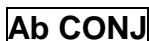
Mikrotiterplate/ Mikrotiterplatte/ plaque de microtitrage/ Piastra di microtitolazione/ Placa de microtitulación/ Placa de Microtitulação/ microtiterplaat/ Mikrotiterplade/ mikrotiterplatta/ microtiterplaat/ Płytko microtiter/ Mikrotiter lap/ Mikrotitračná podložka/ Mikrotitrační podložka/ Микротитърна плака/ Mikrotiiterplaat/ Τρυβλίο μικροπιλοδότησης/ Microplacă/ Mikrotitrská plošča/ Mikrotitruslevy



Reconstitute in/ Rekonstituieren in/ Reconstituer dans/ Ricostituire nel/ Reconstituir en/ Reconstituir em/ reconstituieren in/ Rekonstituér i/ rekonstituera/ Rekonstituować w/ Helyreállítás/ Znovu připravit za/ Znovu pripravít za/ Разтваряне в/ Moodustada uuesti / Ανασυστήστε σε/ Reconstituire în/ Predelava v /rekonstituoi



Sample/ Probe /Echantillon/ campione/ Muestra/ Amostra/ monster/ Prøve/ prov/ Próbká/ Minta/ Vzorka/ Vzorek/ Проба/ Proov/ Δείγμα/ Probă/Vzorec/Näyte



Antibody and Enzyme Conjugate/ Antikörper und Enzym Konjugat/ anticorps conjugué et conjugué enzymatique/ Coniugato di anticorpo ed enzima/ Conjugado de anticuerpos y enzimas/ Conjugado Anticorpo-Enzima/ antilichaam- en enzymconjugaat/ Antistoffer og enzym-konjugat/ antikropps- och enzymkonjugat (antikropp och enzym, konjugat)/ Koniugat antyciał i enzymów/ Antitest és enzim páros/ Protílátkový a enzymatický konjugát/ Protílátkový a enzymatický konjugát/ Антицяло и ензим конюгат/ Antikehad ja ensüümi konjugaat/ Σύμπλοκο αντισώματος-ενζύμου/ Compuși din anticorpi și enzime/ Antitelesa in konjugat encima/ Vasta-aine ja entsyymi konjugaati



Dilute in Buffer X/ Verdünnen in Puffer X/ Diluer dans le tampon X/ Diluire nel tampone X/ Diluir en tampón X/ Diluir no Tampão X/ verdunnen in buffer X/ Fortyndes i buffer X/ späd i buffert X/ Rozcieńczenie w buforze X/ Hígítás X pufferben/ Riedit' v pufrí X/ Ředit v pufru X/ Разреждане в буфер X/ Lahjendada puhvis X/ Αραιώστε σε ρυθμιστικό διάλυμα X/ Diluați în tamponul X/ Razredčiti v pufru X/ laimennetaan x puskuriin



Standard X /Standard X/ Etalon X/ Standard X/ Estándar X/ Standard X/ standaard X/ Standard X/ standard X/ Standard X/ Standard X/ Štandard X/ Standard X/ Стандарт X/ Standard X/ Πρότυπο X/ Standard X/ Standardni X/ Kalibraattori X

Control	KS1/ KS2	Control Serum / Kontrollserum/ Contôle sérique/ Siero di controllo/ Suero de control/ Soro de Controlo/ controleserum/ Kontrollserum/ Kontrollserum/ Serum kontrolne/ Ellenőrző szérum/ Kontrolné sérum/ Kontrolni sérum/ Контролен серум/ Kontrollseerum/ Ορός ελέγχου/Ser de control/ Kontrolni serum/ Kontrolli seerumi
WASHBUF 20x	WP	Washing Buffer Concentrate/ Waschpufferkonzentrat/ Tampon de lavage conc./ Tampone di lavaggio concentrato/ Tampón de lavado concentrado/ Tampão de Lavagem Concentrado/ wasbuffer, geconcentreerd/ Vaskebufferkoncentrat/ Vaskebufferkoncentrat/ tvättbuffertkoncentrat/ Bufor płukania koncentrat/ Mosópufer koncentrátum/ Koncentrát γυμνacieho pufra/ Концентрат на промивен буфер/ Pesupuhvri kontsentraat/ Συμπύκνωμα ρυθμιστικού διαλύματος έκπλυσης/ Concentrat pentru tamponul de spălare/ Koncentrat izpiralnega pufra/ Pesuliuositiiviste
WASHBUF		Washing Buffer / Waschpuffer/ Tampon de lavage/ Tampone di lavaggio/ Tampón de lavado/Tampão de Lavagem/ wasbuffer/ Vaskebuffer/ tvättbuffer/ Bufor płukania/ Mosópufer/ Γυμνáci puffer/ Γυμνáci pufri/ Промивен буфер/ Pesupuhver/ Ρυθμιστικό διάλυμα έκπλυσης/ Tampon pentru spălare /Izpiralni puffer/ Pesuliuos
SUBST TMB	S	Substrate/ Substrat/ Substrat/ Substrato/ Substrato/ Substrato/ substraat/ Substrat/ Substrat/ Substrat/ Szubsztrátum/ Substrát/ Substrát/ Субстрат Substraat/ Υπόστρωμα/ Substrat/ Substrat/ Substraattiliuos
H₂SO₄	SL	Stop Solution/ Stopp Lösung/ Stop Solution/ Soluzione di stop/ Stop Solución/ Solução Stop/ stopoplossing/ Stopopløsning/ Stopplösning/ Stop roztwór/ Megállító oldat/ Roztok na ukončenie/ Roztok pro ukončení/ Стопират разтвор/ Stopp-lahus/ Διάλυμα διακοπής/ Soluție de oprire/ Stop raztopina/ Pysäytysliuos
TAPE		Cover Plate with sealing tape /Platte abkleben/ Recouvrir la microplaque avec bande adhésive/ Coprire la piastra con nastro adesivo/ Cubrir la placa con una cinta adhesiva/ Cobrir a Placa com fita adesiva/ plaatje met tape afdekken/ Afdækningsplade med tape/ maskera platta/ Odkleič plynke/ Tányér leragasztása/ Oblepit' podložku lepiacou páskou/ Olepit podložku lepící páskou/ Плака с лента за запечатване/ Katta plaat isoleerklēerlindīga/ Κολληστε το πλακίδιο με κολλητική ταινία/ Αορεγίτι placa cu o bandă adezivă/ Prelepiti ploščo/ Peitä mikrotitrauslevy oheisella teipillä
MEASURE		Measure plate within 30 min at 450 nm (Referencefilter ≥590nm)/Ausmessung innerhalb von 30 min bei 450 nm (Referenzfilter ≥ 590 nm)./ Mesure lábsorbance en l'éspace de 30 min à450 nm avec ≥590nm longueur d'onde pour référence/Misurazione entro 30 min. a 450 nm (filtro di riferimento ≥ 590 nm)./ Medición de la placa dentro de los siguientes 30 min a 450 nm (filtro de referencia ≥ 590nm)/ Medir a placa dentro de 30 min a 450 nm (Filtro de referência ≥590nm)/ Binnen 30 minuten bij 450 nm meten (referentiefilter ≥ 590 nm)./ Mål plade i løbet af 30 min ved 450 nm (referencefilter ≥590nm)/ Mät inom 30 min vid 450 nm (referensfilter ≥ 590 nm)./ Pomiar w ciągu 30 min przy 450 nm (filtr odniesienia ≥ 590 nm)./ Ki mérés 30 percen belül 450 nm-nél (referenciaszűrő ≥ 590 nm)./ Merať 30 minút pri 450 nm/Měřit 30 minut při 450 nm/ Отчитане в рамките на 30 min при 450 nm (референтен филтър ≥ 590 nm)./ Mõõtmise 30 min jooksul 450 nm korral (võrdlusfilter ≥ 590 nm). Μέτρηση εντός 30 min στα 450 nm (φίλτρο αναφοράς ≥ 590 nm)./ Măsurare în decurs de 30 min la 450 nm (filtru de referință ≥ 590 nm)./ Izmerite ploščico v 30 min pri 450 nm (referenčni filter ≥590nm) / Mittaa 30 minuutin aikana 450 nm:ssä (referenssi suodatin ≥ 590 nm)
Literatur		Literature/ Literatur/ Bibliographie/ Letterario/ Bibliografia/ Literatura documentação/ literatuur/ Litteratur/ litteratur/ Literatura/ Irodalom/ Literatúra/ Literatura/ Литература// Kirjandus/ Βιβλιογραφία/ Bibliografie/ literatura/ Lähdeluettelo
International Test description		International test description/ internationale Testanleitung/ description internationale de test/ Istruzioni per il test internazionali/ Descripción de ensayo internacional/ Descrição internacional do teste/ internationale testbeschrijving/ internationell testbeskrivning/ Opis testu międzynarodowego/ nemzetközi teszt-útmutató/ Medzinárodný návod k testu/ Mezinárodní návod k testu/ rahvusvaheline katse kirjeldus/ Διεθνείς οδηγίες για εργαστηριακές δοκιμές/ instrucțiuni internaționale pentru testare/ mednarodna navodila za preizkus/ Kansainvälinen käyttöohje
End		in all required wells/ in allen benötigten Vertiefungen/ dans tous les godets requis/ in tutti i pozzetti richiesti/ en todos los pozos requeridos/ em todos os tubos necessários/ in alle nodige putjes/ i alle nødvendige brønde/ i alla nödvändiga brunnar/ we wszystkich potrzebnych wgłębieniach/ minden szükséges forrásban/ vo všetkých potrebných miestach/ ve všech potřebných místech/ във всички необходими ямки/ kõigis vajalikes süvendites/ σε όλες τις απαραίτητες κοιλότητες/ în toate cavitățile necesare/ v vseh zahtevanih vdolbinah/kaikkiin tarvittaviin mikrotitrauslevyn syvennyksiin

For in vitro use only!

For Research Use Only!

For professional use only!

CAUTION: Not for human or animal therapeutic or diagnostic use.

Read entire protocol before use!

Symbols	2
Package Insert English	
Table of Content	
Intended Use	5
Introduction	5
Reagents Provided	6
Materials Required but not provided	6
WARNINGS AND PRECAUTIONS	7
Method	8
Specimen	8
Storage of the samples	8
Sample Preparation	8
Technical Recommendations	8
Assay Procedure	10
Calculation of Results	11
Performance Characteristics	12
Literatur / Literature	13
SUMMARY – MEDIAGNOST Mouse Adiponectin ELISA E091-M	14

PACKAGE INSERT ENGLISH

Mediagnost Mouse Adiponectin ELISA E091-M

- is suited for Adiponectin determination in **Mouse Serum** and **Plasma** samples,
- is **fast**: incubation time a total of 2 hours and 30 minutes
- Single Standards with **0.025, 0.075, 0.15, 0.3, 0.65 and 1 ng/ml** native Mouse-Adiponectin are provided in the Kit
- 2 Control Sera are provided for quality control
- uses **high affinity monoclonal antibodies** against mouse Adiponectin
- Microtiter plates are separately breakapart

Intended Use

Measurement of Adiponectin in Mouse serum and plasma samples.

Introduction

Adiponectin was described for the first time in the early 90th of the last century as an endocrine factor produced by adipocytes. Adiponectin is involved in regulation of energy- and fat metabolism. So its concentration in the circulation is said to reflect the risk of atherosclerosis and the degree of insulin resistance. Based on the high incidence of these diseases, adiponectin was and still is object of intensive research regarding the underlying biological mechanisms and regarding its value as biomarker. Beside different cell culture models and studies with human patients, mice and rats are suitable model organisms for basic research and pre-clinical studies.

Therefore we developed and validated this testsystem as a tool for adiponectin measurements in mice usable in research and pre-clinical studies.

Even if the comparability of mice and humans is limited we offer some background information on *human* adiponectin physiology in the following section:

Adiponectin is a 30kDa protein which percentage in serum proteins is 0.01%. It is mainly synthesized by Adipocytes, but also muscle cells and hepatocytes have the ability to synthesize Adiponectin. Until now, IGF-I is the only known natural inductor of the synthesis. It consists of a Collagen-like N-terminal and a globular C-terminal domain [1]. In vivo Adiponectin appears with different oligomers. Beside the trimer and dimer also high molecular multimers exist [1-3]. Up to now two different receptors are known, both receptors are ubiquitarily expressed, though the distribution in the tissues varies. The Adiponectin Receptor 1 (AdipoR1) is especially in muscle- and AdipoR2 in liver tissue synthesized [4].

The significance for the organism is not clear until now. First studies show, that adiponectin correlates negatively with BMI and thus it could have relevance for the energy metabolism for example through the regulation of fatty acid oxidation. Beside the correlation with BMI, Adiponectin level is associated with the Insulin-Resistance [5-7] and so also linked with Type II Diabetes. Adiponectin is associated also with glucose- und lipometabolism [8, 9].

The formerly proposed diagnostic value of the high molecular weight form of adiponectin was not verified using a commercially available testsystem for the determination of HMW adiponectin [10]. Blueher et al. clearly demonstrate that regarding the diagnosis of insulin resistance, measured by whole body glucose uptake below 40 $\mu\text{mol}/\text{kg}\cdot\text{min}$, total adiponectin as determined with the Mediagnost E09, is with an area of 0.92 under the receiver-operating curve, of greater diagnostic value [10].

Furthermore it is involved in inflammatory processes [11-15] and therewith it is of importance for appearance of arteriosclerosis [4, 5, 16] and coronary disease [17, 18], thus the determination of Adiponectin level in plasma could serve to estimate the risk of coronary disease [19, 20]. Beside this Adiponectin influences further physiological processes as for example the angiogenesis [21, 22].

Reagents Provided

1)	MTP	Microtiter plate , ready for use: Microtiter plate with 96 wells, dived up in 12 stripes à 8 wells (separately breakapart), coated with anti-Maus Adiponectin antibody.
2)	CAL	Standards A-F , lyophilised, contain native Mouse Adiponectin. Standard values are between 0.025 - 1 ng/ml (0.025, 0.075, 0.15, 0.3, 0.65 und 1 ng/ml) Adiponectin and have to be reconstituted in 1 ml (each) Dilution Buffer VP . 100 µl per well are used in the assay. If the standards are required for more than one assay process we recommend to store the reconstituted Standards frozen at -20°C. Attention: Standards should be thawed only once – where required please store aliquoted in adequate volumes.
3)	DILU	Dilution buffer VP , 125 ml , ready for use, please use for the reconstitution of Standards A-F, Control Sera KS1 & KS2 and for the serum dilution.
4)	Control	Control Serum KS1 & KS2 , each 250 µl lyophilised: Contains mouse Serum and has to be reconstituted in 250 µl Dilution Buffer VP . The Adiponectin target value concentration and the respective range is given the vial label. The dilution of the Control Sera KS 1&2 should be according to the dilution of the respective samples, the target value concentration should be obtained by multiplication with the respective dilution factor .
5)	Ab CONJ	Antibody-POD-Conjugate AK , 12 ml , ready for use , contains a mixture of biotinylated anti-Adiponectin antibody and HRP (Horseradish Peroxidase)-labelled Streptavidin. Use 100 µl/well in the assay.
6)	WASHBUF 20x	Washing Buffer (WP) , 50 ml , 20 X concentrated solution. Washing Buffer (WP) has to be diluted 1:20 with distilled or demineralised water before use (e.g. add the complete contents of the flask (50 ml) into a graduated flask and fill up with A.dest. to 1000 ml). Attention: After dilution the Washing Buffer is only 4 weeks stable, dilute only according to requirements.
7)	SUBST	Substrate (S) , 12 ml , ready for use, horseradish-peroxidase-(HRP)-substrate, stabilised H ₂ O ₂ Tetramethylbencidine.
8)	H ₂ SO ₄	Stopping Solution (SL) , 12 ml , ready for use, 0.2 M sulphuric acid, Caution acid!
9)		Sealing tape for covering of the microtiter plate, 2 x, adhesive.

Materials Required but not provided

Precision pipettes and multichannel pipettes with disposable plastic tips
 Distilled or deionized water for dilution of the Washing Buffer (WP)
 Vortex-mixer
 Microtiterplate shaker (350 rpm)
 Microtiterplate washer (recommended)
 Micro plate reader ("ELISA-Reader") with filter for 450 and ≥590 nm
 Polyethylen PE/Polypropylen PP tubes for dilution of samples

WARNINGS AND PRECAUTIONS

For research use only. For professional use only.

Before starting the assay, read the instructions completely and carefully. Use the valid version of the package insert provided with the kit. Be sure that everything is understood.

Before use, all kit components should be brought **to room temperature at 20 - 25°C**. Precipitates in buffers should be dissolved before use by thorough mixing and warming.

Do not mix reagents of different lots. Do not use expired reagents.

The microplate contains snap-off strips. Unused wells must be stored at 2 - 8°C in the sealed foil pouch and used in the frame provided.

Caution: This kit contains material of human and/or animal origin.

No known test methods can offer total assurance of the absence of infectious agents; therefore all components and specimens should be treated as potentially infectious.

Following components contain < 0.01% 2-Methyl-4-isothiazolin-3-one solution as preservative : **A-F, AK, VP**

- R34 Irritating to eyes and skin
- R43 Sensibilisation through skin contact possible
- S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice
- S36/37 Wear suitable protective clothing and gloves
- S45 In case of accident or if you feel unwell seek medical advice

Following components contain < 0.01%(w/w) 5-chloro-2-methyl 2H isothiazol-3-one and 2-methyl-2H-isothiazol-3-one as preservative: **A-F, AK, VP, WP**

- R36/38 Irritating to eyes and skin
- R43 Sensibilisation through skin contact possible
- S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice
- S28.1 After contact with skin, wash immediately with plenty of water

Stop solution contains 0.2 M Sulfuric Acid (H₂SO₄)

- R36/38 Irritating to eyes and skin
- S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice
- S28.1 After contact with skin, wash immediately with plenty of water
- S36/37 Wear suitable protective clothing and gloves.

Pipetting of samples and reagents must be done as quickly as possible and in the same sequence for each step. Use separate pipette tips for each sample, control and reagent to avoid cross contamination. Use reservoirs only for single reagents. This especially applies to the substrate reservoirs. Using a reservoir for dispensing a substrate solution that had previously been used for the conjugate solution may turn solution colored. Do not pour reagents back into vials as reagent contamination may occur. Mix the contents of the microplate wells thoroughly to ensure good test results. Do not reuse microwells. Do not let wells dry during assay; add reagents immediately after completing the rinsing steps.

TMB-Substrate (S) contains 3,3',5,5' Tetramethylbenzidine. Store and Incubate in the dark.

- R20/21/R22 Harmful by inhalation, in contact with skin and if swallowed
- R36/37/38 Irritating to eyes, respiratory system and skin
- S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice
- S28.1 After contact with skin, wash immediately with plenty of water
- S36/37 Wear suitable protective clothing and gloves

General first aid procedures:

Skin contact: Wash affected area thoroughly with water. Discard contaminated cloths and shoes.

Eye contact: In case of contact with eyes, rinse immediately with plenty of water at least 15 minutes. In order to assure an effectual rinsing spread the eyelids.

Ingestion: If swallowed, wash out mouth thoroughly with water. Immediately see a physician.

Do not eat, drink or smoke in these areas.

Never pipette the materials with the mouth.

Spilled material must be wiped off immediately and should become disinfected. Clean contaminated areas and equipment with a suitable detergent.

Method

The Mediagnost ELISA for Adiponectin E091-M is a so-called Sandwich-Assay using two specific and high affinity antibodies. The Adiponectin in the samples binds to the first antibody coated on the microtiter plate. In the following step the second specific anti-Adiponectin-Antibody binds in turn to the immobilised Adiponectin. The second antibody is biotinylated and will be applied in a mixture with a Streptavidin-Peroxidase-Enzyme Conjugate. In the closing substrate reaction the turn of the colour will be catalysed quantitatively depending on the Adiponectin-level of the samples.

Specimen

Serum and plasma samples of mice and mouse cell culture medium can be used in this assay.

Influence of Heparin (30 IE/mL), EDTA (6,8 mM) and NaCitrat (0,015 M) on the measurement of Adiponectin has been investigated by recovery experiments. PBS was enriched with recombinant mouse Adiponectin and the above mentioned substances. No significant influence on the recovery of adiponectin was detected.

Hemolytic reactions have to be avoided. The blood has to be allowed to clot and after complete clotting, serum is separated by centrifugation.

Storage of the samples

Storage at RT max. 2 days

Storage at -20°C max. 2 years

More than five freeze/thaw cycles are not possible.

Sample Preparation

Samples have to be diluted in Dilution Buffer (VP).

A sample dilution of 1:10 000 is in general suitable. However, the Adiponectin levels can vary individually significantly, we would therefore recommend to check this and adjust the dilution respectively.

Technical Recommendations

The assay has to be conducted strictly according to the test protocol herein.

Reagents with different lot numbers cannot be mixed. The microtiterplate and reagents are stable until the indicated expiry if stored unopened and protected from sunlight at 2 – 8°C.

The shelf life of the components after opening is not affected, if used appropriately.

Bring all reagents to room temperature (20 - 25°C) before use. Possible precipitations in the buffers have to be resolved before usage by mixing and / or warming.

Incubation at room temperature means: Incubation at 20-25°C

The incubation steps should be performed at mean rotation frequency of a particularly suitable microtiter plate shaker. We are recommending 350 rpm. Due to certain technical differences deviations may occur, in case the rotation frequency must become adjusted. Insufficient shaking may lead to inadequate mixing of the solutions and thereby to low optical densities, high variations and/or false values, excessive shaking may result in high optical densities and/or false values.

Proper washing is of basic importance for a secure, reliable and precise performance of the test. Incomplete washing is common and will adversely affect the test outcome. Possible

consequences may be uncontrolled unspecific variations of measured optical densities, potentially leading to false results calculations of the examined samples. Effects like high background values or high variations may indicate washing problems.

All washing must be performed with the provided washing buffer diluted to usage concentration. Washing volume per washing cycle and well must be 300 µl at least.

The danger of handling with potentially infectious material must be taken into account.

When using an automatic microtiter plate washer, the respective instructions for use must be carefully followed. Device adjustments, e.g. for plate geometry and the provided washing parameters, must be performed. Dispensing and aspirating manifold must not scratch the inside well surface. Provisions must be made that the remaining fluid volume of every aspiration step is minimized. Following the last aspiration step of each washing cycle, this could be controlled, and possible remaining fluid could then be removed, by inverting the plate and repeatedly tapping it dry on non fuzzy absorbent tissue.

Manual washing is an adequate alternative option. Washing Buffer may be dispensed via a multistepper device, a multichannel pipette, or a squirt bottle. The fluid may be removed by dynamically swinging out the microtiter plate over a basin. If aspirating devices are used, care has to be taken that the inside well surface is not scratched. Subsequent to every single washing step, the remaining fluid should be removed by inverting the plate and repeatedly tapping it dry on non fuzzy absorbent tissue.

Standards and Controls

For the reconstitution of the lyophilised components (Standards A - F and Control Sera KS1 &KS2) the kit Dilution Buffer VP has to be used. It is recommended to keep reconstituted reagents at room temperature for 15 minutes and then to mix them thoroughly but gently (no foam!) with a Vortex mixer.

The reconstituted standard and controls can be stored for 2 month at –20°C. Repeated freeze/thaw cycles have to be avoided.

Washing Buffer

The required volume of washing buffer is prepared by 1:20 dilution of the provided 20fold concentrate with deionised water. The diluted Washing Buffer is stable for 4 weeks at 2-8°C. It has to be at room temperature for usage!

Microtiter plate

Unused microtiter plate stripes have to be stored airtight together with the desiccant bag at 2-8°C. The labelled expiry is not influenced in case of proper storage.

Assay Procedure

All determinations (Standards, Control Sera KS1 & KS2 and samples) should be assayed in duplicate. For optimal results, accurate pipetting and adherence to the protocol are recommended.

When performing the assay, the Standards, Control Sera and the samples should be pipetted as fast as possible (e.g., <15 minutes).

All incubations have to be conducted at room temperature (20-25°C)

To avoid distortions due to differences in incubation times, Antibody-POD-Conjugate AK as well as the following Substrate Solution S should be added to the plate in the same order and in the same time interval as the samples. Stop Solution SL should be added to the plate in the same order as the Substrate Solution.

- 1) Add **100 µl Dilution Buffer VP** in the first wells. Subsequently add 100 µl Standard or 100 µl of diluted Control Sera or diluted samples.
- 2) Cover the wells with sealing tape and incubate the plate for **1 hour at room temperature** (shake at 350 rpm)
- 3) After incubation aspirate the contents of the wells and wash the wells **5 times 300 µl Washing Buffer WP** / well. The washing buffer should incubate for at least for 15 seconds/cycle.
- 4) Following the last washing step pipette **100 µl** of the **Antibody-POD-Conjugate AK** in each well.
- 5) Cover the wells with sealing tape and incubate the plate for **1 hour at room temperature** (shake at 350 rpm).
- 6) After incubation wash the wells **5 times** with Washing Buffer as described in step 3
- 7) Pipette **100 µl of the TMB Substrate** Solution in each well.
- 8) Incubate the plate for **30 minutes in the dark at room temperature (20 - 25°C)**.
- 9) Stop the reaction by adding **100 µl of Stopping Solution**.
- 10) Measure the colour reaction within 30 minutes at **450 nm (reference filter ≥590 nm)**.

Calculation of Results

Establishing the Standard Curve

For the evaluation of the assay it is preconditioned that the absorbance values of the blank should be below 0.25, these of standard F should be above 1.0.

Samples, which yield higher absorbance values than Standard F are beyond the standard curve, for reliable determinations these samples should be tested anew with a higher dilution.

Standards are provided in the following concentrations (use the concentration unit as preferred):

Standard	A	B	C	D	E	F
ng/ml	0.025	0.075	0.15	0.3	0.65	1

- 1) Calculate the **mean absorbance** value for the blank from the duplicated determination (well A1/A2).
- 2) Subtract the mean absorbance of the blank from the mean absorbances of all other values.
- 3) Plot the standard concentrations on the x-axis versus the mean value of the absorbance of the standards on the y-axis.
- 4) Recommendation: Calculation of the standard curve should be done by using a computer program. **A higher-grade polynomial, or four parametric logistic (4-PL) curve fit or non-linear regression** are usually suitable for the evaluation (as might be spline or point-to-point alignment in individual cases).
- 5) The Adiponectin concentration of the diluted sample or the diluted control sera KS1&2 in ng/ml (or µg/ml according the chosen unit for the standards) is calculated in this way, the Adiponectin concentration of the **undiluted sample** and of KS1 & KS2 is calculated **by multiplication** with the respective dilution factor.

Performance Characteristics

Standards

The Standards of the ELISA E09 are prepared from **native Adiponectin** in concentrations of **0.025, 0.075, 0.15, 0.3, 0.65 and 1 ng/ml**. The native Adiponectin was calibrated against the recombinant Protein (Manufacturer: R&D Systems, Wiesbaden).

Sensitivity

The analytical sensitivity of the ELISA E091-M yields 0.008 ng/mL (equal to < 0.0008 ng per well; as 2xSD of zero standard in 22fold determination).

Interassay Variability

	Sample 1	Sample 2	Sample 3	Sample 4	Sample 5	Sample 6	Sample 7
Mean	1.12	6.12	7.65	8.04	2.15	7.54	6.87
SD	0.09	1.19	0.60	0.56	0.25	0.55	0.42
VC%	7.66	3.79	7.80	6.98	11.80	7.28	6.08
n	8	8	9	9	9	9	6

Intraassay Variability

	Sample 1	Sample 2
Mean [µg/mL]	7.395	1.712
SD	159	37
VC%	2.15	2.16
n	21	21

Linearity

Linearity				
Dilution	Sample 1	Dilution	Sample 2	Sample 3
1:14000	11.673	1:10000	9.59	
1:28000	10.928	1:12000	11.229	11.886
1:56000	10.67	1:24000	9.856	10.822
1:112000	10.421	1:48000	9.915	10.222
		1:96000	9.684	9.999

recalculated Values in ng/mL

Literature

1. Nakano, Y., et al., *Isolation and characterization of GBP28, a novel gelatin-binding protein purified from human plasma*. J Biochem (Tokyo), 1996. **120**(4): p. 803-12.
2. Pajvani, U.B., et al., *Structure-function studies of the adipocyte-secreted hormone Acrp30/adiponectin. Implications for metabolic regulation and bioactivity*. J Biol Chem, 2003. **278**(11): p. 9073-85.
3. Tsao, T.S., et al., *Role of disulfide bonds in Acrp30/adiponectin structure and signaling specificity. Different oligomers activate different signal transduction pathways*. J Biol Chem, 2003. **278**(50): p. 50810-7.
4. Shimada, K., T. Miyazaki, and H. Daida, *Adiponectin and atherosclerotic disease*. Clin Chim Acta, 2004. **344**(1-2): p. 1-12.
5. Higashiura, K., et al., *Correlations of adiponectin level with insulin resistance and atherosclerosis in Japanese male populations*. Clin Endocrinol (Oxf), 2004. **61**(6): p. 753-9.
6. Spranger, J., et al., *Adiponectin is independently associated with insulin sensitivity in women with polycystic ovary syndrome*. Clin Endocrinol (Oxf), 2004. **61**(6): p. 738-46.
7. Zoico, E., et al., *Adipocytokines, fat distribution, and insulin resistance in elderly men and women*. J Gerontol A Biol Sci Med Sci, 2004. **59**(9): p. M935-9.
8. Ye, J.M., et al., *PPARAlpha /gamma ragaglitazar eliminates fatty liver and enhances insulin action in fat-fed rats in the absence of hepatomegaly*. Am J Physiol Endocrinol Metab, 2003. **284**(3): p. E531-40.
9. Yamauchi, T., et al., *Adiponectin stimulates glucose utilization and fatty-acid oxidation by activating AMP-activated protein kinase*. Nat Med, 2002. **8**(11): p. 1288-95.
10. Blüher (Blueher), M., et al., *Total and high-molecular weight adiponectin in relation to metabolic variables at baseline and in response to an exercise treatment program: comparative evaluation of three assays*. Diabetes Care, 2007. **30**(2): p. 280-5.
11. Delaigle, A.M., et al., *Induction of adiponectin in skeletal muscle by inflammatory cytokines: in vivo and in vitro studies*. Endocrinology, 2004. **145**(12): p. 5589-97.
12. Winzer, C., et al., *Plasma adiponectin, insulin sensitivity, and subclinical inflammation in women with prior gestational diabetes mellitus*. Diabetes Care, 2004. **27**(7): p. 1721-7.
13. Xydakis, A.M., et al., *Adiponectin, inflammation, and the expression of the metabolic syndrome in obese individuals: the impact of rapid weight loss through caloric restriction*. J Clin Endocrinol Metab, 2004. **89**(6): p. 2697-703.
14. Motoshima, H., et al., *Adiponectin suppresses proliferation and superoxide generation and enhances eNOS activity in endothelial cells treated with oxidized LDL*. Biochem Biophys Res Commun, 2004. **315**(2): p. 264-71.
15. Wolf, A.M., et al., *Adiponectin induces the anti-inflammatory cytokines IL-10 and IL-1RA in human leukocytes*. Biochem Biophys Res Commun, 2004. **323**(2): p. 630-5.
16. Okamoto, Y., et al., *Adiponectin reduces atherosclerosis in apolipoprotein E-deficient mice*. Circulation, 2002. **106**(22): p. 2767-70.
17. Schlegel, A., *Adiponectin and risk of coronary heart disease*. Jama, 2004. **292**(1): p. 40; author reply 40.
18. Choi, K.M., et al., *Inflammation, Insulin Resistance, and Glucose Intolerance in Acute Myocardial Infarction Patients without a Previous Diagnosis of Diabetes Mellitus*. J Clin Endocrinol Metab, 2004.
19. Nakamura, Y., et al., *Implications of plasma concentrations of adiponectin in patients with coronary artery disease*. Heart, 2004. **90**(5): p. 528-33.
20. Pischon, T., et al., *Plasma adiponectin levels and risk of myocardial infarction in men*. Jama, 2004. **291**(14): p. 1730-7.
21. Shibata, R., et al., *Adiponectin stimulates angiogenesis in response to tissue ischemia through stimulation of amp-activated protein kinase signaling*. J Biol Chem, 2004. **279**(27): p. 28670-4.
22. Fernandez-Real, J.M., et al., *Adiponectin is associated with vascular function independent of insulin sensitivity*. Diabetes Care, 2004. **27**(3): p. 739-45.

SUMMARY – MEDIAGNOST Mouse Adiponectin ELISA E091-M

Reagent preparation:	Reconstitution:	Dilution
Standards A-F	in 1 ml Dilution Buffer VP	
Control Sera KS1 & KS2	in 250 µl Dilution Buffer VP	1:10 000 with Dilution Buffer VP
Washing Buffer WP		1:20 with Aqua. dest. (e.g., add the complete contents of the flask (50 ml) into a graduated flask and fill with A.dest. to 1000 ml).
Sample Dilution: e.g. 1 :10 000		

Assay Procedure for Double Determination:

Pipette	Reagents	Position
100 µl	Dilution Buffer VP	A1/2
100 µl	Standard A (0.025 ng/ml)	B1/2
100 µl	Standard B (0.075 ng/ml)	C1/2
100 µl	Standard C (0.15 ng/ml)	D1/2
100 µl	Standard D (0.3 ng/ml)	E1/2
100 µl	Standard E (0.65 ng/ml)	F1/2
100 µl	Standard F (1 ng/ml)	G1/2
100 µl	Control Serum KS1	H1/2
100 µl	Control Serum KS2	A3/4
100 µl	Sample dilution	following wells
Cover the wells with the sealing tape.		
Incubation: 1 h at RT, 350 rpm		
5x 300 µl	Aspirate the contents of the wells and wash 5x with 300 µl Wash Buffer WP	each well
100 µl	Antibody-POD-Conjugate AK	each well
Incubation: 1 h at RT, 350 rpm		
5x 300 µl	Aspirate the contents of the wells and wash 5x with 300 µl Wash Buffer WP	each well
100 µl	Substrate Solution S	each well
Incubation: 30 min in the Dark at RT		
100 µl	Stopping Solution SL	each well
Measure the absorbance within 30 min at 450 nm with ≥590 nm as reference wavelength.		



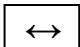
CAL A-F	A -F	Rec in 1 ml VP	
Control	KS1 & KS2	Rec in 250 µl VP	1:10 000 DILU VP
WASHBUF 20x	WP		1:20 DILU A. dest.

SPE	1:10 000 DILU VP
°C 20-25 °C	

100 µl	VP	A1/2
100 µl	CAL A (0.025 ng/ml)	B1/2
100 µl	CAL B (0.075 ng/ml)	C1/2
100 µl	CAL C (0.15 ng/ml)	D1/2
100 µl	CAL D (0.3 ng/ml)	E1/2
100 µl	CAL E (0.65 ng/ml)	F1/2
100 µl	CAL F (1 ng/ml)	G1/2
100 µl	CONTROL KS1 1:10 000 DILU VP	H1/2
100 µl	CONTROL KS2 1:10 000 DILU VP	A3/4
100 µl	SPE 1:10 000 DILU VP	
TAPE		

 1 h  20-25  350 rpm

5x 300 µl	5x WASHBUF WP
100 µl	AbCONJ AK
TAPE	

 1 h  20-25  350 rpm

5x 300 µl	5x WASHBUF WP
100 µl	SUBST TMB S

 0.5 h  20-25 

100 µl	H₂SO₄ SL
MEASURE	