



Product Manual

Adiponectin (human), ELISA kit

Product #: ALX-850-377

USE FOR RESEARCH PURPOSES ONLY

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

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

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

TRADEMARKS AND PATENTS

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

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



Please read entire booklet before proceeding with the assay.



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



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

Intended Use.....	2
Storage, Expiration	2
Introduction	2
Test Principle	3
Precautions	4
Technical Hints	5
Materials Supplied	5
Other Materials Needed.....	6
Preparation of Reagent.....	6
Preparation of Samples	8
Assay Procedure.....	9
Calculations	11
Typical Standard Curve	12
Performance Characteristics.....	13
Definition of the Standard	19
Preliminary Population and Clinical Data	19
Troubleshooting & FAQs.....	22
References.....	22
Explanation of Symbols	27
Assay Procedure Summary	28
Plate Layout.....	29
Contact Information.....	30

INTENDED USE

The Enzo Life Sciences Adiponectin (human), ELISA kit is a sandwich enzyme immunoassay for the quantitative measurement of human adiponectin.

Features

- The total assay time is less than 3 hours
- The kit measures total adiponectin in serum, plasma (EDTA, citrate, heparin), urine, cerebrospinal fluid (CSF)
- Assay format is 96 wells
- Quality Controls are human serum based. No animal sera are used
- Standards are recombinant adiponectin based
- Components of the kit are provided ready to use or concentrated

STORAGE, EXPIRATION

Store the complete kit at 2-8°C. Under these conditions, the kit is stable until the expiration date (see label on the box).

For stability of opened reagents see Performance Characteristics section.

INTRODUCTION

Adiponectin, also referred to as Acrp30, AdipoQ and GBP-28, is a 244 amino acid protein, the product of the *apM1* gene, which is physiologically active and specifically and highly expressed in adipose cells. The protein belongs to the soluble defense collagen superfamily; it has a collagen-like domain structurally homologous with collagen VIII and X and complement factor C1q-like globular domain. Adiponectin forms homotrimers, which are the building blocks for higher order complexes found circulating in serum. Together, these complexes make up approximately 0.01% of total serum protein. Adiponectin receptors AdipoR1 and AdipoR2 have been recently cloned; AdipoR1 is abundantly expressed in skeletal muscle, whereas AdipoR2 is predominantly expressed in the liver. Paradoxically, adipose tissue-expressed adiponectin levels are inversely related to the degree of adiposity. Adiponectin concentrations correlate negatively with glucose, insulin, triglyceride concentrations, liver fat content and body mass index and positively with high-density lipoprotein-cholesterol levels,

hepatic insulin sensitivity and insulin-stimulated glucose disposal. Adiponectin has been shown to increase insulin sensitivity and decrease plasma glucose by increasing tissue fat oxidation.

Clinical studies have shown that low adiponectin levels are associated with insulin resistance and precede the onset of type 2 diabetes. Diabetic patients have low levels of adiponectin and even lower levels of adiponectin were observed in patients with poorly controlled type 2 diabetes and in diabetic patients with coronary heart disease. Hypoadiponectinemia is also closely associated with the metabolic syndrome and with the hypertriglyceridemic waist. Non-alcoholic fatty liver disease is described as part of the metabolic syndrome and levels of adiponectin have inverse association with liver enzymes and fatty liver disease. The key finding is that low adiponectin serum levels predict type 2 diabetes independent of other risk factors.

Adiponectin also inhibits the inflammatory processes of atherosclerosis suppressing the expression of adhesion and cytokine molecules in vascular endothelial cells and macrophages, respectively. This adipokine plays a role as a scaffold of newly formed collagen in myocardial remodelling after ischaemic injury and also stimulates angiogenesis by promoting cross-talk between AMP-activated protein kinase and Akt signalling in endothelial cells. Low serum adiponectin levels are found in patients with coronary artery disease.

Moreover, high circulating levels of adiponectin are associated with decreased risk of myocardial infarction, independent of other factors.

Altogether, monitoring of adiponectin levels and monitoring of processes that affect its production or its receptors are promising targets for prevention and treatment of obesity, insulin resistance, hyperlipidemia and atherosclerosis.

Clinical application and areas of investigation:

- Energy metabolism and body weight regulation
- Metabolic syndrome
- Type 2 diabetes
- Coronary artery disease
- Atherosclerosis

TEST PRINCIPLE

In the Enzo Life Sciences Adiponectin (human), ELISA kit, standards, quality controls and samples are incubated in

microplate wells pre-coated with polyclonal anti-human adiponectin antibody. After 60 minutes incubation and washing, polyclonal anti-human adiponectin antibody, conjugated with horseradish peroxidase (HRP) is added to the wells and incubated for 60 minutes with captured adiponectin. Following another washing step, the remaining HRP conjugate is allowed to react with the substrate solution (TMB). The reaction is stopped by addition of acidic solution and absorbance of the resulting yellow product is measured. The absorbance is proportional to the concentration of adiponectin. A standard curve is constructed by plotting absorbance values against concentrations of standards, and concentrations of unknown samples are determined using this standard curve.

PRECAUTIONS

FOR PROFESSIONAL USE ONLY.



Handle
with care

- Wear gloves and laboratory coats when handling immunodiagnostic materials
- Do not drink, eat or smoke in the areas where immunodiagnostic materials are being handled
- This kit contains components of human origin. These materials were found non-reactive for HBsAg, HCV antibody and for HIV 1/2 antigen and antibody. However, these materials should be handled as potentially infectious, as no test can guarantee the complete absence of infectious agents
- Avoid contact with the acidic Stop Solution and Substrate Solution, which contains hydrogen peroxide and tetramethylbenzidine (TMB). Wear gloves and eye and clothing protection when handling these reagents. Stop and/or Substrate Solutions may cause skin/eyes irritation. In case of contact with the Stop Solution and the Substrate Solution wash skin/eyes thoroughly with water and seek medical attention, when necessary
- The materials must not be pipetted by mouth

TECHNICAL HINTS

- Reagents with different lot numbers should not be mixed
- Use thoroughly clean glassware
- Use deionized (distilled) water, stored in clean containers
- Avoid any contamination among samples and reagents. For this purpose, disposable tips should be used for each sample and reagent
- Substrate Solution should remain colourless until added to the plate. Keep Substrate Solution protected from light
- Stop Solution should remain colourless until added to the plate. The colour developed in the wells will turn from blue to yellow immediately after the addition of the Stop Solution. Wells that are green in colour indicate that the Stop Solution has not mixed thoroughly with the Substrate Solution
- Dispose of consumable materials and unused contents in accordance with applicable national regulatory requirements

MATERIALS SUPPLIED

Kit Components	State	Quantity
Antibody Coated Microtiter Strips	ready to use	96 wells
Conjugate Solution	ready to use	13ml
Set of Standards	ready to use	8 x 1ml
Quality Control HIGH	concentrated	0.1ml
Quality Control LOW	concentrated	0.1ml
Dilution Buffer Conc. (10x)	concentrated	20ml
Wash Solution Conc. (10x)	concentrated	100ml
Substrate Solution	ready to use	13ml
Stop Solution	ready to use	13ml
Product Data Sheet + Certificate of Analysis	-	1 pc

OTHER MATERIALS NEEDED

1. Deionized (distilled) water
2. Test tubes for diluting samples
3. Glassware (graduated cylinder and bottle) for Wash Solution (Dilution Buffer)
4. Precision pipettes to deliver 5 1000 µl with disposable tips
5. Multichannel pipette to deliver 100 µl with disposable tips
6. Absorbent material (e.g. paper towels) for blotting the microtiter plate after washing
7. Vortex mixer
8. Orbital microplate shaker capable of approximately 300 rpm
9. Microplate washer (optional). [Manual washing is possible but not preferable.]
10. Microplate reader with $450 \pm 10\text{nm}$ filter, preferably with reference wavelength 630nm (alternatively another one from the interval 550 - 650nm)
11. Software package facilitating data generation and analysis (optional)

PREPARATION OF REAGENTS

- All reagents need to be brought to room temperature prior to use
- Always prepare only the appropriate quantity of reagents for your test
- Do not use components after the expiration date marked on their label

Assay reagents supplied ready to use:

Antibody Coated Microtiter Strips

Stability and storage: Return the unused strips to the provided aluminium zip-sealed bag with desiccant and seal carefully. Remaining Microtiter Strips are stable 3 months when stored at 2-8°C and protected from the moisture.

Conjugate Solution

Substrate Solution

Stop Solution

Stability and storage: Opened reagents are stable 3 months when stored at 2-8°C.

Human Adiponectin Standards

The Standards are ready to use.

Stability and storage: Opened Standards are stable 3 months when stored at 2-8°C.

Assay reagents supplied concentrated:

Dilution Buffer Conc. (10x)

Dilute only required amount of Dilution Buffer Concentrate. Otherwise dilute all 20ml of Dilution Buffer Concentrate (10x) with 180ml of distilled water to prepare 200ml of Dilution Buffer (1x) for use of all-wells.

Stability and storage: The diluted Dilution Buffer is stable 1 week when stored at 2-8°C. Opened Dilution Buffer Concentrate (10x) is stable 3 months when stored at 2-8°C.

Quality Controls HIGH, LOW

Refer to the Certificate of Analysis for current Quality Control concentration!!!

Dilute Quality Control (HIGH and LOW) 10x with the Dilution Buffer just prior to the assay, e.g. 30µl of QC + 270µl of Dilution Buffer for duplicates. (Quality Controls are supplied diluted 30x). It means the final dilution is 300x and the concentration of Quality Control calculated from the standard curve must be multiplied by a dilution factor of 300.

Mix well (not to foam). Vortex is recommended. Beware of imprecision in pipetting.

Stability and storage: Opened Quality Controls are stable 3 months when stored at 2-8°C.

Do not store the diluted Quality Controls.

Note: Concentration of analyte in Quality Controls need not be anyhow associated with normal and/or pathological concentrations in serum or another body fluid. Quality Controls serve just for control that the kit works in accordance with PDS and CoA and that ELISA test was carried out properly.

Wash Solution Conc. (10x)

Dilute Wash Solution Concentrate (10x) ten-fold in distilled water to prepare a 1x working solution. Example: 100ml of Wash Solution Concentrate (10x) + 900ml of distilled water for use of all 96-wells.

Stability and storage: The diluted Wash Solution is stable 1 month when stored at 2-8°C. Opened Wash Solution Concentrate (10x) is stable 3 months when stored at 2-8°C.

PREPARATION OF SAMPLES

The kit measures adiponectin in serum, plasma (EDTA, citrate, heparin, urine, cerebrospinal fluid (CSF), but also in breast milk.

Samples should be assayed immediately after collection or should be stored at -20°C. Mix thoroughly thawed samples just prior to the assay and avoid repeated freeze/thaw cycles, which may cause erroneous results. Avoid using hemolyzed or lipemic samples.

Serum and plasma samples:

Dilute serum and plasma 300x with the Dilution Buffer prior to the assay in two steps:

Dilution A (10x):

Add 10µl of samples to 90µl of Dilution Buffer. **Mix well** (not to foam).

Dilution B (30x):

Add 10µl of Dilution A into 290µl of Dilution Buffer to prepare final dilution (300x). **Mix well** (not foam).

One step-dilution can be performed (add 5µl of samples to 1495µl of Dilution Buffer). Beware of imprecision in pipetting and mix the samples very thoroughly!

Breast milk, urine and cerebrospinal fluid (CSF) samples:

Dilute samples 3x with Dilution Buffer just prior to the assay, e.g. add 100µl of sample to 200µl of Dilution Buffer for duplicates. **Mix well** (not to foam).

Stability and storage:

Samples should be stored at -20°, or preferably at -70°C for long-term storage. Avoid repeated freeze/thaw cycles.

Stability of milk, urine and CSF samples have not been tested.

Do not store the diluted samples.

See performance characteristics section for stability of serum and plasma samples when stored at 2-8°C, effect of freezing/thawing and effect of sample matrix (serum/plasma) on the concentration of adiponectin.

***Note:** It is recommended to use a precision pipette and a careful technique to perform the dilution in order to get precise results.*

ASSAY PROCEDURE

Adiponectin levels are significantly lower (2-3 orders of magnitude) in breast milk, urine and CSF than in serum and plasma. Therefore, different protocols have to be used.

Protocol (a) for serum and plasma samples:

Sample dilution is 300x

Standard range is 5-100ng/ml (the Standards of 150ng/ml and/or 2ng/ml can be added optionally)

Incubation with Substrate Solution is 10 minutes

Protocol (b) for breast milk, urine and CSF:

Sample dilution 3x

Standard range 1-50ng/ml

Incubation with Substrate Solution is 25-30 minutes

The other assay procedure is same for both ELISA protocols.

1. Pipet 100µl of Standards (5-100ng/ml for serum and plasma samples, 1-50ng/ml for milk, urine and CSF samples), diluted Quality Controls, Dilution Buffer (=Blank) and diluted

samples, preferably in duplicates, into the appropriate wells. See Figure 1 for example of work sheet.

2. Incubate the plate at room temperature (ca. 25°C) for 1 hour, shaking at ca. 300 rpm on an orbital microplate shaker.
3. Wash the wells 3-times with Wash Solution (0.35ml per well). After final wash, invert and tap the plate strongly against paper towel.
4. Add 100µl of Conjugate Solution into each well.
5. Incubate the plate at room temperature (ca. 25°C) for 1 hour, shaking at ca. 300 rpm on an orbital microplate shaker.
6. Wash the wells 3-times with Wash Solution (0.35ml per well). After final wash, invert and tap the plate strongly against paper towel.
7. Add 100µl of Substrate Solution into each well. Avoid exposing the microtiter plate to direct sunlight. Covering the plate with e.g. aluminum foil is recommended.
8. Incubate the plate for 10 minutes (serum and plasma samples) or 25-30 minutes (milk, urine and CSF samples) at room temperature (20-30°C). The incubation time may be extended [up to 20 minutes for serum and plasma samples or up to 50 minutes for milk, urine and CSF samples] if the reaction temperature is below than 20°C. Do not shake the plate during the incubation.
9. Stop the color development by adding 100µl of Stop Solution.
10. Determine the absorbance of each well using a microplate reader set to 450nm, preferably with the reference wavelength set to 630nm (acceptable range: 550 - 650nm). Subtract readings at 630nm (550 - 650nm) from the readings at 450nm. The absorbance should be read within 5 minutes following step 9.

Note: *If some samples and standards have absorbances above the upper limit of your microplate reader, perform a second reading at 405nm. A new standard curve, constructed using the values measured at 405nm, is used to determine adiponectin concentration of off-scale standards and samples. The readings at 405nm should not replace the readings for samples that were “in range” at 450nm.*

Note 2: *Manual washing: Aspirate wells and pipet 0.35ml Wash Solution into each well. Aspirate wells and repeat twice. After final wash, invert and tap the plate strongly against paper towel. Make certain that Wash Solution has been removed entirely.*

	strip 1+2	strip 3+4	strip 5+6	strip 7+8	strip 9+10	strip 11+12
A	Standard 100	QC HIGH	Sample 7	Sample 15	Sample 23	Sample 31
B	Standard 50	QC LOW	Sample 8	Sample 16	Sample 24	Sample 32
C	Standard 20	Sample 1	Sample 9	Sample 17	Sample 25	Sample 33
D	Standard 10	Sample 2	Sample 10	Sample 18	Sample 26	Sample 34
E	Standard 5	Sample 3	Sample 11	Sample 19	Sample 27	Sample 35
F	Standard 2	Sample 4	Sample 12	Sample 20	Sample 28	Sample 36
G	Standard 1	Sample 5	Sample 13	Sample 21	Sample 29	Sample 37
H	Blank	Sample 6	Sample 14	Sample 22	Sample 30	Sample 38

Figure 1: Example of a work sheet.

CALCULATIONS

DETERMINATION OF ADIPONECTIN CONCENTRATIONS

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

1. Calculate the average Net OD for each standard and sample by subtracting the average blank OD from the average OD for each standard and sample.

$$\text{Average Net OD} = \text{Average OD} - \text{Average Blank OD}$$

2. Using linear graph paper, plot the average Net OD for each standard versus adiponectin concentration in each standard. Approximate a straight line through the points.
3. Interpolate the sample concentrations from the standard curve and multiply by the dilution factors for the final sample adiponectin concentrations.

Samples with concentrations outside of the standard curve range will need to be re-analyzed using a different dilution.

The measured concentration of Quality Controls calculated from the standard curve must be multiplied by a dilution factor of 300 and the measured concentration of samples calculated from the standard curve must be multiplied by their respective dilution factor, because Quality Controls and samples have been diluted prior to the assay, e.g. 13.5ng/ml (from standard curve) x 300 (dilution factor for serum and plasma samples) = 4.05 µg/ml

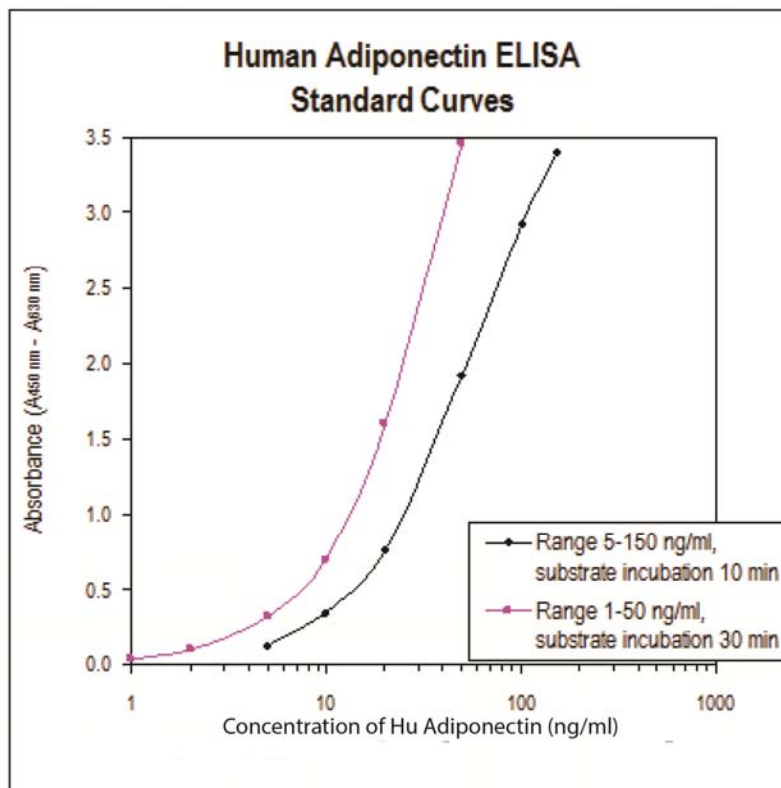


Figure 2: Typical Standard Curves for Adiponectin (human), ELISA kit.

PERFORMANCE CHARACTERISTICS

Typical analytical data of Enzo Life Sciences Adiponectin (human), ELISA kit are presented below.

Sensitivity

Limit of Detection (LOD) (defined as concentration of analyte giving absorbance higher than mean absorbance of blank* plus three standard deviations of the absorbance of blank: $A_{\text{blank}} + 3xSD_{\text{blank}}$) is calculated from the real adiponectin values in wells and is different for two protocols previously described:

For **Protocol (a)** (for serum and plasma samples) recommended sample dilution 300x, calibration range 5–150ng/ml, substrate incubation 10 min: LOD is 0.47ng/ml.

For **Protocol (b)** (for urine and CSF samples) recommended sample dilution 3x, calibration range 1-50ng/ml, substrate incubation 25-30 min: LOD is 0.156ng/ml.

*Dilution Buffer is pipetted into blank wells.

Limit of assay

Results exceeding the calibration range should be repeated with more diluted samples.

The samples with extremely high adiponectin levels can be diluted up to 2 400x. Dilution factor needs to be taken into consideration in calculating the adiponectin concentration.

Specificity

The antibodies used in this ELISA are specific for human adiponectin. The assay recognizes natural and recombinant (full length, mutation-modified trimer-only-forming, and globular domain) human adiponectin.

No cross-reactivity has been observed for human leptin, leptin receptor and resistin at 100ng/ml.

Determination of adiponectin does not interfere with hemoglobin (0.25 mg/ml), bilirubin (85 $\mu\text{mol/l}$) and triglycerides (2.5 mmol/l). Interference over 10% was measured at the higher concentrations.

Sera of several mammalian species were measured in the assay. See results below.

Mammalian serum sample	Observed cross reactivity
Bovine	no
Cat	yes
Dog	yes
Goat	no
Hamster	yes
Horse	no
Monkey	yes
Mouse	no
Pig	no
Rabbit	no
Rat	yes
Sheep	no

Presented results are multiplied by respective dilution factor

For details please contact us at info@enzolifesciences.com.

Precision

Intra-assay (Within-Run) (n=8)

Sample	Mean ($\mu\text{g/ml}$)	SD ($\mu\text{g/ml}$)	CV (%)
1	6.34	0.28	4.4
2	9.41	0.31	3.3

Inter-assay (Run-to-Run) (n=9)

Sample	Mean ($\mu\text{g/ml}$)	SD ($\mu\text{g/ml}$)	CV (%)
1	9.41	0.54	5.8
2	17.74	1.11	6.2

Spiking Recovery

Samples were spiked with different amounts of human adiponectin and assayed.

For **protocol (a)** with **serum** samples:

Sample	Observed ($\mu\text{g/ml}$)	Expected ($\mu\text{g/ml}$)	Recovery O/E (%)
1	4.65	-	-
	21.90	24.34	90.0
	14.40	15.57	92.5
	10.16	10.10	100.6
2	7.79	-	-
	23.87	27.48	86.9
	15.58	18.71	83.3
	11.87	13.24	89.7

For **protocol (b)** with **urine** samples:

Sample	Observed (ng/ml)	Expected (ng/ml)	Recovery O/E (%)
1	15.44	-	-
	76.75	76.30	100.6
	47.53	42.07	113.0
	25.98	25.29	102.7
2	1.05	-	-
	59.91	61.91	96.8
	25.91	27.68	93.6
	12.15	10.90	111.5

For **protocol (b)** with **CSF** samples:

Sample	Observed (ng/ml)	Expected (ng/ml)	Recovery O/E (%)
1	6.53	-	-
	39.46	44.03	89.6
	18.17	21.53	84.4
	13.16	14.03	93.8
2	6.67	-	-
	40.85	44.17	92.5
	19.72	21.67	91.0
	13.61	14.17	96.0

Linearity

Samples were serially diluted with Dilution Buffer and assayed.

For **protocol (a)** with **serum** samples:

Sample	Dilution	Observed ($\mu\text{g/ml}$)	Expected ($\mu\text{g/ml}$)	Recovery O/E (%)
1	-	14.21	-	-
	2x	6.51	7.11	91.6
	4x	4.05	3.55	113.9
	8x	1.73	1.78	97.3
2	-	19.98	-	-
	2x	10.51	9.99	105.2
	4x	5.40	5.00	108.1
	8x	2.35	2.50	94.1

For **protocol (b)** with **urine** samples:

Sample	Dilution	Observed (ng/ml)	Expected (ng/ml)	Recovery O/E (%)
1	-	70.56	-	-
	2x	42.42	35.28	120.2
	4x	21.00	17.64	119.0
	8x	9.18	8.82	104.1
2	-	27.08	-	-
	2x	14.83	13.54	109.5
	4x	7.35	6.77	108.6

For **protocol (b)** with **CSF** samples:

Sample	Dilution	Observed (ng/ml)	Expected (ng/ml)	Recovery O/E (%)
1	-	25.52	-	-
	2x	12.30	12.76	96.4
	4x	5.82	6.38	91.2
2	-	31.97	-	-
	2x	18.39	15.98	115.1
	4x	9.58	7.99	119.9
	8x	4.65	4.00	116.5

Effect of sample matrix

EDTA, citrate and heparin plasmas were compared to respective serum samples from the same 10 healthy/normal individuals. Results are shown below:

Volunteer No.	Serum ($\mu\text{g/ml}$)	Plasma ($\mu\text{g/ml}$)		
		EDTA	Citrate	Heparin
1	9.46	8.49	8.64	10.42
2	7.24	7.9	6.43	6.66
3	6.28	5.99	5.54	6.62
4	9.92	9.86	9.26	8.66
5	19.94	18.95	15.53	19.01
6	18.02	19.58	12.37	17.82
7	8.59	6.76	5.74	17.91
8	18.91	20.16	17.79	20.07
9	18.82	15.34	15.43	17.83
10	8.47	6.64	6.63	8.08
Mean ($\mu\text{g/ml}$)	12.4	12.0	10.3	12.3
Mean Plasma/Serum (%)	-	96.8	83.6	99.6
Coefficient of determination R^2	-	0.96	0.90	0.98

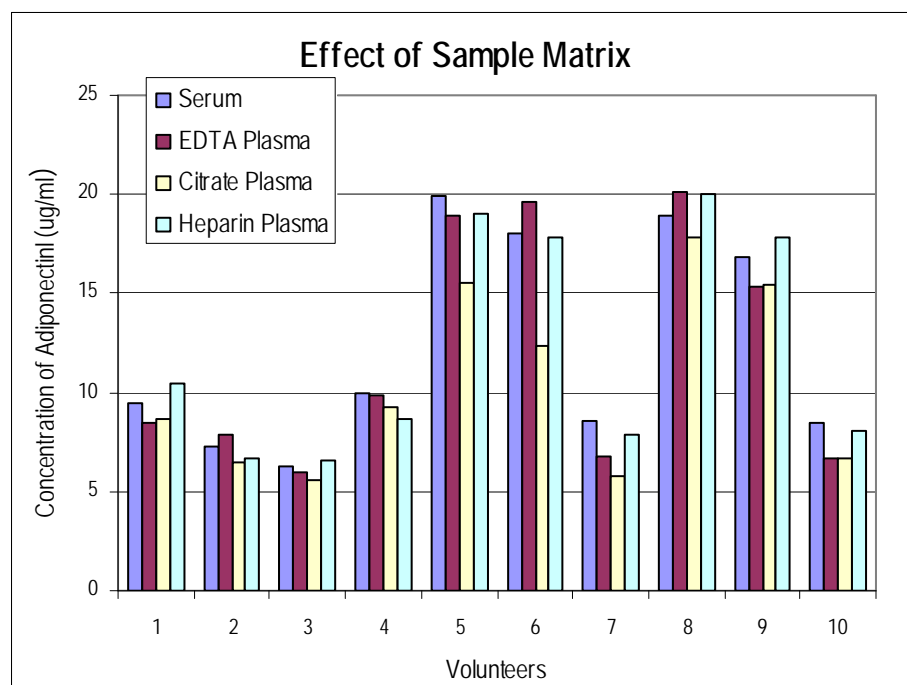


Figure 3: Adiponectin levels measured using Adiponectin (human), ELISA kit from 10 individuals using serum, EDTA, citrate and heparin plasma, respectively.

Stability of samples stored at 2-8°C

Samples should be stored at -20°C. However, no significant decline in concentration of human adiponectin was observed in serum and plasma samples after 7 days when stored at 2-8°C. To avoid microbial contamination, samples were treated with ϵ -aminocaproic acid and sodium azide, resulting in the final concentration of 0.03% and 0.1%, respectively.

Sample	Incubation Temp, Period	Serum ($\mu\text{g/ml}$)	Plasma ($\mu\text{g/ml}$)		
			EDTA	Citrate	Heparin
1	-20°C	3.02	3.13	2.59	3.45
	2-8°C, 1 day	2.95	2.84	2.54	2.95
	2-8°C, 7 days	2.68	2.87	2.5	2.95
2	-20°C	6.77	6.82	6.22	7.06
	2-8°C, 1 day	6.77	6.60	5.83	6.69
	2-8°C, 7 days	7.11	7.06	6.06	7.04
3	-20°C	12.78	12.26	10.81	12.46
	2-8°C, 1 day	13.52	13.29	11.85	13.28
	2-8°C, 7 days	14.05	13.06	12.64	14.12

Effect of Freezing/Thawing

No significant decline was observed in concentration of human adiponectin in serum and plasma samples after repeated (5x) freeze/thaw cycles. However it is recommended to avoid unnecessary repeated freezing/thawing of the samples.

Sample	Number of f/t	Serum ($\mu\text{g/ml}$)	Plasma ($\mu\text{g/ml}$)		
			EDTA	Citrate	Heparin
1	1x	7.97	9.02	7.55	10.28
	3x	8.33	9.06	8.27	9.03
	5x	7.68	9.42	6.77	7.81
2	1x	12.92	14.60	10.87	13.31
	3x	13.34	13.32	12.61	12.78
	5x	12.38	15.11	13.31	15.78
3	1x	11.57	12.23	11.03	14.45
	3x	10.55	14.55	11.03	15.74
	5x	11.97	15.14	10.80	13.88

DEFINITION OF THE STANDARD

The recombinant human adiponectin is used as the Standard. The recombinant human adiponectin is produced in HEK293 cell line and contains 225 amino acid residues of the human adiponectin and 8 extra AA.

PRELIMINARY POPULATION AND CLINICAL DATA

Reference range for serum samples

It is recommended that each laboratory include its own panel of control samples in the assay. Each laboratory should establish its own normal and pathological reference ranges for adiponectin levels with the assay.

Tissue extract

Adiponectin was detected in adipose tissue extracts. Concentrations of 0.2 to 2 μ g/ml were found (total protein concentration 1 mg/ml).

Milk samples

Adiponectin concentrations measured in breast milk samples (n = 18) were in the range of 7 – 40ng/ml.

Urine samples

Adiponectin concentrations were measured in proteinuremic urine samples (n = 10) and non-proteinuremic urine samples (n = 10). Significant differences between the two groups were observed.

<i>Urine samples</i>	<i>Sample ID</i>	<i>Adiponectin (ng/ml)</i>	<i>Mean (ng/ml)</i>	<i>SD (ng/ml)</i>
Proteinuremic samples	4	24.1	59.7 (n=10)	83.9 (n=10)
	19	114.3		
	37	1.1		
	40	3.7		
	41	135.1		
	50	18.5		
	69	5.3		
	83	21.2		
	176	256.8		
	196	16.7		
Non-Proteinuremic samples	51	1.3	3.1 (n=10)	6.5 (n=10)
	52	ND		
	54	19.1		
	59	ND		
	73	ND		
	87	ND		
	103	ND		
	128	ND		
	146	ND		
	163	10.5		

ND - Adiponectin concentrations were below 0.5ng/ml (not detectable).

Cerebrospinal fluid samples

Adiponectin concentrations were measured in serum and CSF samples obtained from the same persons (n = 36).

Sample ID	Adiponectin in CSF (ng/ml)	Adiponectin in serum (ng/ml)
1	17.9	13 810
3	19.8	1 727
4	84.0	9 140
5	17.7	20 080
6	8.1	11 250
11	11.4	6 300
12	200.0	11 700
13	54.0	8 180
14	70.9	15 780
15	4.8	14 870
16	5.3	13 820
23	16.4	10 070
24	3.5	9 180
25	10.8	15 480
27	26.7	20 960
29	29.0	22 900
30	4.4	11 220
31	179.9	32 540
32	13.0	7 580
33	23.6	16 320
34	172.7	8 170
35	14.1	6 940
36	27.3	10 520
41	14.9	18 930
42	5.2	7 660
43	8.1	7 460
46	4.6	12 040
48	32.6	19 620
51	17.8	11 950
52	0.0	7 370
53	30.9	33 200
54	31.2	25 650
55	7.6	13 900
56	10.4	7 530
57	12.2	22 890
58	8.9	13 920
Mean	33.3	14 340
SD	48.3	6 782
n	n = 36	n = 36

TROUBLESHOOTING AND FAQs

Weak signal in all wells

Possible explanations:

- Omission of a reagent or a step
- Improper preparation or storage of a reagent
- Assay performed before reagents were allowed to come to room temperature
- Improper wavelength when reading absorbance

High signal and background in all wells

Possible explanations:

- Improper or inadequate washing
- Overdeveloping; incubation time with Substrate Solution should be decreased before addition of Stop Solution
- Incubation temperature over 30°C

High coefficient of variation (CV)

Possible explanation:

- Improper or inadequate washing
- Improper mixing Standards, Quality Controls or samples

REFERENCES

References to adiponectin:

- Zhu W, Cheng KK, Vanhoutte PM, Lam KS, Xu A. Vascular effects of adiponectin: molecular mechanisms and potential therapeutic intervention. *Clin Sci (Lond)*. 2008 Mar;114(5):361-74.
- Takemura Y, Walsh K, Ouchi N. Adiponectin and cardiovascular inflammatory responses. *Curr Atheroscler Rep*. 2007 Sep;9(3):238-43.
- Kadowaki T, Yamauchi T, Kubota N. The physiological and pathophysiological role of adiponectin and adiponectin receptors in the peripheral tissues and CNS. *FEBS Lett*. 2008 Jan 9;582(1):74-80. Epub 2007 Dec 3.
- Wang ZV, Scherer PE. Adiponectin, cardiovascular function, and hypertension. *Hypertension*. 2008 Jan;51(1):8-14. Epub 2007 Nov 12.

- Behre CJ. Adiponectin, obesity and atherosclerosis. *Scand J Clin Lab Invest.* 2007;67(5):449-58.
- Whitehead JP, Richards AA, Hickman IJ, Macdonald GA, Prins JB. Adiponectin--a key adipokine in the metabolic syndrome. *Diabetes Obes Metab.* 2006 May;8(3):264-80.
- Berner HS, Lyngstadaas SP, Spahr A, Monjo M, Thommesen L, Drevon CaA, Syversen U., Reseland J.E.: Adiponectin and its receptors are expressed in bone-forming cells. *Bone* 2004; 35: 842-849
- Takahashi T, Zhu SJ, Sumino H, Saegusa S, Nakahashi T, Iwai K, Morimoto S, Kanda T: Inhibition of cyclooxygenase-2 enhances myocardial damage in a mouse model of viral myocarditis. *Life Sci* 2005
- Blüher M, Fasshauer M, Kralisch S, Schön MR, Krohn K, Paschke R: Regulation of adiponectin receptor R1 and R2 gene expression in adipocytes of C57BL/6 mice. *Biochem Biophys Res Commun* 2005; 329: 1127-1132
- Pilz S, Maerz W, Weihrauch G, Sargsyan K, Almer G, Nauck M, Boehm BO, Winkelmann BR, Mangge H: Adiponectin serum concentrations in men with coronary artery disease: The Ludwigshafen Risk in Cardiovascular Health (LURIC) study. *Clin Chim Acta* 2005
- Haluzik M, Colombo C, Gavrilova O, Chua S, Wolf N, Chen M, Stannard B, Dietz KR, Le Roith D, Reitmanml. Genetic background (C57BL/6J versus FVB/N) strongly influences the severity of diabetes and insulin resistance in ob/ob mice. *Endocrinology* 2004; 145: 3258-3264
- Ouchi N, Kobayashi H, Kihara S, Kumada M, Sato K, Inoue T, Funahashi T, Walsh K. Adiponectin stimulates angiogenesis by promoting cross-talk between AMP-activated protein kinase and Akt signaling in endothelial cells. *J Biol Chem.* 2004; 279: 1304-1309
- Ishikawa Y, Akasaka Y, Ishii T, Yoda-Murakami M, Choi-Miura NH, Tomita M, Ito K, Zhang L, Akishima Y, Ishihara M, Muramatsu M, Taniyama M. Changes in the distribution pattern of gelatin-binding protein of 28 kDa (adiponectin) in myocardial remodelling after ischaemic injury. *Histopathology.* 2003; 42: 43-52

- Waki H, Yamauchi T, Kamon J, Ito Y, Uchida S, Kita S, Hara K, Hada Y, Vasseur F, Froguel P, Kimura S, Nagai R, Kadowaki T. Impaired multimerization of human adiponectin mutants associated with diabetes. Molecular structure and multimer formation of adiponectin. *J Biol Chem.* 2003; 278: 40352-40363
- Yamauchi T, Kamon J, Ito Y, Tsuchida A, Yokomizo T, Kita S, Sugiyama T, Miyagishi M, Hara K, Tsunoda M, Murakami K, Ohteki T, Uchida S, Takekawa S, Waki H, Tsuno NH, Shibata Y, Terauchi Y, Froguel P, Tobe K, Koyasu S, Taira K, Kitamura T, Shimizu T, Nagai R, Kadowaki T. Cloning of adiponectin receptors that mediate antidiabetic metabolic effects. *Nature.* 2003; 423: 762-769
- Combs TP, Pajvani UB, Berg AH, Lin Y, Jelicks LA, Laplante M, Nawrocki A, Rajala MW, Parlow AF, Cheeseboro L, Ding Y, Russell RG, Lindemann D, Hartley A, Baker GR, Obici S, Deshaies Y, Ludgate ME, Rossetti L, Scherer PE. A transgenic mouse with deletion in the collagenous domain of adiponectin displays elevated circulating adiponectin and improved insulin sensitivity. *Endocrinology* 2003
- Yamauchi T, Kamon J, Waki H, Imai Y, Shimozawa N, Hioki K, Uchida S, Ito Y, Takakuwa K, Matsui J, Takata M, Eto K, Terauchi Y, Komeda K, Tsunoda M, Murakami K, Ohnishi Y, Naitoh T, Yamamura K, Ueyama Y, Froguel P, Kimura S, Nagai R, Kadowaki T. Globular adiponectin protected ob/ob mice from diabetes and ApoE-deficient mice from atherosclerosis. *J Biol Chem.* 2003; 278: 2461-2468
- Pajvani UB, Du X, Combs TP, Berg AH, Rajala MW, Schulthess T, Engel J, Brownlee M, Scherer PE. Structure-function studies of the adipocyte-secreted hormone Acrp30/adiponectin. Implications for metabolic regulation and bioactivity. *J Biol Chem.* 2003; 278: 9073-9085
- Spranger J, Kroke A, Mohlig M, Bergmann MM, Ristow M, Boeing H, Pfeiffer AF. Adiponectin and protection against type 2 diabetes mellitus. *Lancet.* 2003; 361: 226-228
- Kondo H, Shimomura I, Matsukawa Y, Kumada M, Takahashi M, Matsuda M, Ouchi N, Kihara S, Kawamoto T, Sumitsuji S, Funahashi T, Matsuzawa Y. Association of adiponectin mutation with type 2 diabetes: a candidate gene for the insulin resistance syndrome. *Diabetes.* 2002; 51: 2325-2328







- Wang Y, Xu A, Knight C, Xu LY, Cooper GJ. Hydroxylation and glycosylation of the four conserved lysine residues in the collagenous domain of adiponectin. Potential role in the modulation of its insulin-sensitizing activity. *J Biol Chem.* 2002; 277: 19521-19529
- Berg AH, Combs TP, Du X, Brownlee M, Scherer PE. The adipocyte-secreted protein Acrp30 enhances hepatic insulin action. *Nat Med.* 2001; 7: 947-953
- Yamauchi T, Kamon J, Waki H, Terauchi Y, Kubota N, Hara K, Mori Y, Ide T, Murakami K, Tsuboyama – Kasaoka N, Ezaki O, Akanuma Y, Gavrilu O, Vinson C, Reitmanml, Kagechika H, Shudo K, Yoda M, Nakano Y, Tobe K, Nagai R, Kimura S, Tomita M, Froguel P, Kadowaki T. The fat-derived hormone adiponectin reverses insulin resistance associated with both lipoathropy and obesity. *Nat Med.* 2001; 7: 941-946
- Fruebis J, Tsao TS, Javorschi S, Ebbets - Reed D, Erickson MR, Yen FT, Bihain BE, Lodish HF. Proteolytic cleavage product of 30-kDa adipocyte complement-related protein increases fatty acid oxidation in muscle and causes weight loss in mice. *Proc Natl Acad Sci USA.* 2001; 98: 2005-2010
- Das K, Lin Y, Widen E, Zhang Y, Scherer PE. Chromosomal localization, expression pattern, and promoter analysis of the mouse gene encoding adipocyte-specific secretory protein Acrp30. *Biochem Biophys Res Commun.* 2001; 280: 1120-1129
- Arita Y, Kihara S, Ouchi N, Takahashi M, Maeda K, Miyagawa J, Hotta K, Shimomura I, Nakamura T, Miyaoka K, Kuriyama H, Nishida M, Yamashita S, Okubo K, Matsubara K, Muraguchi M, Ohmoto Y, Funahashi T, Matsuzawa Y. Paradoxical decrease of an adipose-specific protein, adiponectin, in obesity. *Biochem Biophys Res Commun.* 1999; 257: 79-83
- Maeda K, Okubo K, Shimomura I, Funahashi T, Matsuzawa Y, Matsubara K. cDNA cloning and expression of a novel adipose specific collagen-like factor, apM1 (adipose most abundant gene transcript 1). *Biochem Biophys Res Commun.* 1996; 221: 286-289
- Kadowaki T, Yamauchi T. Adiponectin and adiponectin receptors. *Endocr Rev.* 2005 May;26(3):439-51. Scherer PE, Williams S, Fogliano M, Baldini G, Lodish HF. A novel serum protein similar to C1q, produced exclusively in adipocytes. *J Biol Chem.* 1995; 270: 26746-26749

References to this product:

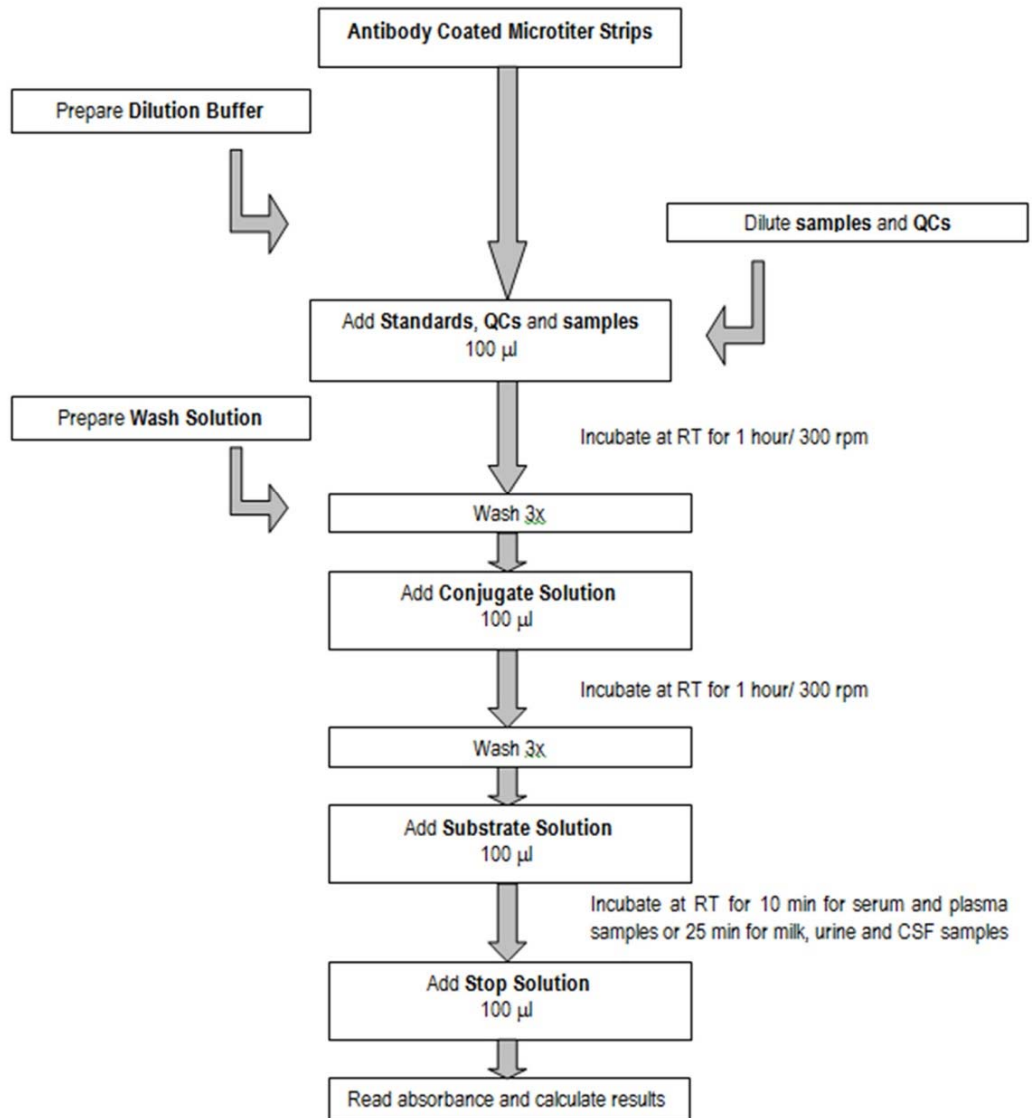
- Risch L, Saely C, Hoefle G, Rein P, Langer P, Gouya G, Marte T, Aczel S, Drexel H: Relationship between glomerular filtration rate and the adipokines adiponectin, resistin and leptin in coronary patients with predominantly normal or mildly impaired renal function. *Clin Chim Acta* 376 (1-2):108-13 (2007). Epub Jul 29 (2006)
- Bronsky J, Karpisek M, Bronska E, Pechova M, Jancikova B, Kotolova H, Stejskal D, Prusa R, Nevoral J: Adiponectin, adipocyte fatty acid binding protein, and epidermal fatty acid binding protein: proteins newly identified in human breast milk. *Clin Chem* 52 (9):1763-70 (2006)
- Tsioufis C, Dimitriadis K, Chatzis D, Vasiliadou C, Tousoulis D, Papademetriou V, Toutouzas P, Stefanadis C, Kalikazaros I. Relation of microalbuminuria to adiponectin and augmented C-reactive protein levels in men with essential hypertension. *Am J Cardiol*. 2005, 96(7), 946-951.
- Karaduman M, Sengul A, Oktenli C, Pekel A, Yesilova Z, Musabak U, Sanisoglu SY, Gunay C, Baysan O, Kocar IH, Tatar H, Ozata M. Tissue levels of adiponectin, tumour necrosis factor-alpha, soluble intercellular adhesion molecule-1 and heart-type fatty acid-binding protein in human coronary atherosclerotic plaques. *Clin Endocrinol (Oxf)*. 2006, 64(2), 196-202.
- Risch L, Guenter H, Saely C, Berchthold S, Weber M, Gouya G, Rein P, Langer P, Marte T, Aczel S, Drexel H: Evaluation of two fully automated novel enzyme-linked immunosorbent assays for the determination of human adiponectin in serum. *Clin Chim Acta* Nov;373(1-2):121-6 (2006) (PubMed)
- Efstathiou SP, Tsioulos DI, Tsiakou AG, Gratsias YE, Pefanis AV, Mountokalakis TD: Plasma adiponectin levels and five-year survival after first-ever ischemic stroke. *Stroke* Sep;36(9):1915-9 (2005) (PubMed)

For more references on this product see our website at www.enzolifesciences.com

EXPLANATION OF SYMBOLS

REF	Catalogue number
Cont.	Content
LOT	Lot number
	See instructions for use
	Biological hazard
	Expiry date
	Storage conditions
	Identification of packaging materials
IVD 	In vitro diagnostic medical device

Assay Procedure Summary





Product Manual

www.enzolifesciences.com

Enabling Discovery in Life Science®

GLOBAL HEADQUARTERS

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

EUROPE/ASIA

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

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