



EAGLE
BIOSCIENCES

Competitive MGO ELISA Kit

Catalog Number: MGO39-K01

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

v. 1.1 (16 OCT 23)

EAGLE BIOSCIENCES, INC.
20A Northwest Blvd., Suite 112, Nashua, NH 03063
Phone: 617-419-2019 Fax: 617-419-1110
WWW.EAGLEBIO.COM



INTENDED USE

The Eagle Biosciences Competitive MGO ELISA Kit is to be used for quantitative determination of MGO in plasma and faeces samples. The Eagle Biosciences Competitive MGO ELISA Kit is for research use only and not for diagnostic or therapeutic procedures.

For further information about this kit, its application, or the procedures in this insert, please contact the Technical Service Team at Eagle Biosciences, Inc at www.EagleBio.com or at 866-411-8023.

ASSAY BACKGROUND

Glycation, or nonenzymatic glycation, is the nonenzymatic reaction of glucose and other reducing sugars with free amino groups of proteins, lipids and nucleic acids. The amino groups of the side chains of arginine and lysine are the primary targets for this type of modification. Over time, the initial glycation products may undergo intramolecular rearrangements and oxidation reactions (glycosidation) and ultimately transform in stable, so-called advanced glycation end products (AGEs). Several compounds, e.g. N ϵ -carboxymethyl-lysine (CML), pentosidine, or methylglyoxal (MGO) derivatives, serve as examples of well-characterized and widely studied AGEs.

AGEs have the potential to interact with a specific receptor (RAGE), a member of the immunoglobulin superfamily, initiating signal pathways that amplify inflammation and oxidative stress, and thereby leading to cellular injury and death. High levels of circulating AGEs are associated with cardiovascular disease, diabetes, chronic kidney disease, and increased mortality. MGO is a highly reactive dicarbonyl metabolite formed during glucose, protein and fatty acid metabolism. Several nonenzymatic as well as enzymatic reactions are involved in MGO formation. MGO can be formed by nonenzymatic elimination of phosphate from triose phosphates, glyceraldehyde-3-phosphate and glycero phosphate, as well as enzymatically from dihydroxyacetone phosphate by MGO synthase. MGO reacts with protein by initial reversible reactions: with arginine and lysine residues forming glycosylamine, with cysteine forming hemithioacetal. The irreversible reaction of MGO with lysine residues of protein forms N ϵ -carboxyethyl-lysine (CEL), N ϵ -carboxymethyl-lysine (CML) and 1,3-di(N-lysino)-4-methyl-imidazolium (MOLD). MGO reacts with arginine residues of protein to form the non-fluorescent products argpyrimidine and hydroxyimidazolone.

MGO-induced AGEs have been linked to the aging process and age-related diseases such as cardiovascular complications of diabetes, neurodegenerative disease and connective tissue disorders. Furthermore, MGO levels are elevated in hyperglycemia and has been implicated in hypertension and atherosclerosis.

The MGO assay can be used to measure MGO-adducts as HSA-MGO arbitrary units. In normal citrate plasma ~200-400 AU/ml can be measured.

ASSAY PRINCIPLE

The competitive MGO ELISA is a ready-to-use solid-phase enzyme-linked immunosorbent assay based on the inhibition principle with a working time of 3½ hours. The efficient format of a plate with twelve disposable 8-well strips allows free choice of batch size for the assay. Samples and standards are pre-incubated with biotinylated tracer antibody in U-shape microtiter plate. Pre-incubated samples and standards with the biotinylated tracer antibodies are incubated on coated strips. Streptavidin-peroxidase conjugate will bind to the biotinylated tracer antibody. Streptavidin-peroxidase will react with the substrate, tetramethylbenzidine (TMB). The enzyme reaction is stopped by the addition of



oxalic acid. The absorbance at 450 nm is measured with a spectrophotometer. A standard curve is obtained by plotting the absorbance (linear) versus the corresponding concentrations of the known standards of the HSA-MGO standards (log). The MGO concentration of samples, which are run concurrently with the standards, can be determined from the standard curve.

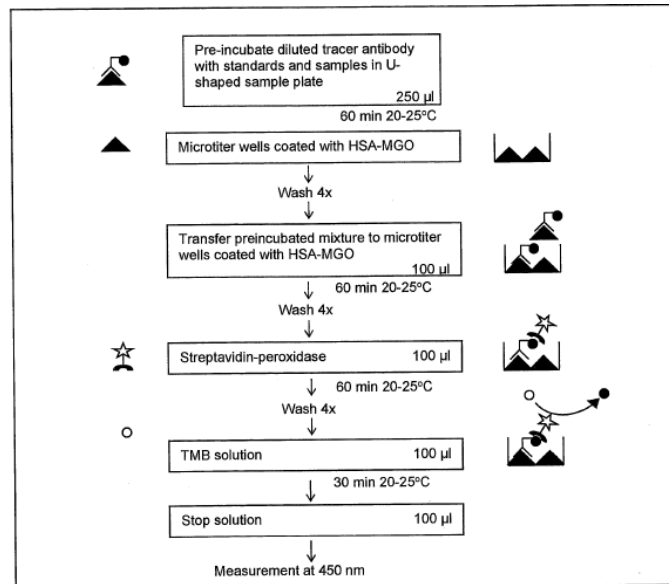
KIT FEATURES

- Working time of 3.5 hours
- Minimum concentration which can be measured is 15.6AU/ml
- Measurable concentration range of 15.6 to 1000 AU/ml.
- Working volume 100µl/well

CROSS REACTIVITY

The ELISA detects MGO containing proteins. Since the assay detects a modified amino acid the assay is useful for proteins of all species.

PROTOCOL OVERVIEW



- The competitive MGO ELISA is a ready-to-use solid-phase enzyme-linked immunosorbent assay based on the inhibition principle with a working time of 3½ hours.
- The efficient format of a plate with twelve disposable 8-well strips allows free choice of batch size for the assay.
- Samples and standards are pre-incubated with biotinylated tracer antibody in U-shape microtiter plate.
- Pre-incubated samples and standards with the biotinylated tracer antibodies are incubated on coated strips.
- Streptavidin-peroxidase conjugate will bind to the biotinylated tracer antibody.
- Streptavidin-peroxidase will react with the substrate, tetramethylbenzidine (TMB).
- The enzyme reaction is stopped by the addition of oxalic acid.



- The absorbance at 450 nm is measured with a spectrophotometer. A standard curve is obtained by plotting the absorbance (linear) versus the corresponding concentrations of the known standards of the HSA-MGO standards (log).
- The MGO concentration of samples, which are run concurrently with the standards, can be determined from the standard curve.

KIT COMPONENTS

Kit Component	Quantity	Color Code
Wash Buffer 20x	1 vial (60ml)	Colorless
Dilution Buffer 10x	1 vial (15 ml)	Green
Standard	2 vials, lyophilized	White
Tracer, biotinylated 100x	2 vials 0.25 ml lyophilized	Blue
Streptavidin-peroxidase 100x	1 tube, 0.25 ml in solution	Brown
TMB Substrate	1 vial (11 ml)	Brown
Stop Solution	1 vial (22 ml)	Red
12 Microtiter strips, pre-coated	1 plate	
U-Shaped Microtiter Plate	1 plate	
Certificate of Analysis	1	
Package Insert	1	

- Upon receipt, store individual components at 2 - 8°C. Do not freeze.
- Do not use components beyond the expiration date printed on the kit label.
- The standard and conjugate in lyophilized form and the streptavidin-peroxidase in concentrated solution are stable until the expiration date indicated on the kit label, if stored at 2 - 8°C.
- The exact amount of standard is indicated on the label of the vial and the Certificate of Analysis.
- The standard is single use. After reconstitution the standard cannot be stored for repeated use.
- The tracer is single use. After reconstitution the conjugate cannot be stored for repeated use.
- The streptavidin-peroxidase can only be stored in concentrated solution and is not stable when stored diluted.
- Upon receipt, foil pouch around the plates should be vacuum-sealed and unpunctured. Any irregularities to aforementioned conditions may influence plate performance in the assay.
- Return unused strips immediately to the foil pouch containing the desiccant pack and reseal along the entire edge of the zip-seal. Quality guaranteed for 1 month if stored at 2 - 8°C.

MATERIALS REQUIRED BUT NOT PROVIDED

- Calibrated micropipettes and disposable tips.
- Distilled or de-ionized water.
- Plate washer: automatic or manual.
- Polypropylene tubes.
- Calibrated ELISA plate reader capable of measuring absorbance at 450 nm.
- Adhesive covers.
- Centrifuge for 1 ml tubes.



WARNING AND PRECAUTIONS

- For research use only, not for diagnostic or therapeutic use.
- This kit should only be used by qualified laboratory staff.
- Do not under any circumstances add sodium azide as preservative to any of the components.
- Do not use kit components beyond the expiration date.
- Do not mix reagents from different kits and lots. The reagents have been standardized as a unit for a given lot. Use only the reagents supplied by manufacturer.
- The assay has been optimized for the indicated standard range. Do not change the standard range.
- Open vials carefully: vials are under vacuum.
- It is advised to spin down streptavidin-peroxidase tubes before use.
- Do not ingest any of the kit components.
- Kit reagents contain 2-chloroacetamide as a preservative. 2-Chloroacetamide is harmful in contact with skin and toxic if swallowed. In case of accident or if you feel unwell, seek medical advice immediately.
- The TMB substrate is light sensitive, keep away from bright light. The solution should be colorless until use.
- The stop solution contains 2% oxalic acid and can cause irritation or burns to respiratory system, skin and eyes. Direct contact with skin and eyes should be strictly avoided. If contact occurs, rinse immediately with plenty of water and seek medical advice.
- Incubation times, incubation temperature and pipetting volumes other than those specified may give erroneous results.
- Do not reuse micro wells or pour reagents back into their bottles once dispensed.
- Handle all biological samples as potentially hazardous and capable of transmitting diseases.
- Use polypropylene tubes for preparation of standard and samples. Do not use polystyrene tubes or sample plates.

SAMPLE PREPARATION

Collection and handling

Plasma

Collect blood using normal aseptic techniques. Blood samples should be kept on ice. If serum is used, separate serum from blood after clotting at room temperature within 1 hour by centrifugation (1500xg at 4°C for 15 min). Transfer the serum to a fresh polypropylene tube.

If plasma is used, separate plasma from blood within 20 minutes after blood sampling by centrifugation (1500xg at 4°C for 15 min). Transfer the plasma to a fresh polypropylene tube.

Faeces

MGO can be measured in faeces if samples are extracted, for example using the following extraction buffer: 0.1 M Tris, 0.15 M NaCl, 1.0 M urea, 10 mM CaCl₂, 5 g/l BSA and 0.25 mM thimerosal (pH 8.0). Add 5 ml extraction buffer to 100 mg sample (giving a dilution factor of 51, assuming the density of faeces to be 1 g/ml). Vortex samples and



filter the samples to remove coarse particles (> 0.6 mm). Shake the filtrate for 20 minutes and centrifuge samples: 10,000xg at 4°C for 20 minutes. Use supernatant for analysis.

Storage

Store samples below -20°C, preferably at -70°C in polypropylene tubes. Storage at -20°C can affect recovery of inhibitory capacity. Use samples within 24 hours after thawing. Avoid multiple freeze-thaw cycles which may cause loss of endoblock LBP activity and give erroneous results. Before performing the assay, samples should be brought to room temperature (18 – 25°C) and mixed gently. Prepare all samples (controls and test samples) prior to starting the assay procedure. Avoid foaming.

DILUTION PROCEDURES

Plasma samples

Samples can be measured accurately undiluted.

Faeces

Samples can be measured accurately undiluted.

Comment regarding recommended sample dilution

The mentioned dilution for samples is a minimum dilution and should be used as a guideline. The recovery of inhibitory compound from an undiluted sample is not 100% and may vary from sample to sample. When testing less diluted samples it is advisable to run recovery experiments to determine the influence of the matrix on the detection of inhibitory compound.

Do not use polystyrene tubes or sample plates for preparation or dilution of the samples.

Guideline for dilution of samples

Please see the table below for recommended sample dilutions. Volumes are based on a total volume of at least 230 µl of diluted sample, which is sufficient for one sample in duplicate in the ELISA. For dilution of samples we recommend to use at least 10 µl of sample.

Dilution	Pre-Dilution	Amount of sample or pre-dilution required	Amount of Dilution buffer required
10x	Not Necessary	25 µl	225 µl
20x	Not necessary	15 µl (sample)	285 µl
50x	Not necessary	10 µl (sample)	490 µl
100x	Not necessary	10 µl (sample)	990 µl
500x	Recommended 10x (see nr.1)	10 µl (pre-dilution)	490 µl
1000x	Recommended 10x (see nr.1)	10 µl (pre-dilution)	990 µl
2000x	Recommended 20x (see nr.2)	10 µl (pre-dilution)	990 µl
5000x	Recommended 50x (see nr.3)	10 µl (pre-dilution)	990 µl

REAGENT PREPARATION

Allow all the reagents to equilibrate to room temperature (20 – 25°C) prior to use. Return to proper storage conditions immediately after use.



Wash buffer

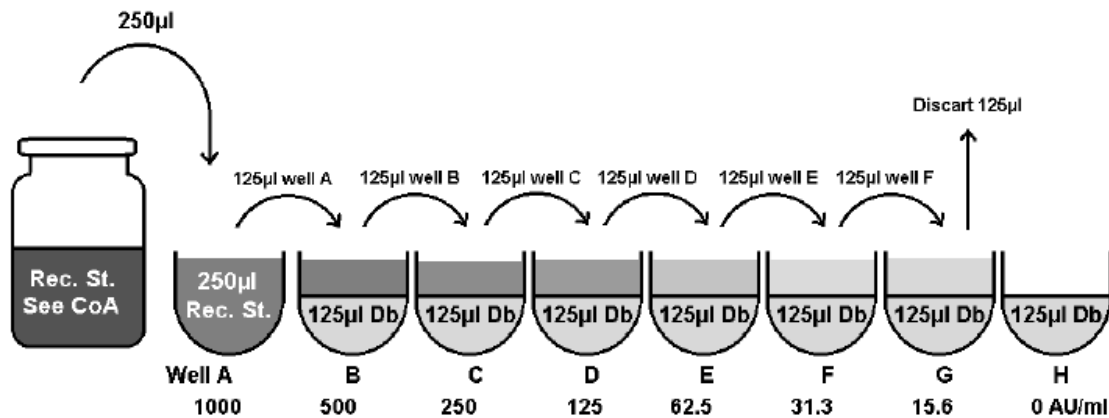
Prepare wash buffer by mixing 60 ml of 20x wash buffer with 1140 ml of distilled or de-ionized water, which is sufficient for 1 x 96 tests. In case less volume is required, prepare the desired volume of wash buffer by diluting 1 part of the 20x wash buffer with 19 parts of distilled or de-ionized water

Dilution buffer

Prepare dilution buffer by mixing 15 ml of the 10x dilution buffer with 135 ml of distilled or de-ionized water, which is sufficient for 1 x 96 tests. In case less volume is required, prepare the desired volume of dilution buffer by diluting 1 part of the 10x dilution buffer with 9 parts of distilled or de-ionized water. Concentrated dilution buffer may contain crystals. In case the crystals do not disappear at room temperature within 1 hour, concentrated dilution buffer can be warmed up to 37°C. Do not shake the solution

Standard solution

The standard is reconstituted by pipetting the amount of dilution buffer mentioned on the CoA in the standard vial. Transfer 250 µl of the reconstituted standard to well A1 of the U-shaped microtiter plate. Prepare a serial dilution of the reconstituted standard with dilution buffer in the U-shaped microtiter plate as shown in the Figure below.



Tracer solution

The tracer is reconstituted by pipetting 0.25 ml distilled or de-ionized endotoxin free water. After reconstitution the tracer cannot be stored for repeated use. Dilute the reconstituted 0.25 ml tracer with 24.75 ml dilution buffer, which is sufficient for 1 x 96 tests. In case less volume is required, prepare the desired volume of tracer by diluting 1 part of the reconstituted tracer with 99 parts of dilution buffer.

Streptavidin-peroxidase solution

It is advised to spin down streptavidin-peroxidase tubes before use. Prepare the streptavidin-peroxidase solution by mixing 0.25 ml of the 100x streptavidin-peroxidase solution with 24.75 ml dilution buffer, which is sufficient for 2 x 96 tests. In case less volume is required, prepare the desired volume of streptavidin-peroxidase solution by diluting 1 part of the 100x streptavidin-peroxidase solution with 99 parts of dilution buffer.

ELISA PROTOCOL

Bring all reagents to room temperature (20 - 25°C) before use.



1. Determine the number of test wells required and fill out the data collection sheet. Return the unused strips of the HSA-MGO coated microtiter plate to the storage bag with desiccant, seal and store at 2 - 8°C.

Please notice: the amount of standards and samples with tracer (total of 250µl) in the U-shape plate is sufficient for a duplicate determination of standards, samples and controls. Therefore, double the amount of wells for the coated microtiter plate are needed.

2. Prepare the standard dilution series in the U-shaped plate. Prepare the samples and add 125 µl to the U-shaped plate.
3. Add 125 µl of the diluted tracer to the 125 µl of standards and samples in the U-shaped plate. Gently mix by pipetting.
4. Cover the tray and tap the tray to eliminate any air bubbles. Be careful not to splash liquid onto the cover.
5. Incubate U-shaped plate for 1 hour at room temperature.
6. Wash the microtiter plate coated with HSA-MGO 4 times with wash buffer using a plate washer or as follows
 - a. Carefully remove cover, avoid splashing.
 - b. Empty the plate by inverting plate and shaking contents out over the sink, keep inverted and tap dry on a thick layer of tissues.
 - c. Add 200 µl of wash buffer to each well, wait 20 seconds, empty the plate as described in 6b.
 - d. Repeat the washing procedure 6b/6c three times.
 - e. Empty the plate and intensively tap on thick layer of tissues. Tap the plate as dry as possible.
7. Transfer 100 µl in duplicate of standard, samples, or controls mixed with tracer from the U-shaped plate into appropriate wells in the coated microtiter plate. Do not touch the side or bottom of the wells.
8. Cover the tray and tap the tray to eliminate any air bubbles. Be careful not to splash liquid onto the cover.
9. Incubate for 1 hour at room temperature.
10. Repeat the wash procedure described in step 6.
11. Add 100 µl of diluted streptavidin-peroxidase to each well, using the same pipetting order as applied in step 7. Do not touch the side or bottom of the wells.
12. Cover the tray and incubate the tray for 1 hour at room temperature.
13. Repeat the wash procedure described in step 6.
14. Add 100 µl of TMB substrate to each well, using the same pipetting order as applied in step 7. Do not touch the side or bottom of the wells.
15. Cover the tray and incubate the tray for 30 minutes at room temperature. It is advised to control the reaction on the plate regularly. In case of strong development the TMB reaction can be stopped sooner. Avoid exposing the



microwell strips to direct sunlight. Covering the plate with aluminium foil is recommended.

16. Stop the reaction by adding 100 μ l of stop solution with the same sequence and timing as used in step 14. Mix solutions in the wells thoroughly by gently swirling the plate. Gently tap the tray to eliminate any air bubbles trapped in the wells.
17. Read the plate within 30 minutes after addition of stop solution at 450 nm using a plate reader, following the instructions provided by the instrument's manufacturer.

In case plate washer is used, please note: use of a plate washer can result in higher background and decrease in sensitivity. We advise validation of the plate washer with the manual procedure. Make sure the plate washer is used as specified for the manual method.

INTERPRETATION OF RESULTS

- Calculate the mean absorbance for each set of duplicate standards, control and samples.
- If individual absorbance values differ by more than 15% from the corresponding mean value, the result is considered suspect and the sample should be retested.
- The mean absorbance of the zero standard should be less than 0.35.
- Create a standard curve using computer software capable of generating a good curve fit. The mean absorbance for each standard concentration is plotted on the vertical (Y) axis versus the corresponding concentration on the horizontal (X) axis (logarithmic scale).
- If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.
- Samples that give a mean absorbance above the absorbance for the highest standard concentration are out of range of the assay. These samples should be retested at a higher dilution.

TECHNICAL HINTS

- User should be trained and familiar with ELISA assays and test procedure.
- If you are not familiar with the ELISA technique it is recommended to perform a pilot assay prior to evaluation of your samples. Perform the assay with a standard curve only following the instructions.
- Improper or insufficient washing at any stage of the procedure will result in either false positive or false negative results. Completely empty wells before dispensing wash buffer, fill with wash buffer as indicated for each cycle and do not allow wells to sit uncovered or dry for extended periods.
- Since exact conditions may vary from assay to assay, a standard curve must be established for every run. Samples should be referred to the standard curve prepared on the same plate.
- Do not mix reagents from different batches, or other reagents and strips. Remainders should not be mixed with contents of freshly opened vials.
- Each time the kit is used, fresh dilutions of standard, sample, tracer, streptavidin-peroxidase and buffers should be made.
- Caps and vials are not interchangeable. Caps should be replaced on the corresponding vials.



- To avoid cross-contaminations, change pipette tips between reagent additions of each standard, between sample additions, and between reagent additions. Also, use separate reservoirs for each reagent.
- The waste disposal should be performed according to your laboratory regulations.

WARRANTY INFORMATION

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.