

OXItek TBARS Assay Kit

THIOBARBITURIC ACID REACTIVE SUBSTANCES

Catalog #: ALX-850-287

(160 tests)

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
Principles of the System	3
Reagents.....	4
Precautions	4
Preparation of Reagents	5
Test Procedure	6
Sample Preparation	7
Calculation and Interpretation of Results	10
References.....	12
Procedural Flow Chart	13
Contact Information.....	14

INTENDED USE

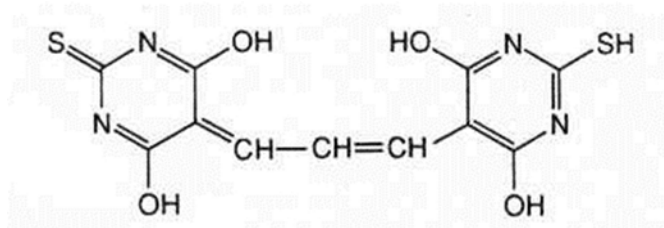
The sensitivity of measuring Thiobarbituric Acid Reactive Substances (TBARS) has made this assay the method of choice for screening and monitoring lipid peroxidation, a major indicator of oxidative stress⁽¹⁻³⁾. This rapid, easy-to-use procedure has been modified by researchers for use with many types of samples including drugs, food products and human and animal biological tissues⁽⁴⁻⁷⁾. The assay has provided important information regarding free radical activity in disease states and has been used for measurement of anti-oxidant activity of several compounds⁽⁸⁻⁹⁾. Although much controversy has appeared in the literature regarding the specificity of TBARS toward compounds other than MDA, it remains the most widely employed assay used to determine lipid peroxidation⁽¹⁰⁾. If lipoprotein fractions are first acid precipitated from the sample, interfering soluble TBARS are minimized. Nonetheless, if TBARS are increased, it is recommended that a more specific assay such as HPLC be performed.

The OXItek TBARS Kit is designed to provide a standardized, reproducible assay with consistent results. Each lot of reagents is quality controlled as a kit, which includes an MDA standard. It is recommended that additional in-house controls are included in each test run.

The OXItek TBARS Assay Kit is for Research Purposes Only.

PRINCIPLE OF THE SYSTEM

Malondialdehyde (MDA) forms a 1:2 adduct with thiobarbituric acid and produces the following:



which can be measured by fluorometry or spectrophotometry. Although this reaction has a much higher sensitivity when measured via fluorometry, protocols for both methods are provided in the Test Procedure section of this insert.

Biological specimens contain a mixture of thiobarbituric acid reactive substances (TBARS), including lipid hydroperoxides and aldehydes, which increase as a result of oxidative stress. TBARS return to normal levels over time, depending upon the presence of anti-oxidants. In practice, TBARS are expressed in terms of malondialdehyde (MDA) equivalents⁽¹¹⁾. In this assay, an MDA standard is used to construct a standard curve against which unknown samples can be plotted.

The OXItek TBARS Assay Kit provides all the necessary reagents to perform 160 tests and is designed for researchers studying oxidative stress and anti-oxidant activity. It is recommended that in-house controls be run with each sample test.

Depending on geographic location, normal plasma and serum TBARS should be <1.5 and <2.0 MDA units respectively. Mean and \pm SD or SE must be established by each laboratory. It is recommended that samples be run in duplicate.

Limitation to Procedure:

1. Hemolyzed, icteric or grossly lipemic plasma samples are not suitable for use in TBARS analysis.
2. Non-lipid TBARS may be present in the sample. It is recommended that a sample with elevated TBARS levels be tested by a more specific test for lipid peroxidation such as HPLC.
3. Normal tissues contain very low levels of free malondialdehyde.

REAGENTS

Materials Supplied:

- ALX-850-287-1 – MDA Diluent (100 ml/kit): Contains sterile deionized water
- ALX-850-287-2 – TBARS Diluent 1 (4 x 50 ml/kit): Contains acetic acid
- ALX-850-287-3 – TBARS Diluent 2 (4 x 50 ml/kit): Contains sodium hydroxide
- ALX-850-287-4 – Thiobabituric Acid (4 vials/kit): Contains 0.53 grams thiobarbituric acid
- ALX-850-287-5 – MDA Standard (20 ml/kit): Contains 100 nmol/ml Malondialdehyde Bis (dimethyl acetal)
- ALX-850-287-6 – SDS Solution (30 ml/kit): Contains sodium dodecyl sulfate
- ALX-850-287-7 – Marbles

Storage:

Store all kit reagents at 2-8°C. The components should be used before the expiration date indicated on the outside of the box.

Materials Required but not Supplied:

- 12 x 75 mm glass test tubes and metal racks
- Disposable gloves
- Adjustable pipettes
- Graduated cylinders and assorted beakers
- Stir plate
- Heat block, incubator or water bath set at 95° C
- Fluorometer or Spectrophotometer

PRECAUTIONS

- To avoid cross contamination, use separate pipette tips for each specimen.
- TBARS Diluent 1 contains acetic acid. Handle and dispose of according to applicable legal and safety guidelines.
- TBARS Diluent 2 contains sodium hydroxide. Handle and dispose of according to applicable legal and safety guidelines.
- If reusing glass tubes, be certain to rinse thoroughly in deionized water.

PREPARATION OF REAGENTS

Note: Prepare fresh for each analysis.

TBA/Buffer Reagent:

Use 1 bottle of TBARS Diluent 1 (50 ml), 1 bottle of TBARS Diluent 2 (50 ml) and 1 vial of Thiobarbituric Acid (TBA). One hundred ml is sufficient for approximately 40 tubes.

Add TBA to a mixing vessel containing half a bottle of TBARS Diluent 1. Rinse the vial with the remaining half of TBARS Diluent 1 and add to the mixing vessel. Add a full bottle of TBARS Diluent 2. Mix until the TBA is completely dissolved. If necessary, very low heat may be used.

MDA Standard For Fluorometer:

Dilute MDA Standard 1:10 by mixing 1.8 ml of MDA Diluent and 0.2 ml of MDA Standard. Mix thoroughly. Prepare a series of 5 standards in MDA Diluent. The dilution scheme for these standards is given below in Table 1.

Table 1: Preparation of MDA Standard For Fluorometer

Standard Number	Concentration of MDA	Volume of MDA Standard	Volume of Diluent
4	4 nmol/ml	400 µl	600 µl
3	3 nmol/ml	300 µl	700 µl
2	2 nmol/ml	200 µl	800 µl
1	1 nmol/ml	100 µl	900 µl
0	0 nmol/ml	0 µl	1000 µl

* The concentration of the 1:10 standard is 10 nmol/ml, and may be used directly as an additional standard when assaying samples expected to have elevated TBARS, such as urines.

MDA Standard For Spectrophotometer:

Use MDA Standard undiluted for 100 nmol/ml standard. Prepare a series of 5 standards in MDA Standard Diluent. The dilution scheme for these standards is given below in Table 2.

Table 2: Preparation of MDA Standard For Spectrophotometer

Standard Number	Concentration of MDA	Volume of 1:10 MDA Standard	Volume of Diluent
4	100 nmol/ml	1000 µl	0 µl
3	50 nmol/ml	500 µl	500 µl
2	25 nmol/ml	250 µl	750 µl
1	12.5 nmol/ml	125 µl	875 µl
0	0 nmol/ml	0 µl	1000 µl



Reagents
require
separate
storage
conditions.

TEST PROCEDURE

Allow all reagents to reach room temperature before use. SDS Solution will take at least one hour if stored at 2-8°C. Heating the SDS Solution at 37°C briefly will redissolve precipitated SDS. SDS Solution can then be stored at room temperature.

Step 1: Collect EDTA plasma (lavender top Vacutainer®). For preparation of other sample types, see sample preparation section of this insert.

Step 2: Label each disposable glass test tube with the standard number or sample identification.

Step 3: Add 100 µl sample or standard to properly labeled tube.

Step 4: Add 100 µl SDS Solution to each tube and swirl to mix.

Step 5: Add 2.5 ml TBA/Buffer Reagent forcefully down the side of each tube.

Step 6: Cover each tube with a glass marble and incubate at 95°C for 60 min.

Step 7: Remove from incubation and cool to room temperature in an ice bath for 10 min.

Step 8: Centrifuge samples at 3000 rpm for 15 min.

Step 9: Remove supernatant from samples for analysis.

Step 10: Fluorescent Analysis: read supernatants with excitation set at 530 nm and emission at 550 nm. It is recommended sensitivity be set at high with a slit width of 5 nm.

Spectrophotometer Analysis: Read absorbance of supernatants at 532 nm.

SAMPLE PREPARATION

Note: Samples should be tested immediately. If serum/plasma samples are not tested immediately, remove from clot, aliquot and store at -70°C.

Platelets:

Collect 5ml specimen of heparinized venous blood (green top Vacutainer®) from patients who have fasted overnight. Remove stopper and let RBCs sediment by gravity. Pipet off platelet-rich plasma (PRP) into clean, plastic centrifuge tube after 1 ml has formed and continue until separation is complete. Centrifuge PRP at 1400 rpm for 5 min, collect supernatant and centrifuge the supernatant at 5000 rpm for 15 min to sediment platelets.

Leukocytes:

Collect blood in EDTA (lavender top Vacutainer®). Allow RBCs to sediment; collect upper "Buffy Coat". Wash leukocytes contained in the Buffy Coat once in an isotonic solution such as saline or Dulbecco's PBS. Resuspend leukocytes in isotonic solution for testing⁽¹⁵⁾.

Erythrocyte Ghosts:

Wash RBC pellet from leukocyte separation with 20 ml of TRIS Buffer, pH 7.4 (6.05 g TRIS, 6.42 g NaCl, 420 ml 0.1M HCl, 580 ml deionized water), centrifuge at 3500 rpm for 10 min, discard supernatant and repeat twice. Add an equal volume of TRIS Buffer to final pellet and incubate a minimum of 4 hours at 4°C. Lysis of erythrocytes is performed on ice with pre-cooled conditions. Add 15 ml lysing solution (301 mg MgSO₄, 372 mg KCl in 500 ml of sterile water) to 0.5 ml of the cell suspension. Immediately add 1 ml of resealing solution (53.7 g KCl, 10.5 g NaCl in 400 ml of deionized water). Keep the suspension on ice for 5 min and then at 37°C for 30 min. Pellet the ghosts by centrifuging for 10 min at 3500 rpm⁽¹²⁾.

Oxidized LDL:

Plasma: Collect fasting heparinized whole blood. Centrifuge at 3500 rpm for 10 min at 5-10°C, carefully remove plasma and place on ice for immediate analysis or freeze several aliquots at -70°C for later analysis. Samples can be safely stored for 1-2 months. Process as described below for serum ⁽³⁾.

Serum: Collect fasting whole blood in a red top vacutainer®. Incubate at room temperature for at least 30 min for clots to form. Centrifuge at 3500 rpm for 10 min. Carefully remove serum and place on ice for immediate analysis or freeze aliquots at -70°C for later analysis. Add 1.5 ml of isotonic saline to 1.5 ml of plasma/serum. Prepare a sodium heparin solution in water containing 40,000 USP units per ml. Prepare a solution of 1.06M manganese chloride in water. A working reagent is prepared with 300 µl sodium heparin solution to 5 ml manganese chloride solution. Scale up as necessary. Add 300 µl of working reagent to the diluted serum/plasma. Vortex and allow the LDL/VLDL to precipitate for 15 min at room temperature. Centrifuge the serum/plasma at 8000 rpm for 15 min. Decant the serum/plasma supernatant and resuspend the pellet in 1.5 ml normal saline or PBS^(3,13). If required, sonicate each sample 5 sec at low setting over ice. Use these samples to start the TBARS assay, substituting water for SDS in the reaction mixture.

Note: When analyzing samples that may be non-fasting, triglycerides measuring above 300-350 mg/dL will yield a turbid supernatant and cannot be measured.

Other Body Fluids:

The TBARS kit is suitable for analysis of CSF, vaginal, synovial, seminal, vitreous, tears, saliva, sperm, pulmonary and gastrointestinal fluids and lavages.

Tissues:

Freeze tissue in liquid nitrogen and immediately crush in a pre-chilled mortar and pestle. Resuspend tissue at 50 mg/ml in normal saline or PBS. Disrupt in a Potter-Elvehjem glass homogenizer. If necessary, sonicate for 15 sec at 40V setting over ice and use uncentrifuged whole homogenate for analysis. Alternatively, homogenize in isotonic media appropriate for sub-cellular fractionation to study TBARS in plasma membranes, nuclear membranes or organelles. Spun supernates may be used for enzyme analyses. Recommend normalizing TBARS values to another constituent such as protein.

Cell Cultures:

Suspend 20 million cells in 1 ml of cell culture medium or buffer of choice such as PBS. Sonicate for 5 second intervals at 40V setting over ice. Use whole homogenates in the assay, being sure to use culture fluid as a sample blank.

Anti-Oxidant Screening:

The TBARS assay may be used for testing anti-oxidants and drugs⁽¹³⁾. Oxidize 2 100 µl aliquots of a sample, such as serum or plasma, with 5mM ferric chloride. Add the anti-oxidant compound to be tested to one. Incubate plasma alone plus the 2 test samples at 37°C for 30 min. Run the TBARS assay and compare to a distilled water control. Calculate percentage of inhibition⁽¹⁴⁾. Drugs may be added directly to plasma samples and compared to plasma alone in the TBARS assay.

CALCULATIONS AND INTERPRETATION OF RESULTS

Typical Standard Curve (Fluorometer):

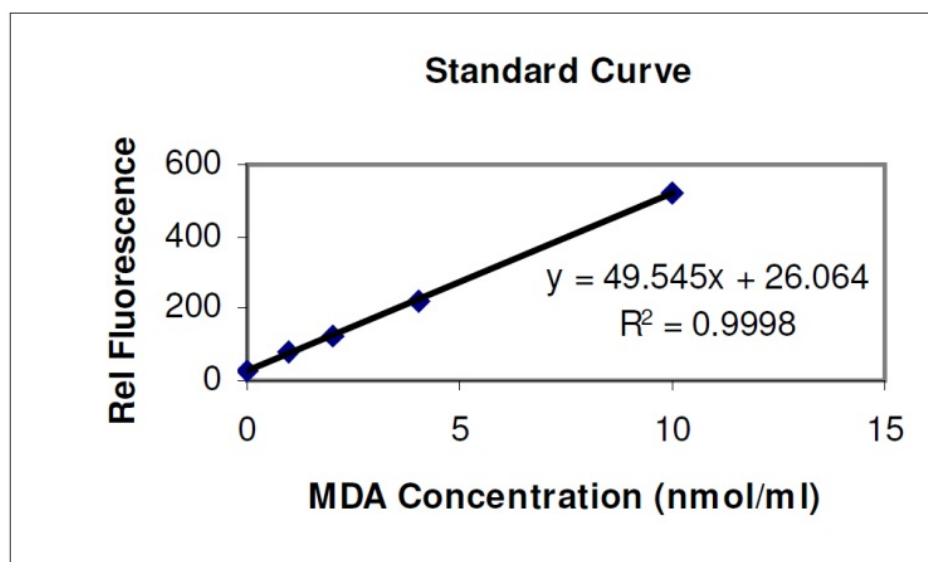
This is an example of a typical standard curve and is not to be used for interpretation of results. Variation may occur in individual laboratories due to pipetting, laboratory and incubator temperatures, etc.

Table 3: Sample Standard Curve

MDA Concentration	Fluorometer Readings
10 nmol/ml	521.15
4 nmol/ml	223.68
2 nmol/ml	126.48
1 nmol/ml	78.69
0 nmol/ml	22.5

For Fluorometer:

Using linear graph paper, plot mean MDA equivalents for each standard used on the X-axis versus the corresponding fluorometer reading on the Y-axis. Determine the concentration of MDA equivalents in nmol/ml in specimens by interpolation from the standard curve. Correct sample values for any other dilutions performed during specimen preparation.



Typical Standard Curve (Spectrophotometer):

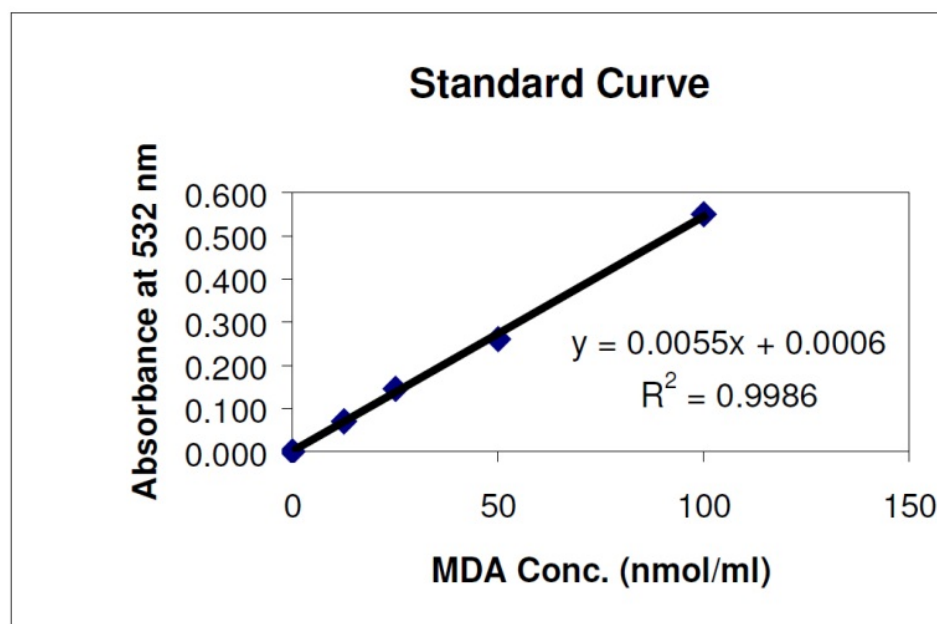
This is an example of a typical standard curve and is not to be used for interpretation of results. Variation may occur in individual laboratories due to pipetting, laboratory and incubator temperatures, etc.

Table 4: Sample Standard Curve

MDA Concentration	Spectrophotometer Readings
100 nmol/ml	0.550
50 nmol/ml	0.260
25 nmol/ml	0.145
12.5 nmol/ml	0.070
0 nmol/ml	0.000

For Spectrophotometer:

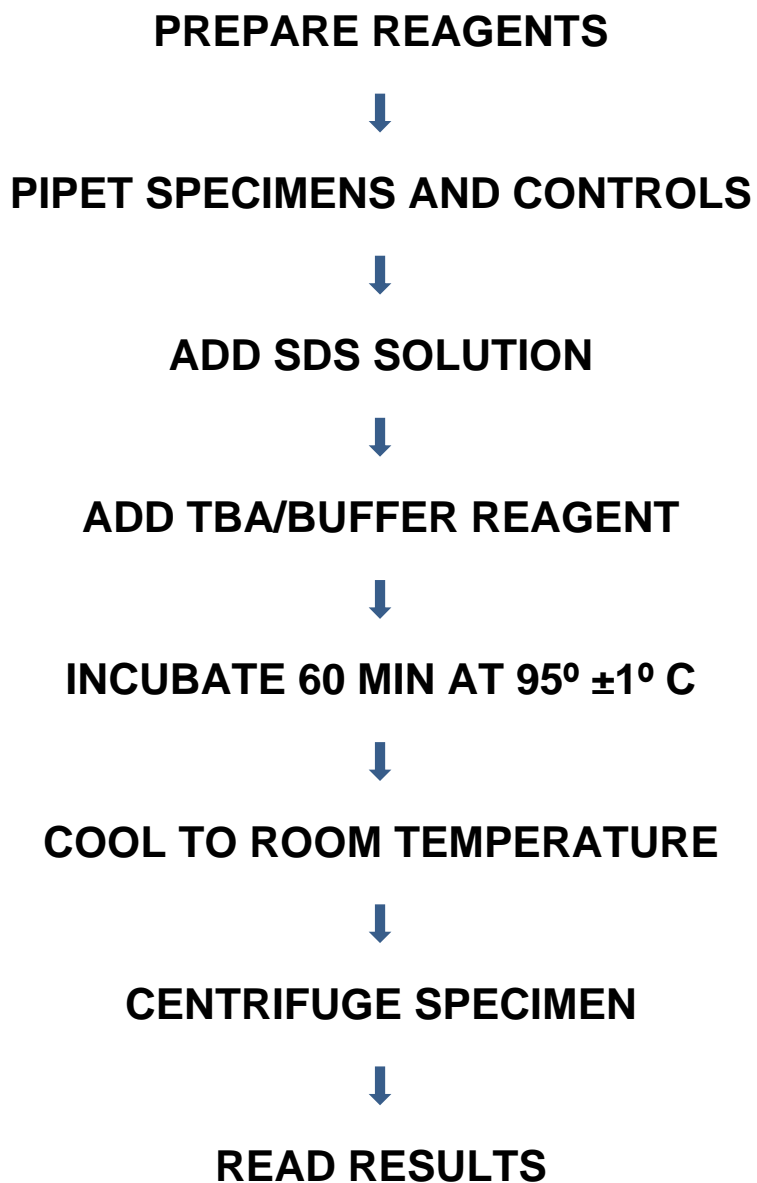
Using linear graph paper, plot mean MDA equivalents for each standard used on the X-axis versus the corresponding spectrophotometer reading on the Y-axis. Determine the concentration of MDA equivalents in nmol/ml in specimens by interpolation from the standard curve. Correct sample values for any other dilutions performed during specimen preparation.



REFERENCES

1. Yagi, K. Simple procedure for specific assay of lipid hydroperoxides in serum or plasma. *Free Radical and Antioxidant Protocols*, 108: 101- 106; 1998
2. Armstrong, D. and Browne, R. The analysis of free radicals, lipid peroxidases, antioxidant enzymes and compounds related to oxidative stress as applied to the clinical chemistry laboratory. *Free Radicals in Diagnostic Medicine*, 366:43-58; 1994
3. Lef'evre G., et.al. Evaluation of lipid peroxidation by measuring thiobarbituric acid reactive substances. *Annals de Biologie Clinique (Paris) May-June*; 56(3): 305-19; 1998
4. Janero, D. Malondialdehyde and thiobarbituric acid-reactivity as diagnostic indices of lipid peroxidation and peroxidative tissue injury. *Free Radical Biology & Medicine*, 9:515-540; 1998
5. Callaway, J.K. et. al. A reliable procedure for comparison of antioxidants in rat brain homogenates. *Journal of Pharmacology Toxicology Methods*, April; 39(3): 155-62; 1998
6. Jentzsh, AM., et. al. Improved analysis of malondialdehyde in human body fluids. *Free Radio Biol Med.*, 20(2): 251-6; 1996
7. Jo, C. et. al. Fluorometric analysis of 2-thiobarbituric acid reactive substances in turkey. *Pout. Sci.*, March; 77(3): 475-80; 1998
8. Villa-Caballero, L. et. al. Oxidative Stress. Should it be measured in the diabetic patient? *Gac Med Mex*, May-June; 136(3): 249-56; 2000
9. Hunnisett A. et. al. Lipoperoxides as an index of free radical activity in bone marrow transplant recipients. Preliminary Observations. *Biol Trace Elem Res*, Jan-March; 47(1-3): 125-32; 1995
10. Armstrong, D. and Browne, R. The analysis of free radicals, lipid peroxidases, antioxidant enzymes and compounds related to oxidative stress as applied to the clinical chemistry laboratory. *Free Radicals in Diagnostic Medicine*, 366:46; 1994
11. Kwon,T. and Watts,B. Malonaldehyde in aqueous solution and its role as a measure of lipid oxidation in foods. *Journal of Food Science*, 29:294-302; 1964
12. Braun,D. and Fromherz,P. Fluorescence interference-contrast microscopy of cell adhesion on oxidized silicon. *Applied Physics A*, 1997
13. Gidez,L. et al. Separation and quantitation of subclasses of human plasma high density lipoproteins by simple precipitation procedure. *Journal of Lipid Research*, 23:1206-1223; 1982
14. Armstrong,D. et al. In vitro screening for antioxidant activity. *Free Radical and Antioxidant Protocols*, 108:315-324; 1998
15. Boyum, A. Separation of leukocytes from blood and bone marrow. *Scandinavian Journal of Clinical Investigation* 21, Supplement 97; 1966.

PROCEDURAL FLOW CHART



The OXltek TBARS Assay Kit is for Research Purposes Only.



Product Manual

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