



## **Glutathione (GSSG/GSH), detection kit**

**Catalog # ADI-900-160**

**Sufficient Reagents for 384 tests with 4 x 96 well-plates**

For use with mammalian cell and tissue extracts, whole blood, plasma, urine, and saliva



## **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.**

## TABLE OF CONTENTS

Introduction.....	4
Principle.....	5
Materials Provided.....	6
Storage.....	6
Other Materials Needed.....	6
Reagent Preparation .....	7
Sample Handling .....	8
Procedural Notes.....	10
Calculation of Results .....	12
Trouble Shooting.....	14
References.....	15
Contact Information .....	16



Check our website for additional protocols, technical notes and FAQs.



Please contact Enzo Life Sciences Technical Support if necessary.

## INTRODUCTION

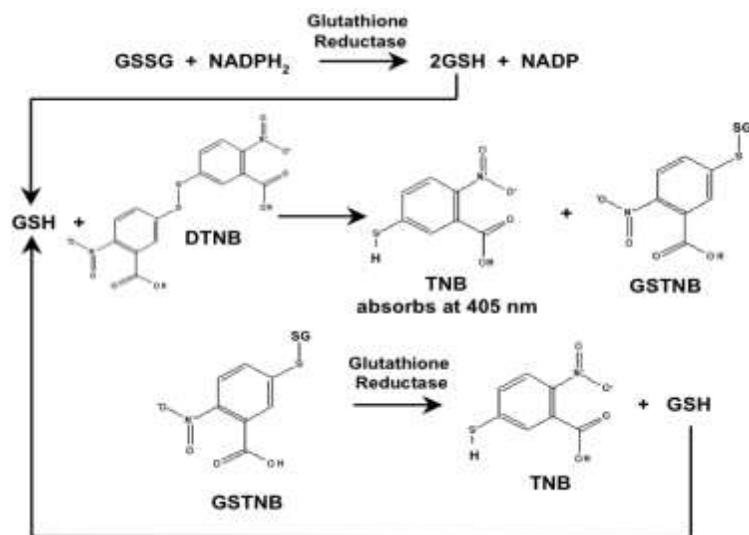
The Glutathione (GSSG/GSH), detection kit is a complete kit for the measurement of total, reduced, and oxidized glutathione concentration in mammalian cell and tissue extracts, whole blood, plasma, saliva, and urine.

Increased oxidative damage of proteins, lipids, and DNA by free radicals is one of the pathogenic mechanisms of diseases such as cancer, atherosclerosis, inflammation, and neurodegenerative disorders<sup>1,2</sup>. Glutathione, a major intracellular non-protein thiol, protects against free radical damage by providing reducing equivalents for several key antioxidant enzymes. In addition Glutathione acts as a scavenger of hydroxyl radicals and nascent oxygen. Elevated Glutathione levels are associated with a reduced rate of illness, higher levels of self-rated health, lower cholesterol, lower body mass index, and lower blood pressure among the elderly<sup>3</sup>. Glutathione provides a primary defense system for the removal of oxidants in the brain. Studies reveal a correlation between low Glutathione levels and damage to neurons that manufacture dopamine, suggesting a link to Parkinson's disease<sup>4</sup>. The concentration of Glutathione ranges from 10  $\mu\text{M}$  to 1 mM in cells and is in the micromolar range in plasma.

The Glutathione Assay utilizes a carefully optimized enzymatic recycling method for the quantification of glutathione. Glutathione Reductase reduces oxidized glutathione (GSSG) to reduced glutathione (GSH). As shown in Figure 1, the sulfhydryl group of GSH reacts with DTNB (5,5'-dithiobis-2-nitrobenzoic acid, Ellman's reagent) to produce a yellow colored 5-thio-2-nitrobenzoic acid (TNB) that absorbs at 405 or 414 nm. The rate of TNB production is directly proportional to the concentration of glutathione in the sample. The measurement of the absorbance of TNB at 405 or 414 nm provides an accurate estimation of glutathione in the sample.

## PRINCIPLE

1. Samples and diluted standards are added to wells of a 96-well plate.
2. Reaction Mix is added to the wells to initiate the reaction.
3. The plate is transferred to a plate reader and absorbance readings are taken at 405 nm or 414 nm every minute for 10 minutes.



## MATERIALS PROVIDED



Do not mix components from different kit lots or use reagents beyond the expiration date of the kit.



The physical, chemical, and toxicological properties of the chemicals and reagents contained in this kit may not yet have been fully investigated. Therefore, we recommend the use of gloves, lab coats, and eye protection while using any of these chemical reagents.

1. **Clear Microtiter Plate Pack, 2 Each, Catalog No. 80-2556**

Two packs of four plates of 96 wells. Clear uncoated solid plates.

2. **25X Assay Buffer, 12 mL, Catalog No. 80-1669**

3. **Reaction Mix, 8 bottles, Catalog No. 80-1667**

One vial is sufficient for 53 wells of a 96 well plate.

4. **GSSG, 2.5 mL (4  $\mu$ M), Catalog No. 80-1666**

5. **Glutathione Reductase, 80  $\mu$ L, Catalog No. 80-1668**

6. **Total Glutathione Assay Layout**

See Figure 2.

## STORAGE



All reagents should be stored at 4°C.



Check our website for additional protocols, technical notes and FAQs.

All components of this kit are stable at 4°C. All kit components are stable at their recommended storage temperatures until the kit's expiration date.

## OTHER MATERIALS NEEDED



For proper performance, use the insert provided with each individual kit received.

1. PBS
2. Distilled water
3. Meta-Phosphoric acid (Sigma-Aldrich Catalog No. 239275)
4. 4-Vinylpyridine (Sigma-Aldrich Catalog No. V3204-5ML)
5. Reagent alcohol
6. Microtubes, 0.5 and 1.5 MI
7. 15 mL conical tubes (adherent and suspension cell preparation)
8. 50 mL conical tubes (tissue preparation)
9. Precision pipettes for volumes between 5-1000  $\mu$ L
10. Multichannel pipettor for volumes between 1 - 50  $\mu$ L and 50  $\mu$ L - 200  $\mu$ L

11. Microplate reader capable of reading at 405 or 414 nm and taking readings every minute for ten minutes and exporting data to an Excel spreadsheet
12. Centrifuge and microfuge for processing samples
13. Sonicator or Homogenizer

## REAGENT PREPARATION

### 1. 1X Assay Buffer

Allow the assay buffer to come to room temperature for a minimum of 30 minutes. Dilute the 25X Assay Buffer to 1X (1:25) with distilled water. The 1X Assay Buffer is used to prepare dilutions of Total Glutathione Standard Curve and to dilute each experimental sample.

### 2. Reaction Mix

Reconstitute one or more bottles of Reaction Mix with 8 mL of distilled water. Periodically swirl the bottle gently over a 15 minute period to dissolve contents. **Immediately before use in the assay, vortex the vial of Glutathione Reductase and add 10  $\mu$ L of the Glutathione Reductase to the bottle of Reaction Mix.** Each bottle of Reaction Mix is sufficient for 53 wells in a 96-well plate, or little more than half a plate. Pool the reconstituted Reaction Mix together into one tube if more than one bottle is used.

### 3. 5% (w/v) meta-Phosphoric acid

Prepare 5% (w/v) meta-Phosphoric acid (Sigma Aldrich Catalog No. 239275, not provided) in distilled water.

### 4. 2M 4-Vinylpyridine

Prepare 2M 4-vinylpyridine solution (Sigma Aldrich Catalog No. V3204-5ML, not provided) by thawing at room temperature, then mixing 108  $\mu$ L 4-vinylpyridine with 392  $\mu$ L ethanol (solution should be prepared and subsequently used only in a chemical fume hood). Use immediately and discard any unused portion. Note: It is recommended that you use 4-vinylpyridine within 1 month of purchase and store at  $-20^{\circ}\text{C}$ . This reagent blocks free thiols present in the reaction, thus eliminating any contribution to the cycling reaction caused by GSH.

### 5. Experimental Samples

After preparing the samples as outlined in the Sample



All solutions must be prepared just before use.



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.



Warning: Prepare and use 2M 4-Vinylpyridine solution in a chemical fume hood.

Handling section that follows, dilute each sample 1:10 with 1X Assay Buffer. Some samples such as whole blood, liver, or red blood cells, may need to be diluted 1:20, 1:40, or more. It may be necessary to make serial dilutions of your extracts to obtain a satisfactory change in absorbance readings with time.

## SAMPLE HANDLING

All samples are treated with 5% (w/v) meta-Phosphoric acid to remove proteins which interfere with the assay.

### Cell Lysate Preparation

1. Detach **adherent cells** by gentle trypsinization. Count the cells and centrifuge at 300 x g for 10 minutes at 4°C. Wash the cells once with cold 1X PBS. Centrifuge **suspension cells** at 300x g for 10 minutes at 4°C. Discard the supernatant. Wash the cells once with cold 1X PBS.
2. Suspend the pellet with 500 µL of cold 5% (w/v) meta-Phosphoric acid per 2-5 x 10<sup>6</sup> cells. Mix thoroughly by repeated pipetting. Homogenize or sonicate the cell suspension and store on ice for 5 minutes.
3. Transfer the suspension to a 1.5 mL tube and centrifuge at 12,000-14,000 x g for 5 minutes at 4°C. Place the supernatant into a clean 1.5 mL tube. Store on ice if you intend to immediately assay for Glutathione, or freeze at -80°C for future use. To thaw, leave samples at room temperature until thawed.

### Tissue Lysate Preparation

1. Remove tissue. Repeatedly wash the tissue with cold isotonic saline (150 mM) or 1x PBS with 0.16 mg/mL heparin to prevent coagulation.
2. Blot tissue on filter paper and weigh.
3. Add ice-cold 5% (w/v) meta-Phosphoric acid (20 mL/g tissue) and homogenize using a cold glass or teflon pestle.
4. Centrifuge the homogenate at 12,000-14,000 x g for 10-15 minutes at 4°C.
5. Collect the clarified supernatant. Store on ice if you intend to immediately assay for Glutathione, or freeze at -80°C for future use. To thaw, leave samples at room temperature until thawed.

## **Erythrocyte Lysate Preparation**

1. Collect blood in Vacutainers containing heparin or sodium citrate as anticoagulant. Centrifuge at 3,000 x g for 10-15 minutes at 4°C.
2. Discard as much of the plasma supernatant as possible. Remove the white buffy coat (leukocytes) on the surface of the erythrocytes.
3. Resuspend the erythrocyte pellet in four volumes of ice-cold 5% (w/v) Meta-Phosphoric. Mix thoroughly and store on ice for 15 minutes.
4. Centrifuge the suspension at 12,000-14,000 x g for 10-15 minutes at 4°C.
5. Collect the clarified supernatant. Store on ice if you intend to immediately assay for Glutathione, or freeze at -80°C for future use. To thaw, leave samples at room temperature until thawed.

## **Whole Blood Lysate Preparation**

1. Collect blood in tubes containing heparin or sodium citrate as anticoagulant.
2. Add four volumes of ice-cold 5% (w/v) meta-Phosphoric acid. Mix thoroughly and store on ice for 15 minutes.
3. Centrifuge at 12,000-14,000 x g for 10-15 minutes at 4°C.
4. Collect the clarified supernatant. Store on ice if you intend to immediately assay for Glutathione, or freeze at -80°C for future use. To thaw, leave samples at room temperature until thawed.

## **Urine, Plasma, and Saliva Preparation**

1. Collect urine, plasma, or saliva and immediately add four volumes of ice-cold 5% (w/v) meta-Phosphoric acid. Mix thoroughly and store on ice for 15 minutes.
2. Centrifuge at 12,000-14,000 x g for 10-15 minutes at 4°C.
3. Collect the clarified supernatant. Store on ice if you intend to immediately assay for Glutathione, or freeze at -80°C for future use. To thaw, leave samples at room temperature until thawed.

## PROCEDURAL NOTES

### Total Glutathione Assay



All standards, controls, and samples should be run in triplicate.



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.



Pipet the reagents to the sides of the wells to avoid possible contamination.

1. Set up the Glutathione standard curve:
  - a) Add 50  $\mu\text{L}$  of 1X Assay Buffer to all the wells in rows A through E, columns 1, 2, and 3 of the microtiter plate (see Figure 2).
  - b) Add 50  $\mu\text{L}$  of the 4  $\mu\text{M}$  GSSG to wells A1, A2, and A3 with a multichannel pipettor. Mix well by pipetting the solution up and down at least ten times.
  - c) Transfer 50  $\mu\text{L}$  from wells A1, A2, and A3 to wells B1, B2, and B3, respectively. Mix well at least 10 times and transfer 50  $\mu\text{L}$  from row B to row C. Continue in this fashion to row D. Mix and discard the last 50  $\mu\text{L}$  from row D. Wells E1, E2, and E3 are set aside as Blank wells. The GSSG content in rows A, B, C, and D, is 100 pmoles/well, 50 pmoles/well, 25 pmoles/well, and 12.5 pmoles/well, respectively.
2. Add 50  $\mu\text{L}$  of your diluted experimental samples to the wells in columns 4 to 12. **Note:** It may be necessary to make serial dilutions of your extracts to obtain a satisfactory change in absorbance readings with time.
3. Prior to the next step, set up the parameters of your plate reader to measure absorbance at 405 nm or 414 nm and to read the required wells. Include a 10 second orbital shake prior to the initial read.
4. Using a multichannel pipettor, add 150  $\mu\text{L}$  of freshly-prepared Reaction Mix to each well.
5. Immediately record the absorbance in the wells at 405 nm or 414 nm using a plate reader at 1 minute intervals over a 10 minute period. **Note:** If you intend to use all the wells on one plate in the assay, it may be necessary to record the absorbance at 2 minute intervals.

## Oxidized Glutathione Assay

1. Add 1  $\mu\text{L}$  of 2 M 4-vinylpyridine per 50  $\mu\text{L}$  of sample and 4  $\mu\text{M}$  GSSG. Incubate for one hour at room temperature (cell lysates should be diluted at least 1:10 prior to 4-vinylpyridine treatment).
2. Serially dilute the 4-vinylpyridine-treated GSSG standard as described above in the total glutathione assay protocol (Step A.1).
3. Serially dilute your 4-vinylpyridine-treated experimental samples as described above in the total glutathione assay protocol (Step A.2).
4. Follow steps 3, 4, and 5 as described above for the total glutathione assay.

	1	2	3	4	5	6	7	8	9	10	11	12
A	GSSG 100 pmole	GSSG 100 pmole	GSSG 100 pmole	Sample 2 1:10	Sample 2 1:10	Sample 2 1:10	Sample 4 1:40	Sample 4 1:40	Sample 4 1:40	Sample 7 1:20	Sample 7 1:20	Sample 7 1:20
B	GSSG 50 pmole	GSSG 50 pmole	GSSG 50 pmole	Sample 2 1:20	Sample 2 1:20	Sample 2 1:20	Sample 5 1:10	Sample 5 1:10	Sample 5 1:10	Sample 7 1:40	Sample 7 1:40	Sample 7 1:40
C	GSSG 25 pmole	GSSG 25 pmole	GSSG 25 pmole	Sample 2 1:40	Sample 2 1:40	Sample 2 1:40	Sample 5 1:20	Sample 5 1:20	Sample 5 1:20	Sample 8 1:10	Sample 8 1:10	Sample 8 1:10
D	GSSG 12.5 pmole	GSSG 12.5 pmole	GSSG 12.5 pmole	Sample 3 1:10	Sample 3 1:10	Sample 3 1:10	Sample 5 1:40	Sample 5 1:40	Sample 5 1:40	Sample 8 1:20	Sample 8 1:20	Sample 5 1:20
E	Back - ground	Back - ground	Back - ground	Sample 3 1:20	Sample 3 1:20	Sample 3 1:20	Sample 6 1:10	Sample 6 1:10	Sample 6 1:10	Sample 8 1:40	Sample 8 1:40	Sample 8 1:40
F	Sample 1 1:10	Sample 1 1:10	Sample 1 1:10	Sample 3 1:40	Sample 3 1:40	Sample 3 1:40	Sample 6 1:20	Sample 6 1:20	Sample 6 1:20	Sample 9 1:10	Sample 9 1:10	Sample 9 1:10
G	Sample 1 1:20	Sample 1 1:20	Sample 1 1:20	Sample 4 1:10	Sample 4 1:10	Sample 4 1:10	Sample 6 1:40	Sample 6 1:40	Sample 6 1:40	Sample 9 1:20	Sample 9 1:20	Sample 9 1:20
H	Sample 1 1:40	Sample 1 1:40	Sample 1 1:40	Sample 4 1:20	Sample 4 1:20	Sample 4 1:20	Sample 7 1:10	Sample 7 1:10	Sample 7 1:10	Sample 9 1:40	Sample 9 1:40	Sample 9 1:40

**Figure 2.** Suggested 96-well plate format for GSSG Standards and samples.

## CALCULATION OF RESULTS

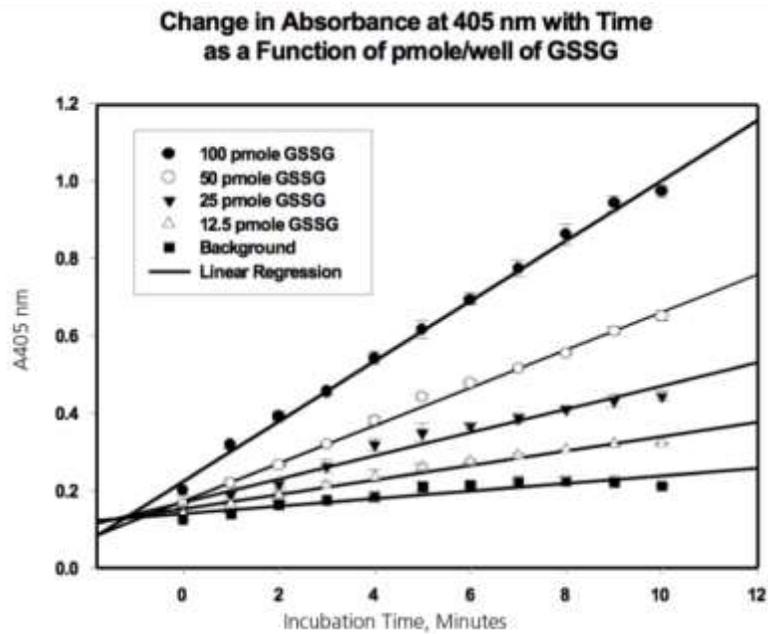
### Determination of Total Glutathione Concentration

1. Take the average of the triplicate absorbance readings for each standard, sample, and Blank at each time point.
2. Plot the average of each standard, sample, and background absorbance (A405 nm) versus incubation time and determine the slope from the linear portion of each curve (Figure 3).
3. Subtract the background slope from the slopes of the standards and the experimental samples.
4. Plot the net slopes of the GSSG standards versus pmoles of Glutathione (Figure 4).
5. Compare the net slopes of the experimental samples with those of the standard curve from Figure 4 to determine the pmoles of GSSG (equivalent to total glutathione) for each experimental sample.

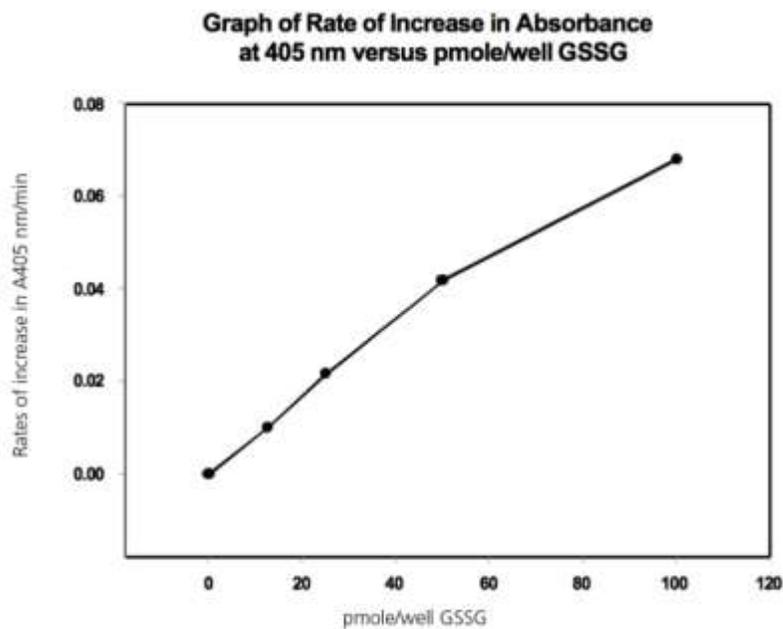
### Determination of Oxidized Glutathione Concentration

1. Follow the procedure described above for generating the Standard GSSG curves (Figures 3 and 4) for the 4-vinylpyridine treated standards.
2. Compare the net slopes of the 4-vinylpyridine-treated experimental samples with those of the 4-vinylpyridine-treated standard curve from Figure 4 to determine the pmoles of oxidized Glutathione for each experimental sample.
3. Subtract the pmole of oxidized glutathione in your sample from the pmole of total glutathione to obtain the pmole of reduced glutathione in your sample.

$$\text{Reduced GSH} = \text{Total glutathione} - \text{Oxidized GSSG}$$



**Figure 3.** Plot of absorbance at 405 nm versus incubation time as a function of pmoles of GSSG/well.



**Figure 4.** Rate of increase in the absorbance at 405 nm as a function of pmole/well GSSG.

## TROUBLE SHOOTING

<b>Problem</b>	<b>Cause</b>	<b>Solution</b>
<b>Erratic within sample values</b>	Inconsistent pipetting technique	Reproducible pipetting is absolutely required for consistent results
	Reaction Mix incompletely solubilized	Periodically vortex the Reaction Mix over a 15 minute period
<b>No color development in standards and samples</b>	Failure to add Glutathione Reductase to Reaction Mix	Vortex Glutathione Reductase briefly and add 10 $\mu$ l to each bottle of reconstituted Reaction Mix
	NADPH is oxidized	Contact Technical Support
<b>No color development in the samples but standards give color</b>	Concentration of glutathione in the samples is below the sensitivity of the assay	Extend incubation time to 30 minutes
<b>Sample absorbance values higher than those of standard curve</b>	Glutathione levels in the sample very high or other thiols are present	Further dilute your samples with 1X assay buffer and reassay

## REFERENCES

1. Onyango IG., and Khan SM. Oxidative stress, mitochondrial dysfunction, and stress signaling in Alzheimer's disease. *Curr Alzheimer Res.* 2006 3:339-349.
2. Medina S., et al. 2003. Antioxidants inhibit the human cortical neuron apoptosis induced by hydrogen peroxide, tumor necrosis factor alpha, dopamine and beta-amyloid peptide 1-42. *Free Radic Res.* 36:1179-84.
3. Julius M., et al. 1994. Glutathione and morbidity in a community-based sample of elderly. *J Clin Epidemiol.* 47: 1021-1026.
4. Mytilineou C, et al. 2002. Glutathione depletion and oxidative stress. *Parkinsonism Relat Disord.* 8:385-387.



# 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](http://www.enzolifesciences.com)**