

ARBOR ASSAYS™
Interactive Assay Solutions™



DetectX[®]

GLUTATHIONE Fluorescent Detection Kit

1 Plate Kit Catalog Number K006-F1

5 Plate Kit Catalog Number K006-F5

Species Independent

Sample Types Validated:

**Whole Blood, Serum, Plasma,
Erythrocytes, Urine, Cell Lysates
and Tissue Samples**

Please read this insert completely prior to using the product.
For research use only. Not for use in diagnostic procedures.

www.ArborAssays.com   

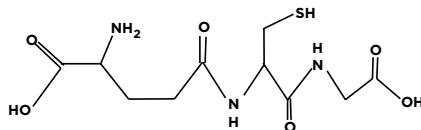
TABLE OF CONTENTS

Background	3
Assay Principle	4
Related Products	4
Supplied Components	5
Storage Instructions	5
Other Materials Required	6
Precautions	6
Sample Types	7
Sample Preparation	7-8
Reagent Preparation	9-10
Assay Protocol	11
Calculation of Results	12
Typical Data	13-15
Validation Data Sensitivity, Linearity, etc.	15-17
Kit Correlation	18
Warranty & Contact Information	19
Plate Layout Sheet	20

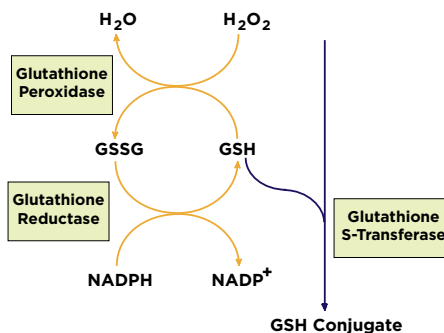


BACKGROUND

Glutathione (L- γ -glutamyl-L-cysteinylglycine; GSH) is the highest concentration non-protein thiol in mammalian cells and is present in concentrations of 0.5 - 10 mM¹. GSH plays a key role in many biological processes, including the synthesis of proteins and DNA, the transport of amino acids, and the protection of cells against oxidation. Harmful hydrogen peroxide cellular levels are minimized by the enzyme glutathione peroxidase (GP) using GSH as a reductant².



The oxidized GSH dimer, GSSG, is formed from GSH and peroxide by the GP reaction (see below). An important role of GSSG in the NF κ B activating signal cascade is suggested by the facts that the potent NF κ B inducer, tetradecanoyl phorbol acetate, increases intracellular GSSG levels and GSSG/GSH ratios³.



Glutathione S-transferases (GST) are an important group of enzymes that catalyze the nucleophilic addition of GSH to electrophiles. They are encoded by 5 gene families; 4 encode cytosolic GST and one encodes the microsomal form of GST. They have been implicated in a number of diseases. In asthma arachidonic acid is converted to unstable leukotriene A₄ (LTA₄). LTA₄ is either hydrated to form LTB₄ or it is conjugated to GSH by a GST, leukotriene C₄ synthase, to form leukotriene C₄, LTC₄ and its derivative LTD₄ are important molecules in bronchial asthma. Leukotriene C₄ synthase is therefore an important therapeutic target. It has also been shown that increased expression of GSTs can lead to drug resistance. Three glutathione adducts of the drug melphalan, used to treat ovarian cancer and multiple myeloma, have been isolated from reactions involving human microsomal GSTs.

1. Meister, A. "On the Discovery of Glutathione." Trends Biochem. Sci. 1988 13(5): 185-188.
2. Meister, A. "The Glutathione-Ascorbic Acid Antioxidant Systems in Animals" J. Biol. Chem. 1994 269:9397-9400.
3. Dröge W, et al., "Functions of Glutathione and Glutathione Disulfide in Immunology and Immunopathology" FASEB J., 1994 8:1131-1138.



ASSAY PRINCIPLE

The DetectX[®] Glutathione kit is designed to quantitatively measure glutathione (GSH), and oxidized glutathione (GSSG) present in a variety of samples. The kit is unique in that both free and oxidized glutathione are detected in the same well in the microtiter plate. No separation or washing is required. Total glutathione is the sum of GSSG plus GSH. Please read the complete kit insert before performing this assay. A GSH standard is provided to generate a standard curve for the assay and all samples should be read off the standard curve. The kit utilizes a proprietary non-fluorescent molecule, ThioStar[®], that will covalently bind to the free thiol group on GSH to yield a highly fluorescent product. After mixing the sample or standard with ThioStar[®] and incubating at room temperature for 15 minutes, the fluorescent product is read at 510 nm in a fluorescent plate reader with excitation at 390 nm. The concentration of the GSH in the sample is calculated, after making a suitable correction for any dilution of the sample, using software available with most fluorescence plate readers.

Free glutathione, GSH, is read first after 15 minutes, followed by addition of a reaction mixture that converts all the oxidized glutathione, GSSG, into free GSH, which then reacts with the excess ThioStar[®] to yield the signal related to Total GSH content. The total concentration of GSH generated in the sample is calculated from the generated signal. We have provided a 96 well plate for measurement but this assay is adaptable for higher density plate formats. The end user should ensure that their HTS black plate is suitable for use with these reagents prior to running samples.

RELATED PRODUCTS

Kits	Catalog No.
Glutathione Colorimetric Detection Kit	K006-H1
Glutathione S-Transferase Fluorescent Activity Kit	K008-F1
Glutathione Reductase Fluorescent Activity Kit	K009-F1

Reagents	Catalog No.
Glutathione Mouse Monoclonal Antibody, 50 µg Mouse IgG2a, Clone L4H raised to glutathione conjugated to KLH Applications: Western blotting, Immunoassay and Immunoprecipitation	A001-50UG
DyLight[®] 488 Glutathione Mouse Monoclonal Antibody, 50 µg Purified monoclonal labeled with a stable FITC like fluorescent dye Applications: Flow cytometry and direct immunofluorescence	A001F-50UG



SUPPLIED COMPONENTS

Black Half Area 96 Well Plate

See: www.arborassays.com/resources/#general-info for plate dimension data.

Kit K006-F1 **or** -F5 1 **or** 5 Each Catalog Number X023-1EA **or** -5EA

Glutathione Standard

Glutathione at 250 μ M in a special stabilizing solution.

Kit K006-F1 **or** -F5 100 μ L **or** 300 μ L Catalog Number C018-100UL **or** -300UL

ThioStar[®] Detection Reagent

ThioStar thiol detection substrate stored in a ziploc pouch with desiccant. Reconstitute with dry DMSO.

Kit K006-F1 2 Plastic vials Catalog Number C021-1EA

Kit K006-F5 4 Glass vials Catalog Number C036-1EA

Dry DMSO

Dry Dimethyl sulfoxide solvent over molecular sieves. May be stored at room temperature.

Kit K006-F1 **or** -F5 4 mL **or** 20 mL Catalog Number X022-4ML **or** -20ML

Assay Buffer Concentrate

A 2X buffer concentrate containing detergents and stabilizers that must be diluted with deionized or distilled water.

Kit K006-F1 **or** -F5 35 mL **or** 200 mL Catalog Number X051-35ML **or** -200ML

NADPH Concentrate

Reduced β -nicotinamide adenine dinucleotide 2'-phosphate (NADPH) as a stable solution.

Kit K006-F1 **or** -F5 300 μ L **or** 1.4 mL Catalog Number X044-300UL **or** -1.4ML

Glutathione Reductase Concentrate

Glutathione Reductase (GR) as a stable solution.

Kit K006-F1 **or** -F5 300 μ L **or** 1.4 mL Catalog Number X048-300UL **or** -1.4ML

Oxidized Glutathione Control

Oxidized Glutathione (GSSG) in a special stabilizing solution. This is an optional control solution to ensure NADPH/GR performance.

Kit K006-F1 **or** -F5 300 μ L Catalog Number C020-300UL

STORAGE INSTRUCTIONS

All components of this kit should be stored at 4°C until the expiration date of the kit. DMSO, when stored at 4°C, will freeze. Can be stored tightly capped at room temperature.



OTHER MATERIALS REQUIRED

Distilled or deionized water

Repeater pipet with disposable tips capable of dispensing 25 μ L.

Aqueous 5-sulfo-salicylic acid dihydrate (SSA) solution at 5% weight/volume (1g of SSA per 20 mL of water) for treating samples to remove protein. We recommend Sigma-Aldrich Catalog Number S2130.

Fluorescence 96 well plate reader capable of reading fluorescent emission at 510 nm, with excitation at 390 nm. Please contact your plate reader manufacturer for suitable filter sets. Set plate parameters for a 96-well Corning Costar 3650 plate. See: www.arborassays.com/resources/#general-info for plate dimension data.

The sensitivity of fluorescent assays is dependant on the capabilities of the plate reader. If your plate reader has adjustable gain you can modify the signals obtained from the assay by increasing or decreasing the gain settings, by changing the aperture settings for monochromator based readers, or by changing the band pass width of the emission and/or excitation filters on some readers. Please review the plate reader manual for details.

Signals expressed by plate readers are Relative Fluorescent Units (RFU) and the values given in the insert were obtained on our plate readers. The RFU numbers you obtain may be different from these, but the assay results should be similar.

Software for converting raw relative fluorescent unit (FLU) readings from the plate reader and carrying out four parameter logistic curve (4PLC) fitting. Contact your plate reader manufacturer for details.

PRECAUTIONS

As with all such products, this kit should only be used by qualified personnel who have had laboratory safety instruction. The complete insert should be read and understood before attempting to use the product.

Sulfosalicylic acid is a strong acid solution and should be treated like any other laboratory acid.

Dimethyl sulfoxide is a powerful aprotic organic solvent that has been shown to enhance the rate of skin absorption of skin-permeable substances. Wear protective gloves when using the solvent especially when it contains dissolved chemicals. *NOTE: DMSO can dissolve certain plastics used in troughs used for holding solutions for multichannel pipets.*

ThioStar[®] Thiol Detection Reagent should be stored at 4°C in the desiccated pouch. Allow desiccated pouch to warm to room temperature prior to opening. ThioStar will react with strong nucleophiles. Buffers containing the preservatives sodium azide, Proclin[™] and Kathon[™] will react with the substrate.

Reconstituted ThioStar in DMSO stored at 4°C in the desiccated pouch. It can be used up to 2 months later. The background on the reconstituted ThioStar will increase slowly over time but the increase will not affect the assay results obtained.



SAMPLE TYPES

GSH is identical across species and we expect this kit may measure GSH from sources other than human. The end user should evaluate recoveries of GSH in samples from other species being tested.

If samples need to be stored after collection, we recommend storing them at -70°C or lower, preferably after being frozen in liquid nitrogen. This assay has been validated for human whole blood, serum, EDTA and heparin plasma, urine, and isolated erythrocytes. Most cell lysates and tissue homogenates should also be compatible. Samples containing visible particulate should be centrifuged prior to using.

All samples will be deproteinized with 5% SSA (see page 6 for preparation), please see sample specific information below for details. This treatment removes any protein thiols present in the samples and also slows oxidation of free GSH.

SAMPLE PREPARATION

All samples must be treated with the SSA solution prepared on page 6. All of the SSA treated centrifuged supernatants must have their SSA concentration brought down to 1% SSA by dilution with Assay Buffer. Further dilutions of the sample, using Sample Diluent (see page 9 for preparation), may be necessary to allow the GSH concentration to be measurement in the assay. Detailed instructions follow.

All samples and standards must be in Sample Diluent before starting the assay.

Use all samples within 2 hours of dilution.

Whole Blood, Serum, EDTA or Heparin Plasma, or Urine

Thoroughly mix sample with an equal volume of cold 5% SSA. Incubate for 10 minutes at 4°C . Centrifuge at 14,000 rpm for 10 minutes at 4°C . Collect the supernatant. If the supernatant contains particulates, re-centrifuge the supernatant for 15 minutes and collect the clarified second supernatant. Samples can be stored in aliquots at $\leq -70^{\circ}\text{C}$ or analyzed immediately. At this point the SSA concentration will be 2.5%.

The supernatant must be diluted 1:2.5 with Assay Buffer by mixing one part with 1.5 parts of Assay Buffer. The SSA concentration will be 1%. The sample will have been diluted 1:5 at this point.

All final dilutions are to be made in Sample Diluent. Treated Whole Blood must be further diluted at least 1:20 for a recommended final dilution of $\geq 1:100$. For Treated Plasma and Treated Urine a final dilution of $\geq 1:5$ is recommended, but further dilutions in Sample Diluent may be necessary.

Tissue Samples

Fresh tissue is washed with ice cold PBS to remove blood then blotted on filter paper before recording wet weight. *NOTE: Samples that have been frozen will contain lysed cells. The PBS wash may contain substantial amounts of GSH and/or GSSG.*

- For Samples Where a Protein Determination is to be Obtained: Homogenize at 10 mg/250 μ L in ice cold 100mM phosphate buffer, pH 7. Centrifuge at 14,000 rpm for 10 minutes at 4°C and remove an aliquot of the supernatant for protein determination. Thoroughly mix a second aliquot of the supernatant with an equal volume of cold 5% SSA. Incubate for 10 minutes at 4°C. Centrifuge at 14,000 rpm for 10 minutes at 4°C to remove precipitated protein. Collect the supernatant. The supernatant must be diluted 1:2.5 with Assay Buffer by mixing one part with 1.5 parts of Assay Buffer. The SSA concentration will be 1%.
- For Samples Not Requiring a Protein Determination: Homogenize at 10 mg/250 μ L in ice cold 5% SSA, incubate at 10 minutes at 4°C, then centrifuge at 14,000 rpm for 10 minutes at 4°C to remove precipitated protein. Collect the supernatant. The supernatant must be diluted 1:5 with Assay Buffer by mixing one part with 4 parts of Assay Buffer. The SSA concentration will be 1%.

Further sample dilutions must be determined by the end-user since it will be dependent upon the tissue type and the amount of tissue used. These dilutions must be made in the prepared Sample Diluent.

Erythrocytes, Red Blood Cells (RBC's)

Collect blood with heparin or EDTA. Centrifuge the sample, remove and discard the plasma and white cell layer. Wash the RBC's 2 times by suspending in 3 volumes of isotonic saline (0.9%), centrifuging at 600 x g for 10 minutes and discarding the saline wash.

After the 2 washes, mix 250 μ L RBC's with 1mL of cold 5% SSA. Incubate for 10 minutes at 4°C. Centrifuge at 14,000 rpm for 10 minutes at 4°C. Collect the supernatant. At this point the SSA concentration will be 4%. The supernatant must be diluted 1:4 with Assay Buffer by mixing one part with 3 parts of Assay Buffer. The SSA concentration will now be 1%. The sample will have been diluted 1:20 at this point. Further dilutions are made in Sample Diluent. *NOTE: Human RBC's require a final dilution of 1:100-1:200 to read within the standard curve.*

Cell Lysates

Washed cell pellets are resuspended at 1-10x10⁶ cells/mL in cold 5% SSA (we used Jurkats at 5x10⁶ cells/mL) and are lysed and deproteinized by vigorous vortexing, freeze/thaw cycling or other suitable disruption method. Incubate cells at 4°C for 10 minutes followed by centrifugation for 10 minutes at 14,000 rpm and 4°C. *NOTE: Samples that have been frozen will contain lysed cells. The PBS wash may contain substantial amounts of GSH and/or GSSG.*

The centrifuged supernatants must be diluted 1:5 with Assay Buffer by mixing one part with 4 parts of Assay Buffer. The SSA concentration will be 1%. The sample will have been diluted 1:5 at this point. Further sample dilutions must be done in Sample Diluent and need to be determined by the end-user since it will be dependent upon the cell type and number of cells used. The recommended final dilution is \geq 1:20.

Use all samples within 2 hours of dilution.



REAGENT PREPARATION

Allow the kit reagents to come to room temperature for 30 minutes. Ensure that all samples have reached room temperature and have been diluted as appropriate prior to running them in the kit.

Assay Buffer

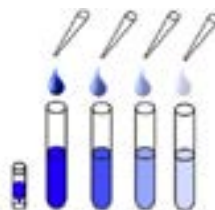
Prepare the Assay Buffer by diluting the supplied Assay Buffer Concentrate with an equal volume of deionized water. Mix thoroughly. Stable at 4°C for 3 months.

Sample Diluent

Prepare the Sample Diluent by diluting one part 5% SSA 1:5 with four parts diluted Assay Buffer and vortex thoroughly. The pH of the Sample Diluent **must** be > 6. Sample Diluent can be stored at 4°C for one month.

Standard Preparation

GSH Standards are prepared by labeling eight test tubes as #1 through #8. Briefly vortex to mix and then spin the vial of standard in a microcentrifuge to ensure contents are at bottom of vial. Pipet 450 μL of Sample Diluent into tube #1 and 250 μL into tubes #2 to #8. Carefully add 50 μL of the Glutathione Standard to tube #1 and vortex completely. Take 250 μL of the GSH solution in tube #1 and add it to tube #2 and vortex completely. Repeat this for tubes #3 through #8. The concentration of GSH in tubes 1 through 8 will be 25, 12.5, 6.25, 3.125, 1.56, 0.781, 0.391 and 0.195 μM .



Use all Standards within 1 hour of preparation.

	Std 1	Std 2	Std 3	Std 4	Std 5	Std 6	Std 7	Std 8
Sample Diluent Vol (μL)	450	250	250	250	250	250	250	250
Addition	Stock	Std 1	Std 2	Std 3	Std 4	Std 5	Std 6	Std 7
Vol of Addition (μL)	50	250	250	250	250	250	250	250
Final Conc (μM)	25	12.5	6.25	3.125	1.56	0.781	0.391	0.195



Control Preparation (Optional)

This optional control solution for ensuring complete conversion of GSSG to GSH can be prepared by adding 5 μL of Oxidized Glutathione Control to 245 μL of Sample Diluent. **Use within 2 hours.**

The Control Preparation ensures that the NADPH and Glutathione Reductase system prepared below in the Reaction Mixture section will adequately reduce GSSG to GSH. If this optional control is run it should yield a value for Total Glutathione of approximately $10 \pm 2 \mu\text{M}$.

ThioStar[®] Detection Reagent

Allow the ziploc bag to warm **completely** to room temperature prior to opening and remove the vial of ThioStar Reagent. Add the volume of DMSO provided to the vial according to the table below. Vortex thoroughly. Store any unused reconstituted Detection Reagent at 4°C in the ziploc pouch with desiccant and use within 2 months.

	Kit K006-F1	Kit K006-F5
Vial Part Number	CO21-1EA, Plastic vial	CO36-1EA, Glass vial
Volume of DMSO to add per vial	1.5 mL	3.5 mL
For # of Wells	Up to 60	Up to 140

Reaction Mixture

Prepare the Reaction Mixture by vortexing the vials of Glutathione Reductase and NADPH Concentrates and then diluting one part each NADPH and Glutathione Reductase Concentrates 1:10 into eight parts Assay Buffer. Vortex thoroughly. See Table for suitable volumes. Store any unused Reaction Mixture at 4°C in an amber vial for no more than 2 days.

Reaction Mix Dilution Table

	1/2 Plate	One Plate
NADPH Concentrate	150 μL	275 μL
Glutathione Reductase Concentrate	150 μL	275 μL
Assay Buffer	1.2 mL	2.2 mL



ASSAY PROTOCOL - FREE AND TOTAL GSH

We recommend that all standards and samples be run in duplicate to allow the end user to accurately determine GSH concentrations.

1. Use the plate layout sheet on the back page to aid in proper sample and standard identification. Set plate parameters for a 96-well Corning Costar 3686 plate. See: www.arborassays.com/resources/#general-info for plate dimension data.
2. Pipet 50 μ L of treated samples, standards or control into wells in the plate.
3. Pipet 50 μ L of Sample Diluent into Zero wells in the plate.
4. Add 25 μ L of the ThioStar Reagent to each well using a repeater pipet.
5. Gently tap the sides of the plate to ensure adequate mixing of the reagents.
6. Incubate at room temperature for 15 minutes.
7. Read the fluorescent signal from each well in a plate reader capable of reading the fluorescent emission at 510 nm with excitation at 370-410 nm. This data will be used to determine Free GSH concentration.
8. Add 25 μ L of the Reaction Mixture to each of the wells using a repeater pipet.
9. Gently tap the sides of the plate to ensure adequate mixing of the reagents.
10. Incubate at room temperature for 15 minutes.
11. Read the fluorescent emission at 510 nm with excitation at 370-410 nm. This data will be used to determine Total GSH concentration.

Total GSH Content Only

Total GSH content can be determined directly by leaving out steps 5, 6 and 7.



CALCULATION OF RESULTS

Average the duplicate FLU readings for each standard and sample. Create a standard curve by reducing the data using the 4PLC fitting routine on the plate reader, after subtracting the mean FLUs for the zero standard. The sample concentrations obtained should be multiplied by the dilution factor to obtain neat sample values.

Or use the online tool from MyAssays to calculate the data:

www.myassays.com/arbor-assays-glutathione-fluorescent-detection-kit.assay



MyAssays

Free glutathione (GSH) concentrations are calculated from the data obtained from step 7 on page 11 utilizing the curve fitting routine supplied with the plate reader.

Total glutathione concentrations of the samples are calculated from the data obtained from step 11 on page 11 utilizing the curve fitting routine supplied with the plate reader. Ensure that the Reaction Mixture is added to all the wells used, including the standard and control wells. The volumes must be the same in the standard, control and samples wells.

Oxidized glutathione (GSSG) concentrations are obtained by subtracting the Free GSH levels from the Total GSH concentrations and dividing by 2. See Below:

$$\text{GSSG} = \frac{(\text{Total GSH} - \text{Free GSH})}{2}$$

2

Total GSH Note:

When Free GSH and Total GSH levels are almost identical, we suggest that you block the free GSH by addition of 2-Vinylpyridine (2VP) to an aliquot of the sample. 2VP is prepared by adding 27 μL of 2-vinylpyridine (such as Sigma Catalog Number 132292) to 98 μL of ethanol. Use immediately and discard remaining unused solutions.

2VP is TOXIC and may cause burns. 2VP solutions should be prepared in a fume hood. Use immediately and discard remaining unused solutions by mixing with copious amounts of water.

To 250 μL of 5% SSA treated samples add 5 μL of the ethanolic solution of 2VP and allow to incubate at room temperature for 1 hour. The 2VP treated samples should then be diluted in Assay Buffer and Sample Diluent according to the dilutions recommended for each sample type on pages 7 and 8 prior to using in the assay.



*The MyAssays logo is a registered trademark of MyAssays Ltd.

WEB INSERT

180117

EXPECT ASSAY ARTISTRY

TYPICAL DATA - FREE GSH

Sample	Mean FLU	Net FLU	GSH Conc. (μM)
Zero	299	0	0
Standard 1	40,945	40,646	25
Standard 2	19,737	19,438	12.5
Standard 3	10,006	9,707	6.25
Standard 4	5,269	4,970	3.125
Standard 5	2,671	2,372	1.56
Standard 6	1,571	1,272	0.781
Standard 7	1,009	710	0.391
Standard 8	639	394	0.195
Sample 1	1,568	1,269	0.77
Sample 2	5,448	5,149	3.41

Always run your own standard curve for calculation of results. Do not use this data.



TYPICAL DATA - TOTAL GSH

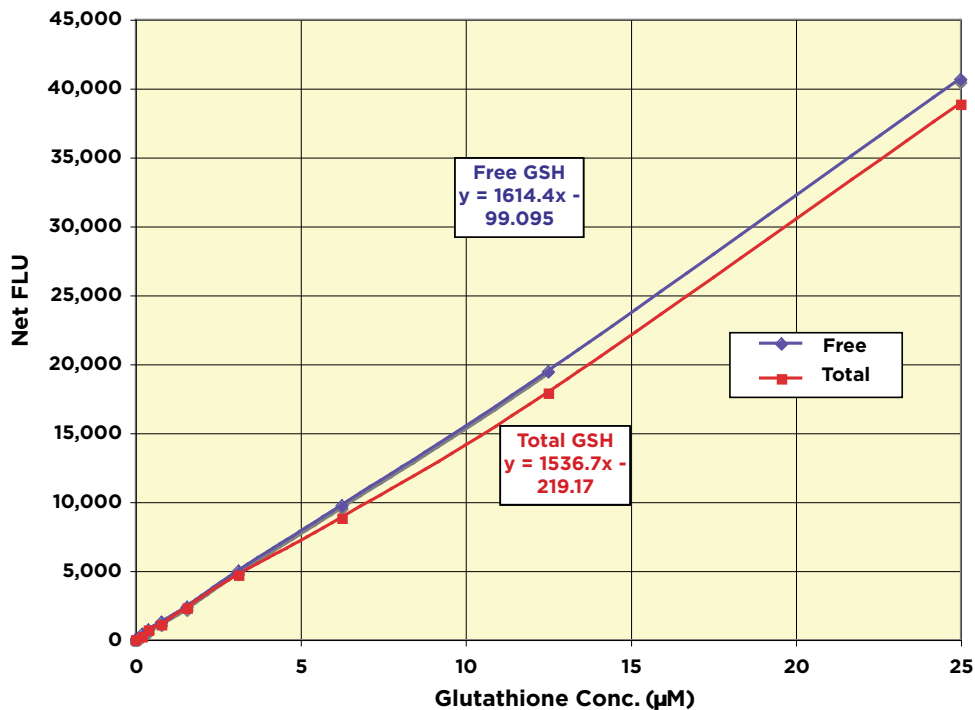
Sample	Mean FLU	Net FLU	GSH Conc. (μM)
Zero	1,127	0	0
Standard 1	39,976	38,849	25
Standard 2	19,034	17,907	12.5
Standard 3	9,958	8,831	6.25
Standard 4	5,814	4,687	3.125
Standard 5	3,429	2,302	1.56
Standard 6	2,209	1,082	0.781
Standard 7	1,816	689	0.391
Standard 8	1,339	212	0.195
Sample 1	4,553	3,426	2.53
Sample 2	10,428	9,301	6.64

Always run your own standard curve for calculation of results. Do not use this data.

See NOTE on Page 12 concerning Free and Total GSH levels in samples.



Typical Standard Curves



Always run your own standard curves for calculation of results. Do not use these data.

See NOTE on Page 12 concerning Free and Total GSH levels in samples.

VALIDATION DATA

Sensitivity and Limit of Detection

Sensitivity was calculated by comparing the FLUs for twenty wells run for each of the zero and standard #8. The detection limit was determined at two (2) standard deviations from the zero along the standard curve.

Sensitivity was determined as 45 nM in the Free GSH and 48 nM in the Total GSH assays.

The Limit of Detection was determined in a similar manner by comparing the FLUs for twenty wells run for each of the zero and a low concentration human serum sample.

The Limit of Detection was determined as 38 nM in the Free GSH and 42 nM in the Total GSH assays.

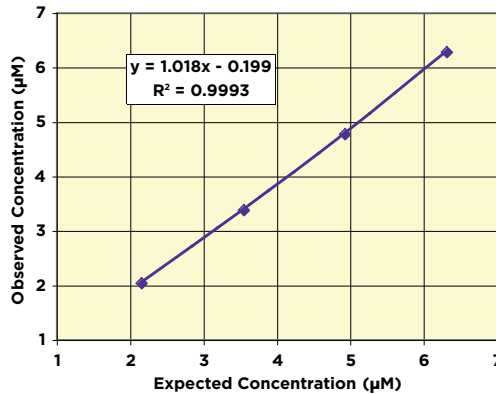


Linearity

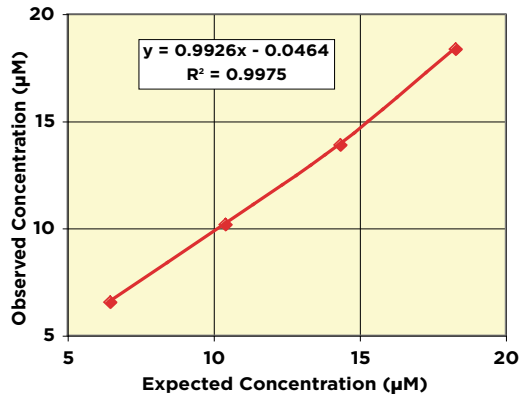
Linearity was determined by taking human RBCs at two different concentrations and mixed in the ratios given below. The measured concentrations were compared to the expected values based on the ratios used.

High RBC Sample	Low RBC Sample	Observed Conc. (µM)		Expected Conc. (µM)		% Recovery	
		Free	Total	Free	Total	Free	Total
80%	20%	6.28	18.35	6.32	18.29	99.3%	100.3%
60%	40%	4.77	13.89	4.93	14.35	96.7%	96.8%
40%	60%	3.38	10.18	3.55	10.41	95.3%	97.8%
20%	80%	2.04	6.55	2.16	6.47	94.5%	101.2%
Mean Recovery						96.5%	99.0%

Free GSH Linearity



Total GSH Linearity



Intra Assay Precision

Two each of SSA treated human urine and whole blood samples were further diluted in 1% SSA Sample Diluent and run in replicates of 20 in an assay. The mean and precision of the calculated GSH concentrations were:

Sample	GSH Conc. (μM)		%CV	
	Free	Total	Free	Total
1	1.27	2.30	4.0	4.7
2	2.00	3.80	3.1	4.7
3	8.33	9.77	4.6	2.7
4	3.89	4.45	3.0	2.3

Inter Assay Precision

Two each of SSA treated human urine and blood samples were further diluted in 1% SSA Sample Diluent and run in duplicates in twenty assays run over multiple days by two operators. The mean and precision of the calculated GSH concentrations were:

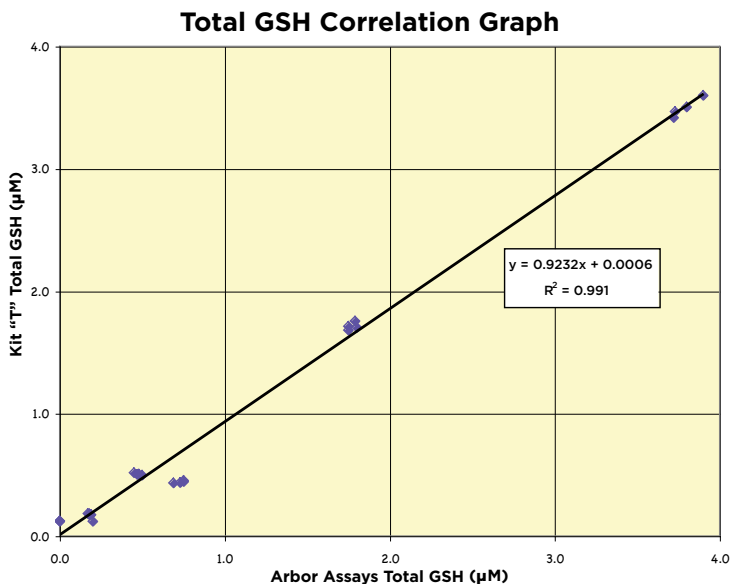
Sample	GSH Conc. (μM)		%CV	
	Free	Total	Free	Total
1	1.30	2.40	8.6	8.3
2	1.83	3.57	14.7	10.0
3	9.38	11.67	6.0	6.0
4	4.89	5.89	7.2	8.0



KIT CORRELATION DATA

We purchased and compared a popular colorimetric total glutathione assay kit (kit "T") that uses Ellman's reagent to detect free glutathione in the sample. Initial experiments used random human urine samples that were processed as described in each kit insert. With kit "T", the values obtained for urine after the recommended treatment with 4 volumes of 5% metaphosphoric acid and subsequent 10 fold dilution with assay buffer put all the values well below the lowest standard. However, the urine samples run in the DetectX[®] kit gave Total GSH values between 0.63 and 4.04 μM .

We also took a Jurkat cell pellet and processed the cells either through the 5% metaphosphoric acid treatment for the kit "T" Ellman's based test or as described on page 9 for the DetectX[®] kit. Cell samples ranged from 25 to 0.78×10^6 cell/mL. Twenty-four samples were run according to manufacturers directions for both kits and the correlation of these samples is shown below.



Many of the cell lysate values for the Ellman's based kit, kit "T", read either below the lowest standard (0.25 μM) or above the highest one (2 μM). This data was calculated via extrapolation from the kinetic method required by kit "T". The lysate values for the DetectX[®] kit were calculated directly from the endpoint standard curve.



LIMITED WARRANTY

Arbor Assays warrants that at the time of shipment this product is free from defects in materials and workmanship. This warranty is in lieu of any other warranty expressed or implied, including but not limited to, any implied warranty of merchantability or fitness for a particular purpose.

We must be notified of any breach of this warranty within 48 hours of receipt of the product. No claim shall be honored if we are not notified within this time period, or if the product has been stored in any way other than outlined in this publication. The sole and exclusive remedy of the customer for any liability based upon this warranty is limited to the replacement of the product, or refund of the invoice price of the goods.

CONTACT INFORMATION

For details concerning this kit or to order any of our products please contact us:

Arbor Assays

1514 Eisenhower Place
Ann Arbor, Michigan 48108 USA

Phone: 734-677-1774

Fax: 734-677-6860

Web: www.ArborAssays.com

E Mail Addresses:

Info@ArborAssays.com

Orders@ArborAssays.com

Technical@ArborAssays.com

Contracts@ArborAssays.com



OFFICIAL SUPPLIER TO ISWE

Arbor Assays and the International Society of Wildlife Endocrinology (ISWE) signed an exclusive agreement for Arbor Assays to supply ISWE members with EIA kits for wildlife conservation research.

DetectX[®], ThioStar[®] and the Arbor Assays logo are all registered trademarks.

www.ArborAssays.com

WEB INSERT

180117 ARBOR ASSAYS



H	G	F	E	D	C	B	A	
								1
								2
								3
								4
								5
								6
								7
								8
								9
								10
								11
								12



Printed on Forest Stewardship Council certified paper

