



ARBOR ASSAYS™
Interactive Assay Solutions™



DetectX[®]

Glutathione S-Transferase Fluorescent Activity Kit

1 Plate Kit Catalog Number K008-F1

Species Independent

Sample Types Validated:

Serum, Plasma, Urine and Cell Lysates

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

www.ArborAssays.com   

K008-F1 WEB 190503

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
Reagent Preparation	8
Assay Protocol	9
Calculation of Results	9
Typical Data	10-11
Validation Data Sensitivity, Linearity, etc.	11-13
Sample Values	14
Interferents and Kinetic Assay	14
Warranty & Contact Information	15
Plate Layout Sheet	16



BACKGROUND

The Glutathione S-Transferase (GST) family of isozymes function to detoxify and neutralize a wide variety of electrophilic molecules by mediating their conjugation with reduced glutathione¹. Human GSTs are encoded by several families, and are expressed in almost all tissues.

Given its pivotal role in ameliorating oxidative stress/damage, GST activity has been repeatedly investigated as a biomarker for arthritis, asthma, COPD, and multiple forms of cancer, as well as an environmental marker³⁻⁷. Examination of GST isoforms and activity in human cancers, tumors and tumor cell lines has revealed the predominance of the acidic pi class. Furthermore, this activity is thought to substantially contribute to the innate or acquired resistance of specific neoplasms to anticancer therapy^{8,9}.

1. Habig, W, et al. "Glutathione S-Transferases: The First Enzymatic Step in Mercapturic Acid Formation" J. Biol. Chem. 1974 249(22):7130-7139.
2. Cook, JA, et al., "Differential Specificity of Monochlorobimane for Isozymes of Human and Rodent Glutathione S-Transferases" Cancer Res. 1991 51:1606-1612.
3. Dalle-Donne, I, et al. "Biomarkers of Oxidative Damage in Human Disease" Clinical Chemistry, 2006 52(4):601-623.
4. Surapneni, KM & VSC Gopan, "Lipid Peroxidation and Antioxidant Status in Patients with Rheumatoid Arthritis" Ind. J. Clin. Biochem. 2008 23(1):41.44.
5. Mohan, SK & O Venkataramana. "Status of Lipid Peroxidation, Glutathione, Ascorbic Acid, Vitamin E and Antioxidant Enzymes in Patients with Osteoarthritis" Ind. J. Med. Sci. 2007 61:9-14.
6. Ferrandina, G., et al., "Glutathione S-Transferase Activity in Epithelial Ovarian Cancer: Association with Response to Chemotherapy and Disease Outcome" Ann. Oncol. 1997 8:343-350.
7. Otitoju, O & Onwarah, INE. "Glutathione S-transferase (GST) Activity as a Biomarker in Ecological Risk Assessment of Pesticide Contaminated Environment" African J. Biotech. 2007 6(12) 1455-1459.
8. Shea, TC, et al., "Identification of an Anionic Form of Glutathione Transferase Present in Many Human Tumors and Human Tumor Cell Lines" Cancer Res. 1988 48:527-533.
9. Shea, TC, et al., "Glutathione Transferase Activity and Isozyme Composition in Primary Human Breast Cancers" Cancer Res. 1990 50:6848-6853.

ASSAY PRINCIPLE

The DetectX® Glutathione S-Transferase Fluorescent Activity Kit is designed to quantitatively measure the activity of GST present in a variety of samples. Please read the complete kit insert before performing this assay. A GST 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 non-fluorescent molecule that is a substrate for the GST enzyme which covalently attaches to glutathione (GSH) to yield a highly fluorescent product. Mixing the sample or standard with the supplied Detection Reagent and GSH and incubating at room temperature for 30 minutes yields a fluorescent product which is read at 460 nm in a fluorescent plate reader with excitation at 390 nm. The activity of the GST in the sample is calculated, after making a suitable correction for any dilution of the sample, using software available with most fluorescence plate readers.

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 Fluorescent Detection Kit	K006-F1/F5
Glutathione Colorimetric Detection Kit	K006-H1
Glutathione Reductase Fluorescent Activity Kit	K009-F1
Urinary Creatinine Detection Kits (2 or 10 Plates)	K002-H1/H5

Reagents	Catalog No.
Glutathione Mouse Monoclonal Antibody, 50 µg Mouse IgG _{2a} , Clone L ₄ H raised to glutathione conjugated to KLH Applications: Western blotting, Immunoassay and Immunoprecipitation	A001-50UG



SUPPLIED COMPONENTS

Black Half Area 96 well Plate

See: www.ArborAssays.com/resources/#general-info for plate dimensions.

One Plate

Catalog Number X023-1EA

Glutathione S-Transferase Standard

Equine Glutathione S-Transferase at 10 U/mL in a special stabilizing solution.

50 μ L

Catalog Number C026-50UL

GST Detection Reagent

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

1 vial

Catalog Number C035-1EA

Dry DMSO

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

2 mL

Catalog Number X022-2ML

Assay Buffer

Phosphate buffer containing proteins and stabilizers.

45 mL

Catalog Number X033-45ML

Glutathione (GSH)

Glutathione supplied as a 20 mM stable solution.

300 μ L

Catalog Number C028-300UL

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. DMSO can be stored tightly capped at room temperature.

OTHER MATERIALS REQUIRED

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

Fluorescence 96 well plate reader capable of reading fluorescent emission at 460 nm, with excitation at 390 nm. Set plate parameters for a 96-well Corning Costar 3686 plate.

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

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.

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.

ACTIVITY STANDARDIZATION

The Glutathione S-Transferase standard used in this kit has been calibrated using an enzymatic method adapted from reference 1.



ARBOR
ASSAYS

SAMPLE TYPES

This assay has been validated for human urine, serum, EDTA, heparin plasma, toadfish liver (*Opsanus tau*) and oyster hemolymph samples. Most cell lysates should also be compatible. Samples that are not clear or that contain visible particulate should be centrifuged prior to using.

GST activity varies across tissues and species, however we expect this kit to measure GST activity from sources other than human. The end user should evaluate recoveries of GST activity in samples from other species being tested.

SAMPLE PREPARATION

Any samples requiring larger dilutions or with GST activities outside the standard curve range should be diluted further with Assay Buffer to obtain readings within the standard curve.

Urine Samples

Samples that are not clear or that contain visible particulate should be centrifuged prior to using. Urine samples should be diluted $\geq 1:2$ in Assay Buffer by adding one part of urine to one part of Assay Buffer. Sample values should be normalized for urinary volume to urinary creatinine levels. Our Urinary Creatinine Detection Kits, K002-H1 and K002-H5, allow simple, safe and accurate determination of urinary creatinine levels.

Serum and Plasma Samples

Samples that are not clear or that contain visible particulate should be centrifuged prior to using. Fresh serum or EDTA and heparin plasma are separated by centrifugation at $600 \times g$ for 10 minutes. Transfer the serum or plasma from the red blood cells into fresh tubes. The serum or plasma may be stored at -80°C or analyzed immediately. Serum or plasma should be diluted with assay Buffer at a dilution of $\geq 1:2$.

Cell Lysates

Washed cell pellets are resuspended at $10\text{-}40 \times 10^6$ cells/mL in Assay Buffer (we used Jurkats at 10×10^6 cells/mL) and are lysed by vigorous vortexing, freeze/thaw cycling or other suitable disruption method. Resulting centrifuged lysate supernatants are measured at appropriate dilutions. The protocol might require adjustment for other cell types. If protein determinations are to be made on the samples, we would recommend using higher number of cells and lysing in your normal PBS-based lysis buffer and determining protein concentration, prior to additional dilutions in Assay Buffer to measure GST activity.

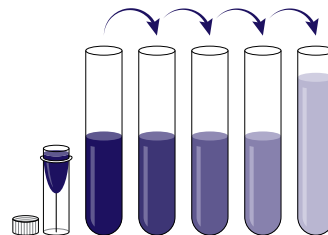
Use all samples within 2 hours of dilution.

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.

Standard Preparation

GST Standards are prepared by labeling test tubes as #1 through #7. Briefly spin vial of standard in a microcentrifuge to ensure contents are at bottom of vial. Pipet 380 μL of Assay Buffer into tube #1 and 200 μL into tubes #2 to #7. Carefully add 20 μL of the Glutathione S-Transferase Standard to tube #1 and vortex completely. Take 200 μL of the GST solution in tube #1 and add it to tube #2 and vortex completely. Repeat these serial dilutions for tubes #3 through #7. The concentration of GST in tubes 1 through 7 will be 500, 250, 125, 62.5, 31.25, 15.61 and 7.81 mU/mL.



Use all Standards within 1 hour of preparation.

	Std 1	Std 2	Std 3	Std 4	Std 5	Std 6	Std 7
Buffer Volume (μL)	380	200	200	200	200	200	200
Addition	Stock	Std 1	Std 2	Std 3	Std 4	Std 5	Std 6
Volume of Addition (μL)	20	200	200	200	200	200	200
Final Conc (mU/mL)	500	250	125	62.5	31.25	15.625	7.81

Detection Reagent

Allow the ziploc pouch to warm **completely** to room temperature prior to opening. Remove the vial of Detection Reagent and add 300 μL of the dry DMSO to the vial. Vortex thoroughly. Store any unused reconstituted Detection Reagent at 4°C in the desiccated pouch for no longer than 2 weeks.

Dilute one part of reconstituted Detection Reagent 1:10 into nine parts of Assay Buffer. 150 μL of Detection Reagent should be diluted with 1.35 mL of Assay Buffer to use half the plate. Discard any excess diluted Detection Reagent.

Glutathione

Dilute one part Glutathione stock provided 1:10 into nine parts of Assay Buffer. 150 μL of Glutathione stock should be diluted with 1.35 mL of Assay Buffer to have enough GSH to be able to read half the plate. Discard any excess diluted glutathione.



ASSAY PROTOCOL

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

1. Use the plate layout sheet on the back page of the insert 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 or standards into duplicate wells in the plate.
3. Pipet 50 μ L of Assay Buffer into duplicate wells as the Zero standard.
4. Add 25 μ L of the Detection Reagent to each well using a repeater pipet.
5. Add 25 μ L of GSH to each well using a repeater pipet.
6. Gently tap the sides of the plate to ensure adequate mixing of the reagents.
7. Incubate at room temperature for 30 minutes.
8. Read the fluorescent signal from each well in a plate reader capable of reading the fluorescent emission at 460 nm with excitation at 370-410 nm. Please contact your plate reader manufacturer for suitable filter sets.
9. Use the plate reader's built-in 4PLC software capabilities to calculate GST activities for each sample.

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 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-s-transferase-fluorescent-kit.assay



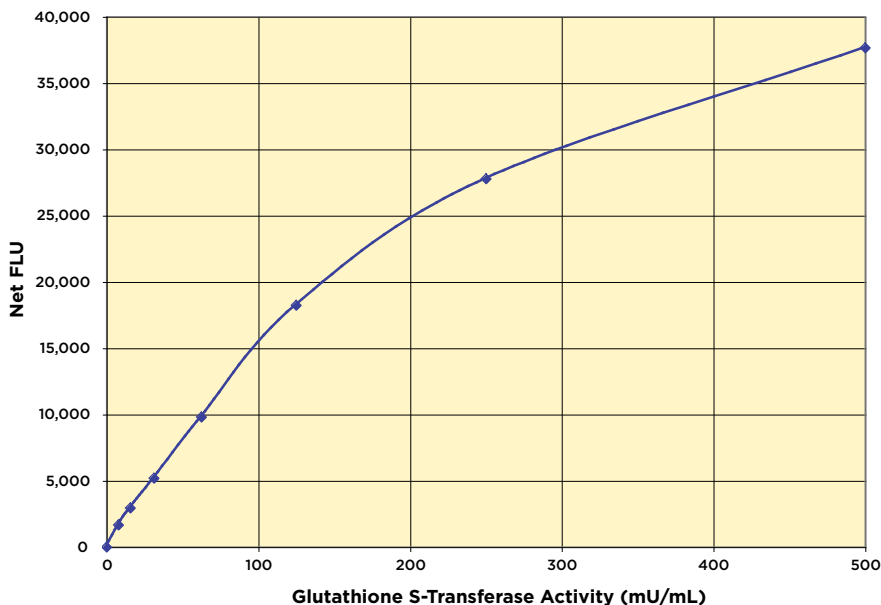
*The MyAssays logo is a registered trademark of MyAssays Ltd.

TYPICAL DATA

Sample	Mean FLU	Net FLU	GST Activity (mU/mL)
Zero	2,749	0	0
Standard 1	40,408	37,659	500
Standard 2	30,529	27,780	250
Standard 3	20,993	18,244	125
Standard 4	12,546	9,797	62.5
Standard 5	7,937	5,188	31.25
Standard 6	5,689	2,940	15.61
Standard 7	4,425	1,676	7.81
Sample 1	7,148	4,399	25.7
Sample 2	5,266	2,517	13.55

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

Typical Standard Curve



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

VALIDATION DATA

Sensitivity and Limit of Detection

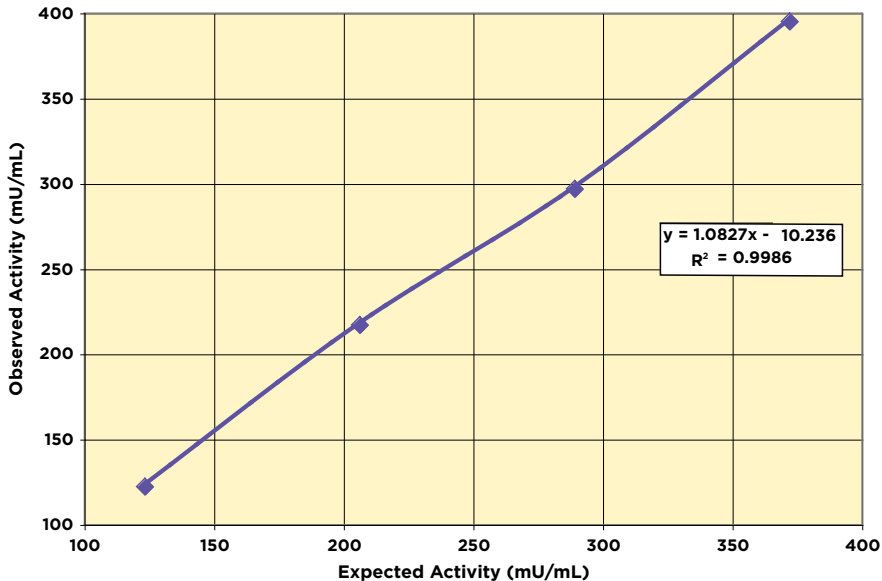
Sensitivity was calculated by comparing the FLUs for twenty wells run for each of the zero and standard #7. The detection limit was determined at two (2) standard deviations from the zero along the standard curve. **Sensitivity was determined as 2.70 mU/mL.**

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 serum sample. **The Limit of Detection was determined as 1.90 mU/mL.**

Linearity

Linearity was determined by taking Jurkat cell lysates at 10×10^6 and 0.8×10^6 cells/mL and mixing in the ratios given below. The measured activities were compared to the expected values based on the ratios used.

Low Cell #	High Cell #	Expected Activity (mU/mL)	Observed Activity (mU/mL)	% Recovery
80%	20%	123.3	122.4	99.3
60%	40%	206.2	217.2	105.3
40%	60%	289.1	297.1	102.7
20%	80%	372.1	395.1	106.2
Mean Recovery				103.4%



Intra Assay Precision

Four serum samples were diluted 1:2 with Assay Buffer and run in replicates of 16 in an assay. The mean and precision of the calculated GST activities were:

Sample	GST Activity (mU/mL)	%CV
1	315.9	4.6
2	221.2	5.6
3	88.2	4.2
4	22.7	6.6

Inter Assay Precision

Four serum samples were diluted 1:2 with Assay Buffer and run in duplicates in twenty assays over multiple days by four operators. The mean and precision of the calculated GST activities were:

Sample	GST Activity (mU/mL)	%CV
1	291.7	12.6
2	218.5	11.0
3	89.6	10.4
4	23.0	15.9

SAMPLE VALUES

Twenty random human serum, heparin and EDTA plasma samples were tested in the assay. Values ranged from 26.8 to 59.8 mU/mL with an average of 39.1 mU/mL. Five random human urine samples were tested in the assay and values were normalized to urinary Creatinine levels utilizing the Arbor Assays Urinary Creatinine Detection kit, K002-H1 or K002-H5. Levels ranged from almost 24 to over 102 mU/mL. When normalized for urine volume using creatinine values, the levels ranged from 2.83 to 20.2 μ U/mg creatinine.

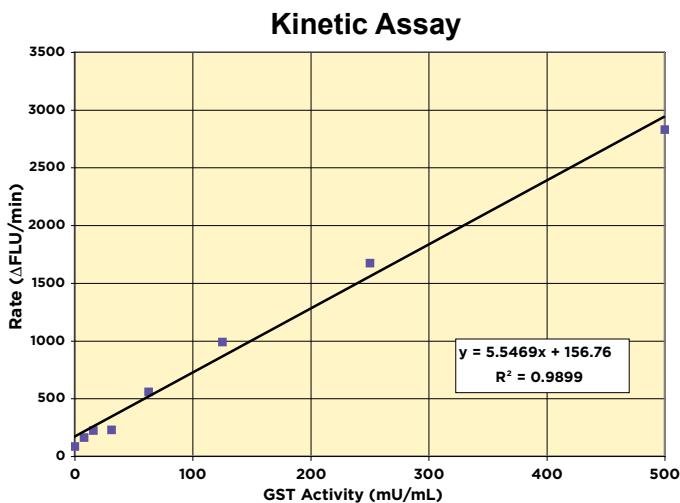
INTERFERENCES

A variety of solvents and detergents were tested as possible interfering substances in the assay. Approximately 10% change was seen in the GST activity in the presence of 1% methanol or DMSO in the sample. Three detergents were also tested: Triton X-100, Tween 20 and SDS. At 0.01% concentration in the sample both SDS and Tween showed no change in activity, whereas Triton showed > 47% decrease at 0.01%.

Bilirubin levels of 2.5 μ g/mL in the sample showed < 5% decrease in GST activity.

END POINT VERSUS KINETIC ACTIVITY

The assay can also be run as a kinetic assay. A human serum sample was read in both an end point and in a kinetic assay. In the end point measurement it had a reading of 12.12 mU/mL and in the kinetic assay a reading of 11.92 mU/mL. A typical standard curve for the kinetic assay is shown below.



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

Email 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

15

K008-F1 WEB 190503



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