

Product Information

Glycine Assay Kit

Catalog Number **MAK261**Storage Temperature -20°C

TECHNICAL BULLETIN

Product Description

Glycine (Gly) is a non-essential amino acid used to help create muscle tissue and convert glucose into energy. The basic functions of glycine include the participation in the synthesis of creatine, glutathione, heme groups, and conjugated bile acids (bile salts). It is also used in the synthesis of DNA and RNA. It is one of the most abundant residues in the triple-helical structure of collagen, which contributes to the elasticity of skin and connective tissue. It acts as a glucogenic amino acid regulating blood sugar levels. Glycine is necessary for central nervous system function. It possesses both inhibitory and excitatory neurotransmitter functions in the brain stem and spinal cord. Recent studies suggest glycine may play a role in tumorigenesis and malignancy.

The Glycine Assay Kit is a simple and reliable high throughput assay that detects the physiological concentration of glycine in a variety of biological fluids. In the assay, glycine is oxidized generating a fluorometric ($\lambda_{\text{ex}} = 535 \text{ nm}/\lambda_{\text{em}} = 587 \text{ nm}$) product proportional to the glycine present. The assay is specific, and other standard and non-standard amino acids do not interfere with the assay. The kit is sensitive to $1 \mu\text{M}$ of glycine.

This kit is suitable for use with plasma, serum, urine, and other biological samples.

Components

The kit is sufficient for 100 assays in 96 well plates.

| | |
|--|--------|
| Glycine Assay Buffer Catalog Number MAK261A | 25 mL |
| Glycine Probe Catalog Number MAK261B | 0.4 mL |
| Glycine Enzyme Mix Catalog Number MAK261C | 1 vL |

Glycine Developer
Catalog Number MAK261D

1 vL

Glycine Standard
Catalog Number MAK261E

1 vL

Reagents and Equipment Required but Not Provided.

- 96 well flat-bottom plate – It is recommended to use black plates with clear bottoms for fluorescence assays
- Spectrophotometric multiwell plate reader
- 10 kDa Molecular Weight Cut-Off (MWCO) Spin Filter
- 10% Glycerol - Molecular Biology Grade (G5516)

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Briefly centrifuge vials before opening. Use ultrapure water for the preparation of reagents. To maintain reagent integrity, avoid repeated freeze/thaw cycles.

Glycine Assay Buffer – Allow buffer to come to room temperature before use. Store at -20°C .

Glycine Probe – Allow probe to come to room temperature before use. Store protected from light at -20°C . Upon thawing, the Glycine Probe is ready-to-use as supplied.

Glycine Enzyme Mix - Dilute $10 \mu\text{L}$ of 100% Glycerol with $90 \mu\text{L}$ of Glycerol Assay Buffer. Vortex for 30 seconds. Reconstitute the Glycine Enzyme Mix with $55 \mu\text{L}$ of Glycine Assay Buffer with 10% Glycerol. Mix well, then aliquot and store at -20°C , protected from light. Keep on ice during use and limit to two freeze/thaw cycles.

Glycine Developer - Reconstitute with 220 μL of Glycine Assay Buffer. Aliquot and store at $-20\text{ }^{\circ}\text{C}$. Keep on ice during use and limit to one freeze/thaw cycle.

Glycine Standard - Reconstitute with 100 μL of water to generate a 100 mM Glycine Standard. Dissolve completely. Store at $-20\text{ }^{\circ}\text{C}$ and use within 2 months of reconstitution.

Storage/Stability

The kit is shipped on wet ice. Storage at $-20\text{ }^{\circ}\text{C}$, protected from light, is recommended.

Procedure

All samples and standards should be run in duplicate.

Glycine Standards for Fluorometric Detection

Dilute 10 μL of the 100 mM Glycine Standard with 990 μL of water to prepare a 1 mM standard solution. Further dilute to 50 μM by adding 50 μL of 1 mM Glycine Standard to 950 μL of water. Add 0, 2, 4, 6, 8, and 10 μL of the 50 μM Glycine standard solution into a 96 well plate, generating 0 (blank), 0.1, 0.2, 0.3, 0.4, and 0.5 nmole/well standards. Add Glycine Assay Buffer to each well to bring the final volume to 50 μL .

Sample Preparation

The fluorometric assay requires 50 μL of sample for each reaction (well).

Homogenize tissue (10–20 mg) or lyse cells (1×10^6) in 100 μL of Glycine Assay Buffer. Centrifuge the samples at $10,000 \times g$ at $2-8\text{ }^{\circ}\text{C}$ for 5 minutes to remove insoluble material. Collect the supernatant.

Centrifuge liquid samples at $10,000 \times g$ at $2-8\text{ }^{\circ}\text{C}$ for 5 minutes.

Add 1–50 μL samples into wells of a 96 well plate. Bring samples to a final volume of 50 μL with Assay Buffer.

Notes: For unknown samples, it is suggested to test several sample volumes to make sure the readings are within the range of the standard curve. Glycine concentration varies over a wide range depending on the sample. For example - human urine: 44–300 μM /mM creatinine; human serum: 126–490 μM glycine; saliva: 10–300 μM glycine. The recommended dilution factor for urine is 50–500 \times , for serum is 16–250 \times , and for saliva is 4–50 \times .

Metabolites found in biological samples can interfere with the assay. Dilute biological samples with Glycine Assay Buffer. If interference is observed in the diluted samples, prepare parallel sample well(s) as sample background control(s) by omitting the Glycine Enzyme Mix.

For samples with high protein content, deproteinize using a 10 kDa MWCO spin filter.

To ensure accurate determination of Glycine in the test samples or for samples having low concentrations of glycine, we recommend spiking samples with a known amount of Glycine Standard (e.g., 0.3 nmole).

Assay Reaction

1. Set up the Master Reaction Mix according to the scheme in Table 1. 50 μL of the Master Reaction Mix is required for each reaction (well).
2. Dilute the Glycine Enzyme Mix 10-fold. (Dilute 2 μL of Glycine Enzyme Mix with 18 μL of Glycine Assay Buffer). 5 μL will be needed for each reaction (well) to be assayed.

Table 1.
Master Reaction Mix

| Reagent | Samples and Positive Control | Background Control Mix |
|------------------------------|------------------------------|------------------------|
| Glycine Assay Buffer | 42 μL | 47 μL |
| Glycine Enzyme Mix (diluted) | 5 μL | – |
| Glycine Developer | 2 μL | 2 μL |
| Glycine Probe | 1 μL | 1 μL |

3. Add 50 μL of the Master Reaction Mix to each sample and positive control well containing the Glycine standard. If using a sample control, add 50 μL of Sample Control Mix to sample control reactions (wells). Mix well using a horizontal shaker or by pipetting.

4. Incubate the plate for 1 hour at 25 °C. Protect the plate from light during the incubation.
5. Measure fluorescence intensity ($\lambda_{\text{ex}} = 535 \text{ nm}$ /
 $\lambda_{\text{em}} = 587 \text{ nm}$).

Results

Calculations

The background is the value obtained for the 0 (assay blank) Glycine Standard. Correct for the background by subtracting the 0 (assay blank) value from all readings. Background values can be significant and must be subtracted from all readings. Subtract the Sample Background Control value from the sample readings.

Use the values obtained from the appropriate Glycine standards to plot a standard curve. The amount of glycine present in the samples may be determined from the standard curve.

Notes: A new standard curve must be set up each time the assay is run.

For spiked samples, correct for interference by using the following equation:

$$\frac{\text{RFU}_{\text{sample (corrected)}}}{(\text{RFU}_{\text{sample+Gly Std (corrected)}}) - (\text{RFU}_{\text{sample (corrected)}})} * \text{Gly spike (nmol)}$$

Concentration of Glycine

$$S_a/S_v = C$$

S_a = Amount of Glycine in the unknown sample (nmole) from standard curve

S_v = Sample volume (μL) added into the wells

C = Concentration of Glycine in sample

Glycine molecular weight: 75.07 g/mole

Sample Calculation

Amount of Glycine (S_a) = 0.384 nmole
(from standard curve)

Sample volume (S_v) = 50.0 μL

Concentration of Glycine in sample

$$0.384 \text{ nmole}/50.0 \mu\text{L} = 0.00768 \text{ nmole}/\mu\text{L}$$

$$0.00768 \text{ nmole}/\mu\text{L} \times 75.07 \text{ ng/nmole} = 0.577 \text{ ng}/\mu\text{L}$$

Troubleshooting Guide

| Problem | Possible Cause | Suggested Solution |
|--|---|--|
| Assay Not Working | Cold assay buffer | Assay Buffer must be at room temperature |
| | Omission of step in procedure | Refer and follow Technical Bulletin precisely |
| | Plate reader at incorrect wavelength | Check filter settings of instrument |
| | Type of 96 well plate used | For fluorescence assays, use black plates with clear bottoms. |
| Samples with erratic readings | Samples prepared in different buffer | Use the Assay Buffer provided or refer to Technical Bulletin for instructions |
| | Cell/Tissue culture samples were incompletely homogenized | Repeat the sample homogenization, increasing the length and extent of homogenization step. |
| | Samples used after multiple freeze-thaw cycles | Aliquot and freeze samples if samples will be used multiple times |
| | Presence of interfering substance in the sample | If possible, dilute sample further |
| | Use of old or inappropriately stored samples | Use fresh samples and store correctly until use |
| Lower/higher readings in samples and standards | Improperly thawed components | Thaw all components completely and mix gently before use |
| | Use of expired kit or improperly stored reagents | Check the expiration date and store the components appropriately |
| | Allowing the reagents to sit for extended times on ice | Prepare fresh Reaction Mix before each use |
| | Incorrect incubation times or temperatures | Refer to Technical Bulletin and verify correct incubation times and temperatures |
| | Incorrect volumes used | Use calibrated pipettes and aliquot correctly |
| Non-linear standard curve | Use of partially thawed components | Thaw and resuspend all components before preparing the reaction mix |
| | Pipetting errors in preparation of standards | Avoid pipetting small volumes |
| | Pipetting errors in the Reaction Mix | Prepare a Reaction Mix whenever possible |
| | Air bubbles formed in well | Pipette gently against the wall of the plate well |
| | Standard stock is at incorrect concentration | Refer to the standard dilution instructions in the Technical Bulletin |
| | Calculation errors | Recheck calculations after referring to Technical Bulletin |
| | Substituting reagents from older kits/lots | Use fresh components from the same kit |
| Unanticipated results | Samples measured at incorrect wavelength | Check the equipment and filter settings |
| | Samples contain interfering substances | If possible, dilute sample further |
| | Sample readings above/below the linear range | Concentrate or dilute samples so readings are in the linear range |

KVG,JAC,MAM 02/18-1