

Product Information

Glycerol 3-Phosphate Colorimetric Assay Kit

Catalog Number **MAK207**

Storage Temperature -20°C

TECHNICAL BULLETIN

Product Description

Glycerol 3-phosphate (G3P) is an important intermediate of carbohydrate and lipid metabolic pathways. It is produced from glycerol by glycerol kinase or from dihydroxyacetone phosphate by glycerol 3-phosphate dehydrogenase. G3P may enter the G3P shuttle to generate NAD^{+} , or may be converted to glyceraldehyde 3-phosphate and enter glycolysis or the lipid biosynthesis pathway.^{1,2}

The Glycerol 3-Phosphate Colorimetric Assay Kit provides a simple assay for measuring G3P in various tissues and cells (ranging from 2–10 nmole/well). G3P is determined by measuring a colorimetric product with absorbance at 450 nm (A_{450}) proportional to the amount of G3P present.

Components

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

G3P Assay Buffer Catalog Number MAK207A	25 mL
G3P Enzyme Mix Catalog Number MAK207B	1 vL
G3P Probe Catalog Number MAK207C	1 vL
G3P Standard Catalog Number MAK207D	1 vL

Reagents and Equipment Required but Not Provided

- 96 well flat-bottom plate – It is recommended to use clear plates for colorimetric assays.
- Spectrophotometric multiwell plate reader

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.

G3P Assay Buffer – Warm buffer to room temperature before use.

G3P Enzyme Mix – Reconstitute with 220 μL of G3P Assay Buffer. Mix well by pipetting. Aliquot and store at -20°C . Keep on ice while in use.

G3P Probe – Reconstitute with 220 μL of water. Mix well by pipetting and store at -20°C . Use within 2 months.

G3P Standard – Reconstitute with 100 μL of water to generate a 100 mM (100 nmole/ μL) G3P Standard Solution. Store at -20°C . Use within 2 months. Keep on ice while in use.

Storage/Stability

The kit is shipped on wet ice and storage at -20°C , protected from light, is recommended.

Procedure

All samples and standards should be run in duplicate.

G3P Standards for Colorimetric Detection

Dilute 10 μL of the 100 mM G3P Standard with 990 μL of water and mix well to make a 1 mM (1 nmole/ μL) G3P Standard Solution. Add 0, 2, 4, 6, 8, and 10 μL of the 1 mM (1 nmole/ μL) G3P Standard Solution into a 96 well plate generating 0 (blank), 2, 4, 6, 8, and 10 nmole/well standards. Add G3P Assay Buffer to each well to bring the volume to 50 μL .

Sample Preparation

Clear liquid samples may be assayed directly.

Tissue samples (10 mg) or cells (1×10^6) can be homogenized in 100 μL of ice cold G3P Assay Buffer. Keep on ice for 10 minutes. Centrifuge the samples at $12,000 \times g$ for 5 minutes to remove insoluble material. Transfer supernatant to a fresh tube.

Add 1–50 μL of the samples into duplicate wells. Bring samples to a final volume of 50 μL using G3P Assay Buffer.

Notes: For unknown samples, it is suggested to test several sample dilutions to ensure the readings are within the linear range of the standard curve.

For samples exhibiting significant background, especially samples containing NADH, include a Sample Blank for each sample by omitting the G3P Enzyme Mix. The Sample Blank readings can then be subtracted from the sample readings.

Assay Reaction

1. Set up Reaction Mixes according to the scheme in Table 1. 50 μL of Reaction Mix is required for each reaction (well).

Table 1.
Reaction Mixes

Reagent	Standards and samples	Sample Blank
G3P Assay Buffer	46 μL	48 μL
G3P Enzyme Mix	2 μL	–
G3P Probe	2 μL	2 μL

2. Add 50 μL of the appropriate Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting.
3. Incubate the plate for 40 minutes at 37 °C.
4. Measure the absorbance (A_{450}) in a microplate reader.

Results

Calculations

Correct for the background by subtracting the measurement obtained for the 0 (blank) G3P standard from that of the standards and samples. Background values can be significant and must be subtracted from all readings. Use the values obtained from the appropriate G3P Standards to plot a standard curve.

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

Subtract the Sample Blank value from the Sample value to obtain the corrected measurement. Using the corrected measurement, determine the amount of G3P (nmole/well) generated by the assay (S_a).

Concentration of G3P

$$C = S_a / S_v$$

where:

S_a = Amount of G3P in unknown sample well (nmole) from standard curve

S_v = Sample volume (μ L) added into the well

C = Concentration of G3P in sample (nmole/ μ L)

Sample Calculation

Amount of G3P (S_a) = 5.841 nmole
(from standard curve)

Sample volume (S_v) = 10 μ L

Concentration of glycerol 3-phosphate in sample:

$$5.841 \text{ nmole} / 10 \mu\text{L} = 0.5841 \text{ nmole}/\mu\text{L}$$

Molecular weight of G3P = 172.1 g/mole

$$0.5841 \text{ nmole}/\mu\text{L} \times 172.1 \text{ ng/nmole} = 100.5 \text{ ng}/\mu\text{L}$$

References

1. Crabtree, B., and Newsholme, E.A., The activities of phosphorylase, hexokinase, phospho-fructokinase, lactate dehydrogenase and the glycerol 3-phosphate dehydrogenases in muscles from vertebrates and invertebrates. *Biochem. J.*, **126**, 49–58 (1972).
2. Larsson, C. et al., The importance of the glycerol 3-phosphate shuttle during aerobic growth of *Saccharomyces cerevisiae*. *Yeast*, **14**, 347–357 (1998).
3. Kunst, L. et al., Altered regulation of lipid biosynthesis in a mutant of *Arabidopsis* deficient in chloroplast Glycerol 3-phosphate acyltransferase activity. *Proc. Natl. Acad. Sci. U. S. A.*, **85**, 4143–4147 (1988).

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 colorimetric assays, use clear plates
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 needed to use 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 Mixes 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 Mixes
	Pipetting errors in preparation of standards	Avoid pipetting small volumes
	Pipetting errors in the Reaction Mix	Prepare Reaction Mixes 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

GD,KVG,LS,MAM 10/14-1