

Gamma Glutamyl Transferase (GGT) Activity Colorimetric Assay Kit

(Catalog #K784-100; 100 reactions; Store kit at -20°C)

I. Introduction:

The Gamma-Glutamyl Transferase (GGT; EC 2.3.2.2) is an enzyme that transfers gamma-glutamyl functional groups. It is found in many tissues, the most notable one being the liver, and has significance in medicine as a diagnostic marker. BioVision's Gamma-Glutamyl Transferase Assay Kit provides a convenient tool for sensitive detection of the GGT in a variety of samples. The GGT in sample will recognize L-γ-Glutamyl-pNA as a specific substrate leading to proportional color development. The activity of GGT can be easily quantified colorimetrically (λ = 418 nm). This assay detects GGT activity as low as 0.5 mIU.

II. Kit Contents:

Components	K784-100	Cap Code	Part Number
GGT Assay Buffer	25 ml	WM	K784-100-1
GGT Substrate	1 Bottle	NM	K784-100-2
GGT Positive Control	1 vial	Green	K784-100-3
pNA Standard (2mM)	400 µl	Yellow	K784-100-4

III. Storage and Handling:

Store the kit at -20°C, protected from light. Allow Assay Buffer to warm to room temperature before use. Briefly centrifuge vials before opening. Read the entire protocol before performing the assay.

IV. Reagent Reconstitution and General Consideration:

GGT Substrate Solution: Add 10 ml assay buffer into substrate bottle and mix well. Take out enough substrate solution (90 µl per assay) for the number of assays to be performed in experiment. Store the rest of the GGT Substrate Solution into -20°C quickly. **Note:** The GGT Substrate solution is unstable at room temperature (can be hydrolyzed by itself) which increases the assay background.

GGT Positive Control: Reconstitute with 100 µl diH₂O. Pipette up and down several times to completely dissolve the pellet into solution (**Don't vortex**). Aliquot enough GGT Positive Control (10 µl per assay) for the number of assays to be performed in each experiment and aliquot and freeze the rest immediately at -20°C for future use. The GGT Positive Control is stable for up to 1 month at -20°C after reconstitution or freeze-thaw cycles (< 5 times). Keep the GGT Positive Control on ice during the preparation.

V. GGT Activity Assay Protocol:

1. pNA Standard Curve: (Warm for 1-2 min at 37°C to completely melt DMSO):

Add 0, 4, 8, 12, 16, 20 µl of the 2 mM pNA standard solution into a 96-well plate in duplicate to generate 0, 8, 16, 24, 32, 40 nmol/well standard. Adjust the final volume to 100 µl with GGT Assay Buffer.

2. Sample Preparations:

Tissues (10 mg) or cells (1×10⁶) can be homogenized in the 200 µl GGT Assay Buffer then centrifuged (13,000 x g, 10 min.) to remove insoluble material. Serum samples (10 µl) can be directly added into each well. Prepare test samples to 10 µl/well with GGT Assay Buffer in a 96-well plate. We suggest testing several doses of your sample to make sure the readings are within the linear range of the standard curve.

3. Reaction Mix:

Add 90 µl GGT Substrate Mix into each well containing the test samples and positive controls. Mix well. **Do not add to pNA Standards.**

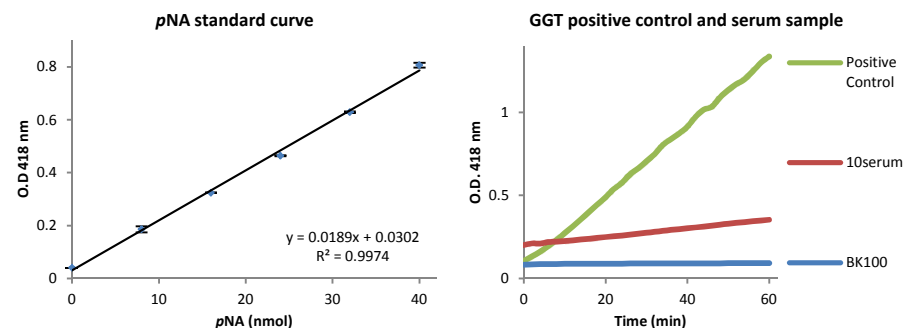
4. Measurement: For pNA Standard Curve, measure OD at 418 nm in a microplate reader. For the samples and positive controls, incubate the mix for 3 min at 37°C, then measure OD at 418 nm in a microplate reader (A₀), incubate for another 30 min to 2 hr at 37°C to measure OD at 418 nm again (A₁); incubation times will depend on the GGT activity in the samples. We recommend measuring the OD in a kinetic method (preferably every 3- 5 min) and choose the period of linear range which falls within pNA Standard Curve to calculate the GGT activity of the samples.

5. Calculation: Plot the pNA standard Curve, then calculate the GGT activity of the test samples: ΔOD = A₁ - A₀, apply the ΔOD to the pNA standard curve to get B nmol of pNA generated by GGT in the given time.

$$\text{GGT Activity} = \frac{B}{T \times V} \times \text{Sample Dilution Factor} = \text{nmol/min/ml} = \text{mU/ml}$$

Where: **B** is the pNA amount from standard Curve (in nmol)
T is the time incubated (in min)
V is the sample volume added into the reaction well (in ml)

Unit Definition: One unit GGT will generate 1.0 µmol of pNA per min at 37°C.
Note: One pNA unit ≈ 1.5 IU.



RELATED PRODUCTS:

- ADP/ATP Ratio Assay Kit
- Glucose Assay Kit
- Uric Acid Assay Kit
- Creatine Assay Kit
- Ammonia Assay Kit
- Triglyceride Assay Kit
- Nitric Oxide Assay Kit
- GGT Activity Fluorometric Assay Kit
- Ethanol Assay Kit

- Ascorbic Acid Quantification Kit
- Fatty Acid Assay Kit
- Pyruvate Assay Kit
- Creatinine Assay Kit
- Free Glycerol Assay Kit
- Total Antioxidant Capacity (TAC) Assay Kit
- Glutamate Kit
- Formate Assay Kit
- Cholesterol Assay Kit

FOR RESEARCH USE ONLY! Not to be used on humans.

GENERAL TROUBLESHOOTING GUIDE:

Problems	Cause	Solution
Assay not working	<ul style="list-style-type: none"> • Use of ice-cold assay buffer • Omission of a step in the protocol • Plate read at incorrect wavelength • Use of a different 96-well plate 	<ul style="list-style-type: none"> • Assay buffer must be at room temperature • Refer and follow the data sheet precisely • Check the wavelength in the data sheet and the filter settings of the instrument • Fluorescence: Black plates (clear bottoms) ; Luminescence: White plates ; Colorimeters: Clear plates
Samples with erratic readings	<ul style="list-style-type: none"> • Use of an incompatible sample type • Samples prepared in a different buffer • Cell/ tissue samples were not completely homogenized • Samples used after multiple free-thaw cycles • Presence of interfering substance in the sample • Use of old or inappropriately stored samples 	<ul style="list-style-type: none"> • Refer data sheet for details about incompatible samples • Use the assay buffer provided in the kit or refer data sheet for instructions • Use Dounce homogenizer (increase the number of strokes); observe for lysis under microscope • Aliquot and freeze samples if needed to use multiple times • Troubleshoot if needed • Use fresh samples or store at correct temperatures until use
Lower/ Higher readings in Samples and Standards	<ul style="list-style-type: none"> • Improperly thawed components • Use of expired kit or improperly stored reagents • Allowing the reagents to sit for extended times on ice • Incorrect incubation times or temperatures • Incorrect volumes used 	<ul style="list-style-type: none"> • Thaw all components completely and mix gently before use • Always check the expiry date and store the components appropriately • Always thaw and prepare fresh reaction mix before use • Refer datasheet & verify correct incubation times and temperatures • Use calibrated pipettes and aliquot correctly
Readings do not follow a linear pattern for Standard curve	<ul style="list-style-type: none"> • Use of partially thawed components • Pipetting errors in the standard • Pipetting errors in the reaction mix • Air bubbles formed in well • Standard stock is at an incorrect concentration • Calculation errors • Substituting reagents from older kits/ lots 	<ul style="list-style-type: none"> • Thaw and resuspend all components before preparing the reaction mix • Avoid pipetting small volumes • Prepare a master reaction mix whenever possible • Pipette gently against the wall of the tubes • Always refer the dilutions in the data sheet • Recheck calculations after referring the data sheet • Use fresh components from the same kit
Unanticipated results	<ul style="list-style-type: none"> • Measured at incorrect wavelength • Samples contain interfering substances • Use of incompatible sample type • Sample readings above/below the linear range 	<ul style="list-style-type: none"> • Check the equipment and the filter setting • Troubleshoot if it interferes with the kit • Refer data sheet to check if sample is compatible with the kit or optimization is needed • Concentrate/ Dilute sample so as to be in the linear range
<p>Note: The most probable list of causes is under each problem section. Causes/ Solutions may overlap with other problems.</p>		