

Glucose Oxidase Activity Assay Kit

PromoKine

Convenient kit for sensitive colorimetric/fluorometric detection of glucose oxidase activity in a variety of samples

Instruction Manual

| Catalog Number | PK-CA577-K788 | | | | | | | | | | | | | | | | | | | | | |
|---|---|--------------------|--------------|--------------------|------------------|------------------|------------------|---------------|-----------------|-----------------|-----------------------------|-----------------|-----------------|---------------|------------------|-------|---|--------|--------|------------------------------------|--------|--------|
| Description | <p>The glucose oxidase enzyme (GOx) (EC 1.1.3.4) is an oxidoreductase commonly found in a wide variety of microorganisms that catalyzes the oxidation of glucose to hydrogen peroxide and D-glucono-δ-lactone. GOx aids in breaking the sugar down into its metabolites.</p> <p>PromoKine's Glucose Oxidase Activity Assay Kit provides a convenient tool for sensitive detection of the GOx in a variety of samples. Glucose oxidase in samples recognizes D-glucose as a specific substrate leading to proportional color development. The activity of GOx can be easily quantified colorimetrically ($\lambda = 570$ nm) or fluorometrically (Ex/Em = 535/585 nm). GOx assay detects glucose oxidase activity as low as 0.01mU.</p> | | | | | | | | | | | | | | | | | | | | | |
| Quantity | 100 assays | | | | | | | | | | | | | | | | | | | | | |
| Kit Components | <table border="1"><thead><tr><th>Components</th><th>Quantity</th><th>Color Code</th></tr></thead><tbody><tr><td>GOx Assay Buffer</td><td>25 ml</td><td>WM</td></tr><tr><td>OxiRed Probe</td><td>0.2 ml</td><td>Red</td></tr><tr><td>GOx Developer (lyophilized)</td><td>1 vial</td><td>Green</td></tr><tr><td>GOx Substrate</td><td>1 ml</td><td>Blue</td></tr><tr><td>H₂O₂ Standard (0.88 M)</td><td>0.1 ml</td><td>Yellow</td></tr><tr><td>GOx Positive Control (lyophilized)</td><td>1 vial</td><td>Purple</td></tr></tbody></table> | Components | Quantity | Color Code | GOx Assay Buffer | 25 ml | WM | OxiRed Probe | 0.2 ml | Red | GOx Developer (lyophilized) | 1 vial | Green | GOx Substrate | 1 ml | Blue | H ₂ O ₂ Standard (0.88 M) | 0.1 ml | Yellow | GOx Positive Control (lyophilized) | 1 vial | Purple |
| Components | Quantity | Color Code | | | | | | | | | | | | | | | | | | | | |
| GOx Assay Buffer | 25 ml | WM | | | | | | | | | | | | | | | | | | | | |
| OxiRed Probe | 0.2 ml | Red | | | | | | | | | | | | | | | | | | | | |
| GOx Developer (lyophilized) | 1 vial | Green | | | | | | | | | | | | | | | | | | | | |
| GOx Substrate | 1 ml | Blue | | | | | | | | | | | | | | | | | | | | |
| H ₂ O ₂ Standard (0.88 M) | 0.1 ml | Yellow | | | | | | | | | | | | | | | | | | | | |
| GOx Positive Control (lyophilized) | 1 vial | Purple | | | | | | | | | | | | | | | | | | | | |
| Storage and Reagents Preparation | <p>Store kit at -20°C, protect from light. Warm Assay Buffer to room temperature before use. Briefly centrifuge all small vials prior to opening. Read entire protocol before performing assay.</p> <p>GOx Developer and GOx Positive Control: Reconstitute with 220 μl Assay Buffer. Pipette up and down several times to completely dissolve the pellet (Don't vortex). Aliquot and freeze at -20°C. Stable for up to 2 months at -20°C after reconstitution or freeze-thaw cycles (< 5 times). Keep GOx Positive Control on ice while in use.</p> | | | | | | | | | | | | | | | | | | | | | |
| Assay Protocol | <p>1. H₂O₂ Standard Curve: Add 10 μl 0.88 M H₂O₂ Standard to 870 μl dH₂O to make 10 mM H₂O₂ Standard. Dilute 10 mM H₂O₂ Standard further to 1:10 with Assay Buffer to make a 1 mM H₂O₂ Standard. Add 0, 2, 4, 6, 8, 10 μl of the diluted 1 mM H₂O₂ Standard into a series of wells of 96-well plate to generate 0, 2, 4, 6, 8, 10 nmol/well H₂O₂ Standard.</p> <p><u>For the fluorometric assay,</u> dilute 1 mM H₂O₂ Standard 1:10 with Assay Buffer to make 100 μM H₂O₂ standard. Add 0, 2, 4, 6, 8, 10 μl of the diluted 100 μM H₂O₂ standard into a series of wells of 96-well plate to generate 0, 0.2, 0.4, 0.6, 0.8, 1.0 nmol/well H₂O₂ Standard. Adjust the final volume to 50 μl with Assay Buffer.</p> <p>2. Sample Preparations: Homogenize cells (1×10^6) with 100-200 μl Assay Buffer. Centrifuge at 13,000 g for 10 minutes to remove the insoluble material. 5-50 μl serum samples can be directly diluted in the Assay Buffer. Add 1-50 μl sample per well, adjust final volume to 50 μl with Assay Buffer. For samples having high background, prepare a parallel sample well as the background control. Note: For unknown samples, we suggest testing several doses to ensure the readings are within the standard curve range.</p> <p>3. Positive Control: Add 2-10 μl of Positive Control into the desired well(s) and adjust final volume to 50 μl with Assay Buffer.</p> <p>4. Reaction Mix: Mix enough reagents for the number of assays to be performed. For each well, prepare 50 μl Reaction Mix containing:</p> <table border="1"><thead><tr><th></th><th>Reaction Mix</th><th>Background Control</th></tr></thead><tbody><tr><td>GOx Assay Buffer</td><td>36 μl</td><td>46 μl</td></tr><tr><td>GOx Developer</td><td>2 μl</td><td>2 μl</td></tr><tr><td>OxiRed Probe</td><td>2 μl</td><td>2 μl</td></tr><tr><td>GOx Substrate</td><td>10 μl</td><td>-----</td></tr></tbody></table> <p>Add 50 μl of the reaction mix to each well containing samples, Positive Control, and standards. Mix well. Note: The fluorometric assay is ~ 10-fold more sensitive than the colorimetric assay so dilute the probe 10 times in Assay Buffer & use the same volume (2 μl). Background control mix is recommended for samples having high background.</p> | | Reaction Mix | Background Control | GOx Assay Buffer | 36 μl | 46 μl | GOx Developer | 2 μl | 2 μl | OxiRed Probe | 2 μl | 2 μl | GOx Substrate | 10 μl | ----- | | | | | | |
| | Reaction Mix | Background Control | | | | | | | | | | | | | | | | | | | | |
| GOx Assay Buffer | 36 μl | 46 μl | | | | | | | | | | | | | | | | | | | | |
| GOx Developer | 2 μl | 2 μl | | | | | | | | | | | | | | | | | | | | |
| OxiRed Probe | 2 μl | 2 μl | | | | | | | | | | | | | | | | | | | | |
| GOx Substrate | 10 μl | ----- | | | | | | | | | | | | | | | | | | | | |

5. Measurement: Incubate the plate for 5 minutes at 37°C and measure OD at 570 nm or fluorescence at Ex/Em = 535/585 nm (A1). Incubate for another 15 minutes to 2 hours at 37°C and again measure (A2). Note: Incubation time depends on the glucose oxidase activity in the samples. We recommend measuring in a kinetic method (preferably every 1–2 minutes) and choose the period of linear range to calculate the glucose oxidase activity of the samples. If the absorbance exceeds 0.7 OD at about ~15 minutes, dilute the sample and rerun the assay. The H₂O₂ Standard curve can read in end point mode (i.e. at the end of incubation time).

6. Calculation: Subtract the 0 standard reading from all readings. Plot H₂O₂ Standard Curve. Calculate the glucose oxidase activity of the test sample: $\Delta OD = A2 - A1$. Apply the ΔOD to the H₂O₂ Standard Curve to get B nmol of H₂O₂ generated by Glucose Oxidase during the reaction time ($\Delta T = T2 - T1$).

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

Where: B is the H₂O₂ amount from Standard Curve (nmol).

ΔT is the time incubated (min).

V is the sample volume added into the reaction well (ml).

Unit Definition: One unit of GOx is the amount of enzyme that generates 1.0 μmol of H₂O₂ per min at 37°C.

Intended Use

For in vitro research use only. Not for diagnostic or therapeutic procedures.

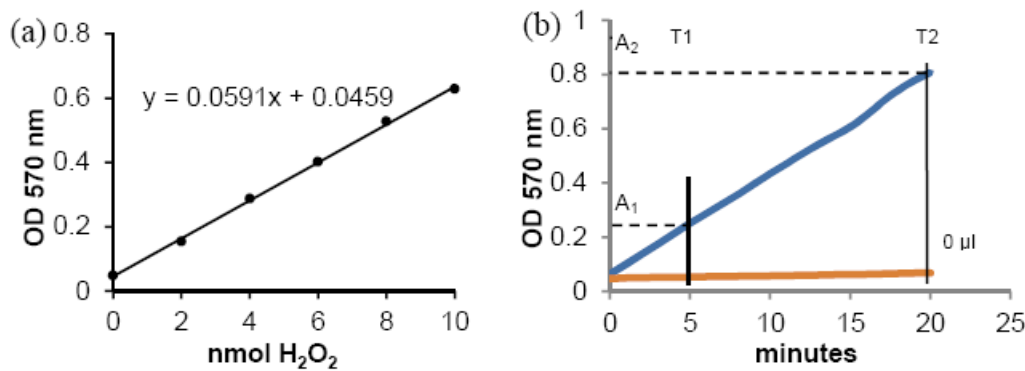


Figure: H₂O₂ Standard Curve (a). Glucose oxidase activity in sample (b). Assays were performed following the kit protocol.

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: info@promokine.info
www.promokine.info

USA/Canada

Phone: 1 – 866 – 251 – 2860 (toll free)
Fax: 1 – 866 – 827 – 9219 (toll free)

Deutschland

Telefon: 0800 – 776 66 23 (gebührenfrei)
Fax: 0800 – 100 83 06 (gebührenfrei)

France

Téléphone: 0800 90 93 32 (ligne verte)
Téléfax: 0800 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 – 96 03 33 (toll free)
Fax: 0800 – 169 85 54 (toll free)

Other Countries

Phone: +49 6221 – 649 34 0
Fax: +49 6221 – 649 34 40