

Glucose Oxidase Activity Assay Kit (Colorimetric)

ab219924

[3 References](#) [1 Image](#)

Overview

Product name	Glucose Oxidase Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Cell Lysate
Assay type	Enzyme activity (quantitative)
Assay time	0h 20m
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Glucose Oxidase Activity Assay Kit ab219924 uses a simple colorimetric assay to measure glucose oxidase (GOx) activity in cell and tissue extracts, and biological fluids. The assay can be performed in a convenient 96-well or 384-well plate format and is easily adapted to automation without a separation or wash step.</p> <p>The glucose oxidase assay protocol uses a red probe to monitor glucose oxidase activity, leading to an increase in absorbance that can be easily detected at OD575 nm.</p> <p>Glucose oxidase assay protocol summary:</p> <ul style="list-style-type: none">- add samples and standards to wells- add glucose oxidase reaction mix- analyze every 2-3 mins with a microplate reader in kinetic mode for at least 10-30 mins at 37°C
Notes	Glucose oxidase (notatin, GOx, EC 1.1.3.4) is a dimeric enzyme that catalyzes oxidation of beta-D-glucose into hydrogen peroxide and D-glucono-1,5-lactone, which is hydrolyzed to gluconic acid.
Platform	Microplate reader

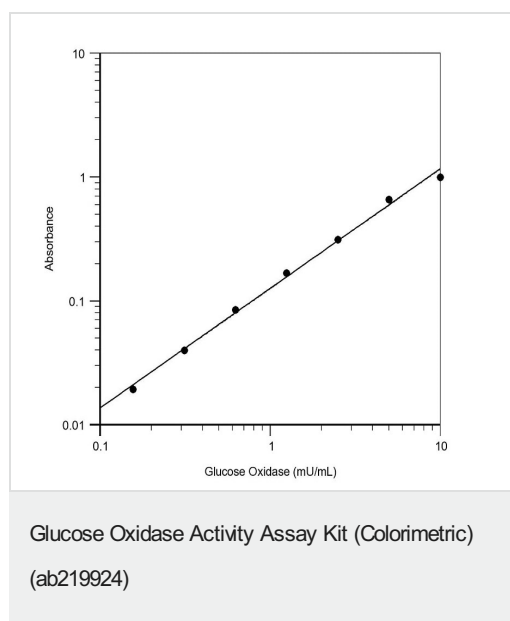
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	500 tests
Assay Buffer	1 x 50ml

Components	500 tests
DMSO	1 x 200µl
Glucose	1 vial
Glucose Oxidase (100 U)	1 vial
Horseradish Peroxidase (HRP)	1 vial
Red Dye (light-sensitive)	1 vial

Images



Glucose Oxidase Activity Assay Kit (Colorimetric) (ab219924).
Typical glucose oxidase dose response curve. Absorbance was measured on a 96-well clear bottom plate using a SpectraMax microplate reader with path check on.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors