

## Product datasheet

# Glucose Oxidase Assay Kit (Fluorometric) ab138884

[1 Image](#)

### Overview

---

<b>Product name</b>	Glucose Oxidase Assay Kit (Fluorometric)
<b>Detection method</b>	Colorimetric/Fluorometric
<b>Sample type</b>	Urine, Plasma, Other biological fluids
<b>Assay type</b>	Enzyme activity
<b>Sensitivity</b>	0.05 mU/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	<p>Glucose Oxidase Assay Kit ab138884 uses a quick and sensitive fluorometric assay for the measurement of glucose oxidase activity. It can be performed in a 96-well or 384-well microtiter plate format and is readily adapted to automation without a separation step.</p> <p>The glucose oxidase assay protocol uses a red dye which enables a dual recordable mode. The fluorescent signal can be easily read by either a fluorescence microplate reader at Ex/Em = 540/590 nm or an absorbance microplate reader at 576 nm.</p> <p>Glucose oxidase assay protocol summary:</p> <ul style="list-style-type: none"><li>- add standards and samples to wells</li><li>- add reaction mix to wells and incubate for 10-30 min at 37°C</li><li>- analyze with a microplate reader</li></ul>
<b>Notes</b>	<p>Glucose oxidase is a dimeric protein that catalyzes the oxidation of beta-D-glucose into hydrogen peroxide and D-glucono-1,5-lactone, which is hydrolyzed to gluconic acid. It is widely used for the determination of glucose in body fluids and in removing residual glucose and oxygen from beverages, food and other agricultural products. Furthermore, Glucose oxidase is commonly used in biosensors to detect glucose.</p>
<b>Platform</b>	Microplate reader

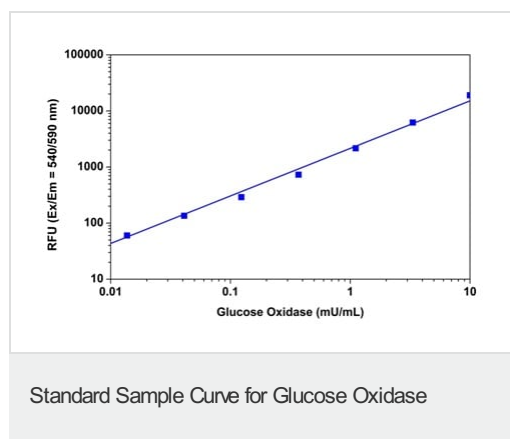
### Properties

---

<b>Storage instructions</b>	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	500 tests
AbRed Indicator	1 vial
Assay Buffer	1 x 50ml
DMSO	1 x 200µl
Glucose oxidase	1 vial
Glucose standard	1 vial
Horseradish Peroxidase	1 vial

## Images



Glucose oxidase dose response was measured with ab138884 in a solid black 96-well plate using a microplate reader. As low as 0.05 mU/ ml glucose oxidase was detected with 30 minutes incubation (n=3).

**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