

# Glucose Dehydrogenase Assay Kit (Colorimetric)

## ab102532

[1 References](#) [2 Images](#)

Overview

|                  |  |
|------------------|--|
| Product name     | Glucose Dehydrogenase Assay Kit (Colorimetric)   |
| Detection method | Colorimetric   |
| Sample type      | Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell culture media   |
| Assay type       | Enzyme activity  |
| Sensitivity      | > 0.01 mU/well   |
| Assay time       | 2h 00m   |
| Product overview | Abcam's Glucose Dehydrogenase Assay Kit (Colorimetric) provides a convenient tool for sensitive detection of the GDH in a variety of samples. The GDH present in sample will recognize D-glucose as a specific substrate leading to a proportional color development. The activity of GDH can be easily quantified colorimetrically ( $\lambda = 450\text{ nm}$ ). Visit our <a href="#">FAQs page</a> for tips and troubleshooting. |
| Notes            | <p>This product is manufactured by BioVision, an Abcam company and was previously called K786 Glucose Dehydrogenase Activity Colorimetric Assay Kit. K786-100 is the same size as the 100 test size of ab102532.</p> <p>This assay detects GDH activity as low as 0.01 mU.</p> <p>Developer (Lyophilized) is light sensitive and should be protected from light during storage and preparation.</p>                                  |
| Platform         | Microplate reader  |

Properties

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

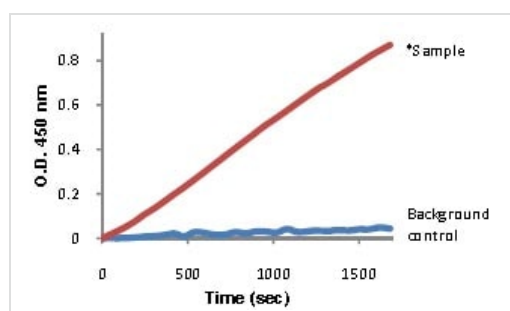
| Components              | 100 tests |
|-------------------------|-----------|
| Developer (Lyophilized) | 1 vial    |
| GDH Assay Buffer        | 1 x 25ml  |
|                         |           |

| Components                         | 100 tests |
|------------------------------------|-----------|
| GDH Positive Control (Lyophilized) | 1 vial    |
| Glucose                            | 1 x 1ml   |
| NADH Standard (Lyophilized)        | 1 vial    |

## Relevance

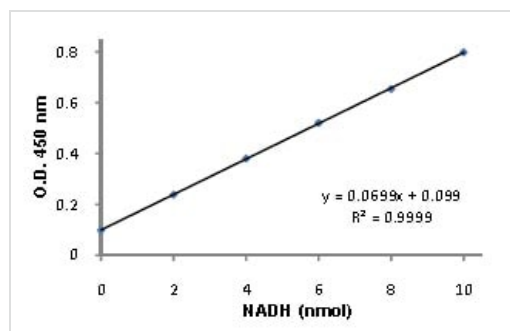
Glucose dehydrogenase (NAD<sup>+</sup>) (EC 1.1.1.118) is an enzyme that catalyzes the chemical reaction: D-glucose + NAD<sup>+</sup> = D-glucono-1,5-lactone + NADH + H<sup>+</sup>. This enzyme belongs to the family of oxidoreductases, specifically those acting on the CH-OH group of donor with NAD<sup>+</sup> or NADP<sup>+</sup> as acceptor.

## Images



Example of Bovine Liver Sample test curves obtained using ab102532

Functional Studies - Glucose Dehydrogenase  
Activity Detection Kit (ab102532)



Example of NADH standard curve obtained using ab102532

Functional Studies - Glucose Dehydrogenase  
Activity Detection Kit (ab102532)

**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