

# Glucose Dehydrogenase Assay Kit (Colorimetric) ab102532

## 1 References [画像数 2](#)

### 製品の概要

製品名	Glucose Dehydrogenase Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell culture media
アッセイタイプ	Enzyme activity
検出感度	> 0.01 mU/well
全工程の試験時間	2h 00m
製品の概要	<p>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 (<math>\lambda = 450 \text{ nm}</math>).</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
特記事項	<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>
試験プラットフォーム	Microplate reader

### 製品の特性

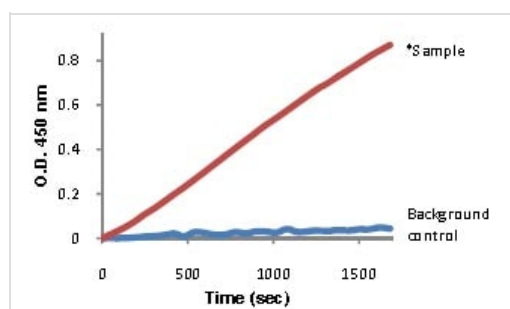
保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Developer (Lyophilized)	1 vial
GDH Assay Buffer	1 x 25ml
GDH Positive Control (Lyophilized)	1 vial

内容	100 tests
Glucose	1 x 1ml
NADH Standard (Lyophilized)	1 vial

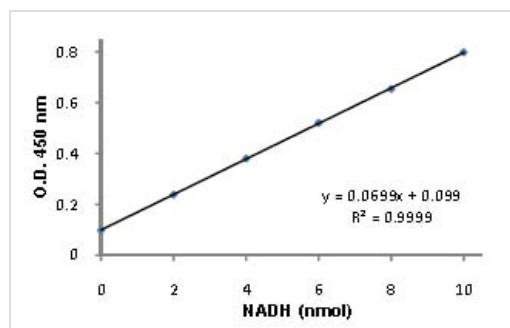
**関連性** Glucose dehydrogenase (NAD<sup>+</sup>) (EC 1.1.1.118) is an enzyme that catalyzes the chemical reaction: D-glucose + NAD<sup>+</sup> = D-glucono-1,5-lactone + NADH + H<sup>+</sup>. This enzyme belongs to the family of oxidoreductases, specifically those acting on the CH-OH group of donor with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

## 画像



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.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

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