

Glucose Dehydrogenase Assay Kit (Colorimetric) ab102532

1 References 2 图像

概述	
产品名称	Glucose Dehydrogenase Assay试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell culture media
检测类型	Enzyme activity
灵敏度	> 0.01 mU/well
检测时间	2h 00m
产品概述	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 FAQs page for tips and troubleshooting.
说明	<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>
平台	Microplate reader
性能	

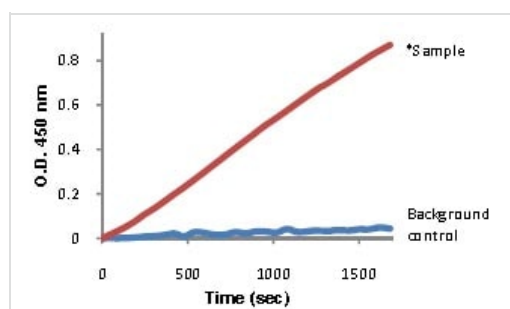
存放说明Store at -20°C. Please refer to protocols.

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

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

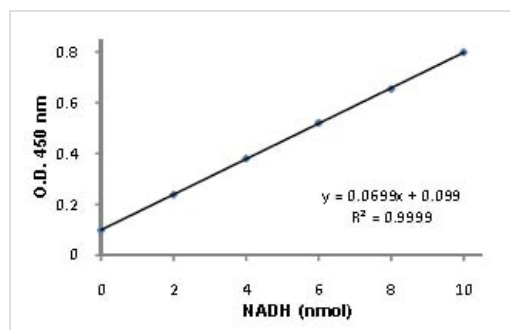
相关性 Glucose dehydrogenase (NAD⁺) (EC 1.1.1.118) is an enzyme that catalyzes the chemical reaction: D-glucose + NAD⁺ = D-glucono-1,5-lactone + NADH + H⁺. This enzyme belongs to the family of oxidoreductases, specifically those acting on the CH-OH group of donor with NAD⁺ or NADP⁺ as acceptor.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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