

Glutamate Dehydrogenase Activity Assay Kit ab102527

[22 References](#) [4 图像](#)

概述

产品名称	Glutamate Dehydrogenase Activity Assay试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
检测类型	Enzyme activity
灵敏度	> 0.01 mU/well
产品概述	<p>Glutamate Dehydrogenase Activity Assay Kit (ab102527) provides a convenient tool for sensitive detection of GDH in a variety of samples. GDH in sample will consume glutamate as a specific substrate and generate NADH stoichiometrically, resulting in a proportional color development. The GDH activity is easily quantified colorimetrically ($\lambda = 450 \text{ nm}$). This assay detects GDH activity as low as 0.01mU in serum or tissue and cell extracts.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K729 Glutamate Dehydrogenase Activity Colorimetric Assay Kit. K729-100 is the same size as the 100 test size of ab102527.</p> <p>Glutamate dehydrogenase (GDH) is an enzyme that converts glutamate to α-Ketoglutarate, and vice versa. It represents a key link between catabolic and metabolic pathways, and is therefore ubiquitous in eukaryotes.</p>
平台	Microplate reader

性能

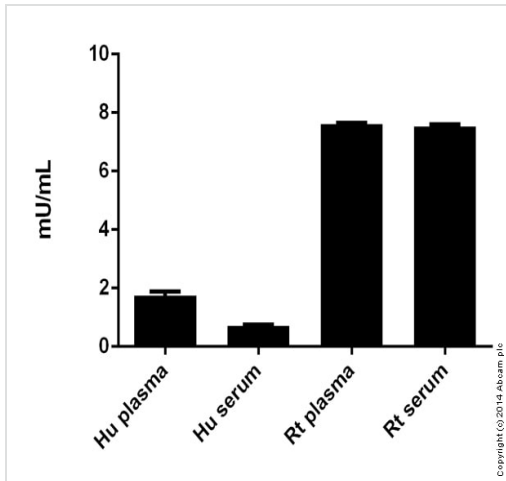
存放说明 Store at -20°C. Please refer to protocols.

组件	100 tests
Developer Solution III	1 vial
GDH Assay Buffer	1 x 25ml
GDH Positive Control	1 vial
Glutamate Solution	1 x 1ml

组件	100 tests
NADH Standard I	1 vial

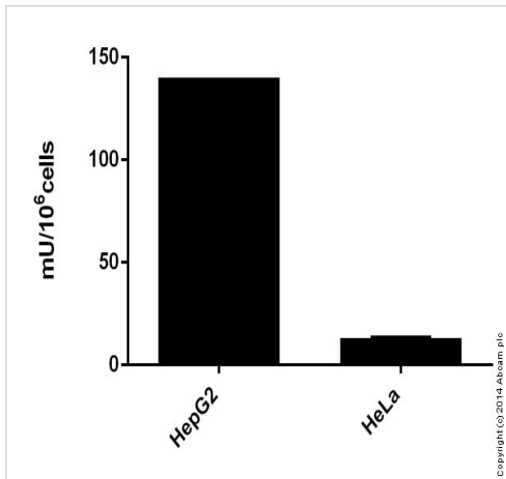
相关性 Glutamate dehydrogenase (GDH) is an enzyme that converts glutamate to alpha-Ketoglutarate, and vice versa. It represents a key link between catabolic and metabolic pathways, and is therefore ubiquitous in eukaryotes.

图片



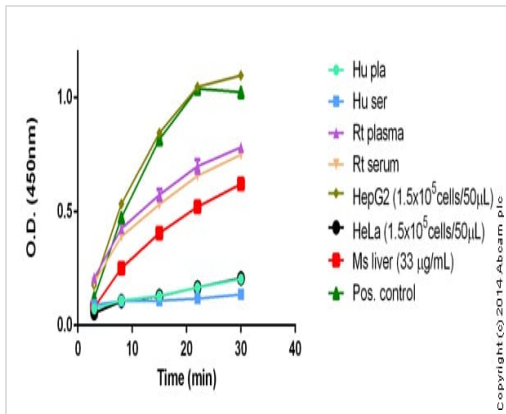
GDH measured showing activity (mU) per mL of tested sample

Functional Studies - Glutamate dehydrogenase
Detection Kit (ab102527)



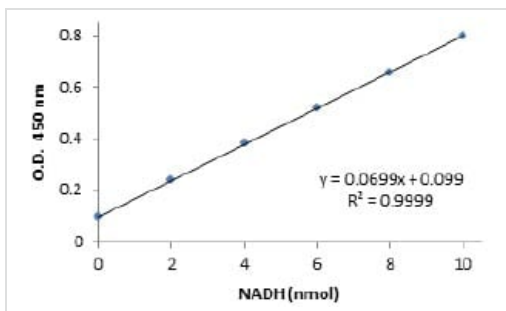
GDH measured in cell lysates showing activity (mU) per 10⁶ cells tested

Functional Studies - Glutamate Dehydrogenase
Detection Kit (ab102527)



GDH measured in various samples showing O.D.'s in time course

Functional Studies - Glutamate Dehydrogenase
Detection Kit (ab102527)



NADH Standard Curve

Functional Studies - Glutamate Dehydrogenase
Detection Kit (ab102527)

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