

Glutamate Dehydrogenase Activity Assay Kit ab102527

[22 References](#) [4 Images](#)

Overview

Product name	Glutamate Dehydrogenase Activity Assay Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Enzyme activity
Sensitivity	> 0.01 mU/well
Product overview	<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>
Notes	<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>
Platform	Microplate reader

Properties

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

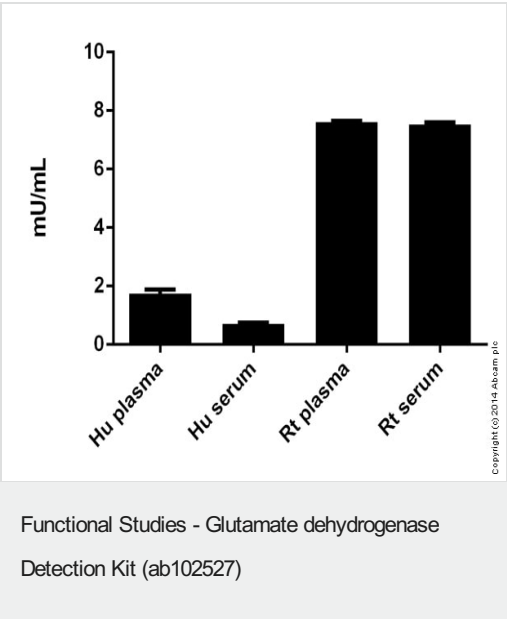
Components	100 tests
Developer Solution III	1 vial
GDH Assay Buffer	1 x 25ml
GDH Positive Control	1 vial
Glutamate Solution	1 x 1ml

Components	100 tests
NADH Standard I	1 vial

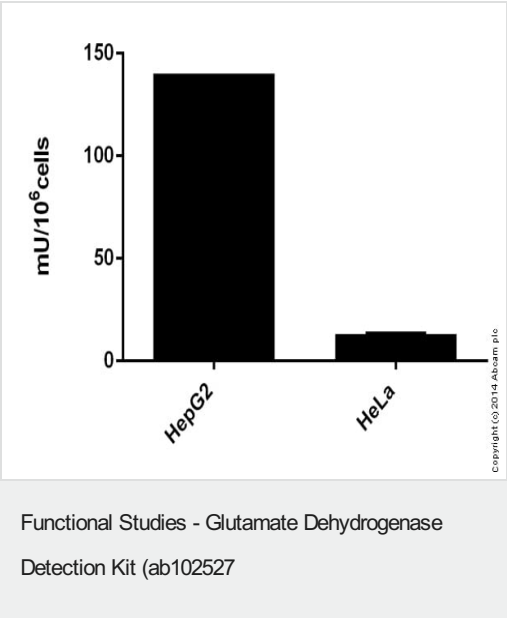
Relevance

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.

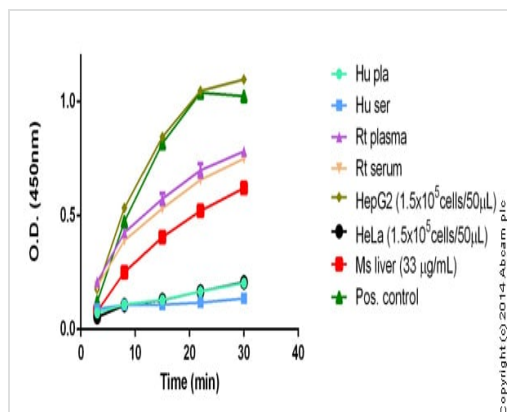
Images



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

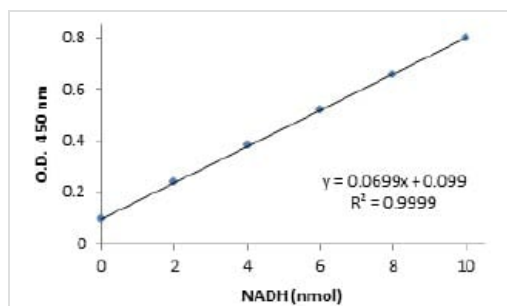


GDH measured in cell lysates showing activity (mU) per 10⁶ cells tested



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.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