

Glutamate Dehydrogenase Activity Assay Kit ab102527

22 References 画像数 4

製品の概要

製品名	Glutamate Dehydrogenase Activity Assay Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Enzyme activity
検出感度	> 0.01 mU/well
製品の概要	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. Visit our FAQs page for tips and troubleshooting.
特記事項	<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>
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.

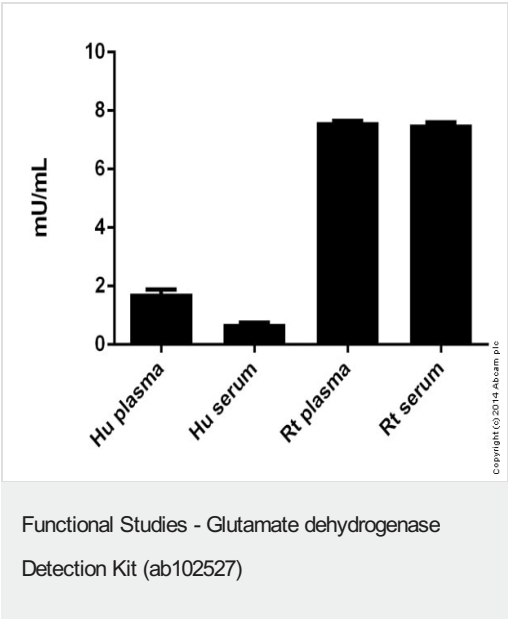
内容	100 tests
Developer Solution III	1 vial
GDH Assay Buffer	1 x 25ml
GDH Positive Control	1 vial
Glutamate Solution	1 x 1ml

内容	100 tests
NADH Standard I	1 vial

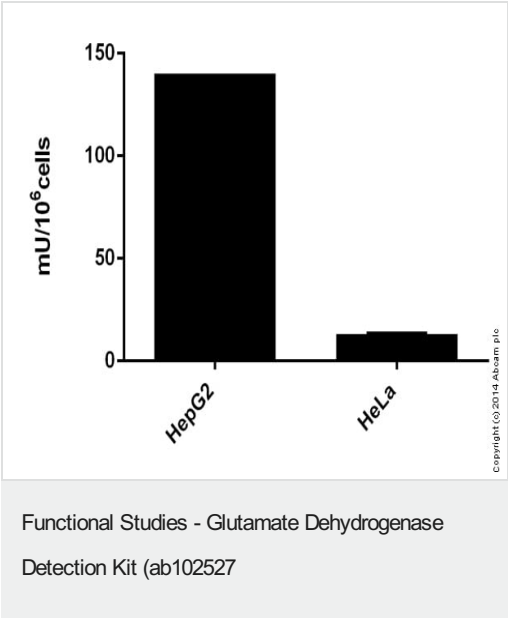
関連性

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.

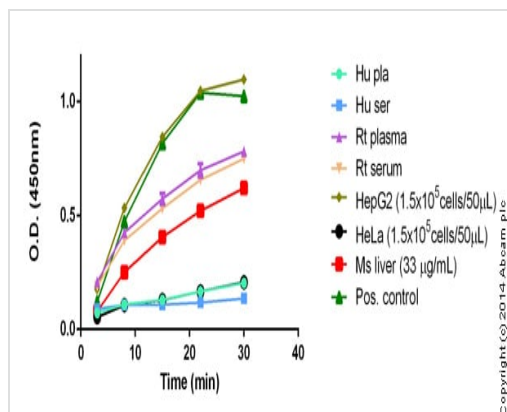
画像



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

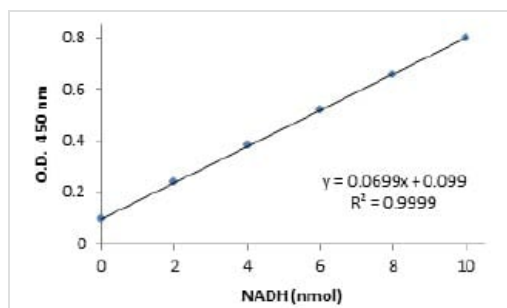


GDH measured in cell lysates showing activity (mU) per 10⁶ cells tested



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.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