

Glucose Assay Kit ab272532

1 References 画像数 1

製品の概要

製品名	Glucose Assay Kit
検出方法	Colorimetric
サンプルの種類	Serum, Plasma, Other biological fluids, Food samples
アッセイタイプ	Quantitative
検出範囲	39 μ M - 16600 μ M
製品の概要	<p>Glucose Assay Kit (ab272532) is a simple, direct and automation-ready procedures for measuring glucose concentrations find wide applications in research and drug discovery. This assay is designed to measure glucose directly in serum or plasma without any pretreatment. The improved otoluidine method utilizes a specific color reaction with glucose. The absorbance at 630nm is directly proportional to glucose concentration in the sample.</p> <p>Sensitive and accurate: Use as little as 5 μL samples. Linear detection range 0.7 mg/dL (39 μM) to 300 mg/dL (16.6 mM) glucose in 96-well plate.</p> <p>Simple and convenient: The procedure involves addition of a single working reagent and incubation for 8 min in a boiling water bath. Improved reagent stability. The optimized formulation has greatly enhanced the reagent and signal stability.</p> <p>Low interference in biological samples: No pretreatments are needed. Assays can be directly performed on serum and plasma samples.</p>
アプリケーション	適用あり: Functional Studies

製品の特性

保存方法 Please refer to protocols.

内容	100 tests
Reagent	1 x 50ml
Standard	1 x 1ml

関連性

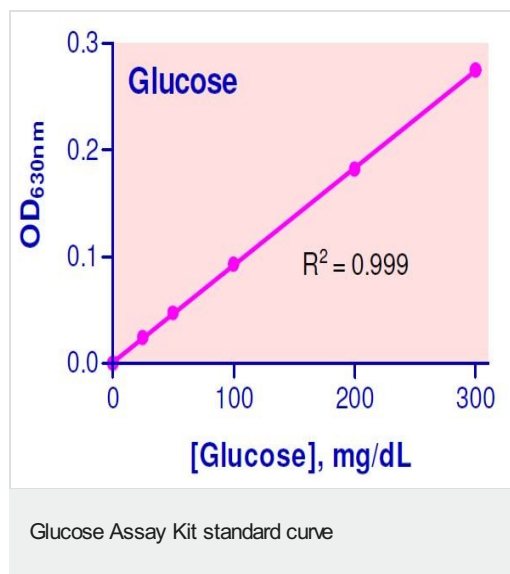
Glucose (C₆H₁₂O₆; FW: 180.16) is a ubiquitous energy source in most organisms, from bacteria to humans. The breakdown of carbohydrates produces mono- and disaccharides, most of which is glucose. Through glycolysis and TCA (citric acid cycle), glucose is oxidized to eventually form CO₂ and water, generating the universal energy molecule ATP. Glucose is a primary source of energy for the brain and a critical component in the production of proteins and in lipid metabolism and therefore measurement of glucose level is a key diagnostic parameter for many metabolic disorders.

アプリケーション

The Abpromise guarantee **Abpromise保証は、次のテスト済みアプリケーションにおけるab272532の使用に適用されます**
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご確認ください。

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

画像



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