

Product datasheet

Glucose Uptake Assay Kit (Colorimetric) ab136955

★★★★★ 1 Abreviews 82 References 3 Images

Overview

Product name	Glucose Uptake Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Sensitivity	<= 0.01 nmol/well
Assay time	3h 00m
Species reactivity	Reacts with: Mammals, Other species
Product overview	Glucose Uptake Assay Kit (Colorimetric) (ab136955) is a highly sensitive and easy to use non-radioactive assay kit which can detect glucose uptake as low as 10 pmol/well in a variety of cell types.

2-deoxyglucose (2-DG) is used in glucose uptake assay protocols because of its structural similarity to glucose. 2-DG is taken up by glucose transporters and metabolized to 2-DG-6-phosphate (2-DG6P). 2-DG6P cannot be further metabolized, and thus accumulates within cells. The accumulated 2-DG6P is directly proportional to 2-DG (or glucose) uptake by cells. In this assay, the 2-DG6P is oxidized to generate NADPH, the level of which can be determined by an enzymatic recycling amplification reaction.

Glucose uptake assay protocol summary:

- prepare cells with suitable glucose starvation / uptake stimulation depending on experimental set-up
- add 2-DG to cells and incubate for 20 mins at 37°C
- wash cells with PBS to remove exogenous 2-DG
- lyse cells with extraction buffer and repeated pipetting
- freeze/thaw lysates and heat at 85°C for 40 min
- cool on ice for 5 min
- add neutralizing buffer, spin and transfer supernatant to new tubes
- add supernatants and standards to wells
- add reaction mix A and incubate for 1 hr at 37°C
- add extraction buffer and heat to 90°C for 40 min
- cool on ice for 5 min and add neutralizing buffer
- add reaction mix B
- analyze every 2-3 mins on microplate reader in kinetic mode at 37°C

Notes

This product is manufactured by BioVision, an Abcam company and was previously called K676 Glucose Uptake Colorimetric Assay Kit. K676-100 is the same size as the 100 test size of ab136955.

Review our [Metabolism Assay Guide](#) to learn about assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also about how to assay metabolic function in live cells using your plate reader.

Platform

Microplate reader

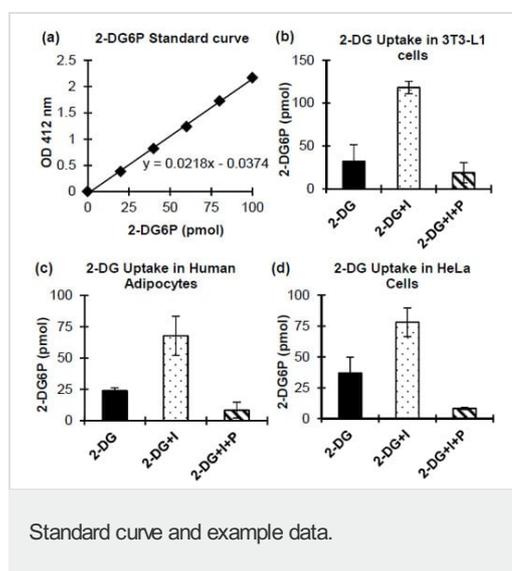
Properties

Storage instructions

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

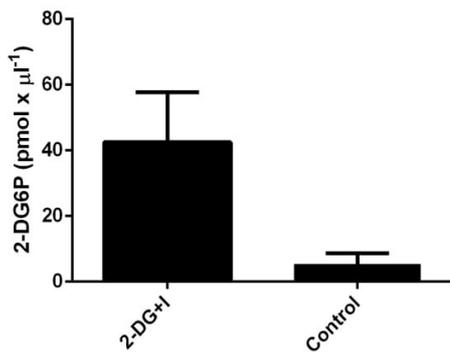
Components	Identifier	100 tests
2-Deoxyglucose	Purple	1 x 1ml
2-DG6P Standard (Lyophilized)	Yellow	1 vial
Assay Buffer	WM	1 x 25ml
Enzyme Mix (Lyophilized)	Orange	1 vial
Extraction Buffer	NM	1 x 17ml
Glutathione Reductase (Lyophilized)	Green	2 vials
Neutralizing Buffer	Clear	1 x 2.5ml
Recycling Mix(Lyophilized)	Blue	1 vial
Substrate	Red	2 vials

Images



2-DG6P Standard curve (a) and 2-DG uptake in 3T3-L1 cells (b), Human adipocytes (c) and HeLa cells (d) respectively. I=Insulin; P=Phloretin.

2-DG Uptake in 3T3-L1 Adipocytes (1500 cells/well)



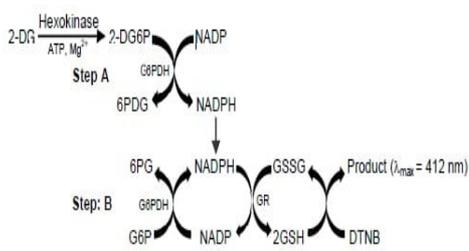
Functional Studies - Glucose Uptake Assay Kit
(ab136955)

Glucose uptake in 3T3-L1 adipocytes stimulated with insulin (I).
3T3-L1 adipocytes were differentiated using:

Dexamethasone [ab120743](#) (1mM, 1:1000)

IBMX [ab120840](#) (11.5 mg/mL, 1:100)

Insulin [ab123768](#) (1 mg/mL, 1:1000)



Assay Procedure

Step A: 2-DG oxidation to generate NADPH; Step B: NADPH recycling amplification Reaction.

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