

Product datasheet

2-NBDG Glucose Uptake Assay Kit ab235976

[7 References](#) [1 Image](#)

Overview

Product name	2-NBDG Glucose Uptake Assay Kit
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Product overview	2-NBDG Glucose Uptake Assay Kit ab235976 provides a convenient tool for studying cellular glucose uptake. The kit uses 2-NBDG, a fluorescently-labeled deoxyglucose analog, as a probe for the detection of glucose taken up by cultured cells.

The 2-NBDG assay kit is suitable for analysis with flow cytometer, fluorescent microscope or microplate reader.

2-NBDG protocol summary:

- treat cells with experimental treatment in glucose-free medium
- add 2-NBDG and incubate for 10 min
- remove cell culture supernatant and wash cells with cell-based assay buffer
- add propidium iodide at this point to exclude dead cells if analyzing with flow cytometry
- add cell-based assay buffer and analyze immediately

2-NDG can be detected with filters designed to detect FITC (Ex/Em 485/535 nm).

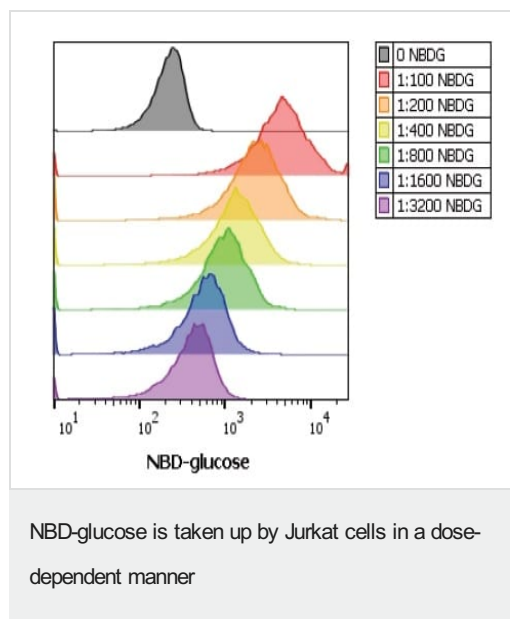
Platform Microplate reader, Fluor. microscope, Flow cyt.

Properties

Storage instructions Please refer to protocols.

Components	1 x 96 tests	1 x 96 tests
Cell-Based Assay Apigenin	1 x 100µl	1 x 100µl
Cell-Based Assay Buffer Tablet	1 tablet	1 tablet
Cell-Based Assay NBD Glucose	1 x 500µl	1 x 500µl
Cell-Based Propidium Iodide Solution	1 x 250µl	1 x 250µl

Images



Jurkat (Human T cell leukemia cell line from peripheral blood) cells were equilibrated in glucose-free medium for two hours prior to being treated with the indicated dilution of NBD-glucose for 10 minutes at 37°C. Cells were washed and data were collected on a MACSQuant cytometer. NBD-glucose fluorescence within the live population is shown.

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