

Glycolysis Assay [Extracellular acidification] ab197244

[26 References](#) [3 图像](#)

概述

产品名称	Glycolysis Assay [Extracellular acidification]
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based (quantitative)
检测时间	1h 30m
产品概述	Glycolysis Assay [Extracellular Acidification] (ab197244) is an easy mix-and-measure, 96 well fluorescence plate reader-based approach for the analysis of extracellular acidification (ECA/ECAR). As lactate production is the main contributor to this acidification, ab197244 is a convenient and informative measure of cellular glycolytic flux. Such measurements offer an important insight into the central role played by altered glycolytic activity in a wide array of physiological and pathophysiological processes, including cellular adaptation to hypoxia and ischemia, and the development and progression of tumorigenesis.

The performance of Glycolysis Assay facilitates sensitive robust microtiter-plate based measurements, thereby overcoming many of the problems associated with the more cumbersome potentiometric pH approach. Rates of extracellular acidification are calculated from changes in fluorescence signal over time and, as the measurement is non-destructive and fully reversible (pH-sensitive reagent is not consumed), measurement of time-courses and multiple drug treatments are possible.

说明 Learn more about the full range of [assays to measure glycolysis, oxygen consumption, fatty acid oxidation and metabolic flux in live cells](#).

Or review the full [metabolism assay guide](#) for other assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress.

平台 Microplate reader

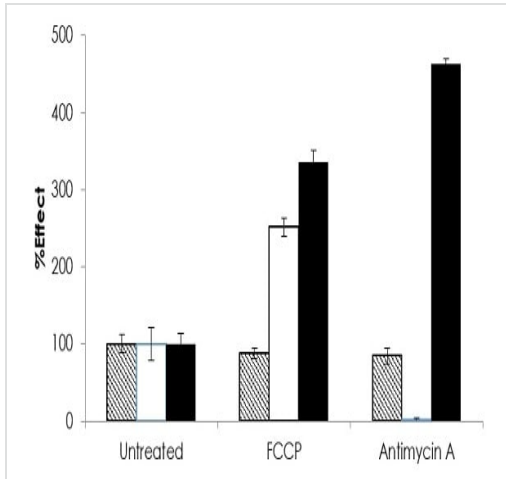
性能

存放说明 Store at +4°C. Please refer to protocols.

组件	96 tests	4 x 96 tests
Glycolysis Assay Reagent (lyophilized)	1 vial	4 vials

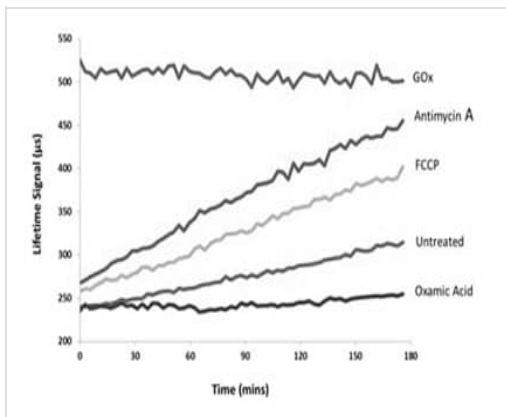
组件	96 tests	4 x 96 tests
Respiration Buffer	1 tablet	4 tablets

图片



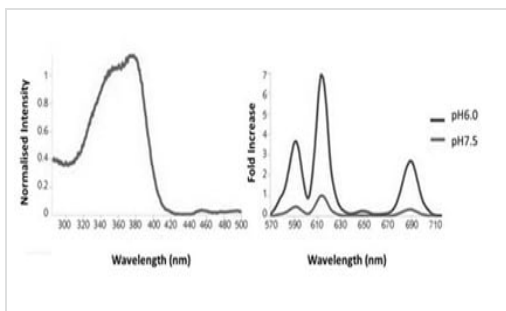
Simultaneous quantification of mitochondrial respiration and glycolytic flux

Cellular Energy Flux for HepG2 cells (seeded at 65,000 per well), treated with a combination of drug compounds modulating the ETC (Antimycin A [1 μ M] and FCCP [2.5 μ M]), shown as a percentage relative to untreated control cells. Comparative measurements were taken with Extracellular Oxygen Consumption Assay ([ab197243](#)) (white column) and Glycolysis Assay [Extracellular acidification] ([ab197244](#)) (black column) show the shift between mitochondrial respiration and glycolysis and the cellular control of energy (ATP; measured 1h post-treatment using Luminescent ATP Detection Assay kit ([ab113849](#)) (striped column)).



Typical lifetime profile

Typical Lifetime profile of Glycolysis Assay for adherent cells, treated with typical control compounds, including Oxamic acid recommended as a negative control. The effect of Glucose Oxidase as a positive signal control is illustrated schematically.



Excitation and emission spectra

Excitation and Emission spectra of Glycolysis Assay. Left panel shows normalized excitation (Ex 340 – 410nm; Peak 360-380nm). Right panel shows emission maxima (Em 590, 615 and 690nm) fold increase between pH6.0 and pH7.5.

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors