

Glycolysis Stress Test Companion Assay ab222945

3 图像

概述

产品名称	Glycolysis Stress Test Companion Assay
检测方法	Fluorescent
样品类型	Adherent cells
检测类型	Cell-based (quantitative)
产品概述	Glycolysis Stress Test Companion Assay (ab222945) allows glycolytic characterization at user-defined glucose concentrations in live cells when used in combination with our <u>Glycolysis Assay [Extracellular Acidification] (ab197244)</u> .

The assay uses 2-Deoxyglucose (2-DG) and oligomycin to facilitate an in-depth analysis of cellular glycolytic flux: 2-DG blocks glycolysis through competitive hexokinase inhibition allowing quantification of non-glycolytic extracellular acidification (ECA); oligomycin inhibits ATP synthase, therefore preventing aerobic ATP generation and forcing cells to increase glycolytic flux to meet ATP demand. The levels of such compensatory glycolysis achievable under test conditions can reveal underlying glycolytic perturbation not evident under basal conditions and facilitates assessment of ATP-demand limited glycolysis capability. Glucose oxidase is included in the kit to provide a convenient cell-free ECA assay signal control.

说明	Learn more about the full range of <u>assays to measure glycolysis, oxygen consumption, fatty acid oxidation and metabolic flux</u> in live cells. Or review the full <u>metabolism assay guide</u> for other assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress.
----	---

平台	Microplate reader
----	-------------------

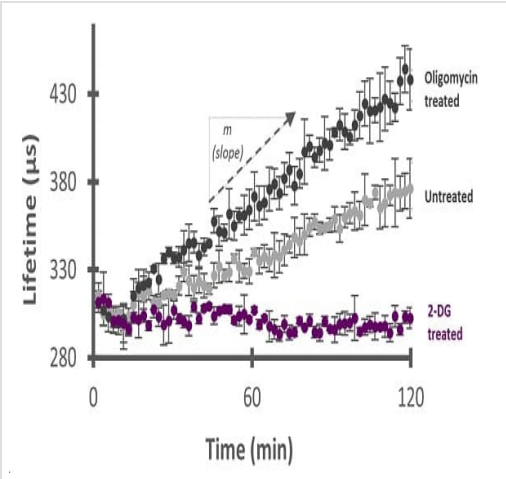
性能

存放说明	Please refer to protocols.
------	----------------------------

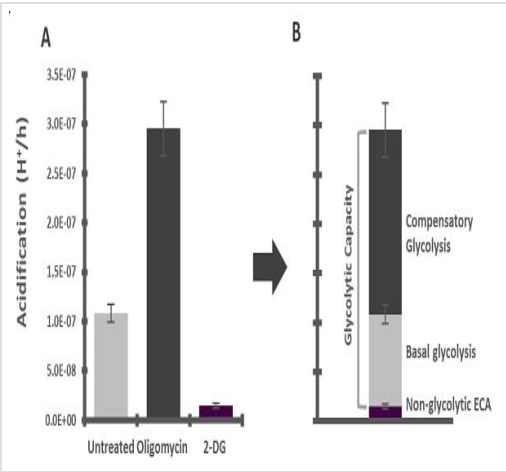
组件	55 tests
2-DG	1 x 600µl
Glucose	1 vial

组件	55 tests
Glucose Oxidase Signal Control	1 vial
Glucose-free Respiration Buffer	1 tablet
Oligomycin	1 vial

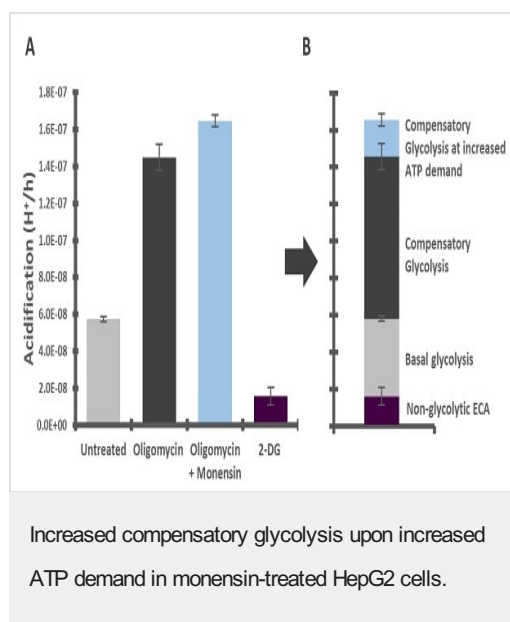
图片



Extracellular acidification of A549 cells treated with 2-DG (purple) and oligomycin (black).



Glycolytic characterization of A549 cells.



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