

Extracellular Oxygen Consumption Assay ab197243

73 References **3 图像**

概述

产品名称	Extracellular Oxygen Consumption Assay
检测方法	Fluorescent
样品类型	Tissue, Adherent cells, Suspension cells, Purified mitochondria
检测类型	Cell-based
检测时间	1h 30m
产品概述	<p>Extracellular Oxygen Consumption Assay Kit ab197243 is a mix-and-read, 96-well fluorescence plate reader assay for the real-time kinetic analysis of extracellular oxygen consumption rates (OCR). The oxygen consumption rate is a measure of the cellular respiration rate, and of mitochondrial function.</p> <p>The assay is optimized for isolated mitochondria and cell cultures, and can be used with tissues, enzyme preparations, and small organisms.</p> <p>The fluorescent dye used in this assay kit is quenched by oxygen. The dye excites at 360-380 nm (max 380) and emits at 630-680 nm (max 650). It is also available separately as ab197242.</p> <p>In the assay, an oil layer is added on top of the assay medium to limit diffusion of oxygen into the assay medium. As mitochondrial respiration depletes the oxygen within the assay medium, quenching of the fluorescent dye is reduced, and the fluorescence signal increases proportionately.</p> <p>The reaction is non-destructive and fully reversible (the oxygen sensitive dye is not consumed) enabling assay time courses and drug treatments.</p>
说明	<p>Learn more about the full range of assays to measure glycolysis, oxygen consumption, fatty acid oxidation and metabolic flux in live cells.</p> <p>Or review the full metabolism assay guide for other assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress.</p>
平台	Microplate reader

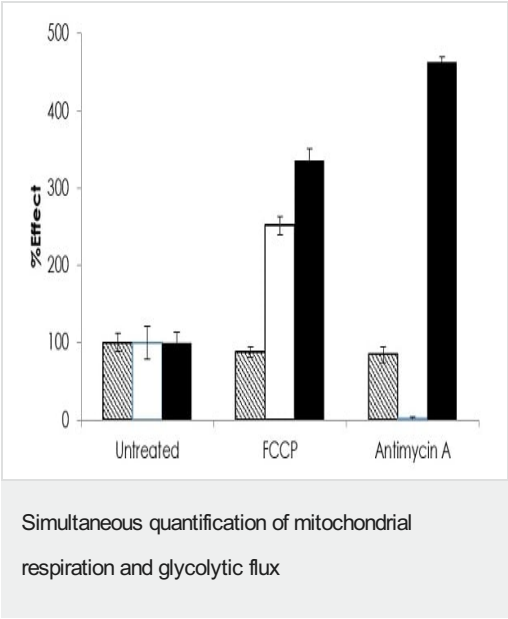
性能

存放说明

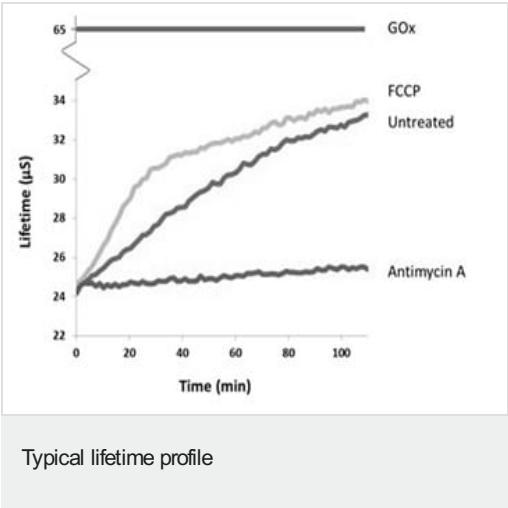
Store at +4°C. Please refer to protocols.

组件	96 tests	4 x 96 tests
Extracellular O2 Consumption Reagent	1 vial	4 vials
High Sensitivity Oil	1 x 15ml	4 x 15ml

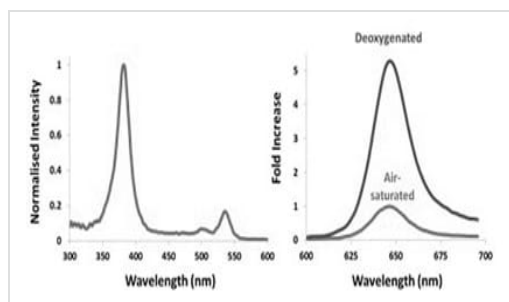
图片



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 of Extracellular O₂ Consumption Assay for adherent cells, treated with different ETC compounds, including Antimycin A (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 Extracellular O₂ Consumption Reagent. Left panel shows normalized excitation (Ex = 360-400nm; Peak 380nm). Right panel shows emission (Em = 630 - 680nm; Peak 650nm) in oxygenated and deoxygenated conditions.

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