

Intracellular Oxygen Concentration Assay ab197245

3 References 画像数 6

製品の概要

製品名	Intracellular Oxygen Concentration Assay
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Quantitative
全工程の試験時間	1h 30m
種交差性	交差種: Human

交差が予測される動物種: Mammals 

製品の概要

Intracellular Oxygen Concentration Assay (ab197245) is an easy mix and measure 96 well fluorescence plate reader based approach for the analysis of intracellular oxygen concentration at the cell monolayer. The assay is based on the ability of oxygen to quench the excited state of the oxygen-sensitive probe. The probe is taken up via nonspecific energy dependent endocytosis and, after washing, the cells are monitored on a fluorescence plate reader (dual-read TR-F required for full oxygen quantitation). The probe phosphorescence is quenched by intracellular oxygen in a non-chemical, reversible manner allowing the measurement of average intracellular O₂ levels and facilitating real-time monitoring of relative changes in cellular oxygen consumption. The probe signal increases with a reduction in intracellular oxygen and decreases with an increase in intracellular oxygen. The probe is excitable at 360-400 or 535 nm and emits at 630-680 nm. Optimal filter combinations are Ex/Em = 340/642 nm.

The flexible plate reader format, allows multiparametric or multiplex combination with a range of other reagents and it is suitable for HTP automation.

特記事項

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.

試験プラットフォーム

Microplate reader

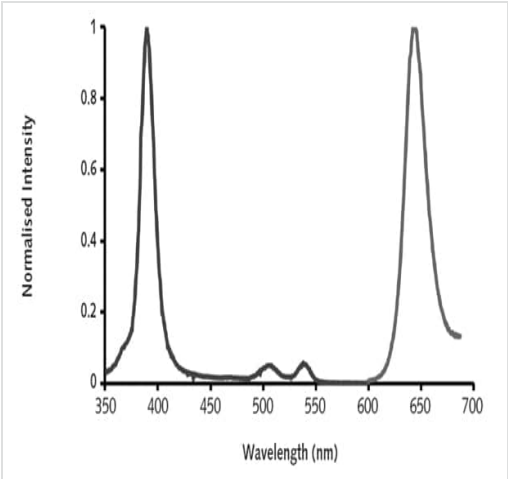
製品の特性

保存方法

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

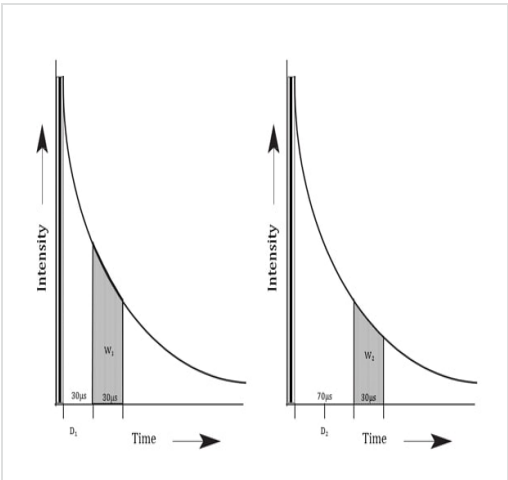
内容	96 tests	4 x 96 tests
Intracellular O2 probe	1 vial	4 vials

画像

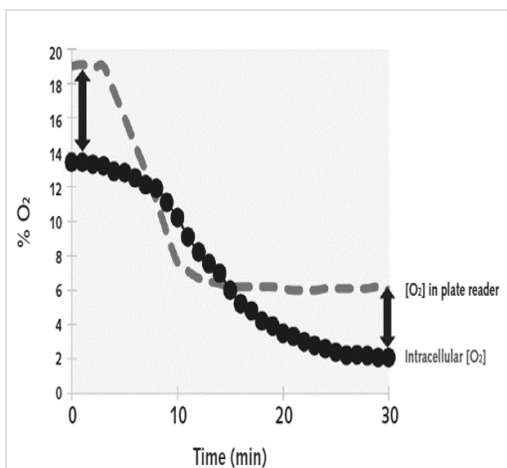


Excitation and Emission spectra of Intracellular O₂ probe, showing normalized excitation (Ex 360-400nm; Peak 380nm) and emission (Em 630-670nm; Peak 650nm).

Excitation and emission spectra of O₂ probe

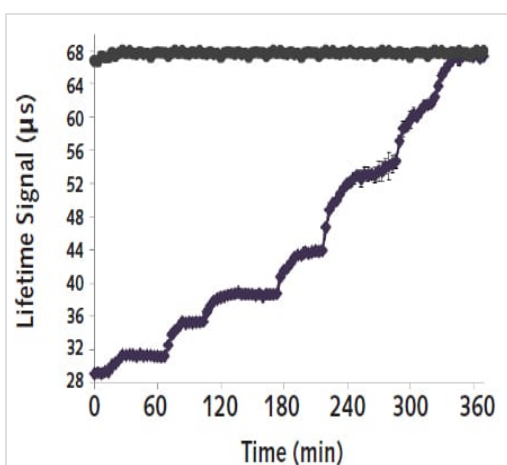


Illustrating dual read TR-F measurement.



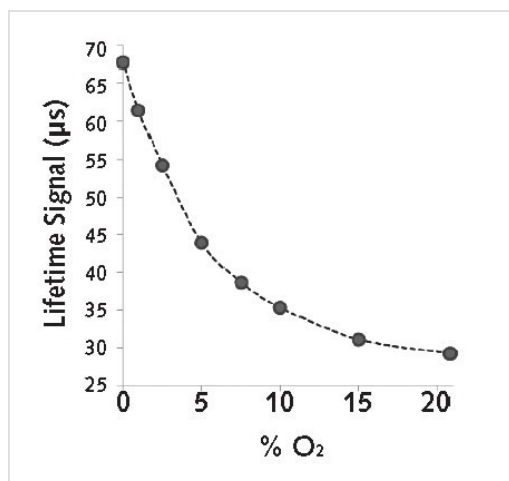
HEK293T cell oxygenation

HEK293T cell oxygenation. HEK293T cells were cultured in 2D and measured at ambient oxygen. Intracellular O₂ levels were ~ 14%. Reducing instrument O₂ to 6% caused cellular oxygenation to drop to ~2%. Assay performed using a CLARIOstart equipped with an ACU module (BMG Labtech).



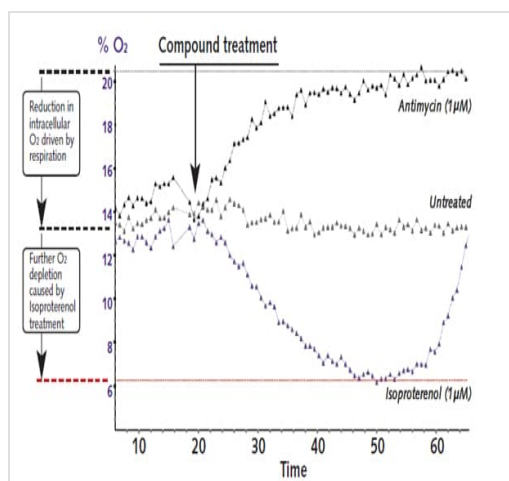
Sample callibration data.

Sample Calibration Data. Intra O₂ probe Lifetime profiles measured at decreasing [O₂] with parallel glucose oxidase treatment to achieve 0% O₂.



Relationship between probe lifetime (τ) and applied $[O_2]$. Applying a first order exponential fit generates a calibration function of $O_2\% = A1 \times \text{Exp}(-\tau / t1)$. Example: $O_2\% = 659.3 \times \text{Exp}(-\tau / 8.475)$.

Relationship between probe lifetime (τ) and applied $[O_2]$.



Measuring the impact of cell metabolism on iPS-derived cardiomyocyte oxygenation. During measurement, cells are treated with antimycin (ETC inhibitor) and isoproterenol (β -adrenoreceptor agonist).

Measuring the impact of cell metabolism on iPS-derived cardiomyocyte oxygenation.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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