

Cell Viability Assay Kit (Fluorometric - Dual Green/Red) ab112121

画像数 1

製品の概要

製品名	Cell Viability Assay Kit (Fluorometric - Dual Green/Red)
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Quantitative
種交差性	交差種: Mammals, Other species
製品の概要	<p>Abcam's Cell Viability assay kits are a set of tools for monitoring cell viability and cellular functions. ab112121 uses two non-fluorescent indicators: The CellGreen fluorescent dye for viable cells and a cell-impermeable DNA-binding dye for the cells with compromised membranes. The non-fluorescent Green Indicator Dye can easily permeate intact live cells and is hydrolyzed by intracellular esterase to generate the strongly fluorescent hydrophilic CellGreen fluorescent dye which is well-retained in the cell cytoplasm. The esterase activity is proportional to the number of viable cells. The DNA-binding dye is quite polar and impermeable for viable cells that have intact membranes. It becomes fluorescent only upon binding to the DNA of dead cells. Cells grown in black-wall plates can be stained and quantified in less than two hours.</p> <p>ab112121 is more robust and accurate than the other viability assays. It can be readily adapted for a wide variety of fluorescence platforms such as microplate assays, fluorescence microscope, and flow cytometry. The kit provides all the essential components with an optimized assay protocol. It is suitable for both proliferating and non-proliferating cells (either suspension or adherent cells). ab112121 comes with sufficient reagents to perform either 200 assays (96-well format) or 800 assays (a 384-well format).</p>
特記事項	<p>ab112121 should be stored desiccated. Assay kit comes with sufficient reagents to perform either 200 assays (96-well format) or 800 assays (a 384-well format).</p> <p>Related assays</p> <p>Review the cell health assay guide to learn about kits to perform a cell viability assay, cytotoxicity assay and cell proliferation assay.</p>
試験プラットフォーム	Microplate reader, Fluor. microscope, Flow cyt.

製品の特性

保存方法

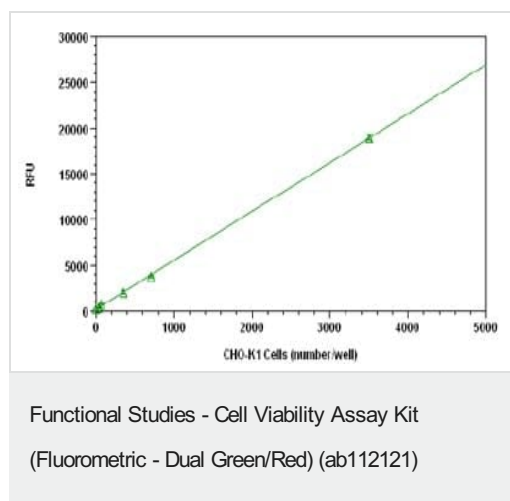
Store at -20°C. Please refer to protocols.

内容	2 x 96 tests
Assay Buffer	1 x 20ml
CellGreen fluorescent dye	2 vials
DMSO	1 x 100µl
Propidium Iodide	1 x 40µl

関連性

Cell viability is a determination of living or dead cells, based on a total cell population. Cell viability assess healthy cells in a sample, with no distinction between dividing or quiescent cells. An increase in cell viability indicates cell growth, while a decrease in viability can generally be interpreted as the result of either toxic effects of compounds/agents or suboptimal culture conditions.

画像



CHO-K1 cell number response was measured with ab112121. CHO-K1 cells at 0 to 5,000 cells/well/100 µL were seeded overnight in a black wall/clear bottom 96-well plate. The cells were incubated with 100 µL/well of Green dye-loading solution for 1 hour at 37 °C. The fluorescence intensity was measured at Ex/Em = 490/ 525 nm. The fluorescence intensity was linear ($R^2 = 1$) to the cell number as indicated. The detection limit was 30 cells/well (n=6).

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