

Cell Viability Assay Kit (Fluorometric - Green) ab112122

1 References [1 图像](#)

概述

产品名称	Cell Viability Assay试剂盒(Fluorometric - Green)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Quantitative
种属反应性	与反应: Mammals, Other species
产品概述	<p>There are a variety of parameters that can be used to monitor cell viability. The green fluorescent dye used in the kit is a hydrophobic compound. It easily permeates intact live cells and gets enhanced fluorescence upon entering into live cells. The hydrolysis of the non-fluorescent substrate by intracellular esterases generates a strongly green fluorescent hydrophilic product that is well-retained in the cell cytoplasm. The green fluorophore generated by the non-fluorescent substrate used in the kit has the spectral properties of fluorescein at Ex/Em = ~490 nm/520 nm. When well excited with the Argon Laser at 488 nm, the fluorophore emits intense green fluorescence at ~520 nm.</p> <p>ab112122 provides all the essential components with an optimized cell-labeling protocol for fluorescence microplate assays. It can also be used with a fluorescence microscope equipped with a FITC filter set.</p> <p>ab112122 provides an effective tool of labeling cells for fluorescence microplate and microscopic investigations of cellular functions. It is useful for a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity. ab112122 is suitable for proliferating and non-proliferating cells.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	Review the cell health assay guide to learn about more kits to perform a cell viability assay , cytotoxicity assay and cell proliferation assay .
平台	Microplate reader, Fluor. microscope, Flow cyt.

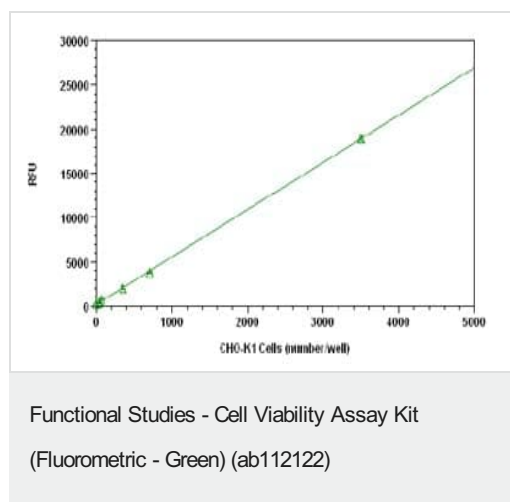
性能

存放说明 Store at -20°C. Please refer to protocols.

组件	5 x 96 tests
Assay Buffer	1 x 50ml
CellGreen fluorescent dye	5 vials
DMSO	1 x 200µl

相关性 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.

图片



CHO-K1 cell number response was measured with ab112122. 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.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