

Cell Viability Assay Kit (Fluorometric - Blue) ab112120

1 Image

Overview

Product name	Cell Viability Assay Kit (Fluorometric - Blue)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Quantitative
Assay time	2h 00m
Species reactivity	Reacts with: Mammals, Other species
Product overview	Abcam Cell Viability Assay Kits are a set of tools for monitoring cell viability and cellular functions. There are a variety of parameters that can be used to monitor cell viability.

ab112120 uses our proprietary cell viability dye whose fluorescence is strongly enhanced upon entering into live cells. The dye is a hydrophobic compound that easily permeates intact live cells. The weakly fluorescent Blue probe is hydrolyzed by intracellular esterase to generate a strongly fluorescent hydrophilic product that is well-retained in the cell cytoplasm. The esterase activity is proportional to the number of viable cells, and thus directly related to the fluorescence intensity of the product generated from the esterase-catalyzed hydrolysis of the fluorogenic substrate. Cells grown in black wall/clear bottom plates can be stained and quantified in less than two hours.

ab112120 is more robust than tetrazolium salt based assays. It can be readily adapted for many different types of fluorescence platforms such as microplate assays, fluorescence microscope, and flow cytometry.

ab112120 is useful for a variety of studies, including cell adhesion, chemotaxis, multidrug resistance, cell viability, apoptosis and cytotoxicity. It provides all the essential components with an optimized cell-labeling protocol and can be used for both suspension and adherent cells.

Visit our [FAQs page](#) for tips and troubleshooting.

Notes Review the [cell health assay guide](#) to learn about more kits to perform a [cell viability assay](#), [cytotoxicity assay](#) and [cell proliferation assay](#).

Platform Microplate reader, Fluor. microscope, Flow cyt.

Properties

Storage instructions

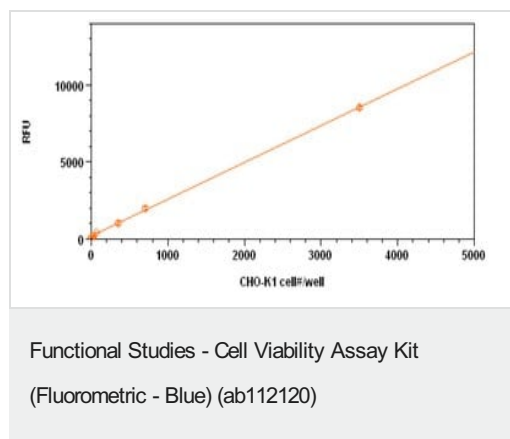
Store at -20°C. Please refer to protocols.

Components	5 x 96 tests
Assay Buffer	1 x 50ml
Cell Blue dye	5 vials
DMSO	1 x 200µl

Relevance

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.

Images



CHO-K1 cell number response was measured using ab112120. 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 blue indicator solution for 1 hour at room temperature. The fluorescence intensity was measured at Ex/Em = 360/450 nm. The fluorescence intensity was linear ($R^2 = 1$) to the cell number as indicated. The detection limit was 70 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.com/abpromise> or contact our technical team.

Terms and conditions

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

