

Cell Viability Assay Kit (Fluorometric - Blue Ex 405 nm) ab176748

[1 References](#) [画像数 2](#)

製品の概要

製品名	Cell Viability Assay Kit (Fluorometric - Blue Ex 405 nm)
検出方法	Fluorescent
サンプルの種類	Suspension cells
アッセイタイプ	Cell-based (quantitative)
種交差性	交差種: Mammals, Other species
製品の概要	<p>Cell Viability Assay Kit (Fluorometric - Blue Ex 405 nm) (ab176748) uses a proprietary cell viability dye whose fluorescence is strongly enhanced upon entering into live cells.</p> <p>The dye is a hydrophobic compound that easily permeates intact live cells. The weakly fluorescent CytoCalcein Violet 450, AM is hydrolyzed by intracellular esterase to generate a strongly fluorescent hydrophilic product that is well-retained in the cell cytoplasm.</p> <p>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.</p> <p>Cells grown in black wall/clear bottom plates can be stained and quantified in less than two hours.</p> <p>The assay is more robust than tetrazolium salt or Alamar Blue® based ones. It can be readily adapted for many different types of fluorescence platforms such as microplate assays, fluorescence microscope, and flow cytometry.</p>
特記事項	Review the cell health assay guide to learn about more kits to perform a cell viability assay , cytotoxicity assay and cell proliferation assay .
試験プラットフォーム	Microplate reader, Fluor. microscope, Flow cyt.

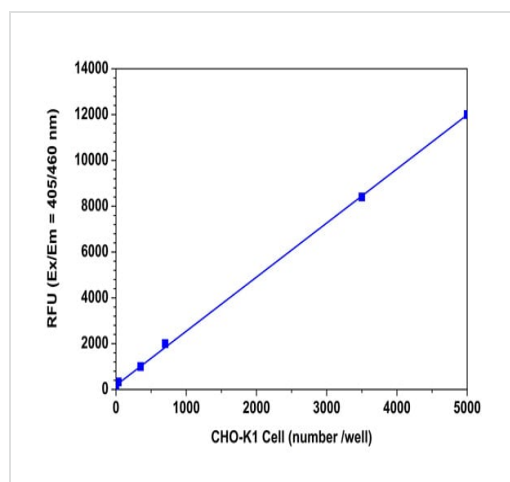
製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	500 tests
Assay Buffer	1 x 50ml
CytoCalcein Violet 450, AM	5 vials
DMSO	1 x 200 μ l

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.

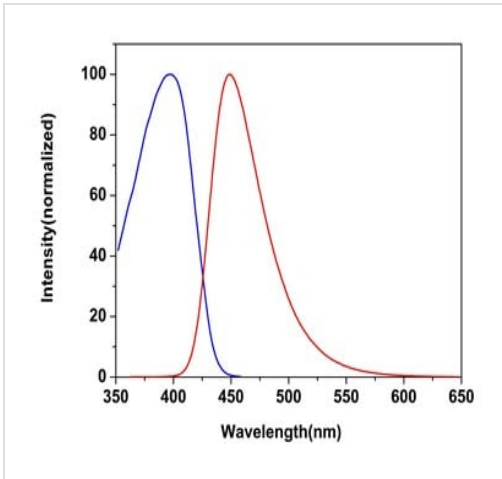
画像



CHO-K1 cell number measured with Cell Viability Assay Kit (Fluorometric-Blue Ex 405nm)

CHO-K1 cell number response was measured with Cell Viability Assay Kit (Fluorometric - Blue Ex 405 nm) (ab176748). CHO-K1 cells were seeded overnight at 0 - 5,000 cells/well/100 μ L in a black wall/clear bottom 96-well plate. The cells were incubated with 100 μ L/well of CytoCalcein Violet 450 dye-loading solution for 1 hour at room temperature. The fluorescence intensity was measured at Ex/Em = 405/460 nm. The fluorescence intensity was linear ($R^2 = 1$) to the cell number as indicated.

The detection limit was 70 cells/well (n=6).



Excitation and Emission Spectra for Cell Viability
Assay Kit (Fluorometric - Blue Ex 405 nm)
(ab176748)

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