

# Resazurin Assay Kit - Cell Cytotoxicity (Fluorometric) ab112119

[3 References](#) [画像数 1](#)

### 製品の概要

製品名	Resazurin Assay Kit - Cell Cytotoxicity (Fluorometric)
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 00m
種交差性	交差種: Mammals, Other species
製品の概要	

#### Recommended alternative kit

We recommend that you purchase [Resazurin Assay Kit ab129732](#) instead of this kit.

Resazurin Assay Kit ab112119 (Cell Cytotoxicity, Fluorometric) ab112119 provides a fast, simple, accurate and homogeneous assay for the fluorometric detection of viable cells.

The resazurin assay protocol is based on the reduction of oxidized non-fluorescent blue resazurin to a red fluorescent dye (resorufin) by the mitochondrial respiratory chain in live cells. The amount of resorufin produced is directly proportional to the number of living cells. Resorufin has Ex/Em of 530-560 / 590 nm.

The resazurin assay is very commonly used to quantify the numbers of live cells, and monitor cell viability / cytotoxicity.

ab112119 is more sensitive for cell proliferation and cytotoxicity than other assays such as MTT. The kit components are stable and have minimal cytotoxicity, thus a longer incubation time (such as 24-48 hours) is possible if required.

Resazurin assay protocol summary:

- add assay solution to cells in 96 or 384 well plates and shake gently for 30 sec
- incubate for 1-24 hr at 37°C
- analyze with microplate reader

### 特記事項

ab112119 should be stored desiccated

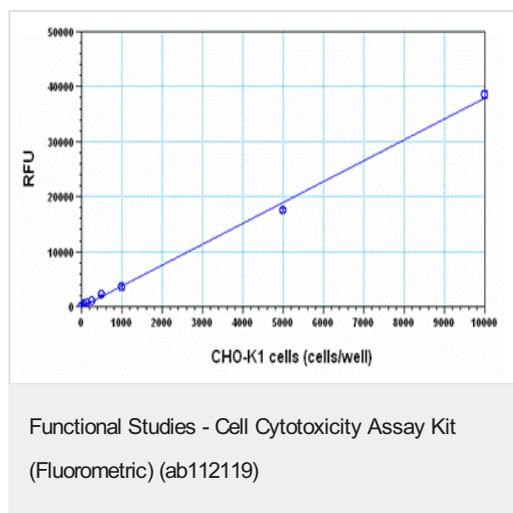
製品の特性

保存方法

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

内容	1000 tests
Assay Solution	1 x 20ml

画像



CHO-K1 cell number response was measured with ab112119. CHO-K1 cells at 0 to 10,000 cells/well/100 µL were seeded overnight in a black wall/clear bottom 96-well plate. The cells were incubated with 20 µL/well of Assay Solution for 3 hours at 37 °C. The absorbance intensity was measured at 570 nm and 605 nm. The fluorescence intensity was linear ( $R^2 = 0.998$ ) to the cell number as indicated. The detection limit from the blank controls was 60 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