

## Product datasheet

# Cell Cytotoxicity Assay Kit (Colorimetric) ab112118

★★★★★ [1 Abreviews](#) [23 References](#) [2 Images](#)

### Overview

---

<b>Product name</b>	Cell Cytotoxicity Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Quantitative
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	<p>Monitoring cell cytotoxicity is one of the most essential tasks for studying cellular functions. There are a variety of parameters that can be used. ab112118 uses a proprietary water-soluble dye that changes its absorption spectra upon cellular reduction. The absorption ratio change is directly proportional to the number of living cells. The characteristics of its high sensitivity, non-radioactivity and no-wash method make ab112118 suitable for high throughput screening of cell proliferation or cytotoxicity against a variety of compounds.</p> <p>ab112118 does not require pre-mixing of components and has higher sensitivity compared to tetrazolium based colorimetric assays (such as MTT and XTT). It comes with reagents sufficient to run 1000 assays. The kit components are quite stable with minimal cytotoxicity, thus longer incubation times (such as 24-48 hours) are possible if required.</p> <p>ab112118 is robust and convenient to use. It can be readily adapted for a wide variety of instrument platforms. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p> <p>Cytotoxicity assay protocol summary:</p> <ul style="list-style-type: none"><li>- add assay solution to cells and shake for 30 s</li><li>- incubate for 1-24 hrs</li><li>- analyze with microplate reader</li></ul>
<b>Notes</b>	<p>ab112118 should be stored desiccated.</p> <p>As low as 300 cells can be accurately quantified using ab112118.</p>
<b>Platform</b>	Microplate reader

### Properties

---

## Storage instructions

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

Components	Identifier	1000 tests
Assay Solution	Component A	1 x 20ml

## Images

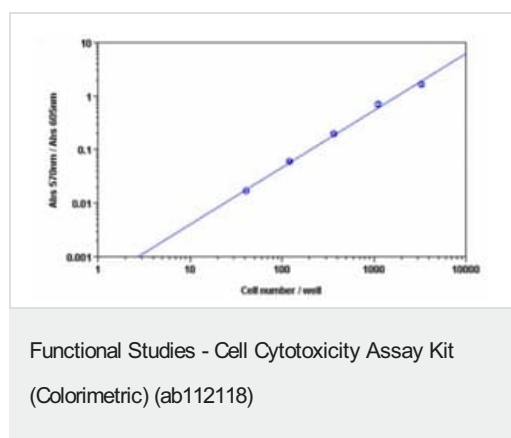
% Cell viability		
	HeLa	NIH3T3
1 $\mu$ M	28.51%	56.03%
10 $\mu$ M	12.78%	43.51%
50 $\mu$ M	10.52%	34.02%

Copyright (c) 2014 Abcam plc

Functional Studies - Cell Cytotoxicity Assay Kit  
(Colorimetric) (ab112118)

•HeLa: day before experiment cells seeded in 96 wp @ 1e4 cells/well and treated with 1  $\mu$ M, 10  $\mu$ M and 50  $\mu$ M Camptothecin (incubated for 24h); on the day of experiment cells incubated with Assay Solution for 2 hours.

•NIH3T3: day before experiment cells seeded in 96 wp @ 1e4 cells/well and treated with 1  $\mu$ M, 10  $\mu$ M and 50  $\mu$ M Camptothecin (incubated for 24h); on the day of experiment cells incubated with Assay Solution for 2 hours.



CHO-K1 cell number response was measured with ab112118. CHO-K1 cells at 0 to 10,000 cells/well/100  $\mu$ L were seeded overnight in a black wall/clear bottom 96-well plate. The cells were incubated with 20  $\mu$ L/well of Assay Solution for 3 hours at 37 °C. The absorbance intensity was measured at 570 nm and 605 nm. The ratio of OD570/OD605 is proportional to the number of cells as indicated.

**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