

# Cell Viability Assay Kit (Fluorometric - Blue) ab102501

[3 References](#) [2 图像](#)

### 概述

产品名称	Cell Viability Assay试剂盒(Fluorometric - Blue)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Quantitative
灵敏度	> 100 cells/well
范围	100 cells/well - 100000 cells/well
实验步骤	One step assay
种属反应性	与反应: Mammals
产品概述	<p>This Cell Proliferation Assay Kit (Fluorometric-Blue) ab102501 (previously known as VisionBlue Quick Cell Viability Fluorometric Assay K303) provides by far the most sensitive and easiest means for quantifying cell proliferation and viability. Simply add the single reagent to cell culture, incubate and read the fluorescence intensity. The assay utilizes the redox dye resazurin which is not fluorescent, but upon reduction by viable metabolically active cells, the dye becomes highly fluorescent (Ex = 530-570 nm; Em = 590-620 nm). Therefore, viable metabolically active cells can be easily measured.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K303 VisionBlue™ Quick Cell Viability Fluorometric Assay Kit. K303-2500 is the same size as the 2500 test size of ab102501.</p> <p>Key Features:</p> <ul style="list-style-type: none"><li>Simple procedure: Just Add-Incubate-Read. No washing, No solubilization.</li><li>Highly Sensitive: Detect as few as 100 cells.</li><li>Longer Stability: Reagent is stable at 4C for at least a year.</li><li>Safe: Non-radioactive, non-toxic, and cells can be further used for other experiments.</li><li>Wide applications: Used in studying a variety of growth stimulations or inhibitions (e.g., by growth factors, cytokine, nutrients, anticancer drugs, apoptosis inducers/inhibitors, toxicityinducing chemicals, etc.).</li></ul>
平台	Microplate

### 性能

**存放说明**

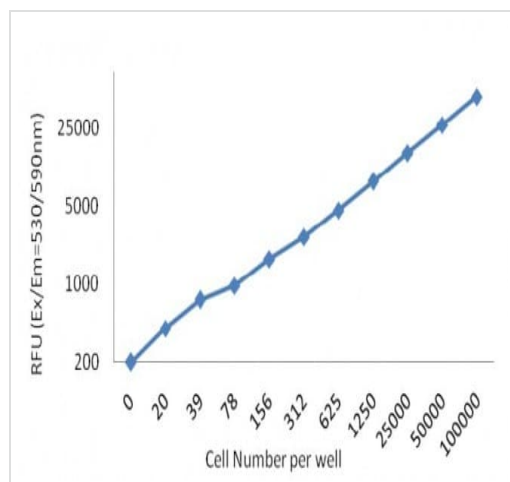
Store at -20°C. Please refer to protocols.

组件	500 tests	2500 tests
Cell Viability Dye	1 x 5ml	1 x 25ml

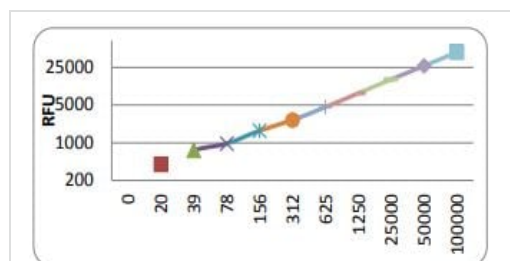
**相关性**

Cell proliferation is the multiplication or reproduction of cells, as a result of cell growth and cell division, resulting in the expansion of a cell population.

**图片**



Cell Viability Assay Kit (Fluorometric - Blue)



Cell Viability Assay Kit (Fluorometric - Blue)

**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