

# WST-8 Assay Kit (Cell Proliferation) ab65475

[20 References](#) [4 图像](#)

### 概述

|      |   |
|------|---|
| 产品名称 | WST-8 Assay试剂盒(Cell Proliferation)  |
| 检测方法 | Colorimetric  |
| 样品类型 | Adherent cells, Suspension cells  |
| 检测类型 | Quantitative  |
| 检测时间 | 4h 00m  |
| 产品概述 | <p>WST-8 Assay Kit ab65475 (previously called WST-1 Assay Kit) is an alternative version of <a href="#">ab65473</a>, our most popular WST-1 Assay Kit (which is also supplied as a ready-to-use reagent <a href="#">ab155902</a>). This kit uses a different WST-8 analog, and also contains a stop solution.</p> |

The WST-8 assay provides by far the easiest and most sensitive means for performing a quantitative cell proliferation assay, cell viability assay, or cytotoxicity assay in mammalian cells.

The WST-8 assay protocol is based on the cleavage of the tetrazolium salt to formazan by cellular mitochondrial dehydrogenase. The amount of the dye generated by activity of dehydrogenase is directly proportional to the number of living cells. The formazan dye produced by viable cells can be quantified by microplate reader by measuring the absorbance of the dye solution at 440 nm.

The WST-8 assay protocol is very simple:

- add the WST-8 assay reagent to the cell culture media and incubate for between 0.5 and 4 hrs
- shake the plate to mix the contents
- analyze the amount of formazan dye produced by measuring the absorbance at 440 nm

This kit contains an optional stop solution to stop the production of formazan dye.

### 说明

This product is manufactured by BioVision, an Abcam company and was previously called K302 Quick Cell Proliferation colorimetric Assay Kit Plus. K302-2500 is the same size as the 2500 test size of ab65475.

The WST-8 assay is add-and-read, and requires no washing, harvesting, or solubilization steps. It is faster, stable, and more sensitive than MTT, XTT, MTS based assays. The assay correlates well with the [3H]-thymidine incorporation assay.

Review our [cell health assays guide](#) to learn more about our other [cell viability](#), [cytotoxicity](#) and [cell proliferation](#) assay kits.

### 平台

Microplate reader

## 性能

### 存放说明

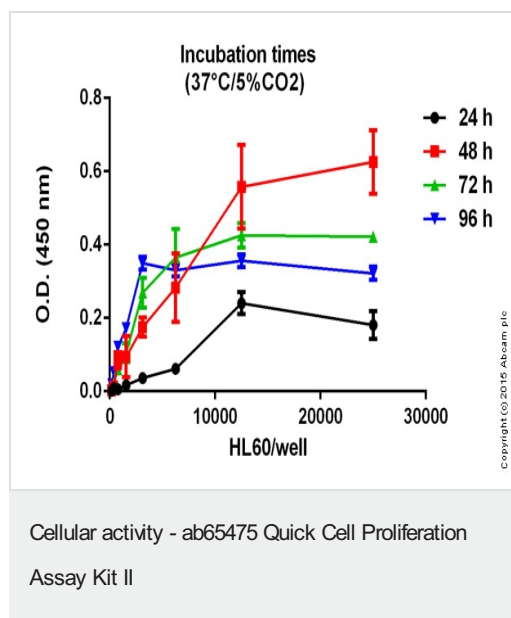
Store at -20°C. Please refer to protocols.

| 组件                       | 500 tests | 2500 tests |
|--------------------------|-----------|------------|
| Electrocoupling Solution | 1 x 5ml   | 1 x 25ml   |
| Stop Solution IV         | 1 x 5ml   | 1 x 25ml   |
| WST Reagent I            | 1 vial    | 5 vials    |

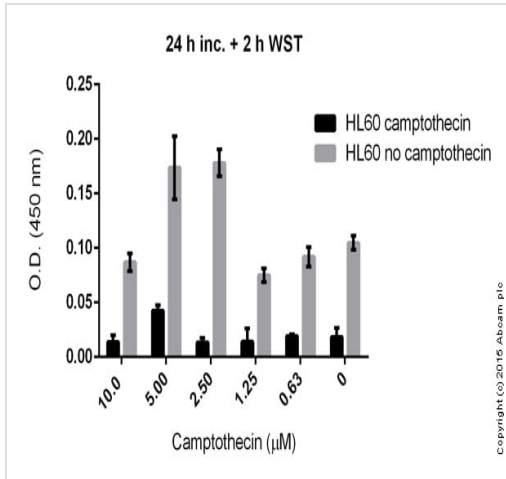
### 相关性

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

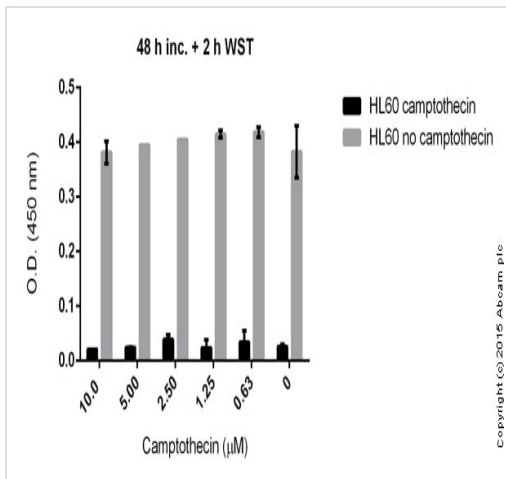


Mitochondrial dehydrogenases checked in various cell concentrations at different incubation times (2h WST for each condition)



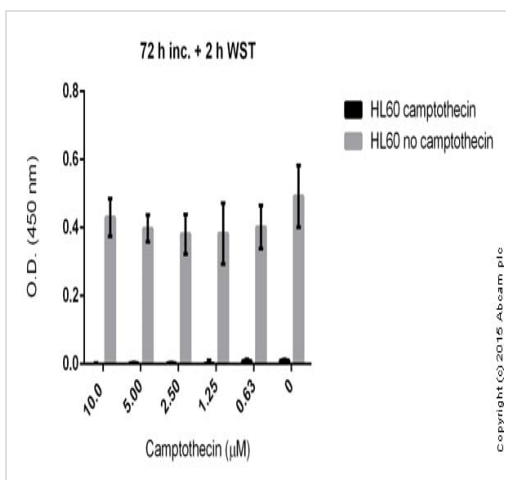
HL60 cell proliferation measured after 24h incubation (37°C/5%CO<sub>2</sub>) and 2h WST at various camptothecin concentrations

Cellular activity - ab65475 Quick Cell Proliferation Assay Kit II



HL60 cell proliferation measured after 48h incubation (37°C/5%CO<sub>2</sub>) and 2h WST at various camptothecin concentrations

Cellular activity - ab65475 Quick Cell Proliferation Assay Kit II



HL60 cell proliferation measured after 72h incubation (37°C/5%CO<sub>2</sub>) and 2h WST at various camptothecin concentrations

Cellular activity - ab65475 Quick Cell Proliferation Assay Kit II

**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