

Product datasheet

WST-1 Assay Kit (Cell Proliferation) ab65475

5 References 4 Images

Overview

Product name	WST-1 Assay Kit (Cell Proliferation)
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells
Assay type	Quantitative
Assay time	4h 00m
Product overview	WST-1 Assay Kit ab65475 is an alternative version of ab65473 , our most popular WST-1 Assay Kit (which is also supplied as a ready-to-use reagent ab155902). This kit uses a different WST-1 analog, and also contains a stop solution.

The WST-1 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-1 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-1 assay protocol is very simple:

- add the WST-1 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.

Notes This kit was previously called Quick Cell Proliferation Assay Kit II, and WST-1 Cell Proliferation Assay Kit.

The WST-1 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.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Microplate reader

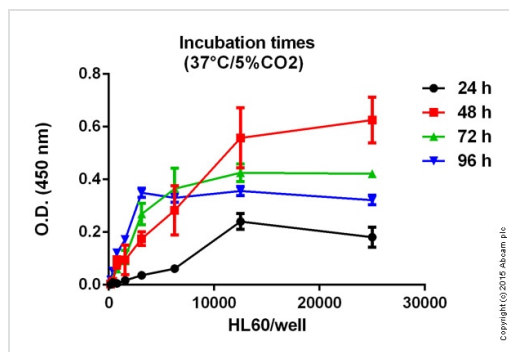
Properties

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

Components	500 tests	2500 tests
Electro Coupling Solution (ECS)	1 x 5ml	1 x 25ml
Stop Solution	1 x 5ml	1 x 25ml
WST Reagent (lyophilized)	1 vial	1 vial

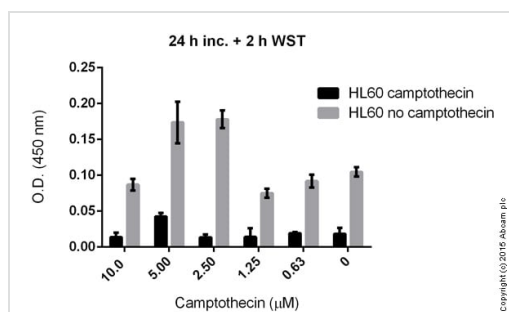
Relevance Cell proliferation is the multiplication or reproduction of cells, as a result of cell growth and cell division, resulting in the expansion of a cell population.

Images



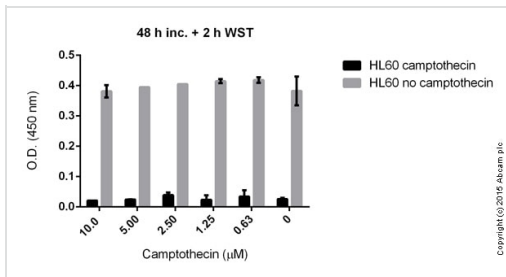
Mitochondrial dehydrogenases checked in various cell concentrations at different incubation times (2h WST for each condition)

Cellular activity - ab65475 Quick Cell Proliferation Assay Kit II



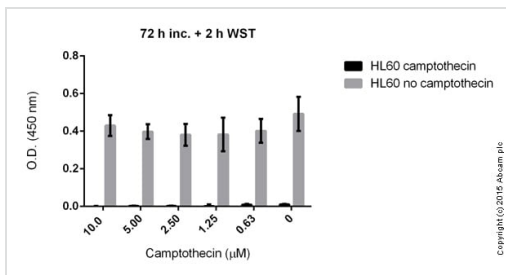
HL60 cell proliferation measured after 24h incubation (37°C/5%CO2) and 2h WST at various camptothecin concentrations

Cellular activity - ab65475 Quick Cell Proliferation Assay Kit II



HL60 cell proliferation measured after 48h incubation (37°C/5%CO₂) 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%CO₂) 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.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