

Product datasheet

WST-1 Assay Reagent - Cell Proliferation (ready to use) ab155902

★★★★★ [3 Abreviews](#) [35 References](#)

Overview

Product name	WST-1 Assay Reagent - Cell Proliferation (ready to use)
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells
Assay time	0h 30m
Product overview	<p>WST-1 Assay Reagent ab155902 provides a simple, accurate and ready-to-use assay to measure cell proliferation, cell viability and cytotoxicity in mammalian cells.</p> <p>The WST-1 assay protocol is based on the cleavage of the tetrazolium salt WST-1 to formazan by cellular mitochondrial dehydrogenases. The larger the number of viable cells, the higher the activity of the mitochondrial dehydrogenases, and in turn the greater the amount of formazan dye formed.</p> <p>The WST-1 assay protocol is very simple:</p> <ul style="list-style-type: none"> - 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 <p>We also offer complete WST-1 assay kits: WST-1 assay kit ab65473 which is the full kit version of this reagent, and WST-1 assay kit ab65475 which uses an alternative WST-1 analog.</p>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K304 Ready-to-use Cell Proliferation Colorimetric Reagent, WST-1. K304-2500 is the same size as the 2500 test size of ab155902.</p> <p>WST-1 assays are non-radioactive, rapid and more sensitive than MTT, XTT, or MTS-based assays. The entire assay can be performed in the same microtiter plate and does not require extra steps like washing, harvesting and cell solubilization.</p> <p>Review our cell health assays guide to learn more about our other cell viability, cytotoxicity and cell proliferation assay kits.</p>
Platform	Microplate reader

Properties

Storage instructions

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

Components	2500 tests
WST-1 Reagent I	1 x 2500 tests

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.

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