

WST-1 Assay Kit (Cell Proliferation) ab65473

39 References

製品の概要

製品名	WST-1 Assay Kit (Cell Proliferation)
検出方法	Colorimetric
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Quantitative
全工程の試験時間	4h 00m
製品の概要	WST-1 Cell Proliferation Assay Kit (ab65473) 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. It is also available in a ready-to-use reagent format: ab155902 .

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 is just add-and-read, requiring no washing, no harvesting, and no 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.

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

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K301 Quick Cell Proliferation Colorimetric Assay Kit. K301-500 is the same size as the 500 test size of ab65473.</p> <p>Review our cell health assays guide to learn more about our other cell viability, cytotoxicity and cell proliferation assay kits.</p>
------	--

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

保存方法

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

内容	500 tests	2500 tests
Electrocoupling Solution	1 x 5ml	1 x 25ml
WST Reagent IV	1 vial	1 vial

関連性

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.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors