

Product datasheet

Cell Counting Kit 8 (WST-8 / CCK8) ab228554

10 References 2 Images

Overview

Product name	Cell Counting Kit 8 (WST-8 / CCK8)
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells
Product overview	<p>Cell Counting Kit 8 (WST-8 / CCK8) (ab228554) provides a convenient and robust way of performing a cell viability assay. The kit uses a water-soluble tetrazolium salt to quantify the number of live cells by producing an orange formazan dye upon bio-reduction in the presence of an electron carrier.</p> <p>WST-8 / CCK8 solution is added directly to the test cells with no pre-mixing of components required.</p> <p>WST-8 / CCK8 tetrazolium salt is reduced by cellular dehydrogenases to an orange formazan product that is soluble in tissue culture medium. The amount of formazan produced is directly proportional to the number of living cells and is measured by absorbance at 460 nm.</p> <p>The excellent stability and little cytotoxicity of the WST-8 / CCK8 solution make the kit useful for assays that require long incubation (such as 24 to 48 hours).</p> <p>The detection sensitivity is higher than with other tetrazolium salt-based assays such as MTT, XTT or MTS etc.</p>
Notes	Review the cell health assay guide to learn about other kits to perform a cell viability assay , cytotoxicity assay or cell proliferation assay .
Platform	Microplate reader

Properties

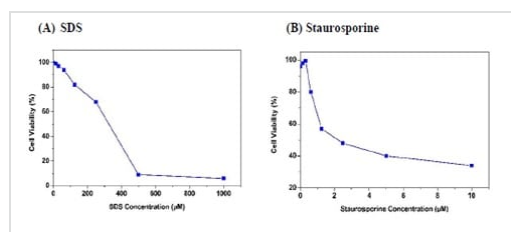
Storage instructions Store at +4°C. Please refer to protocols.

Components	1000 tests
WST-8 Solution	1 x 10ml

Relevance

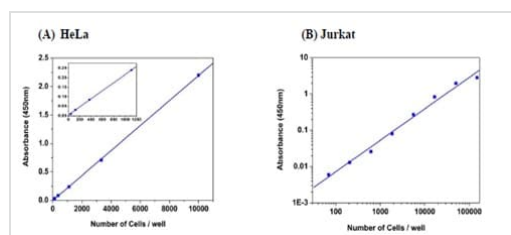
Cell proliferation is the multiplication or reproduction of cells, as a result of cell growth and cell division, resulting in the expansion of a cell population.

Images



Cytotoxicity tests of HeLa cells.

Cytotoxicity tests of HeLa cells in response to (A) SDS and (B) Staurosporine treatment were measured with The cell Counting Kit 8 (WST-8) (ab228554). HeLa cells at 10,000 cells/well/100 µL were seeded overnight in a black wall/clear bottom 96-well plate. Cells were treated with serially diluted SDS for 2 hours or Staurosporine for 4 hours. The absorbance was measured at 460 nm using a plate reader.



Cell Counting Assay.

Cell number was determined with the cell Counting Kit 8 (WST-8) (ab228554). (A) HeLa cells at 0 to 10,000 cells/well/100 µL, and (B) Jurkat cells at 0 to 100,000 cells/well/100 µL were added in a clear bottom 96-well plate. The absorbance was measured at 460 nm using a plate reader.

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