

CELL COUNTING KIT-8



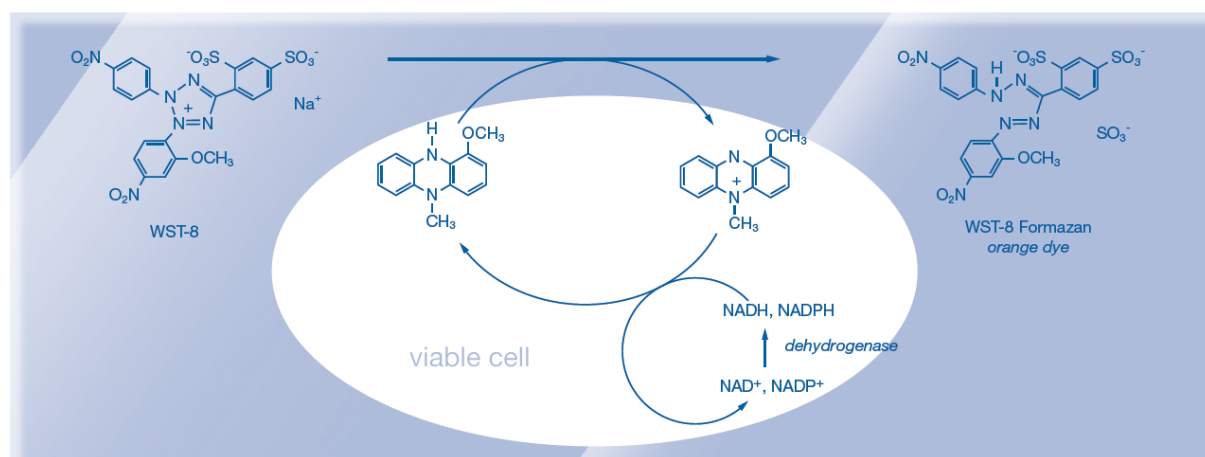
[Code#: CK04-11] 1,000 tests
[Code#: CK04-13] 3,000 tests
[Code#: CK04-20] 10,000 tests

PRODUCT DESCRIPTION

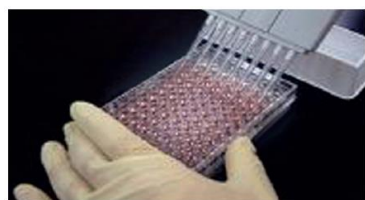
Cell Counting Kit-8 (CCK-8) is a sensitive colorimetric reagent used for the determination of cell viability in cell proliferation and cytotoxicity assays. Dojindo's highly water-soluble tetrazolium salt, Wst-8, is reduced by dehydrogenase activities in cells to give an orange-colored formazan dye, it is soluble in the tissue culture media. The detection sensitivity of CCK-8 is higher than that of other tetrazolium salts such as MTT, XTT or MTS.

PRINCIPLE

The amount of the WST-8 Formazan dyes generated by the activities of dehydrogenases in cells, are directly proportional to the number of living cells.



PROCEDURE



Add CCK-8 solution

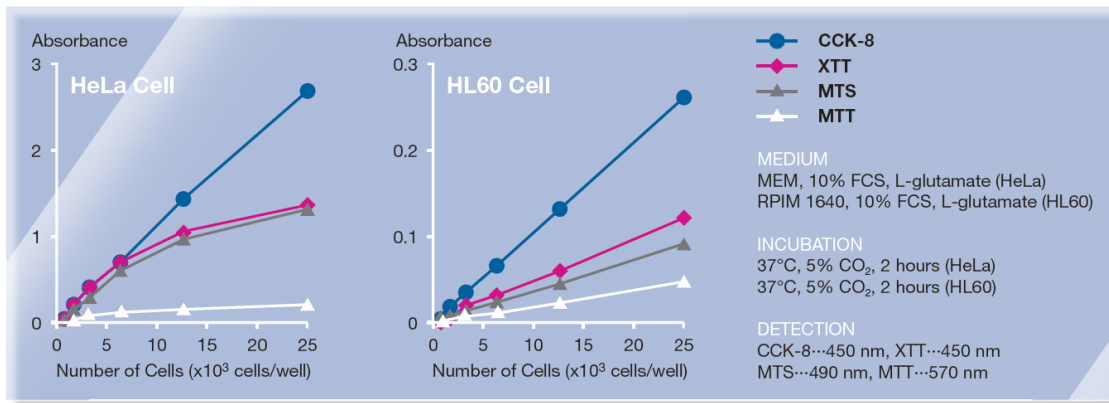


Incubate for 1-4 hours



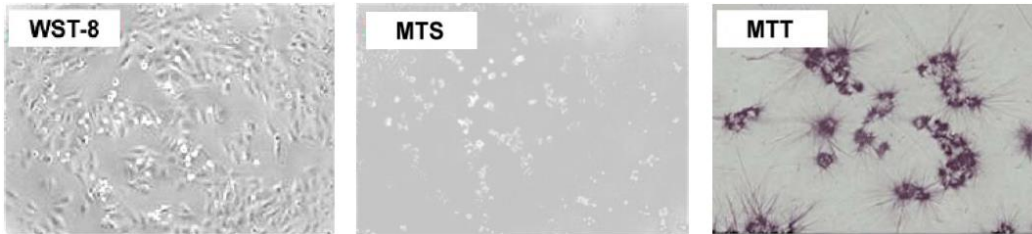
Measure O.D. at 450 nm

SENSITIVITY



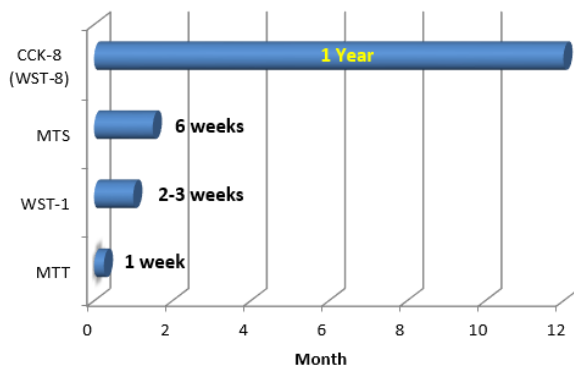
NO TOXICITY TO CELL

As seen in the photos, even after 24 hours of incubation with CCK-8, cells continue to proliferate. Other cell viability assay kits are toxic to cells, and most of the cells are killed during the first 3 hours of incubation.



LONGER STABILITY OF REAGENT

With stability up to a year at 5°C, you do not have to worry about preparing your cell proliferation reagents right before your experiment. Cell Counting Kit-8 is even stable for 6 months at room temperature. There is no one out there who can survive the heat like us.



DOJINDO STATE OF THE ART
LIFE SCIENCE TECHNOLOGIES

EUROPEAN HEADQUARTERS
DOJINDO EU GMBH

Leopoldstr. 254, 80807 Munich, Germany
Phone +49 89 3540-4806
Fax +49 89 3540-4806
email info@dojindo.eu.com

www.dojindo.eu.com

DISTRIBUTED BY