

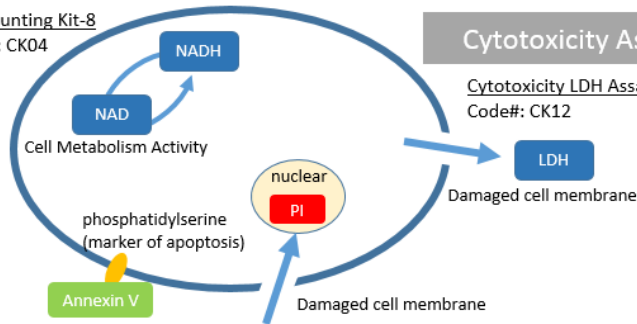
Dojindo Cell Analysis Products

Cell Counting Kit-8 [Code#: CK04]
 Cytotoxicity LDH Assay Kit-WST [Code#: CK12]
 Annexin V, FITC Apoptosis Detection Kit [Code#: AD10]



Cell Proliferatin / Viability Assay

Cell Counting Kit-8
 Code#: CK04



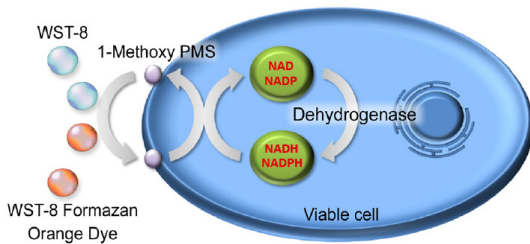
Cytotoxicity Assay

Cytotoxicity LDH Assay Kit-WST
 Code#: CK12

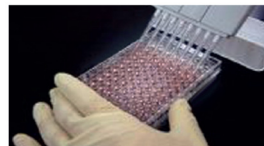
Apoptosis Detection

Annexin V, FITC Apoptosis Detection Kit
 Code#: AD10

CELL COUNTING KIT-8



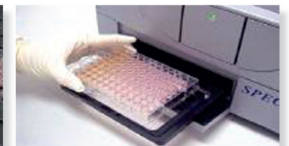
Cell Counting Kit-8 (CCK-8) allows sensitive colorimetric assays for the determination of cell viability in cell proliferation. The detection sensitivity of CCK-8 is higher than the other tetrazolium salts such as MTT, XTT, or MTS. Unlike MTT, XTT, MTS and WST-1, CCK-8 is stable up to one year in a refrigerator (0-5°C).



Add CCK-8 solution



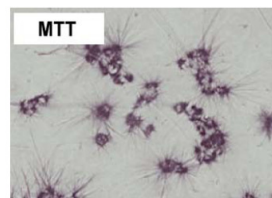
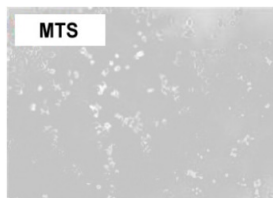
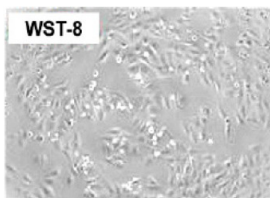
Incubate for 1-4 hours



Measure O.D. at 450 nm

LOW TOXICITY TO CELL

As seen in the photos, even after 24 hours of incubation with CCK-8 (WST-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.



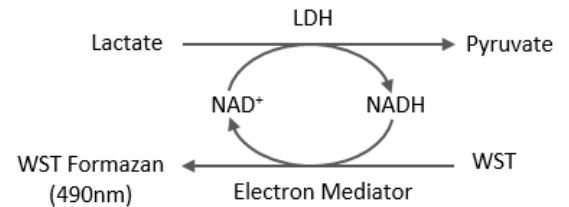
KIT CONTENTS

Product code	CK04-11 (1000 tests)	CK04-13 (3000 tests)
Cell Counting Kit-8	2 x 5 ml	6 x 5 ml



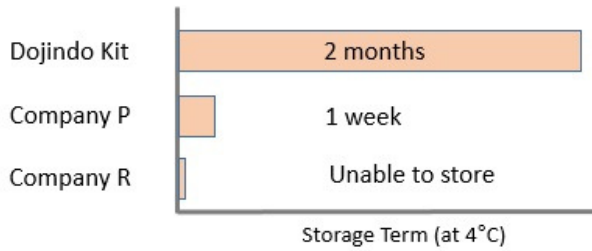
Cytotoxicity LDH Assay Kit-WST

Cytotoxicity LDH Assay Kit-WST is a kit for determination of cytotoxicity by measuring a lactate dehydrogenase (LDH) activity released from damaged cells. LDH is a stable cytoplasmic enzyme presented in all types of cells and released into the cell culture medium through damaged plasma membrane. Cytotoxicity LDH Assay kit-WST can be used to measure the released LDH according to the following scheme.



STABILITY DATA

Unlike competitive products, the reconstituted working solution can be stored for 2 months at refrigerator.



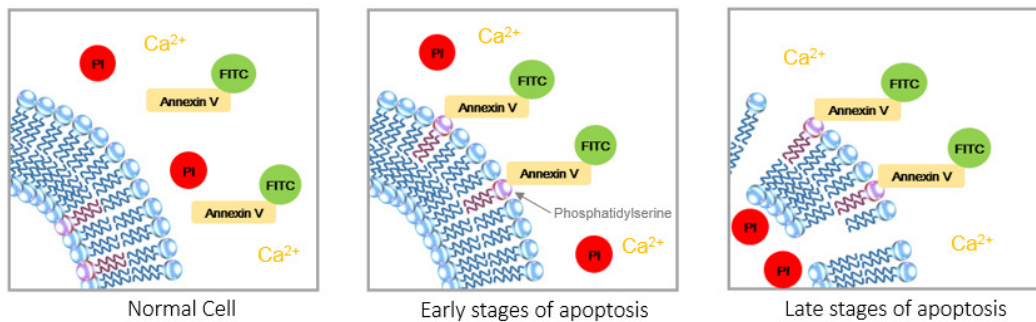
KIT CONTENTS

Product code	CK12 (100 tests)
Dye Mixture	1 tube
Assay Buffer	11 ml x 1 tube
Lysis Solution	1.1 ml x 1 tube
Stop Solution	5.5 ml x 1 tube



Annexin V, FITC Apoptosis Detection Kit

Annexin V stained cells are used to indicate cell membrane changes that occur in the early stage of apoptosis. By using fluorescent-labeled Annexin V, the apoptotic cells can be identified by flow cytometry or fluorescence microscopy.



KIT CONTENTS

Product code	AD10-10 (50 tests)	AD10-12 (2 x 50 tests)
Annexin V, FITC Conjugate	50 assay x 1 bottle	50 assay x 2 bottles
PI Solution	50 assay x 1 bottle	50 assay x 2 bottles
Annexin V Binding Buffer	50 assay x 1 bottle	50 assay x 2 bottles



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