

Cell Counting Kit-8

Simple, fast and accurate assay to measure cell proliferation and cytotoxicity

The Cell Counting Kit-8 is a colorimetric assay kit used to measure cell proliferation and cytotoxicity.

It is a ready-to-use solution that does not require radioisotopes and correlates with the [³H]-thymidine incorporation assay. It can be added directly to the cell media for fast, high-throughput screening without a solubilization process obtaining highly reproducible and accurate results. CCK-8 has shown to achieve higher sensitivity and stability than MTT, MTS or WST-1.

Citations: 108

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-039-0100	100 tests
ALX-850-039-KI01	500 tests
ALX-850-039-KI02	5x500 tests

Manuals, SDS & CofA

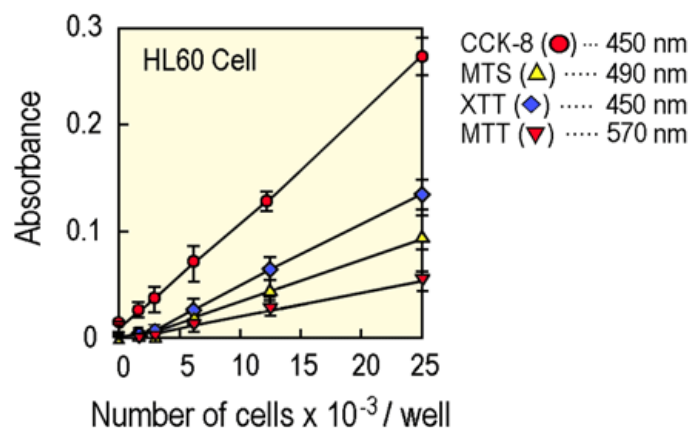
[View Online »](#)

- One-step, ready-to-use solution with no radioisotopes
- High sensitivity that correlates with the [³H]-thymidine incorporation assay
- High-throughput screening without a solubilization step
- More sensitive and stable than MTT, MTS or WST-1

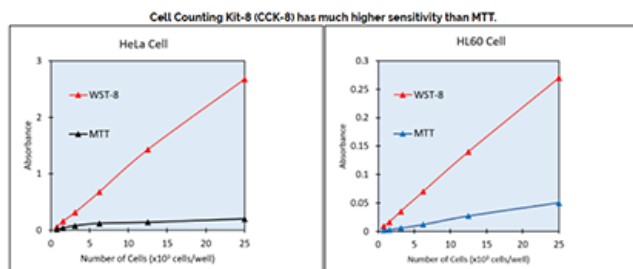
Simple, Fast, and Accurate Method for Measuring Cell Proliferation

Reagent	MTT	MTS WST-1	CCK-8
Preparation	Thaw Reagent Dissolve MTT	Thaw Reagent Dissolve MTS/WST-1	Thaw Reagent
Procedure	Add Reagent Measure Abs.	Add Reagent Measure Abs.	Add Reagent Measure Abs.
Handling Time	40 minutes	30 minutes	15 minutes

Cell Counting kit-8 performance

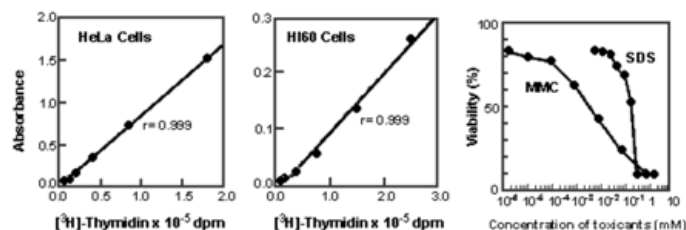


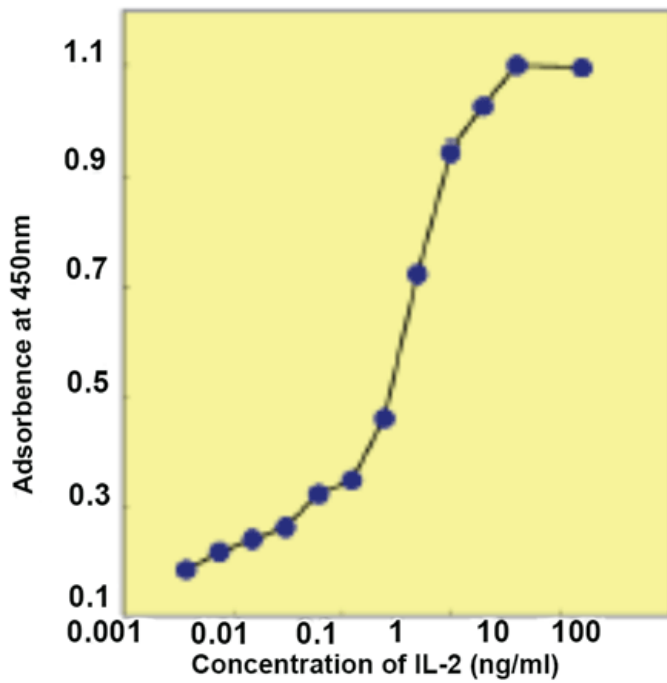
CTLL-2 cells were incubated with various concentrations of IL-2 for 72 hours. CCK-8 solution was added to each well and the absorbance at 450nm was measured. IL-2 exposure resulted in an increase absorbance which correlates to an increase in cell proliferation. CCK-8 shows greatest sensitivity.



Medium: MEM, 10% FCS, L-glutamate (HeLa)
 RPMI 1640, 10% FCS, L-glutamate (HL60)
 Incubation: 37°C, 5%CO₂, 2 hours (HeLa)
 37°C, 5%CO₂, 2 hours (HL60)
 Detection: CCK-8 450 nm, MTT – 570 nm

CCK-8 sensitivity for HeLa and HL60 cells are more sensitive than MTT





CTLL-2 cells were incubated with various concentrations of IL-2 for 72 hours. CCK-8 solution was added to each well and the absorbance at 450nm was measured. IL-2 exposure resulted in an increase absorbance which correlates to an increase in cell proliferation.

Handling & Storage

Handling Protect from light. Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CCKI-8

Application Colorimetric detection

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents WST-8 solution, 1-methoxy-PMS

Quantity
100 tests: 1 mL
500 tests: 5 mL
2500 tests: 5 x 5 mL

Technical Info / Product Notes **Principle:** Employs the tetrazolium salt WST-8 (2-(2-methoxy-4-nitrophenyl)-3-(4-nitrophenyl)-5-(2,4-disulfophenyl)-2*H*-tetrazolium . monosodium salt), that produces a highly water soluble formazan dye upon biochemical reduction in the presence of an electron carrier, 1-methoxy-PMS. The amount of the yellow colored formazan dye generated by dehydrogenases in cells is directly proportional to the number of viable cells in a culture medium.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)