

Cell Transformation Assay Kit (Fluorometric)

08/16

(Catalog # K922-100; 100 assays; Store at -20°C)

I. Introduction:

Transformed cells can proliferate without attaching to surface. Anchorage-independent cell growth is the hallmark of cell transformation. The Soft-Agar Assay is a traditional method for screening cell transformation *in vitro*. However, this method is lengthy (3-4 weeks incubation), laborious (counting colonies) and inconsistent (due to subjective counting). BioVision's Cell Transformation Assay is faster, stable, and more sensitive than the traditional soft-agar assay. The kit uses a quantitative dye that binds to nucleic acid and generates green fluorescence. This one-step method is non-radioactive and simple (just add-and-read, and does not require tedious labor such as counting colonies). The assay is high-throughput adaptable and has wide linear range from 50-60,000 cells. The entire assay can be finished within 7-8 days.

II. Applications:

- Measurement of cell transformation in response to carcinogens, oncogenes, etc.
- Assessments of chemicals that induce or inhibit cell transformation

III. Sample Type:

- Adherent or suspension cells

IV. Kit Contents:

Components	K922-100	Cap Code	Part Number
Agarose Powder	240 mg	NM	K922-100-1
DMEM Solution (10X)	2 X 1.5 ml	Clear	K922-100-2
Staining Solution	1 ml	Brown	K922-100-3
Agarose Solubilization Solution	5 ml	NM	K922-100-4
Quantitative Dye (200X)	0.1 ml	Red	K922-100-5

V. User Supplied Reagents and Equipment:

- 96-well clear tissue culture plate and 96-well white plate
- Sterile dH₂O, PBS, and FBS
- Microscope
- Multi-well spectrophotometer (ELISA reader)

VI. Storage Conditions and Reagent Preparation:

Store kit at -20°C, protected from light. Briefly centrifuge small vials prior to opening. Read entire protocol before performing the assay. *Prepare reagents and perform assays under sterile conditions* (i.e. tissue culture hood/biosafety cabinet).

- **Agarose Powder:** To make a 1.2% agarose solution, add 20 ml of sterile dH₂O into the Agarose Powder bottle. Open the bottle cap, slightly, and heat the bottle on a heat block until the Agarose Powder is completely dissolved (~ 100 °C; 30-40 min is recommended). Gently shake the bottle to solubilize the agarose. Transfer the bottle to a 37°C water bath and keep it for 30 min. to equilibrate to temperature. *Unused 1.2% agarose solution can be stored at 4°C under sterile conditions.*

Note: Keep the Agarose solution in a 37 °C water bath throughout cell-seeding process to prevent solidification of the agarose solution.

- **DMEM Solution (10X):** Dilute 10X DMEM in sterile dH₂O to 1X DMEM containing 10% FBS (1X DMEM/10% FBS). For example, dilute 100 µl of DMEM Solution (10X) into 900 µl dH₂O with 100 µl of FBS. Make as much as needed. Store at 4°C. Before using, warm to 37°C in a water bath.
- **Quantitative Dye (200X):** To make 1X Quantitative Dye Solution, dilute 200X Quantitative Dye with 1X PBS. For example: add 10 µl of 200X Quantitative Dye into 1.9 ml 1X PBS, mix well. *Discard the unused 1X Quantitative Dye Solution and always prepare fresh dilution.*

VII. Cell Transformation Assay Protocol:

1. Sample Preparation:

- a. **Preparation of Base Agarose Layer:** Prepare 75 µl/well base agarose mix. For each well, mix:

1.2% agarose solution	37.5 µl
DMEM Solution (10X)	7.5 µl
100% FBS	7.5 µl
dH ₂ O	22.5 µl

Prepare enough Base Agarose mix for the number of experiments to be performed. Mix well. Add 75 µl of base agarose mix into desired wells in a 96-well clear bottom tissue culture plate. Keep the plate at 4°C for 15 min. to solidify the agarose.

Note: Prior to adding the top agarose layer with cells, warm the plate at room temperature by keeping in tissue culture hood for 10 min.

- b. **Preparation of Top Agarose Layer with Cells:** Prepare a stock solution of cells (1-5 X10⁶ cells/ml) in 1X DMEM/10% FBS medium. Calculate and adjust the desired cell concentration (see note a) based on the number of cells per well for the assay. Prepare 75 µl/well top agarose-cell mix as follows:

1.2% agarose solution	25.0 μ l
DMEM Solution (10X)	5.5 μ l
FBS	5.5 μ l
Cells in 1X DMEM/10% FBS	20 μ l
dH ₂ O	19 μ l

Make as much as needed. Mix by pipetting. Add 75 μ l of agarose-cell mix into each well of the 96-well clear bottom tissue culture plate already containing the solidified base agarose layer. Keep the plate at 4°C for 10 min. to solidify the top agarose-cell mix. Bring the plate to room temperature by keeping it in the tissue culture hood for 10 min. Add total of 100 μ l of 1X DMEM/10% FBS medium with or without test compound into each well and incubate at 37°C for 6-8 days.

Notes:

- *Assay has linear range from 50 to 60,000 cells, depending on the cell type used in the experiment. Adjust the cell numbers to avoid over-seeding.
 - Prepare parallel well(s) as blank control (no cells) with same amount of culture medium and reagents for the reagent background reading.
 - During the process of plating the Base Agarose Layer and Top Agarose Layer with cells, keep 1.2% agarose solution, DMEM solution (10X), sterile dH₂O, and FBS in a 37°C water bath to equilibrate the temperature and to prevent solidification of agarose in case of 1.2% agarose layers.
 - Multi-channel pipette can be used for plating Base Agarose Layer. Add agarose-cell mix carefully to avoid bubbles in both base and top agarose layers.
 - Colony Visualization (Optional): Add 10 μ l Staining Solution into each well and incubate for 60 min. at 37°C incubator with 5% CO₂. Colonies formed by transformed cells can be visualized and imaged under the microscope.
- Cell-Dose Curve:** On day 0: Prepare a cell-dose curve by using the stock made in step 1.b (1-5 X10⁶ cells/ml in 1X DMEM/10% FBS medium). Arrange eight serial dilutions (2-fold) in separate 1.5 ml centrifuge tubes with 1X DMEM/10% FBS medium (150 μ l). Add 50 μ l of Agarose Solubilization Solution to each tube, mix and incubate cells for 15 min at RT. Transfer 20 μ l of the each mixture into 96-well white plate. Add 80 μ l 1X Quantitative Dye Solution to each well, protect from light and shake for 10 min on a shaker. Measure the fluorescence (RFU) of the blank and diluted cell solutions using a microtiter plate reader at Ex/Em= 480/530 nm. Subtract all readings from blank, and plot the Cell-Dose Curve.
 - Measurement:** On day 6-8 (at the end of the desired incubation time, step 1.b): Carefully remove the medium above the top agarose layer by pipetting. Add 50 μ l of Agarose Solubilization Solution into each well and incubate at 37 °C incubator for 1 hr. to solubilize the agarose. Transfer 20 μ l of Solubilized Agarose-cell mix into a 96-well white plate and add 80 μ l of 1X Quantitative Dye Solution. Protect from light and gently shake for 15 min. at room temperature. Measure fluorescence (Ex/Em= 480/530 nm).
 - Calculation:** Subtract 0 Standard reading from all readings. Plot the Cell-Dose Curve (number of cells vs RFU_{480/530}). Apply the sample readings (Δ RFU) to the Cell-Dose Curve (Cell number/well: 1/10th of the original stock solutions, step 2) to get the number of transformed cells (B). The total number of transformed cells can be calculated using the equation RFU_{480/530} = slope*cells +b.

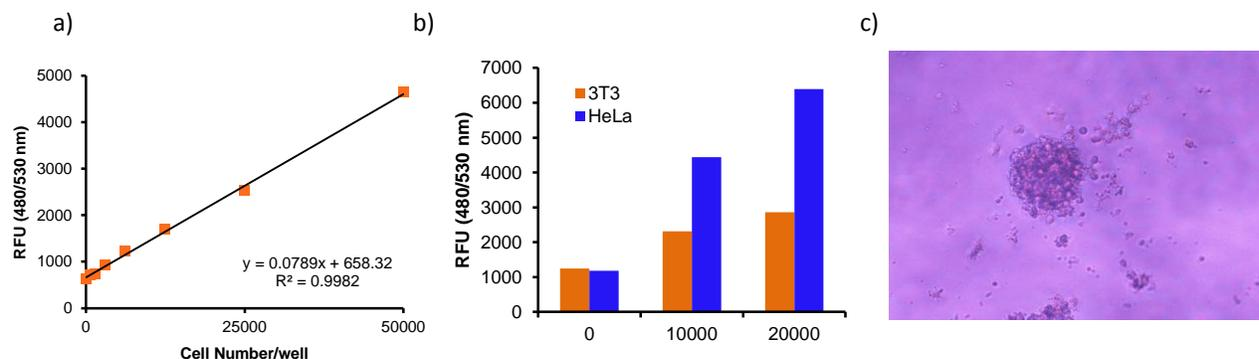


Figure: a) HeLa Cell Dose Curve; b) Anchorage-independent growth of 3T3 and HeLa cells. 3T3 and HeLa cells were serially diluted and seeded in agarose gel. Cells were solubilized and detected by the Quantitative Dye. **c) Image of a HeLa Colony.** HeLa cells were cultured for 7 days according to the kit protocol.

VIII. RELATED PRODUCTS:

Quick Cell Proliferation Colorimetric Assay Kit (K301)
 BrdU Cell Proliferation Assay Kit (K306)
 MTS Cell Proliferation Colorimetric Assay Kit (K300)
 Ready-to-use Cell Proliferation Reagent, WST-1 (K304)
 EZCell™ Cell Cycle Analysis Kit (K920)
 Annexin V Apoptosis Kits (K101-K104)
 ApoSENSOR™ ATP Cell Viability Bioluminescence Assay Kit (K254)
 EZViable™ Calcein AM Cell Viability Assay Kit (Fluorometric) (K305)

Quick Cell Proliferation Colorimetric Assay Kit Plus (K302)
 EZCell™ Cell Migration/Chemotaxis Assay Kits (K906-K912)
 Live-Dead Cell Staining Kit (K501)
 VisionBlue™ Quick Cell Viability Fluorometric Assay Kit (K303)
 Senescence Detection Kit (K320)
 EZSolution™ 7-Aminoactinomycin D (2727)
 StayBrite™ Highly Stable ATP Bioluminescence Assay Kit(K791)
 ADP Colorimetric Assay Kit II (K356)

FOR RESEARCH USE ONLY!



FOR RESEARCH USE ONLY! Not to be used on humans.