

Product datasheet

Cell Transformation Assay Kit (Fluorometric) ab235699

1 Image

Overview

Product name	Cell Transformation Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Species reactivity	

Predicted to work with: Mammals 

Product overview

Transformed cells can proliferate without attaching to a 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). The Cell Transformation Assay Kit (Fluorimetric) (ab235699) 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.

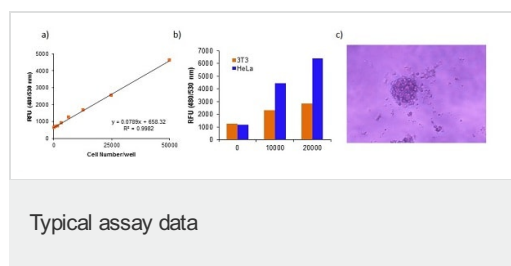
Platform	Microplate reader
-----------------	-------------------

Properties

Storage instructions	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	100 tests
Agarose Powder	1 x 240mg
Agarose Solubilization Solution	1 x 5ml
DMEM Solution (10X)	2 x 1.5ml
Quantitative Dye (200X)	1 x 0.1ml
Staining Solution	1 x 1ml

Images



(a) HeLa (Human epithelial cell line from cervix adenocarcinoma) cell-dose curve; (b) Equal numbers of NIH/3T3 (Mouse embryo fibroblast cell line) and HeLa cells were serially diluted and seeded in agarose gel. Cells were solubilized and detected with Quantitative Dye; (c) Image of HeLa cell colonies. HeLa cells were cultured for 7 days according to the kit protocol.

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