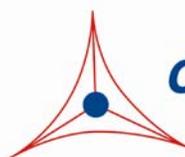

Product Manual

CytoSelect™ 96- Well Cell Transformation Assay (Soft Agar Colony Formation)

Catalog Number

CBA- 130	96 assays
CBA- 130- 5	5 x 96 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

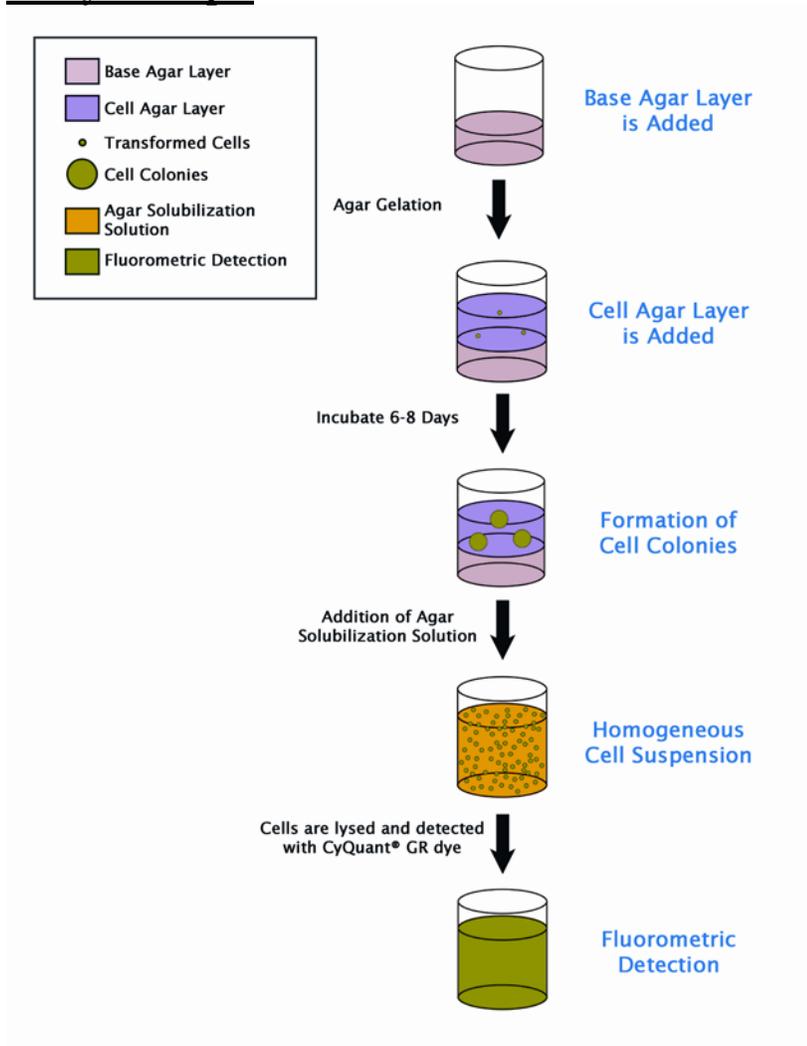
Neoplastic transformation occurs via a series of genetic and epigenetic alterations that yield a cell population that is capable of proliferating independently of both external and internal signals that normally restrain growth. For example, transformed cells show reduced requirements for extracellular growth promoting factors, are not restricted by cell-cell contact, and are often immortal. Anchorage-independent growth is one of the hallmarks of transformation, which is considered the most accurate and stringent in vitro assay for detecting malignant transformation of cells.

Traditionally, the soft agar colony formation assay is a common method to monitor anchorage-independent growth, which measures proliferation in a semisolid culture media after 3-4 weeks by manual counting of colonies. Standard soft agar assays are usually performed in 100-mm or 60 mm dishes, where cells are allowed to grow inside a semisolid culture media for 3-4 weeks before sizable colonies appear. This method is quite cumbersome, time-consuming, and difficult when testing a large number of samples. Additionally, the manual counting of colonies is highly subjective; with varying colony sizes, it's difficult to determine meaningful results.

Cell Biolabs CytoSelect™ 96-well Cell Transformation Assay does **not** involve subjective manual counting of colonies or require a 3-4 week incubation period. Instead cells are incubated only 6-8 days in a semisolid agar media before being solubilized, lysed and detected by the patented CyQuant® GR Dye in a fluorescence plate reader (see Assay Principle below). This format provides a quantitative, high-throughput method to accurately measure cell transformation. Additionally, the short incubation time (6-8 days) makes it possible to assay cells transiently transfected with oncogenes or siRNA.

The CytoSelect™ 96-well Cell Transformation Kit provides a robust system for screening oncogenes and cell transformation inhibitors. Each kit provides sufficient quantities to perform 96 tests in a microtiter plate.

Assay Principle



Related Products

1. CBA-106: CytoSelect™ 96-Well Cell Migration Assay (8µm, Fluorometric)
2. CBA-106-C: CytoSelect™ 96-Well Cell Migration and Invasion Assay (8µm, Fluorometric)
3. CBA-110: CytoSelect™ 24-Well Cell Invasion Assay (Basement Membrane, Colorimetric)
4. CBA-112: CytoSelect™ 96-Well Cell Invasion Assay (Basement Membrane, Fluorometric)
5. CBA-135: CytoSelect™ 96-Well Cell Transformation Assay (Cell Recovery, Colorimetric)
6. CBA-140: CytoSelect™ 96-Well Cell Transformation Assay (Cell Recovery, Fluorometric)
7. CBA-145: CytoSelect™ 384-Well Cell Transformation Assay
8. CBA-150: CytoSelect™ In Vitro Tumor Sensitivity Assay
9. CBA-155: CytoSelect™ Clonogenic Tumor Cell Isolation Kit
10. CBA-320: CytoSelect™ 96-Well Hematopoietic Colony Forming Cell Assay

Kit Components

1. CytoSelect™ Agar Powder (Part No. 113001): One 1.2 g bottle
2. 5X DMEM Solution (Part No. 113002): Two 1.5 mL sterile tubes
3. Agar Solubilization Solution (Part No. 113003): One 6 mL amber bottle
4. 8X Lysis Buffer (Part No. 113004): One 3 mL bottle
5. CyQuant GR Dye (Part No. 10103): One 25 µL tube

Materials Not Supplied

1. Cells and Culture Medium
2. 1X PBS
3. 37°C Incubator, 5% CO₂ Atmosphere
4. Light Microscope
5. 96-well Fluorometer
6. Microwave or Heating Block
7. Water bath
8. (Optional) Positive Control cells such as NIH 3T3 (Ras G12V)

Storage

Store all components at 4°C.

Preparation of Reagents

- 1.2% Agar Solution: Place 1.2 g of Agar Powder in a sterile bottle, add 100 mL of sterile cell culture grade water. Microwave or boil until agar is completely dissolved.
- 2X DMEM/20% FBS Medium: In a sterile tube, dilute the provided 5X DMEM in sterile cell culture grade water to 2X containing 20% FBS. For example, to prepare a 5 mL solution, add 2 mL of 5X DMEM, 1 mL of FBS and 2 mL of sterile cell culture grade water. Sterile filter the 2X media to 0.2 µm.

Note: You may substitute your own medium in place of the DMEM we provide, but ensure that it is at a 2X concentration.

- CyQuant Working Solution: Immediately before use, prepare sufficient amount of the CyQuant Working Solution by diluting the CyQuant GR Dye 1:400 with 1X PBS. For example, add 10 µL to 4 mL of 1X PBS. Use the solution immediately; do not store the CyQuant Working Solution.

Assay Protocol (must be under sterile conditions)

I. Preparation of Base Agar Layer

1. Melt 1.2% Agar Solution in a microwave and cool to 37°C in a water bath.
2. Warm 2X DMEM/20% FBS medium to 37°C in a water bath. Allow at least 30 minutes for the temperature to equilibrate.
3. Mix equal volumes of 1.2% Agar Solution and 2X DMEM/20% FBS medium in a sterile, pre-warmed tube by inverting several times. Immediately transfer 50 µL of the mixture to each well of a 96-well sterile flat-bottom microplate. Gently tap the plate a few times to allow the agar solution to evenly cover the wells.

Notes:

- *Work quickly with the agar solution to avoid gelation. Also, try to avoid adding air bubbles to the well.*
 - *To avoid fast and uneven evaporation that leads to aberrant results, we suggest not using the wells on the plate edge, or filling the edge wells with medium to reduce evaporation.*
4. Transfer the plate to 4°C for 30 minutes to allow the base agar layer to solidify.
 5. Prior to adding the Cell Agar Layer (Section II), allow the plate to warm up for 15 minutes at 37°C.

II. Preparation of Cell Agar Layer (samples should be assayed in triplicate)

1. Melt 1.2% Agar Solution in a microwave and cool to 37°C in a water bath.
2. Warm 2X DMEM/20% FBS medium to 37°C in a water bath. Allow at least 30 minutes for the temperature to equilibrate.
3. Harvest and resuspend cells in culture medium at $0.4 - 4 \times 10^5$ cells/mL, keep the cell suspension warm in a 37°C water bath.
4. Mix equal volumes of 1.2% Agar Solution, 2X DMEM/20% FBS media, and cell suspension (1:1:1) in a sterile, pre-warmed tube by inverting several times. Immediately transfer 75 µL of the mixture to each well of the 96-well flat-bottom microplate already containing the solidified base agar layer (25 µL of cell suspension containing 1000-10000 cells/well will be seeded).

Note: Work quickly with the agar solution to avoid gelation, but gently pipette as not to disrupt the base layer integrity. Also, try to avoid adding air bubbles to the well. Always include negative control wells that contain no cells in the cell agar layer.

5. Transfer the plate to 4°C for 15 minutes to allow the cell agar layer to solidify.

III. Quantitation of Anchorage-Independent Growth

1. Add 100 µL of culture medium containing cell growth activator(s) or inhibitor(s) to each well.
2. Incubate the cells for 6-8 days at 37°C and 5% CO₂. Examine the cell colony formation under a light microscope.

3. Remove culture medium by inverting the plate and blotting on paper towel. Gently tap several times.
4. Add 50 μL of Agar Solubilization Solution to each well of the 96-well plate. Incubate for 1 hr at 37°C.
5. Pipette each well 5-10 times to ensure complete agar solubilization.
6. Add 25 μL of 8X Lysis Buffer to each well. Pipette each well 5-10 times to ensure a homogeneous mixture.
7. Incubate the plate at room temperature for 15 minutes.
8. Transfer 10 μL of the mixture to a 96-well plate suitable for fluorescence measurement.
9. Add 90 μL of the CyQuant Working Solution to each well. Incubate 10 minutes at room temperature.
10. Read the plate in a 96-well fluorometer using a 485/520 nm filter set.

Cell Dose Curve (optional)

1. Harvest and resuspend cells in culture medium at $1 - 5 \times 10^6$ cells/mL.
2. Prepare a serial of 2-fold dilution with culture medium, including a medium blank.
3. Transfer 125 μL of each cell dilution to a microfuge tube. Add 50 μL of Agar Solubilization Solution and 25 μL of 8X Lysis Buffer to each tube. Vortex each tube to ensure a homogeneous mixture. Incubate the tubes at room temperature for 15 minutes.
4. Transfer 10 μL of the mixture to a 96-well plate suitable for fluorescence measurement.
5. Add 90 μL of the CyQuant Working Solution to each well. Incubate 10 minutes at room temperature.
6. Read the plate in a 96-well fluorometer using a 485/520 nm filter set.

Example of Results

The following figures demonstrate typical results with the CytoSelect™ 96-well Cell Transformation Assay Kit. Fluorescence measurement was performed on SpectraMax Gemini XS Fluorometer (Molecular Devices) with a 485/538 nm filter set and 530 nm cutoff. One should use the data below for reference only. This data should not be used to interpret actual results.

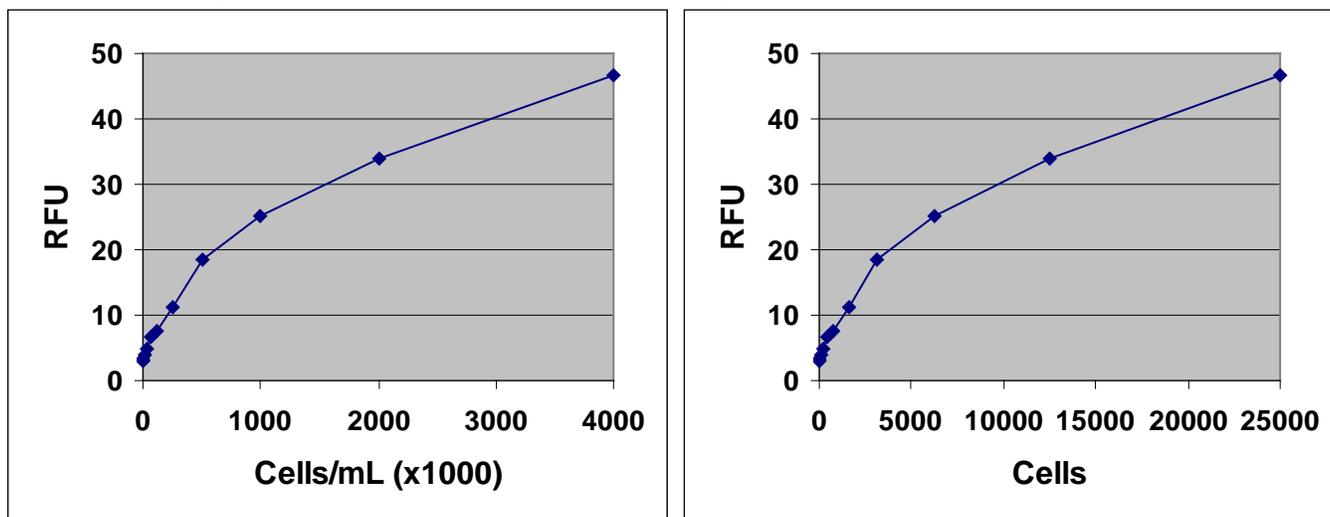


Figure 1. HeLa Cell Dose Curve. Cervical carcinoma HeLa cells were resuspended at 4×10^6 cells/mL and titrated 1:2 in culture medium, followed by addition of Agar Solubilization Solution, Lysis Buffer, and Cyquant® GR Dye detection (as described in the Cell Dose Section). Results were shown by cell concentration or by actual cell number in CyQuant Detection.

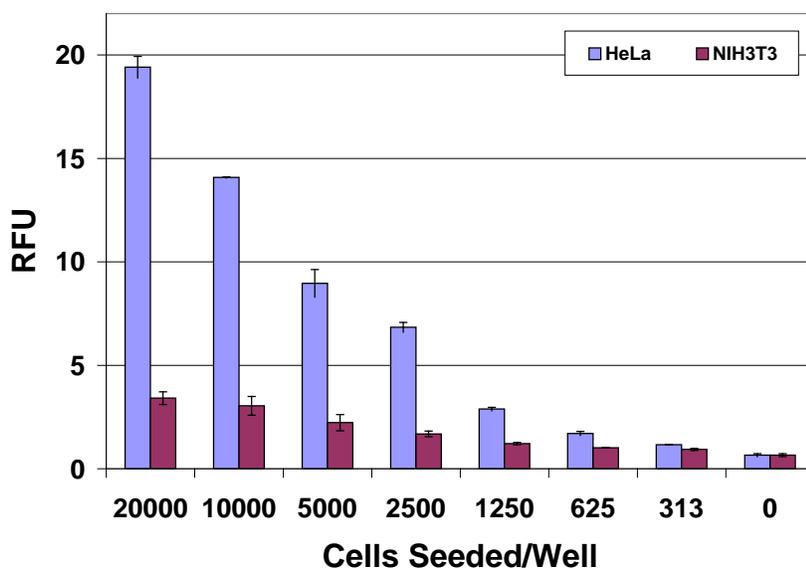


Figure 2. Anchorage-Independent Growth of HeLa Cells. HeLa cells were seeded at various concentrations and cultured for 6 days. HeLa cell transformation is determined according to the assay protocol.

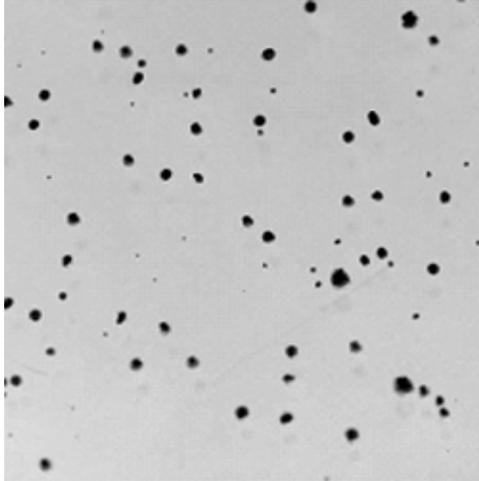


Figure 3. HeLa Colony Formation. HeLa cells were cultured for 14 days according to the assay protocol. Colonies were visualized by 0.1% p-iodonitro tetrazolium violet (INT) staining.

Calculation of Anchorage-Independent Growth

1. Compare RFU values with the Cell Dose Curve and extrapolate the cell concentration in soft agar.
2. Calculate the Total Transformed Cell Number/Well
Total Transformed Cells/Well = cells/mL in soft agar x 0.125 mL/well

For example: If you extrapolate your RFU value from your cell dose curve and determine you have 500,000 cells/mL in your soft agar sample.

Total Transformed Cells/Well = 500,000 cells/mL x 0.125 mL/well = 62,500 cells/well

References

1. Shin SI, Freedman VH, Risser R, and Pollack R. (1975) *Proc Natl Acad Sci U S A.* 72:4435-9.
2. Hahn WC, Counter CM, Lundberg AS, Beijersbergen RL, Brooks MW and Weinberg RA. (1999) *Nature* 400:464-8.

Recent Product Citations

1. Martinez-Cruzado, L. et al. (2017). Trabectedin and Camptothecin Synergistically Eliminate Cancer Stem Cells in Cell-of-Origin Sarcoma Models. *Neoplasia* **19(6)**:460-470.
2. Lee, B. et al. (2017). The long noncoding RNA SPRIGHTLY acts as an intranuclear organizing hub for pre-mRNA molecules. *Sci. Adv.* **3(5)**:e1602505.
3. Manzanares, M.A. et al. (2017). Transforming Growth Factors α and β Are Essential for Modeling Cholangiocarcinoma Desmoplasia and Progression in a Three-Dimensional Organotypic Culture Model. *Am J Pathol.* pii: S0002-9440(17)30236-5. doi: 10.1016/j.ajpath.2017.01.013.
4. Andrio, E. et al. (2017). Identification of cancer-associated missense mutations in hacc1 that impair cell growth control and Rac1 ubiquitylation. *Sci Rep.* **7**:44779. doi: 10.1038/srep44779.

5. Illuzzi, J. L., et al. (2017). Tumor-associated APE1 variant exhibits reduced complementation efficiency but does not promote cancer cell phenotypes. *Environ. Mol. Mutagen.* **58**(2):84-98. doi:10.1002/em.22074
6. Otaka Y, et al. (2017). STXBP4 drives tumor growth and is associated with poor prognosis through PDGF Receptor signaling in lung squamous cell carcinoma. *Clin Cancer Res.* 1815.2016. doi: 10.1158/1078-0432.CCR-16-1815.
7. Thompson, JK, et al. (2017). Asbestos-Induced Mesothelial to Fibroblastic Transition Is Modulated by the Inflammasome. *Am J Pathol.* **187**(3):665-678. doi: 10.1016/j.ajpath.2016.11.008.
8. Wang R, et al. (2017). Basic studies on epigenetic carcinogenesis of low-dose exposure to 1-trichloromethyl-1,2,3,4-tetrahydro- β -carboline (TaClo) in vitro. *PLoS One.* **12**(2):e0172243. doi: 10.1371/journal.pone.0172243.
9. Bell, J.B. et al. (2016). MNK inhibition disrupts mesenchymal glioma stem cells and prolongs survival in a mouse model of glioblastoma. *Mol. Cancer Res.* **14**:984-993.
10. Sheyn, D. et al. (2016). Human induced pluripotent stem cells differentiate into functional mesenchymal stem cells and repair bone defects. *Stem Cells Trans. Med.* **5**:1447-1460.
11. Lu, M. et al. (2016). The high expression of long non-coding RNA PANDAR indicates a poor prognosis for colorectal cancer and promotes metastasis by EMT pathway. *J Cancer Res Clin Oncol.* doi:10.1007/s00432-016-2252-y.
12. Qu, Y. et al. (2016). Axitinib blocks Wnt/ β -catenin signaling and directs asymmetric cell division in cancer. *Proc Natl Acad Sci U S A.* **113**:9339-9344.
13. Ciampi, R. et al. (2016). Classical point mutations of RET, BRAF and RAS oncogenes are not shared in papillary and medullary thyroid cancer occurring simultaneously in the same gland. *J Endocrinol Invest.* doi:10.1007/s40618-016-0526-5.
14. Bai, D. et al. (2016). The ATPase hCINAP regulates 18S rRNA processing and is essential for embryogenesis and tumour growth. *Nat Commun.* doi:10.1038/ncomms12310.
15. Hua, G. et al. (2016). The four and a half LIM domains 2 (FHL2) regulates ovarian granulosa cell tumor progression via controlling AKT1 transcription. *Cell Death Dis.* doi:10.1038/cddis.2016.207.
16. Wolfenson, H. et al. (2016). Tropomyosin controls sarcomere-like contractions for rigidity sensing and suppressing growth on soft matrices. *Nat Cell Biol.* **18**:33-42.
17. Martinez-Cruzado, L. et al. (2016). Aldh1 expression and activity increase during tumor evolution in sarcoma cancer stem cell populations. *Sci Rep.* doi:10.1038/srep27878.
18. Gong, J. et al. (2016). Downregulation of STAT3/NF- κ B potentiates gemcitabine activity in pancreatic cancer cells. *Mol Carcinog.* doi:10.1002/mc.22503.
19. Nakanishi, Y. et al. (2016). Activating mutations in PIK3CB confer resistance to PI3K inhibition and define a novel oncogenic role for p110 β . *Cancer Res.* doi:10.1158/0008-5472.
20. Tsukamoto, Y. et al. (2015). Expression of DDX27 contributes to colony-forming ability of gastric cancer cells and correlates with poor prognosis in gastric cancer. *Am J Cancer Res.* **5**:2998.
21. Bunda, S. et al. (2015). Inhibition of SHP2-mediated dephosphorylation of Ras suppresses oncogenesis. *Nat Commun.* doi:10.1038/ncomms9859.
22. He, C. et al. (2015). The Hippo/YAP pathway interacts with EGFR signaling and HPV oncoproteins to regulate cervical cancer progression. *EMBO Mol Med.* **7**:1426-1449.
23. Ercan, D. et al. (2015). EGFR mutations and resistance to irreversible pyrimidine based EGFR inhibitors. *Clin Cancer Res.* **21**:3913-3923.
24. Mayr, C. et al. (2015). 3-Deazaneplanocin A may directly target putative cancer stem cells in biliary tract cancer. *Anticancer Res.* **35**:4697-4705.

25. He, C. et al. (2015). The Hippo/YAP pathway interacts with EGFR signaling and HPV oncoproteins to regulate cervical cancer progression. *EMBO Mol Med*. doi:10.15252/emmm.201404976.
26. Hua, G. et al. (2015). YAP induces high-grade serous carcinoma in fallopian tube secretory epithelial cells. *Oncogene*. doi:10.1038/onc.2015.288.
27. Ukaji, T. et al. (2015). Inhibition of IGF-1-mediated cellular migration and invasion by migracin A in ovarian clear cell carcinoma cells. *PLoS One*. **10**:e0137663.
28. Fujita, T. et al. (2015). Identification and characterization of CXCR4-positive gastric cancer stem cells. *PLoS One*. **10**:e0130808.
29. Bon, H. et al. (2015). Salt-inducible Kinase 2 regulates mitotic progression and transcription in prostate cancer. *Mol Cancer Res*. **13**:620-635.

Please see the complete list of product citations: <http://www.cellbiolabs.com/96-well-cell-transformation-assays-standard-soft-agar>.

License Information

This product is provided under an intellectual property license from Life Technologies Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased product and components of the product only in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The sale of this product is expressly conditioned on the buyer not using the product or its components, or any materials made using the product or its components, in any activity to generate revenue, which may include, but is not limited to use of the product or its components: (i) in manufacturing; (ii) to provide a service, information, or data in return for payment; (iii) for therapeutic, diagnostic or prophylactic purposes; or (iv) for resale, regardless of whether they are resold for use in research. For information on purchasing a license to this product for purposes other than as described above, contact Life Technologies Corporation, 5791 Van Allen Way, Carlsbad CA 92008 USA or outlicensing@lifetech.com.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2004-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.