Product Manual

CytoSelect[™] 96-Well In Vitro Tumor Sensitivity Assay (Soft Agar Colony Formation)

Catalog Number

CBA-150	96 assays	
CBA-150-5	5 x 96 assays	

FOR RESEARCH USE ONLY Not for use in diagnostic procedures

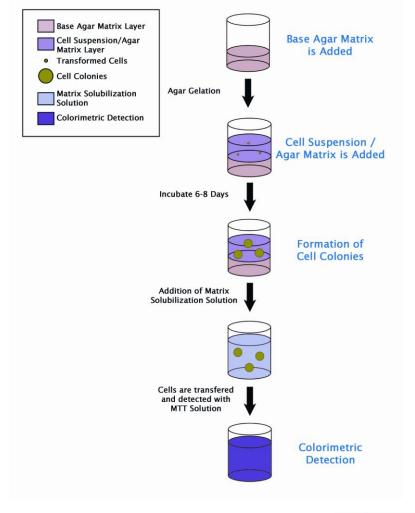


Introduction

Tumor sensitivity assays are intended to help predict the sensitivity of various tumors to chemotherapeutic agents, with the intent of identifying the most effective treatment with the fewest side effects. With this information, physicians can devise tailor made chemotherapy regiments and eliminate ineffective drugs, sparing patients of unnecessary toxicity. Ideally, an in vitro tumor sensitivity assay must be reliable, sensitive, and resemble the 3-D, in vivo environment (such as culturing in collagen gel or soft agar).

Traditionally, the soft agar colony formation assay is a common method to monitor anchorageindependent growth, which measures proliferation in a semisolid culture media after 3-4 weeks by manual counting of colonies. Cell Biolabs' CytoSelect[™] 96-well In Vitro Tumor Sensitivity 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 proprietary semisolid agar media before being solubilized, transferred and detected by the provided MTT Solution in a microtiter plate reader (see Assay Principle below).

The CytoSelect[™] 96-well In Vitro Tumor Sensitivity Assay provides a stringent, anchorageindependent model for chemosensitivity testing and potential anticancer drug screening. Each kit provides sufficient quantities to perform 96 tests in a microtiter plate.





Related Products

- 1. CBA-130: CytoSelect[™] 96-Well Cell Transformation Assay (Soft Agar Colony Formation)
- 2. CBA-135: CytoSelectTM 96-Well Cell Transformation Assay (Cell Recovery, Colorimetric)
- 3. CBA-140: CytoSelect[™] 96-Well Cell Transformation Assay (Cell Recovery, Fluorometric)
- 4. CBA-320: CytoSelect[™] 96-Well Hematopoietic Colony Forming Cell Assay

Kit Components (shipped at room temperature)

- 1. <u>10X CytoSelectTM Agar Matrix Solution</u> (Part No. 114001): One 10 mL sterile bottle
- 2. CytoSelectTM Matrix Diluent (Part No. 114002): One 4 mL sterile bottle
- 3. <u>5X DMEM Medium</u> (Part No. 113005): One 5 mL bottle
- 4. <u>1X Matrix Solubilization Buffer</u> (Part No. 115001): One 20 mL sterile bottle
- 5. Detergent Solution (Part No. 113501): One 10 mL bottle
- 6. <u>MTT Solution</u> (Part No. 113502): One 1 mL tube

Materials Not Supplied

- 1. Tumor Cells (cancer cell line or cells prepared from solid tumor)
- 2. Anticancer Agents (e.g., Taxol, 5-Fluorouracil, anticancer mAb or siRNA)
- 3. 37°C Incubator, 5% CO₂ Atmosphere
- 4. Light Microscope
- 5. 96-well Microtiter Plate Reader
- 6. 37°C and boiling water baths

<u>Storage</u>

Store all components at 4°C.

Preparation of Reagents

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.

• 10X CytoSelect[™] Agar Matrix Solution: Heat the Agar Matrix Solution bottle to 90-95°C in a water bath for 30 minutes, or until agar matrix liquefies (microwaving is optional). Transfer the bottle to a 37°C water bath for 20 minutes and maintain until needed.



Assay Protocol (must be under sterile conditions)

I. Preparation of Base Agar Matrix Layer

- 1. Heat the 10X CytoSelect[™] Agar Matrix Solution to 90-95°C in a water bath for 30 minutes, or until agar matrix liquefies (microwaving is optional). Transfer the bottle to a 37°C water bath for 20 minutes and maintain until needed.
- 2. Warm the 2X DMEM/20% FBS medium (see Preparation of Reagents section) to 37°C in a water bath. Allow at least 30 minutes for the temperature to equilibrate.
- 3. According to Table 1 (below), prepare the desired volume of Base Agar Matrix Layer in the following sequence:
 - a. In a sterile tube, add the appropriate volume of 2X DMEM/20% FBS medium.
 - b. Next, add the corresponding volume of sterile water. Mix well.
 - c. Finally, add the corresponding volume of 10X CytoSelect[™] Agar Matrix Solution. Mix well.

Note: The 10X CytoSelectTM Agar Matrix Solution is slightly viscous; care should be taken in accurately pipetting the appropriate volume.

2X DMEM/20%	Sterile Water	10X	Total Volume of	# of Tests in 96-
FBS Medium	(mL)	CytoSelect TM	Base Agar Matrix	well Plate (50
(mL)		Agar Matrix	Layer (mL)	μL/test)
		Solution (mL)		
2.5	2	0.5	5	100
1.25	1	0.25	2.5	50
0.5	0.4	0.1	1	20

Table 2. Preparation of Base Agar Matrix Layer

- 4. After mixing, maintain the Base Agar Matrix Layer at 37°C to avoid gelation.
- 5. Dispense 50 μL of Base Agar Matrix Layer into each well of a 96-well sterile flat-bottom microplate (samples should be assayed in triplicate). Gently tap the plate a few times to ensure the Base Agar Matrix Layer evenly covers the wells.

Notes:

- Work quickly with the layer 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.
- 6. Transfer the plate to 4°C for 30 minutes to allow the Base Agar Matrix Layer to solidify.
- 7. Prior to adding the Cell Suspension/Agar Matrix Layer (Section II), allow the plate to warm to room temperature for 30 minutes.



II. Addition of Cell Suspension/Agar Matrix Layer (under sterile conditions)

- 1. Heat the 10X CytoSelect[™] Agar Matrix Solution to 90-95°C in a water bath for 30 minutes, or until agar matrix liquefies (microwaving is optional). Transfer the bottle to a 37°C water bath for 20 minutes and maintain until needed.
- Warm the 2X DMEM/20% FBS medium (see Preparation of Reagents section) and CytoSelect[™] Matrix Diluent 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 $0.1 1 \ge 10^6$ cells/mL. Keep the cell suspension warm in a 37°C water bath.
- 4. According to Table 2 (below), prepare the desired volume of Cell Suspension/Agar Matrix Layer in the following sequence:
 - a. In a sterile tube, add the appropriate volume of 2X DMEM/20% FBS medium.
 - b. Next, add the corresponding volume of CytoSelectTM Matrix Diluent. Mix well.
 - c. Next, add the corresponding volume of 10X CytoSelect[™] Agar Matrix Solution. Mix well.
 - d. Finally, add the corresponding volume of cell suspension. Mix well.

Note: The CytoSelectTM Matrix Diluent and 10X CytoSelectTM Agar Matrix Solution are slightly viscous; care should be taken in accurately pipetting the appropriate volumes.

2X DMEM/20% FBS Medium (mL)	CytoSelect™ Matrix Diluent (mL)	10X CytoSelect [™] Agar Matrix Solution (mL)	Cell Suspension (mL)	Total Volume of Cell Suspension/ Agar Matrix Layer (mL)	# of Tests in 96-well Plate (75 µL/test)
3.5	2.75	0.75	0.5	7.5	100
1.75	1.375	0.375	0.25	3.75	50
0.875	0.688	0.188	0.125	1.875	25

 Table 3. Preparation of Cell Suspension/Agar Matrix Layer

- 5. After mixing, incubate the Cell Suspension/Agar Matrix Layer at room temperature for 5 minutes.
- Immediately dispense 75 μL of Cell Suspension/Agar Matrix Layer into each well of the 96well plate, already containing the Base Agar Matrix Layer (Section I). *Notes:*
 - Work quickly with the layer 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 Suspension/Agar Matrix Layer.
- 7. Transfer the plate to 4°C for 20 minutes to allow the Cell Suspension/Agar Matrix Layer to solidify.
- 8. Allow the plate to warm to room temperature for 30 minutes.



- 9. Add 50 μ L of culture medium containing anticancer agents (e.g. Taxol, 5-Fluorouracil, mAb, etc.) to each well.
- 10. Incubate the cells for 6-8 days at 37°C and 5% CO₂. Examine the colony formation under a light microscope.

III. Quantitation of Anchorage-Independent Growth

- 1. Add 125 μ L of the 1X Matrix Solubilization Buffer to each well.
- 2. Pipette the entire volume of the well 10-12 times to mix thoroughly and solubilize the agar matrix completely.
- 3. Transfer 100 μ L of the mixture to a 96-well microtiter plate.
- 4. Add 10 μ L of MTT Solution to each well. Pipette each well 7-10 times to ensure a homogeneous mixture.
- 5. Incubate the plate for 2-4 hours at 37°C and 5% CO₂. *Note: Under the microscope, a purple precipitate should be visible within the cells.*
- 6. Add $100 \,\mu\text{L}$ of Detergent Solution to each well.
- 7. Incubate the plate in the dark for 2-4 hours at room temperature.
- 8. Pipette each well 7-10 times to ensure a homogeneous mixture.
- 9. Measure the absorbance at 570 nm in a 96-well microtiter plate reader.

<u>Cell Dose Curve (optional)</u>

- 1. Heat the 10X CytoSelect[™] Agar Matrix Solution to 90-95°C in a water bath for 30 minutes, or until agar matrix liquefies (microwaving is optional). Transfer the bottle to a 37°C water bath for 20 minutes and maintain until needed.
- 2. Warm the 2X DMEM/20% FBS medium (see Preparation of Reagents section) and CytoSelect[™] Matrix Diluent 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 5 10×10^6 cells/mL.
- 4. Prepare a serial 2-fold dilution in culture medium, including a blank without cells.
- 5. Transfer 50 μ L of each dilution to a 96-well plate.
- 6. According to Table 3 (below), prepare the desired volume of Cell Dose Curve Solution in the following sequence:
 - a. In a sterile tube, add the appropriate volume of 2X DMEM/20% FBS medium.
 - b. Next, add the corresponding volume of sterile water. Mix well.
 - c. Next, add the corresponding volume of CytoSelectTM Matrix Diluent. Mix well.
 - d. Finally, add the corresponding volume of 10X CytoSelect[™] Agar Matrix Solution. Mix well.



Note: The CytoSelectTM Matrix Diluent and 10X CytoSelectTM Agar Matrix Solution are slightly viscous; care should be taken in accurately pipetting the appropriate volumes.

2X	Sterile Water	CytoSelect TM	10X	Total Volume of
DMEM/20%	(mL)	Matrix Diluent	CytoSelect [™]	Cell Dose Curve
FBS Medium		(mL)	Agar Matrix	Solution (mL)
(mL)			Solution (mL)	
1.25	0.45	0.55	0.25	2.5
0.625	0.225	0.275	0.125	1.25

Table 4. Preparation of Cell Dose Curve Solution

- 7. Immediately dispense $125 \ \mu$ L of Cell Dose Curve Solution into the wells of the 96-well plate, already containing the cell serial dilution (from step 5).
- 8. Add 125 μ L of 1X Matrix Solubilization Buffer to each well. Pipette each well 10-12 times to mix thoroughly.
- 9. Transfer 100 μ L of the mixture to a 96-well microtiter plate.
- 10. Add 10 μ L of MTT Solution to each well. Pipette each well 7-10 times to ensure a homogeneous mixture.
- 11. Incubate the plate for 2-4 hours at 37°C and 5% CO₂. *Note: Under the microscope, a purple precipitate should be visible within the cells.*
- 12. Add 100 µL of Detergent Solution to each well.
- 13. Incubate the plate in the dark for 2-4 hours at room temperature.
- 14. Pipette each well 7-10 times to ensure a homogeneous mixture.
- 15. Measure the absorbance at 570 nm in a 96-well microtiter plate reader.

Example of Results

The following figures demonstrate typical results with the CytoSelect[™] Cell Transformation Assay Kit. One should use the data below for reference only. This data should not be used to interpret actual results.



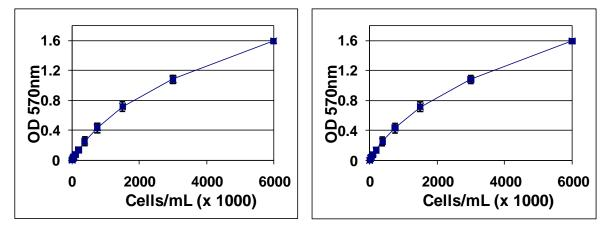


Figure 1. HeLa Cell Dose Curve. Cervical carcinoma HeLa cells were resuspended at $6 \ge 10^6$ cells/mL and titrated 1:2 in culture medium, followed by addition of Cell Dose Curve Solution, Matrix Solubilization Solution, MTT Solution, and Detergent Solution (as described in the Cell Dose Section). Results are shown by cell concentration or by actual cell number in MTT Detection.

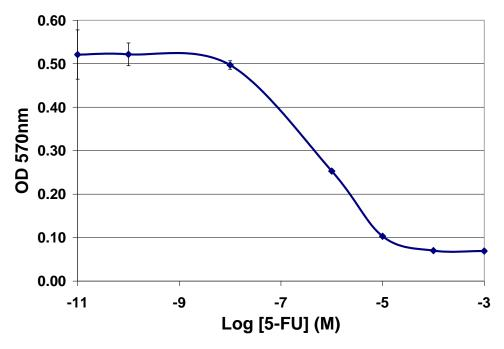


Figure 2. Inhibition of Hela Cell Transformation by 5-Fluorouracil. HeLa cells were seeded at 5000 cells/well and cultured 7 days at various 5-FU concentrations. Cell transformation was determined according to the assay protocol. IC50 value of 5-Fluorouracil on HeLa cell anchorage-independent growth was determined to be $\sim 1 \mu M$.



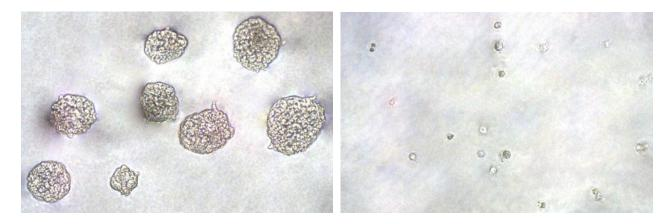


Figure 3. Inhibition of HeLa Cell Anchorage-Independent Growth by Taxol. HeLa cells were cultured for 7 days in the absence (left) or presence (right) of 1 nM Taxol according to the assay protocol.

Calculation of Anchorage-Independent Growth

- 1. Compare OD_{570nm} values with the Cell Dose Curve and extrapolate the cell concentration.
- 2. Calculate the Total Transformed Cell Number/Well Total Transformed Cells/Well = cells/mL x 0.050 mL/well

For example: If you extrapolate your OD_{570nm} value from your cell dose curve and determine you have 500,000 cells/mL in your sample. **Total Transformed Cells/Well** = 500,000 cells/mL x 0.050 mL/well = 25,000 cells/well

References

- 1. Shin SI, Freedman VH, Risser R, and Pollack R. (1975) Proc Natl Acad Sci USA. 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. Nakamura D. (2023). The evaluation of tumorigenicity and characterization of colonies in a soft agar colony formation assay using polymerase chain reaction. *Sci Rep.* **13**(1):5405. doi: 10.1038/s41598-023-32442-6.
- Chandrasekaran, B. et al. (2023). Antiandrogen-Equipped Histone Deacetylase Inhibitors Selectively Inhibit Androgen Receptor (AR) and AR-Splice Variant (AR-SV) in Castration-Resistant Prostate Cancer (CRPC). *Cancers (Basel)*. 15(6):1769. doi: 10.3390/cancers15061769.
- 3. Nakachi, S. et al. (2022). Impact of anti-diabetic sodium-glucose cotransporter 2 inhibitors on tumor growth of intractable hematological malignancy in humans. *Biomed Pharmacother*. doi: 10.1016/j.biopha.2022.112864.
- 4. Katayama, Y. et al. (2022). Heterogeneity among tumors with acquired resistance to EGFR tyrosine kinase inhibitors harboring EGFR-T790M mutation in non-small cell lung cancer cells. *Cancer Med.* doi: 10.1002/cam4.4504.
- 5. Xie, X. et al. (2021). Babao Dan is a robust anti-tumor agent via inhibiting wnt/β-catenin activation and cancer cell stemness. *J Ethnopharmacol*. doi: 10.1016/j.jep.2021.114449.



- Dehghanian, S.Z. et al. (2021). ABT-751 Induces Multiple Anticancer Effects in Urinary Bladder Urothelial Carcinoma-Derived Cells: Highlighting the Induction of Cytostasis through the Inhibition of SKP2 at Both Transcriptional and Post-Translational Levels. *Int J Mol Sci.* 22(2):945. doi: 10.3390/ijms22020945.
- Fujimura, T. et al. (2020). Enhanced antitumor effect of alectinib in combination with cyclindependent kinase 4/6 inhibitor against RET-fusion-positive non-small cell lung cancer cells. *Cancer Biol Ther.* doi: 10.1080/15384047.2020.1806643.
- 8. Chandrasekaran, B. et al. (2020). Chronic exposure to cadmium induces a malignant transformation of benign prostate epithelial cells. *Oncogenesis*. **9**(2):23. doi: 10.1038/s41389-020-0202-7.
- Tyagi, A. et al. (2019). Combination of androgen receptor inhibitor and cisplatin, an effective treatment strategy for urothelial carcinoma of the bladder. *Urol Oncol.* pii: S1078-1439(19)30101-2. doi: 10.1016/j.urolonc.2019.03.008.
- Wei, R.J. et al. (2019). Inhibition of the formation of autophagosome but not autolysosome augments ABT-751-induced apoptosis in TP53-deficient Hep-3B cells. *J Cell Physiol*. 234(6):9551-9563. doi: 10.1002/jcp.27643.
- 11. Wang, L. et al. (2018). Effects of ebv-miR-BART7 on tumorigenicity, metastasis, and TRAIL sensitivity of non-small cell lung cancer. *J Cell Biochem*. doi: 10.1002/jcb.28289.
- 12. von Frowein, J. et al. (2018). MiR-492 regulates metastatic properties of hepatoblastoma via CD44. *Liver Int.* **38**(7):1280-1291. doi: 10.1111/liv.13687.
- 13. Pal, D. et al. (2018). Suppression of Notch1 and AKT mediated epithelial to mesenchymal transition by Verrucarin J in metastatic colon cancer. *Cell Death Dis*. **9**(8):798. doi: 10.1038/s41419-018-0810-8.
- Chandrasekaran, B. et al. (2017). Molecular insights: Suppression of EGFR and AKT activation by a small molecule in non-small cell lung cancer. *Genes Cancer*. 8(9-10):713-724. doi: 10.18632/genesandcancer.154.
- 15. Pal, D. et al. (2017). Inhibition of autophagy prevents cadmium-induced prostate carcinogenesis. *Br. J. Cancer* **117(1)**:56-64.
- 16. Kato, K. et al. (2017). Opposite effects of tumor protein D (TPD) 52 and TPD54 on oral squamous cell carcinoma cells. *Int J Oncol.* doi: 10.3892/ijo.2017.3929
- 17. Wei, R. J. et al. (2016). A microtubule inhibitor, ABT-751, induces autophagy and delays apoptosis in Huh-7 cells. *Toxicol Appl Pharmacol*. doi:10.1016/j.taap.2016.09.021.
- 18. Mukudai, Y. et al. (2016). Methanol and butanol extracts of Paeonia lutea leaves repress metastasis of squamous cell carcinoma. *Evid Based Complement Alternat Med.* doi:10.1155/2016/6087213.
- 19. Damodaran, C. et al. (2016). miR-301a expression: A prognostic marker for prostate cancer. *Urol Oncol.* doi: 10.1016/j.urolonc.2016.03.009.
- 20. Zheng, Y. et al. (2016). Glioma-derived platelet-derived growth factor-BB recruits oligodendrocyte progenitor cells via platelet-derived growth factor receptor-α and remodels cancer stroma. *Am J Pathol.* doi: 10.1016/j.ajpath.2015.12.020.
- 21. Joshi, P. et al. (2015). MicroRNA-148a reduces tumorigenesis and increases TRAIL-induced apoptosis in NSCLC. *Proc Natl Acad Sci U S A*. **112**:8650-8655.
- 22. Meador, C. B. et al. (2015). Optimizing the sequence of anti-EGFR-targeted therapy in EGFRmutant lung cancer. *Mol Cancer Ther.* **14**:542-552.
- 23. Peng, Y. T. et al. (2015). Upregulation of cyclin-dependent kinase inhibitors CDKN1B and CDKN1C in hepatocellular carcinoma-derived cells via goniothalamin-mediated protein stabilization and epigenetic modifications. *Toxicol Rep.* doi: 10.1016/j.toxrep.2015.01.010.



- 24. Akl, M. R. et al. (2015). Araguspongine C induces autophagic death in breast cancer cells through suppression of c-Met and HER2 receptor tyrosine kinase signaling. *Mar Drugs.* **13**:288-311.
- 25. Suman, S. et al. (2014). The pro-apoptotic role of autophagy in breast cancer. *Bri J Cancer*. **111**:309-317.
- 26. Suman, S. et al. (2014). Activation of AKT signaling promotes epithelial–mesenchymal transition and tumor growth in colorectal cancer cells. *Mol Carcinog.* **53**:E151-E160.
- 27. Bard-Chapeau, E. et al. (2013). EVI1 oncoprotein interacts with a large and complex network of proteins and integrates signals through protein phosphorylation. *PNAS*. **110**:E2885-E2894.
- 28. Takezawa. K. et al. (2012). HER2 amplification: a potential mechanism of acquired resistance to EGFR inhibition in EGFR-mutant lung cancers that lack the second-site EGFRT790M mutation. *Cancer Discovery*. **2**: 922-933.
- 29. Li, C. et al. (2012). The root bark of Paeonia moutan is a potential anticancer agent in human oral squamous cell carcinoma cells. *Anticancer Res.* **32**:2625-2630.
- 30. Itamochi, H. et al. (2011). Inhibiting the mTOR pathway synergistically enhances cytotoxicity in ovarian cancer cells induced by etoposide through upregulation of c-Jun. *Clin. Cancer Res.* **17**:4742-4750.

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. 5628 Copley Drive San Diego, CA 92111 Worldwide: +1 858-271-6500 USA Toll-Free: 1-888-CBL-0505 E-mail: <u>tech@cellbiolabs.com</u> www.cellbiolabs.com

©2006-2024: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.

