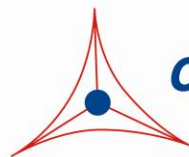

Product Manual

96-Well Cellular Senescence Assay Kit (SA- β -gal Activity, Fluorometric Format)

Catalog Number

CBA-231	120 assays
CBA-231-5	5 x 120 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Normal primary cells proliferate in culture for a limited number of population doublings prior to undergoing terminal growth arrest and acquiring a senescent phenotype. This finite life span correlates with the age of the organism and with the life expectancy of the species from which the cells were obtained; such that the older the age or the shorter the life span, the less the ability of the cells to undergo population doubling. Senescent cells are characterized by an irreversible G₁ growth arrest involving the repression of genes that drive cell cycle progression and the upregulation of cell cycle inhibitors like p16^{INK4a}, p53, and its transcriptional target, p21^{CIP1}. They are resistant to mitogen-induced proliferation, and assume a characteristic enlarged, flattened morphology. Research into the pathways that positively regulate senescence and ways cells bypass senescence is therefore critical in understanding carcinogenesis. Normal cells have several mechanisms in place to protect against uncontrolled proliferation and tumorigenesis.

Senescent cells show common biochemical markers such as expression of an acidic senescence-associated β -galactosidase (SA- β -Gal) activity. While senescence has been characterized primarily in cultured cells, there is also evidence that it occurs *in vivo*. Cells expressing markers of senescence such as SA- β -Gal have been identified in normal tissues.

The 96-well Cellular Senescence Assay Kit provides an easy-to-use and efficient method to determine the cellular senescence by measuring SA- β -Gal activity using a fluorometric substrate. This quantitative assay uses cell lysate for both SA- β -galactosidase activity determination and normalization of samples containing different cell numbers. Each kit provides sufficient quantities to perform up to 120 assays in a 96-well plate.

Related Products

1. CBA-080: CytoSelect™ 24-Well Anoikis Assay
2. CBA-081: CytoSelect™ 96-Well Anoikis Assay
3. CBA-230: Cellular Senescence Assay Kit (SA- β -gal Staining)
4. CBA-232: Quantitative Cellular Senescence Assay (SA β -Gal)
5. CBA-240: CytoSelect™ Cell Viability and Cytotoxicity Assay
6. AKR-100: β Galactosidase Staining Kit

Kit Components

1. 2X Cell Lysis Buffer (Part No. 123101): One 10 mL bottle
2. 2X Reaction Buffer (Part No. 123102): One 10 mL bottle
3. SA- β -Gal Substrate (20X) (Part No. 123103): One 300 μ L amber tube
4. Stop Solution (Part No. 123104): One 25 mL bottle

Materials Not Supplied

1. Senescent cells or tissue samples
2. 37°C Incubator
3. β -mercaptoethanol
4. 96-well plate suitable for a fluorescence plate reader
5. 96-well Fluorometer
6. Protein Assay Reagents

Storage

Store SA- β -gal substrate solution protected from light at -20°C. Store all other components at room temperature.

Preparation of Reagents

- 1X Cell Lysis Buffer: Prepare a 1X Cell Lysis Buffer by diluting the provided 2X stock 1:2 in ddH₂O. Store the diluted solution at room temperature for up to six months. Immediately before use, add proper amount of proteinase inhibitors such as PMSF.
- 2X Assay Buffer: Immediately before use, add β -mercaptoethanol to 2X Reaction Buffer at a final concentration of 10 mM and dilute 20X SA- β -Gal Substrate to 1X with 2X Reaction Buffer containing 10 mM β -mercaptoethanol. Don't store 2X Assay Buffer.

Reagent	96-well	24-well	6-well	10 cm
1X Cell Lysis Buffer	100 μ L	400 μ L	1000 μ L	1500 μ L

Assay Protocol

1. Aspirate the medium from the senescent cells.
2. Wash the cells once with 200 μ L of cold 1X PBS and aspirate the wash.
3. Add 100 μ L of cold 1X Cell Lysis Buffer (see the table above for the required amount of 1X Cell Lysis Buffer of other plate formats). Incubate at 4°C for 5 minutes. Transfer the whole lysate to a microcentrifuge tube and centrifuge 10 minutes at 4°C. Collect supernatant as cell lysate.
4. (optional) Determine the total protein concentration of each cell lysate sample by protein assay such as Pierce's BCA protein Assay.
5. Transfer 50 μ L of the cell lysate to a 96-well plate. Add 50 μ L of freshly prepared 2X Assay Buffer. Incubate the wells at 37°C protected from light for 1- 3 hr.
6. Remove 50 μ L of the reaction mixture to a 96-well plate suitable for fluorescence measurement. Stop the reaction by adding 200 μ L of Stop solution.
7. Read fluorescence with a fluorescence plate reader at 360 nm (Excitation) / 465 nm (Emission).

Example of Results

The following figures demonstrate typical with the 96-well Cellular Senescence Assay Kit. Fluorescence measurement was performed on SpectraMax Gemini XS Fluorometer (Molecular Devices) with a 355 nm/460 nm filter set. One should use the data below for reference only. This data should not be used to interpret actual results.

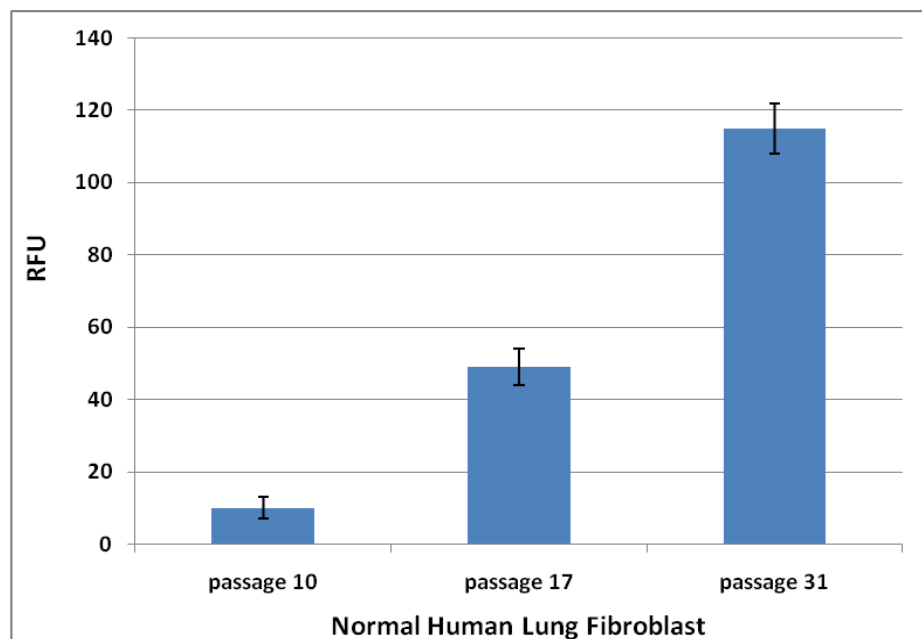


Figure 1: SA-β-Gal activity in Senescent Human Lung Fibroblast HFL-1 Cells. Normal HFL-1 cells with different passage numbers were lysed. Lysates were allowed to incubate with SA-β-Gal Substrate for 1 hr at 37°C. SA-β-Gal activities were measured as described in the Assay Protocol.

References

1. Current Protocols in Molecular Biology, John Wiley & Sons Press.
2. Campisi, J. (2000) *In Vivo* 14, 183-188.
3. Dimri, G. P., X. Lee, G. Basile, M. Acosta, G. Scott, C. Roskelley, E. E. Medrano, M. Linskens, I. Rubelj, O. Pereira-Smith, M. Peacocke, and J. Campisi. (1995) *Proc. Natl. Acad. Sci. USA* 92:9363-9367.

Recent Product Citations

1. Park, J.E., et al. (2017). Standardized *Kaempferia parviflora* Extract Inhibits Intrinsic Aging Process in Human Dermal Fibroblasts and Hairless Mice by Inhibiting Cellular Senescence and Mitochondrial Dysfunction. *Evid Based Complement Alternat Med.* 2017:6861085. doi: 10.1155/2017/6861085.
2. Zhang, X. et al. (2016). Cryptosporidium parvum infection attenuates the ex vivo propagation of murine intestinal enteroids. *PHY2* 4:e13060.

3. Platas, J. et al. (2016). Anti-senescence and anti-inflammatory effects of the C-terminal moiety of PTHrP peptides in OA osteoblasts. *J Gerontol A Biol Sci Med Sci*. doi:10.1093/gerona/glw100.
4. Sato, K. et al. (2016). Exosomes in liver pathology. *J Hepatol*. doi:10.1016/j.jhep.2016.03.004.
5. Kim, D. S. et al. (2016). Cell culture density affects the proliferation activity of human adipose tissue stem cells. *Cell Biochem Funct*. doi:10.1002/cbf.3158.
6. Hu, W. et al. (2015). Mechanistic investigation of bone marrow suppression associated with palbociclib and its differentiation from cytotoxic chemotherapies. *Clin Cancer Res*. doi:10.1158/1078-0432.CCR-15-1421.
7. Chang, Z. et al. (2015). Ascorbic acid provides protection for human chondrocytes against oxidative stress. *Mol Med Rep*. doi:10.3892/mmr.2015.4231.
8. Hirano, T. et al. (2015). A novel interaction between FLICE-associated huge protein (FLASH) and E2A regulates cell proliferation and cellular senescence via tumor necrosis factor (TNF)-alpha-p21WAF1/CIP1 axis. *PLoS One*. **10**:e0133205.
9. Liao, C. K. et al. (2015). Depletion of B cell CLL/lymphoma 11B gene expression represses glioma cell growth. *Mol Neurobiol*. doi:10.1007/s12035-015-9231-1.
10. Shimizu, R. et al. (2015). Cholangiocyte senescence caused by lysophosphatidylcholine as a potential implication in carcinogenesis. *J Hepatobiliary Pancreat Sci*. doi: 10.1002/jhbp.256.
11. Klinkhammer, B. M. et al. (2014). Mesenchymal stem cells from rats with chronic kidney disease exhibit premature senescence and loss of regenerative potential. *PLoS One*. **9**:e92115-e92115.
12. Li, Y. et al. (2014). Inhibition of APE1/Ref-1 redox activity rescues human retinal pigment epithelial cells from oxidative stress and reduces choroidal neovascularization. *Redox Biol*. **2**:485-494.
13. Qi, H. et al. (2014). COH-203, a novel microtubule inhibitor, exhibits potent anti-tumor activity via p53-dependent senescence in hepatocellular carcinoma. *Biochem Biophys Res Commun*. **455**:262-268.
14. Lu, J. et al. (2014). SIRT1 counteracted the activation of STAT3 and NF-κB to repress the gastric cancer growth. *Int J Clin Exp Med*. **7**:5050-5058.

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

©2007-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.