



Cellular Senescence Assay Kit

KAA002

**FOR RESEARCH USE ONLY
Not for use in diagnostic procedures.**

USA & Canada

Phone: +1(800) 437-7500 • Fax: +1 (909) 676-9209 • Europe +44 (0) 23 8026 2233

Australia +61 3 9839 2000 • Germany +49-6192-207300 • • ISO Registered worldwide •

www.chemicon.com • custserv@chemicon.com • techserv@chemicon.com

Introduction

Cellular senescence is one of the most fundamental aspects of cell behavior, and is thought to play a critical role in regulating cellular lifespan both *in vitro* and *in vivo* [1-3]. Primary somatic cells grown *in vitro* do not proliferate indefinitely. Instead, after a period of rapid proliferation, cell division rate slows, and ultimately ceases altogether, with the cells becoming unresponsive to mitogenic stimuli. This process is termed cellular senescence, and senescent cells have a well-defined accompanying phenotype - increased cell size, distinctive flat morphology, accumulated lipofuscin granules, wide changes in gene expression, and activity of senescence-associated β -galactosidase (SA- β -gal) [2,3].

It is generally believed that cellular senescence reflects some of the changes that occur during the aging of organisms, and senescent cells have been detected *in vivo* at sites of age-related pathology, such as benign hyperplastic prostate [4] and atherosclerotic lesions [5]. Recent studies have also provided convincing demonstrations of cellular senescence occurring *in vivo* in response to internally- and externally-induced stress signals [6,7]. In all of these studies, senescence was characterized by the appearance of senescence-associated β -galactosidase (SA- β -gal) activity, in common with the senescent phenotype *in vitro*.

Cellular senescence has become an increasingly important target in the development of novel therapeutics. Emerging data implicates senescence bypass in the development of cancer and suggests that senescence may represent a tumor suppressor mechanism. The demonstration that tumor cells can be induced to undergo replicative senescence following the introduction of negative cell-cycle regulators, anti-telomerase peptides, or drug treatment suggests that induction of senescence can be exploited as a basis for cancer therapy [8,9].

For Research Use Only; Not for use in diagnostic procedures

Test Principle

As outlined above, a classic characteristic of the senescent phenotype is the induction of senescence-associated β -galactosidase (SA- β -gal) activity. SA- β -gal is present only in senescent cells, not in presenescent, quiescent, or proliferating cells. Chemicon's Cellular Senescence Assay Kit provides all the reagents required to efficiently detect SA- β -gal activity at pH 6.0 in cultured cells and tissue sections. In this assay, SA- β -gal catalyzes the hydrolysis of X-gal, which results in the accumulation of a distinctive blue color in senescent cells. Each kit provides sufficient quantities of reagents to perform at least 50 assays in 35 mm wells.

Kit Components

1. 100X Fixing Solution: (Part No. 2004755) One 1.5 mL vial
2. 10X Staining Solution A: (Part No. 2004756) One 15 mL bottle
3. 10X Staining Solution B: (Part No. 2004754) One 15 mL bottle
4. X-gal Solution: (Part No. 2004752) Two 1.5 mL vials

Materials Not Supplied

1. 1X PBS (Phosphate Buffered Saline)
2. 37°C incubator
3. Phase contrast or light microscope
4. Polypropylene tubes
5. Cells / Tissue samples
6. 70% glycerol (optional)

Precautions

- Please refer to the Material Safety Data Sheet at www.chemicon.com for any necessary precautions.

Storage

Store X-gal solution protected from light at -20°C, and other kit components at 4°C. All components supplied are stable for 1 year.

Assay Instructions

The following protocol is designed for one 35mm well of a 6-well plate.

Set-up:

1. Prepare a 1X PBS solution (not provided).
2. Dilute the 100X Fixing Solution with PBS to make a 1X working solution. 1mL fixing solution /well is sufficient (*e.g.* Add 10 μ L fixing solution to 990 μ L PBS).
3. Prepare SA- β -gal Detection Solution by mixing Staining Solutions A and B, X-Gal and PBS as indicated in Table 1. Prepare fresh for each use. See Table 1 for suggested volumes. Mix well before use.

Table 1. Preparation of SA- β -gal Detection Solution

Reagents	1 well (35 mm)	5 wells (35 mm)	10 wells (35 mm)
Staining Solution A (10X)	200 μ L	1 mL	2 mL
Staining Solution B (10X)	200 μ L	1 mL	2 mL
X-Gal	50 μ L	250 μ L	500 μ L
PBS	1.55 mL	7.75 mL	15.5 mL
Total	2 mL	10 mL	20 mL

Procedure:

1. Aspirate the growth medium from the cells.
2. Wash the cells once with 2 mL 1X PBS and aspirate the wash.
3. Add 1 mL 1X Fixing Solution per well. Incubate at room temperature for 10 - 15 minutes.
4. Remove the Fixing Solution and wash the cells twice with 2 mL 1X PBS. Aspirate after each wash.
5. Add 2 mL of freshly prepared 1X SA- β -gal Detection Solution.
6. Incubate the cells at 37°C, without CO₂ and protected from the light, for at least 4 hours. Samples may safely be left overnight.
7. Remove the SA- β -gal Detection Solution. Wash the stained cells twice with 2mL PBS. Aspirate after each wash.
8. Count the blue stained cells under phase contrast, or light microscopy.
9. For long term storage, overlay the stained cells with 70% glycerol – diluted in 1X PBS. Store at 4 - 8°C.

Sample Results



References

1. Ball AJ & Levine F. Telomere-independent cellular senescence in human fetal cardiomyocytes. *Aging Cell*. 2005 4:21-30.
2. Dimri GP *et al*. A biomarker that identifies senescent human cells in culture and in aging skin in vivo. *Proc. Natl Acad. Sci. USA* 1995 92:9363–9367.
3. Duncan EL *et al*. Senescence and immortalization of human cells. *Biogerontology* 2000 1:103–121.
4. Choi J *et al*. Expression of senescence-associated beta-galactosidase in enlarged prostates from men with benign prostatic hyperplasia. *Urology* 2000 56:160–166.
5. Vasile E *et al*. Differential expression of thymosin beta-10 by early passage and senescent vascular endothelium is modulated by VPF/VEGF: evidence for senescent endothelial cells in vivo at sites of atherosclerosis. *FASEB J* 2003 15:458–466.
6. Satyanarayana A *et al*. Telomere shortening impairs organ regeneration by inhibiting cell cycle re-entry of a subpopulation of cells. *EMBO J*. 2003 22:4003–4013.
7. Schmitt CA *et al*. A senescence program controlled by p53 and p16INK4a contributes to the outcome of cancer therapy. *Cell* 2002 109:335–346.
8. Marcotte R & Wang E. Replicative senescence revisited. *J Gerontol A Biol Sci Med Sci*. 2002 57:B257-69.
9. Dimri GP. What has senescence got to do with cancer? *Cancer Cell* 2005 7:505-512.

Warranty

EMD Millipore Corporation (“EMD Millipore”) warrants its products will meet their applicable published specifications when used in accordance with their applicable instructions for a period of one year from shipment of the products. **EMD MILLIPORE MAKES NO OTHER WARRANTY, EXPRESSED OR IMPLIED. THERE IS NO WARRANTY OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE.** The warranty provided herein and the data, specifications and descriptions of EMD Millipore products appearing in EMD Millipore’s published catalogues and product literature may not be altered except by express written agreement signed by an officer of EMD Millipore. Representations, oral or written, which are inconsistent with this warranty or such publications are not authorized and if given, should not be relied upon.

In the event of a breach of the foregoing warranty, EMD Millipore Corporation’s sole obligation shall be to repair or replace, at its option, the applicable product or part thereof, provided the customer notifies EMD Millipore Corporation promptly of any such breach. If after exercising reasonable efforts, EMD Millipore Corporation is unable to repair or replace the product or part, then EDM Millipore shall refund to the Company all monies paid for such applicable Product. **EMD MILLIPORE CORPORATION SHALL NOT BE LIABLE FOR CONSEQUENTIAL, INCIDENTAL, SPECIAL OR ANY OTHER DAMAGES RESULTING FROM ECONOMIC LOSS OR PROPERTY DAMAGE SUSTAINED BY ANY COMPANY CUSTOMER FROM THE USE OF ITS PRODUCTS.**

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

(c) 2009 - 2016: Merck KGaA, Darmstadt. All rights reserved. No part of these works may be reproduced in any form without permission in writing

KAA002

August 2016
Version 3: 4002519

