

# Cellular senescence activity assay

## Highly reproducible SA-β-gal activity kit

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.

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 G<sub>1</sub> growth arrest involving the repression of genes that drive cell cycle progression and the upregulation of cell cycle inhibitors like p16<sup>INK4a</sup>, p53, and its transcriptional target, p21<sup>CIP1</sup>. 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.

Citations: 27

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-KIT129-0120

120 assays

## Manuals, SDS & CofA

[View Online »](#)

- Quantify SA-β-gal activity in a fluorescence plate reader
- Easy-to-use protocol with results in less than 90 minutes
- Fully compatible with all cell types

## Handling & Storage

**Handling** Store SA- $\beta$ -gal substrate solution protected from light at -20°C. Store all other components at room temperature.

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Fluorescent detection, HTS

**Application Notes** For the determination of cellular senescence by measuring SA- $\beta$ -gal activity using a fluorometric substrate.

**Contents**

- 2X Cell lysis buffer
- 2X Reaction buffer
- SA- $\beta$ -gal substrate (20X)
- Stop solution

**UniProt ID** P16278

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)