Cellular senescence activity assay

Highly reproducible SA-\(\beta\)-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_1 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.

Citations: 25

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-β-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-β-gal activity using a fluorometric substrate.

Contents 2X Cell lysis buffer

2X Reaction buffer

SA-β-gal substrate (20X)

Stop solution

UniProt ID P16278

Last modified: May 29, 2024

