

Senescence Staining Assay

To detect to detect β -Gal activity in senescent cells.

Cellular Senescence is the phenomenon wherein cells stop dividing. It can be induced by both internal (oxidative stress, DNA damage, etc.) and external (toxins, UV light, etc.) factors Senescence is a state of arrested growth in which cells remain viable but stop dividing. Senescent cells display increase of cell size, senescence-associated expression of β -galactosidase, p53 and P16 activity, along with altered patterns of gene expression. AkrisBio's Senescence Detection Kit is designed to detect β -Gal activity. β -Gal is present primarily in senescent cells but can give a false positive with cells that normally express β -gal such as mature macrophages.

Ordering Information

[Order Online »](#)

AKR-COS-0102	1Kit
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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