

Cellular Senescence Detection Kit (SA- β - Gal Staining)

Our Cellular Senescence Staining Kit provides an efficient method to visualize Senescence Associated (SA) β -galactosidase. SA- β -Gal catalyzes the hydrolysis of X-gal, which produces a blue color in senescent cells. Visualize results with a standard light microscope.

Ordering Information

[Order Online »](#)

CBI-CBA-230	50 assays
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Stain cells for senescence-associated β -galactosidase
- Assay performed in a standard 35 mm culture dish

Handling & Storage

Use/Stability Store X-gal solution protected from light at -20°C. Store all other components at 4°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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