

Product datasheet

Senescence Detection Kit ab65351

★★★★★ 1 Abreviews 76 References 1 Image

Overview

Product name	Senescence Detection Kit
Detection method	Colorimetric
Sample type	Tissue, Adherent cells
Assay type	Enzyme activity
Assay time	1h 10m
Product overview	Senescence Detection Kit (ab65351) is designed to histochemically detect SA-beta-Gal activity in cultured cells and tissue sections, a known characteristic of senescent cells. The SA-beta-Gal is present only in senescent cells and is not found in presenescent, quiescent or immortal cells.

See Senescence Assay Kit [ab228562](#) to detect beta galactosidase activity in senescent cells by flow cytometry.

Senescence assay protocol summary:

- wash cells / sections with PBS
- fix with fixative solution for 10 min
- wash with PBS
- incubate in staining solution mix for 1 hr
- analyze staining with a microscope

Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K320 Senescence Detection Kit. K320-250 is the same size as the 250 test size of ab65351.</p> <p>Senescence is thought to be a tumor suppressive mechanism and an underlying cause of aging. Senescence represents an arrested state in which the cells remain viable, but not stimulated to divide by serum or passage in culture. Senescent cells display increase of cell size, senescence-associated expression of beta-galactosidase (SA-beta-Gal) activity, and altered patterns of gene expression.</p>
--------------	---

Properties

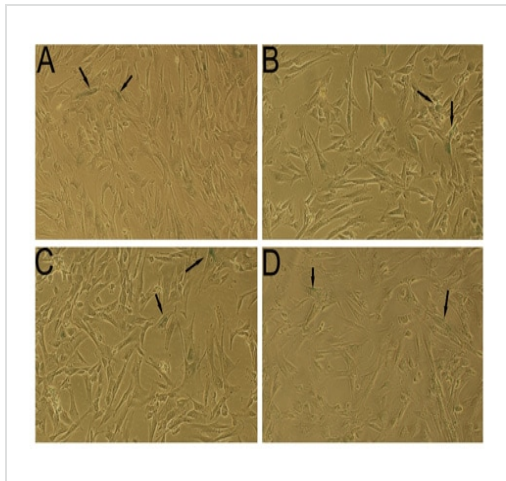
Storage instructions	Store at -20°C. Please refer to protocols.
Storage buffer	Preservative: None

Components	Identifier	250 tests	5000 tests
100X Staining Supplement	Red	1 x 1.5ml	20 x 1.5ml
1X Fixative Solution	NM	1 x 125ml	20 x 125ml
1X Staining Solution	WM	1 x 125ml	20 x 125ml
X-Gal (150 mg, lyophilized)	Green	1 vial	20 vials

Relevance

Cellular senescence is a growth-arrest program by which normal diploid cells lose the ability to divide, and it plays a critical role in regulating lifespan both in vivo and in vitro. Cellular senescence occurs as reflection of organism aging and in response to internal and external stress signals.

Images



Senescence-associated beta-galactosidase staining at the 2nd(A), 4th(B), 8th(C), 12th passage *in vitro* expansion. Cells were plated at a density of 10,000 cells/cm² for 24h before staining. Five representative images (100x) were taken from diverse areas of cell culture, using phase-contrast microscopy to assess the number of positive cells.

Image obtained from Angelucci S et al; Proteome Sci, 2010 Mar 26;8:18

Human Wharton's jelly cells - Senescence Detection

Kit (ab65351)

Angelucci S et al., Proteome Sci, 8, 18, 2010
 Reproduced under the Creative Commons license
<http://creativecommons.org/licenses/by/2.0>

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors