

beta Galactosidase Staining Kit ab102534

★★★★★ [1 Abreviews](#) [15 References](#) [1 图像](#)

概述

产品名称	beta Galactosidase Staining 试剂盒
样品类型	Adherent cells, Suspension cells
检测类型	Direct
产品概述	<p>The LacZ gene from <i>E. coli</i> is one of the most commonly used reporter genes for testing the efficiency of expression vector mediated gene transfer and for studying the regulation of promoters of genes. The LacZ gene encodes the enzyme β-galactosidase, which is very stable, resistant to proteolytic degradation, can utilize a variety of substrates and can be easily assayed <i>in situ</i>.</p> <p>Abcam's beta Galactosidase Detection Kit (ab102534) utilizes X-gal as the substrate for β-galactosidase.</p>

Visit our [FAQs page](#) for tips and troubleshooting.

说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K802 Beta-Galactosidase Staining Kit. K802-250 is the same size as the 250 test size of ab102534.</p> <p>This kit has been optimized for transfected cells. If you want to detect senescence, we recommend using Senescence Detection Kit (ab65351).</p>
----	---

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	250 tests
Fixative Solution III	1 x 125ml
Staining Solution II	1 x 125ml
100X Staining Supplement	1 x 1.5ml
X-Gal	1 x 150mg

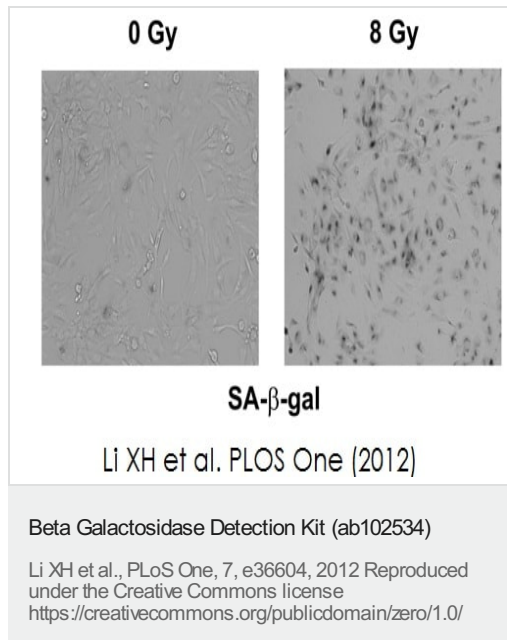
相关性 Beta galactosidase is a hydrolase enzyme that cleaves beta-linked terminal galactosyl residues

from gangliosides, glycoproteins, and glycosaminoglycans. Beta galactosidase is an essential enzyme in the human body. Deficiencies in the protein can result in galactosialidosis or Morquio B syndrome. Senescent cells display senescence-associated expression of beta galactosidase activity.

细胞定位

Isoform 1: Lysosome. Isoform 2: Cytoplasm, perinuclear region. Note=Localized to the perinuclear area of the cytoplasm but not to lysosomes.

图片



Gamma-radiation induced senescence in hFOB cells. The number of senescence-associated-β-galactosidase positive cells (*left panel*) clearly increases 72h after 8 Gy γ irradiation. Representative data was from one of a total of 3 independent experiments. *Image obtained from Li XH et al; PLOS One; 2012;7(5):e36604*

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.cn/abpromise> or contact our technical team.

Terms and conditions

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