

In-Cell ELISA (ICE) Support Pack ab111542

4 References [1 图像](#)

概述

产品名称	In-Cell ELISA (ICE) Support Pack
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based (quantitative)
实验步骤	Multiple steps standard assay
产品概述	<p>ab111542 is for use with suspension, apoptotic/detaching cells and adherent cell lines. The pack contains five 96-well microplates, buffers and protocol to perform ICE with Abcam's ICE-validated antibodies. For an ICE assay, it is necessary to purchase both primary antibody(ies) and labeled secondary antibody(ies). Antibodies are sold separately, allowing customizing the target(s) of interest, method of detection and multiplexing.</p>

In-Cell ELISA uses quantitative immunocytochemistry to measure protein levels or post-translational modifications within cells. The cells are fixed to the bottom of a coated 96-well plate (provided). Targets of interest are detected by primary antibodies, which are in turn quantified with labeled secondary antibody(ies). Abcam offers highly-specific, well-characterized primary antibodies, IRDye®- or HRP-labeled anti-mouse and anti-rabbit secondary antibodies, as well as IRDye®-labeled isotype specific anti-mouse antibodies. By combining antibodies of different species or isotype and appropriate IR-labeled secondary antibodies, two color multiplexing can be achieved in the 800/700 channels. IR imaging and quantification is performed using a LI-COR® Odyssey® or Aeries® system. HRP-labeled complexes are developed and quantified colorimetrically using spectrophotometer.

Two protocols are available when using this kit. Protocol (1) suspension cells and cells likely to detach under experimental conditions (for example, adherent cells undergoing apoptosis readily detach from a culture plate). Protocol (2) - A second protocol is available for normal adherent cells. These protocols can be used with any of Abcam's ICE-validated antibodies. Specific scientific information, background and working concentration for each antibody are detailed in each antibody's corresponding datasheet.

说明

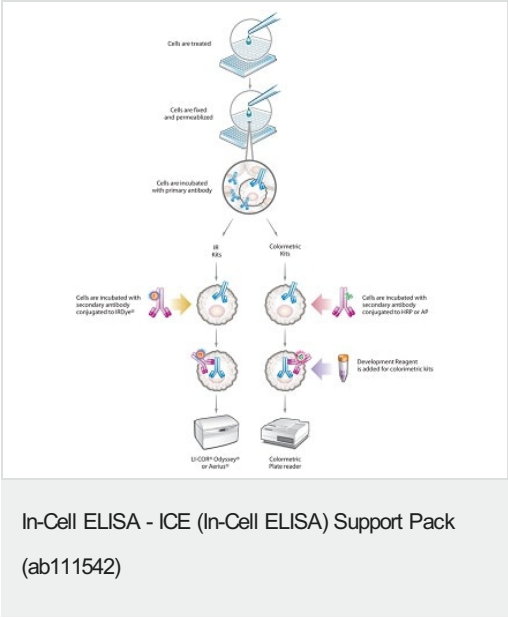
Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

性能

存放说明Store at +4°C. Please refer to protocols.

组件	5 x 96 tests
100X Triton X-100	1 x 1.25ml
10X Blocking Solution	1 x 40ml
10X Phosphate Buffered Saline (PBS)	1 x 250ml
400X Tween-20	1 x 4ml
96-well Assay Plate	5 units
Janus Green Stain	1 x 30ml
Plate Seals	5 units

图片



Cells are grown to ~80% confluency in a 96- or 384-well plate, a drug/other treatment is applied to stimulate a cellular response. The cells are then fixed and permeabilized, effectively "freezing" them. Primary antibodies are then added which bind to their intended targets within the mitochondria or other subcellular compartment. After incubation, the unbound primary antibodies are washed away and secondary antibodies are added. These secondaries are conjugated to either IRDyes® or to an enzyme label (HRP or AP) for the colorimetric versions of the assays. Unbound secondaries are washed away, reaction buffer is added for the colorimetric assays, and the signal is read on a suitable instrument for the kit type.

» [In-cell ELISA diagram](#) in PDF format

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