

Phosphotyrosine In-Cell ELISA Kit ab126431

1 图像

概述

产品名称	Phosphotyrosine In-Cell ELISA试剂盒
检测方法	Colorimetric
样品类型	Adherent cells
检测类型	Cell-based (qualitative)
检测时间	5h 10m
实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse, Rat, Human
产品概述	<p>ab126431 is a very rapid, convenient and sensitive assay kit that can monitor the activation or function of important biological pathways in cells. It can be used for measuring the relative amount of Phosphotyrosine and screening the effects of various treatments, inhibitors (such as siRNA or chemicals), or activators in cultured human, mouse and rat cell lines. By determining Phosphotyrosine protein in your experimental model system, you can verify pathway activation in your cell lines without spending excess time and effort in preparing cell lysate and performing an analysis of Western Blot.</p> <p>In the Cell-Based Phosphotyrosine ELISA kit, cells are seeded into a 96 well tissue culture plate. The cells are fixed after various treatments, inhibitors or activators. After blocking, HRP-Anti-Phosphotyrosine is pipetted into the wells and incubated. The wells are washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of protein. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
平台	Microplate

性能

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

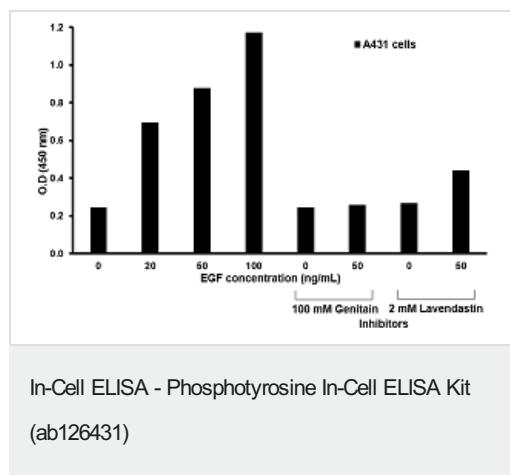
组件	1 x 96 tests
HRP Conjugated Anti-Phosphotyrosine Concentrate (2000X)	1 x 7µl
Blocking Buffer Concentrate (5X)	1 x 20ml

组件	1 x 96 tests
Fixing Solution	1 x 30ml
Quenching Buffer Concentrate (30x)	1 x 2ml
Stop Solution	1 x 14ml
TMB One-Step Substrate Reagent	1 x 12ml
Uncoated 96-well Microplate	1 unit
Wash Buffer A Concentrate (20X)	1 x 30ml
Wash Buffer B Concentrate (20X)	1 x 30ml

相关性

The phosphorylation of specific tyrosine residues has been shown to be a primary mechanism of signal transduction during normal mitogenesis, cell cycle progression and oncogenic transformation, its role in other areas such as differentiation and gap junction communication, is a matter of active and ongoing research. Antibodies that specifically recognize phosphorylated tyrosine residues have proved to be invaluable to the study of tyrosine phosphorylated proteins and the biochemical pathways in which they function.

图片



A431 cells were treated for 30 min with 50 μ L of 100 mM Genistein or 2 mM Lavendustin in appropriate wells at room temperature prior to EGF stimulation. Added 50 μ L different concentrations of rhEGF (0, 20 or 100 ng/ml in serum free DMEM) to appropriate wells. Then incubated for 10 min at 37°C.

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