

画像数 3

醫藥用外劇物

| | |
|------------|---|
| 製品名 | Human STAT5 (Tyr694) In-Cell ELISA Kit |
| 検出方法 | Colorimetric |
| サンプルの種類 | Adherent cells |
| アッセイタイプ | Semi-quantitative |
| 全工程の試験時間 | 5h 10m |
| ステップ | Multiple steps standard assay |
| 種交差性 | 交差種: Human |
| 製品の概要 | <p>ab126429 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 STAT5 (Tyr694) phosphorylation and screening the effects of various treatments, inhibitors (such as siRNA or chemicals), or activators in cultured human cell lines. By determining STAT5 protein phosphorylation 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 In-Cell STAT5 (Tyr694) ELISA kit, cells are seeded into a 96 well tissue culture plate. The cells are fixed after various treatments, inhibitors or activators. After blocking, Anti Phospho-STAT5 (Tyr694) or Anti-STAT5 (primary antibody) is pipetted into the wells and incubated. The wells are washed, and HRP-conjugated anti-mouse IgG (secondary antibody) is added to the wells. The wells are washed again, 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> |
| 試験プラットフォーム | Microplate |

保存方法 Store at -20°C. Please refer to protocols.

| 内容 | 1 x 96 tests |
|---|--------------|
| HRP-conjugated Anti-Mouse IgG Concentrate | 1 x 10µl |

| 内容 | 1 x 96 tests |
|--|--------------|
| Blocking Buffer Concentrate (5X) | 1 x 20ml |
| Fixing Solution | 1 x 30ml |
| Uncoated 96-well Microplate | 1 unit |
| Mouse Anti-Phospho-STAT5 (Tyr694) Concentrate (Item G) | 1 x 7μl |
| Mouse Anti-STAT5 Concentrate (Item H) | 1 x 10μl |
| Quenching Buffer Concentrate (30x) | 1 x 2ml |
| Stop Solution | 1 x 14ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |
| Wash Buffer A Concentrate (20X) | 1 x 30ml |
| Wash Buffer B Concentrate (20X) | 1 x 30ml |

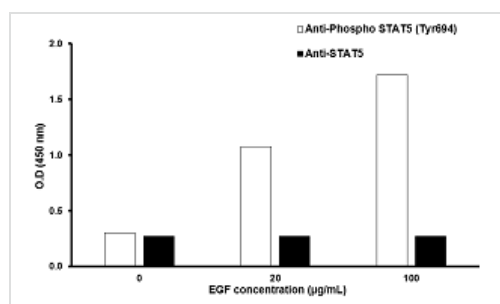
機能 Carries out a dual function: signal transduction and activation of transcription. Binds to the GAS element and activates PRL-induced transcription.

配列類似性 Belongs to the transcription factor STAT family.
Contains 1 SH2 domain.

翻訳後修飾 Tyrosine phosphorylated in response to IL-2, IL-3, IL-7, IL-15, GM-CSF, growth hormone, prolactin, erythropoietin and thrombopoietin. Tyrosine phosphorylation is required for DNA-binding activity and dimerization. Serine phosphorylation is also required for maximal transcriptional activity.

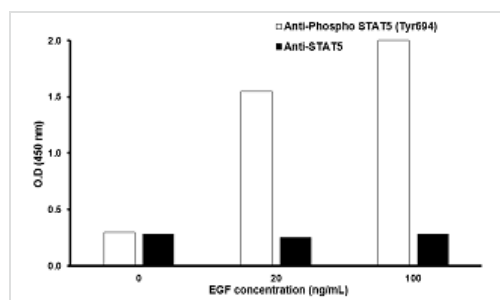
細胞内局在 Cytoplasm. Nucleus. Translocated into the nucleus in response to phosphorylation.

画像



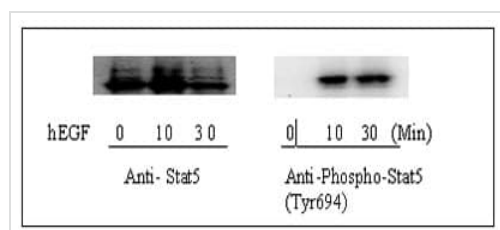
In-Cell ELISA - STAT5 (Tyr694) Human In-Cell ELISA Kit (ab126429)

A431 cells were stimulated by different concentrations of EGF for 10 minutes at 37°C.



In-Cell ELISA - STAT5 (Tyr694) Human In-Cell ELISA Kit (ab126429)

A431 cells were stimulated by different concentrations of EGF for 30 minutes at 37°C.



Western blot - STAT5 (Tyr694) Human In-Cell ELISA Kit (ab126429)

Western blot analysis of extracts from 100 ng/ml hEGF treated A431 cells. Phospho-STAT5 (Tyr694) and STAT5 antibodies were used in both detection assays.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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