

Rapid ELISA-Based Measurement of Protein Phosphorylation Using RayBio® Phosphorylation ELISA Kits

Valerie Sloane Jones, Ph.D.*¹, Ruochun Huang, Ph.D.¹, Ruo-Pan Huang, M.D., Ph.D.^{1,2}

¹RayBiotech, Inc., 3607 Parkway Lane, Suite 100, Norcross, GA 30092, Phone: 770-729-2992; ²RayBiotech, Inc., Guangzhou 510600 China.
*To whom correspondence should be addressed.

Abstract

RayBio® Phosphorylation ELISA is a rapid, convenient and sensitive sandwich ELISA for the in vitro measurement of key signaling pathway phosphoproteins in cell or tissue extracts. Over 20 different kits are available for well-studied pathway molecules such as EGFR, Akt, Erk1/2, p38 alpha, Mek1, STAT1, STAT3, and Met. RayBio Phosphorylation ELISA also features site-specific phospho-antibodies which can detect a single phosphorylated residue (for example, Tyr1086 of EGFR). This method is an improvement on traditional labor-intensive immunoblotting protocols that require 2-3 days of processing time.

Immunoblotting is frequently used for immunodetection of phosphoproteins, in some cases following immunoprecipitation of the cellular extract. The signals from each probe can then be compared and phospho-specific fractions visually assessed or quantified by densitometry. However, the time and effort required in these protocols has driven many researchers to seek out ELISA-based methods, which offer a faster, higher throughput alternative.

In this article, we present a study in which the phosphorylation of two proteins, STAT1 (Tyr701) and Mek1 (Ser217/221), is detected in response to various agonists. Data from the RayBio Phosphorylation ELISA kits will be compared to traditional immunoblotting.

Background

The detection of the phosphorylation of a specific protein is usually accomplished by immunological methods, in which the researcher can measure the fraction of phosphorylated target protein relative to total levels of expressed target protein in the cell. By comparing the phosphorylated fractions in cellular extracts under various experimental conditions, one can identify to what degree and what condition activates or blocks an important signaling pathway. This approach requires the use of a phospho-specific antibody in parallel with a "loading control" to assess relative levels of total cellular protein. Typically, the loading control may be incorporated by using either the corresponding "pan" antibody (recognizing the target protein in both its phosphorylated and non-phosphorylated states), or alternatively an antibody against a "housekeeping" protein such as β -actin or GAPDH. The signals from each antibody can then be compared between samples. This methodology has largely replaced the classical, laborious techniques of directly radiolabeling phosphoproteins with ³²P-orthophosphate, followed by SDS-PAGE gel separation and subsequent autoradiographic analysis.

Assay Format, Procedure and Application

RayBio Phosphorylation ELISA kits contain antibodies for both phosphorylated and total protein (loading control) detection in parallel. Each kit includes a positive control, which is a cell lysate specially prepared to produce a strong phospho-positive signal. The positive control is intended to be used as a reference sample rather than an actual calibration standard, and its inclusion in the assay protocol is optional.

In a single microplate, up to 24 different samples may be simultaneously tested in duplicate (including positive controls, if used), producing both phosphorylated and total protein signal outputs. The protocol involves pipetting fresh or thawed lysates into an antibody-coated 96-well microplate, followed by washes and subsequent incubation with anti-phosphorylated antibody or anti-pan antibody. After further washing, an HRP-conjugated secondary antibody is added. Colorimetric detection of captured protein is carried out using TMB substrate solution where blue color intensity develops in proportion to the amount of phosphorylated protein or pan protein bound.

Some researchers may be interested in screening for only phosphorylated protein signals rather than total and phosphorylated in parallel. For this application, each Phosphorylation ELISA kit is also available in “phospho-only” format, containing only a phospho-specific antibody pair. This format allows the researcher to screen more conditions (48 duplicates per plate) for a lower cost compared to the “phospho plus total” readout format.

The assay procedure has minimal hands-on time and hundreds of samples can be assayed by a single technician in one day. Thus, the RayBio Phosphorylation ELISA represents a more rapid and efficient alternative to multi-step phospho-protein detection by immunoblot analysis, which requires a minimum of two days processing time. Below, we present two examples of analyses using the Phosphorylation ELISA kits.

Site-Specific STAT1 Activation

STAT is a transcription factor implicated in programming gene expression for biological events as diverse as embryonic development, programmed cell death, organogenesis, innate immunity, adaptive immunity and cell growth regulation¹. In the study presented here, cultured A431 cells (an epidermoid carcinoma cell line) were treated with recombinant human epidermal growth factor (EGF) to induce EGFR-mediated signaling cascades and JAK-STAT signaling. Cell lysates were analyzed by both RayBio Phospho-STAT1 (Tyr701) ELISA (Figure 1A) and traditional immunoblot (Figure 1B). After EGF treatment, a robust and rapid phosphorylation of STAT1 at Tyr701 was observed in parallel with total STAT1 levels by both detection methods (Figure 1). We note that immunoblots were incubated with anti-phospho-STAT1 (Tyr701) and anti-pan STAT1 (the same antibodies used in the RayBio Phosphorylation ELISA kit).

To compare the sensitivity of the ELISA kit to traditional Western blot analysis, EGF-treated A431 cell lysates were serially diluted and assayed by each approach. Both methods exhibited similar sensitivity, with lowest detectable limits well below 1 µg/ml of total lysate protein (Figure 2), concluding that the ELISA kit is equally sensitive as traditional immunoblot, but can be completed inside 6 hours.

Site-Specific Mek1 Activation

Mek1 is a protein kinase that functions in the mitogen-activated protein kinase (MAPK) signaling cascade, and is involved in the regulation of normal cell proliferation, survival and differentiation². In this study, HeLa cells were treated with tetradecanoylphorbol-13-acetate (TPA), a potent activator of protein kinase-C (PKC) pathways. Cell lysates were analyzed using both RayBio[®] Phospho-Mek1 (Ser217/221) ELISA, and by immunoblot. After TPA treatment, a robust and rapid phosphorylation of Mek1 was detected in parallel with total Mek1 levels by both methods, concluding that the Phospho-Mek1 (Ser217/221) ELISA has similar sensitivities to the immunoblot method (Figure 3). Immunoblots were incubated with anti-phospho-Mek1 (Ser217/221) or anti-pan Mek1 (the same antibodies used in the RayBio Phosphorylation ELISA kit).

Summary

RayBio[®] Phosphorylation ELISA kits are a rapid, sensitive, high-throughput, cost-effective and convenient method for monitoring the activation of important biological pathways in cells or tissues. Unlike immunoblots, which require several days processing to obtain parallel signals for phosphorylated and total target protein, the Raybiotech's ELISA method achieves comparable sensitivity within 6 hours processing time. Up to 24 different experimental conditions can be simultaneously screened in duplicate on one plate with phosphorylated and total (pan) protein readouts assayed. In this format, several hundred samples can be assayed by a single technician in one day.

References

1. Horvath, CM. STAT proteins and transcriptional responses to extracellular signals. *Trends Biochem Sci* 25(10), 496-502 (2000).
2. Roberts PJ, Der CJ. Targeting the Raf-MEK-ERK mitogen-activated protein kinase cascade for the treatment of cancer. *Oncogene* 26, 3291–3310 (2007).

To Order or Receive Technical Assistance

In the U.S. and Canada, call 1-888-494-8555 or contact us at info@raybiotech.com.

In other countries worldwide, please visit: <http://www.raybiotech.com/distributors.html>

Connect With Us

Join RayBiotech on Facebook and Twitter for the latest updates, news, products, innovations and promotions!



<https://www.facebook.com/pages/RayBiotech/149958741713501?v=wall>



<https://twitter.com/RayBiotech>

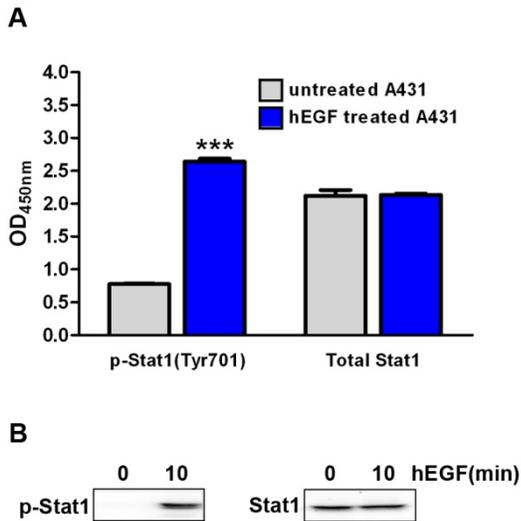


Figure 1: Detection of relative Phospho-STAT1 levels in A431 Cells following stimulation with Epidermal Growth Factor (EGF). A431 cells were treated with 100 ng/ml recombinant human EGF for 10 minutes. (A) Analysis of cell lysates using Human/Mouse Phospho-STAT1 (Y701) and Total STAT1 ELISA (cat # PEL-STAT1-Y701-002). (B) Immunoblot analysis of lysates using antibodies against phospho-STAT1 and STAT1 (the same antibodies used in the ELISA kit). ***=P<0.001, n=3.

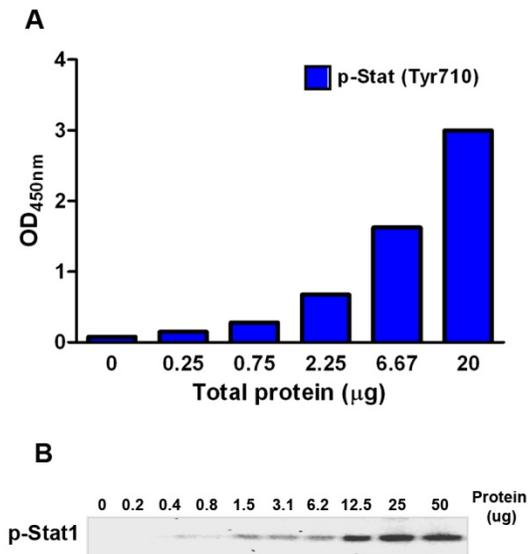


Figure 2: Comparative sensitivity of the detection of Phospho-STAT1 levels in A431 Cells following stimulation with Epidermal Growth Factor (EGF) by RayBio Phospho-STAT1 (Tyr701) ELISA and by traditional immunoblot techniques. A431 cells were treated with 100 ng/mL recombinant human EGF for 20 minutes to induce phosphorylation of STAT1. Serial dilutions of lysates were analyzed by (A) RayBio Human/Mouse Phospho-STAT1 (Y701) and Total STAT1 ELISA (cat # PEL-STAT1-Y701-002) and (B) by immunoblot. Immunoblots were incubated with anti-phospho-STAT1 (the same antibody used in the ELISA kit).

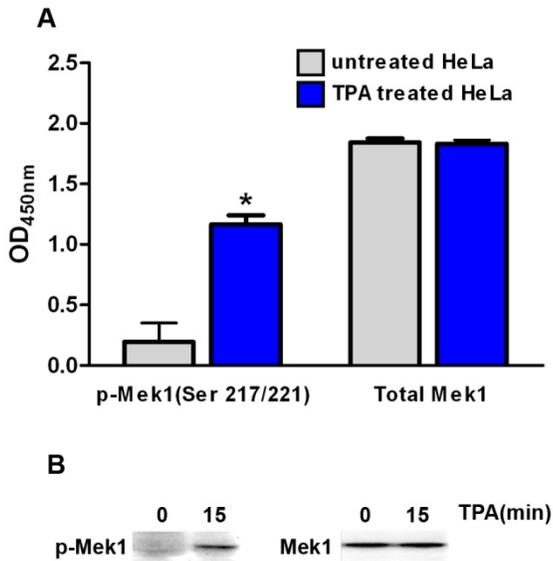


Figure 3: Detection of relative Phospho-Mek levels in HeLa Cells following stimulation with TPA. HeLa cells were treated with TPA for 15 minutes. Cell lysates were analyzed using (A) Human/Mouse/Rat Phospho-Mek (S217/S221) & Total Mek ELISA (cat # PEL-MEK-S217-002) and (B) by immunoblot. Immunoblots were incubated with anti-Mek1 or anti-phospho-Mek1 (the same antibodies used in the ELISA kit). *= $P < 0.05$, $n = 3$.