

Development and Validation of a Novel EGFR Receptor-Neutralizing (D1D4J) Monoclonal Antibody

Krystyna Zuberek Hincman, Christopher A. Manning, Michael Nelson, Michael Lewis,

Roy Scialdone, Jaime Darce, Stephen Lutz, Evans John Burford, Herbert Haack, Matthew Silver

Cell Signaling Technology, Inc., 3 Trask Lane, Danvers MA 01923

Abstract Number: 4595

INTRODUCTION

- The epidermal growth factor receptor (EGFR) is a cell-surface receptor for members of the EGF family. EGFR dimerizes upon ligand binding, which leads to autophosphorylation, downstream signaling and ultimately receptor internalization.
- In cancer, EGFR mutations and amplification have been linked to cell proliferation and survival.
- There are two FDA approved neutralizing monoclonal antibodies: cetuximab (BMS; mouse/human chimeric) and panitumumab (Amgen; fully humanized). However, these antibodies are often unavailable to researchers studying this important pathway.
- Various immunization strategies have been attempted at Cell Signaling Technology, Inc (CST) previously to generate an extracellular EGFR Rabbit mAb with limited success. A specifically designed whole cell immunization protocol described lead to suitable monoclonal candidates.
- Here we describe the development and validation of a new rabbit monoclonal EGFR-neutralizing antibody (clone D1D4J).

CONCLUSIONS

- D1D4J is a highly specific EGFR monoclonal antibody that recognizes the extracellular domain of human and mouse EGFR, and is the only commercially available extracellular Rabbit mAb on the market.
- D1D4J demonstrated strong and specific performance in Western blot (WB), immunofluorescence (IF) (live and fixed), flow cytometry (Flow), and immunoprecipitation (IP) assays.
- D1D4J also shows strong neutralizing ability as determined by multiple assays (WB and IF).
- D1D4J is a useful research tool to support further preclinical development of EGFR targeted therapies.
- EGF Receptor (D1D4J) XP[®] Rabbit mAb (Neutralizing) #54359 was validated using cell-based models and IF imaging. D1D4J binds EGFR and neutralizes EGF-induced activation and internalization. Compared to other Research Use Only (RUO) neutralizing antibodies on the market, D1D4J shows superior performance and is a useful research tool to study EGFR signaling.

Methods

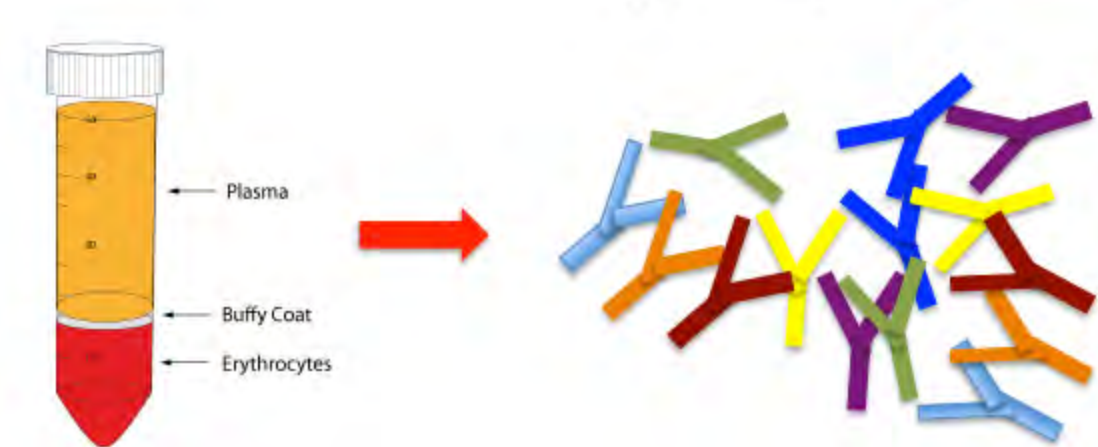
Figure 1A: Whole Cell Immunization

Rabbits were immunized with mammalian cells expressing full length human EGFR protein



Figure 1B: Serum Analysis and Rabbit Selection

Rabbit serum tested for EGFR reactivity by flow cytometry



Cell Systems Used for Analysis	Over-expression or Endogenous	EGFR Expression Level
293 (Mock)	transfected mock	-
293 O/E hEGF Receptor	transfected	++++
A549	endogenous	+
A-431	endogenous	++
Jurkat	endogenous	-
CAMA-1	endogenous	-

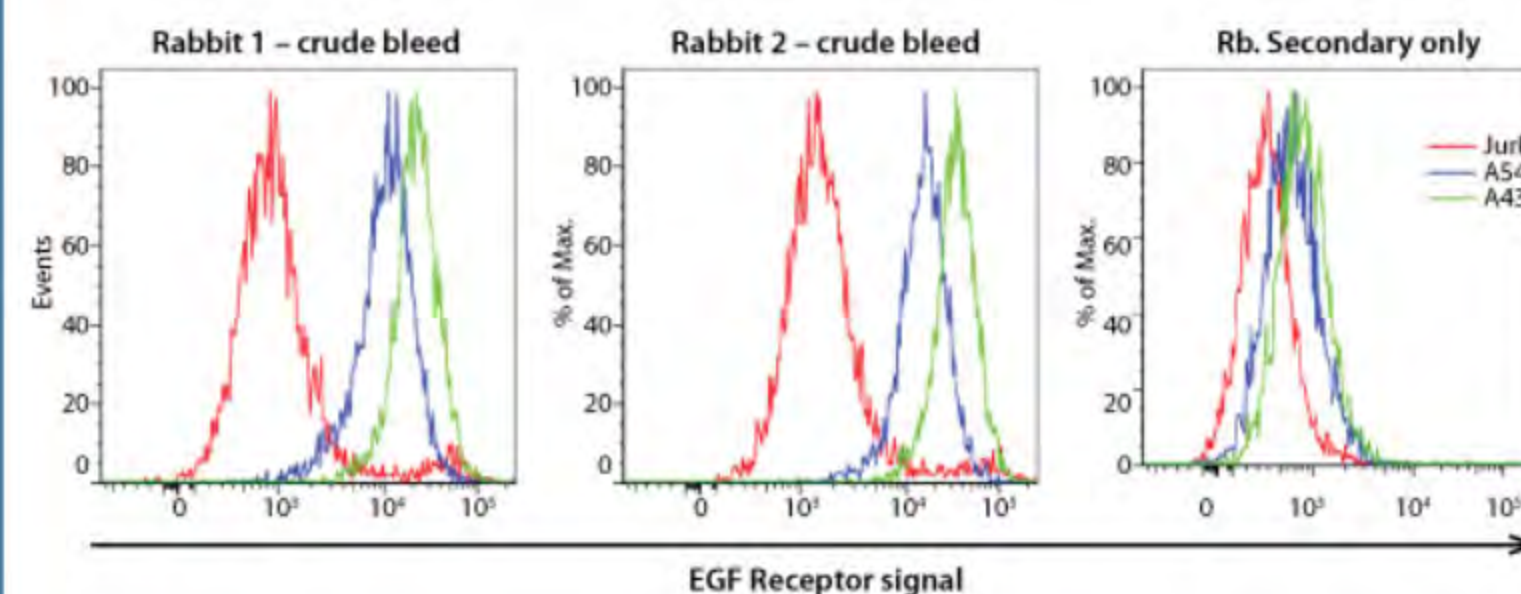
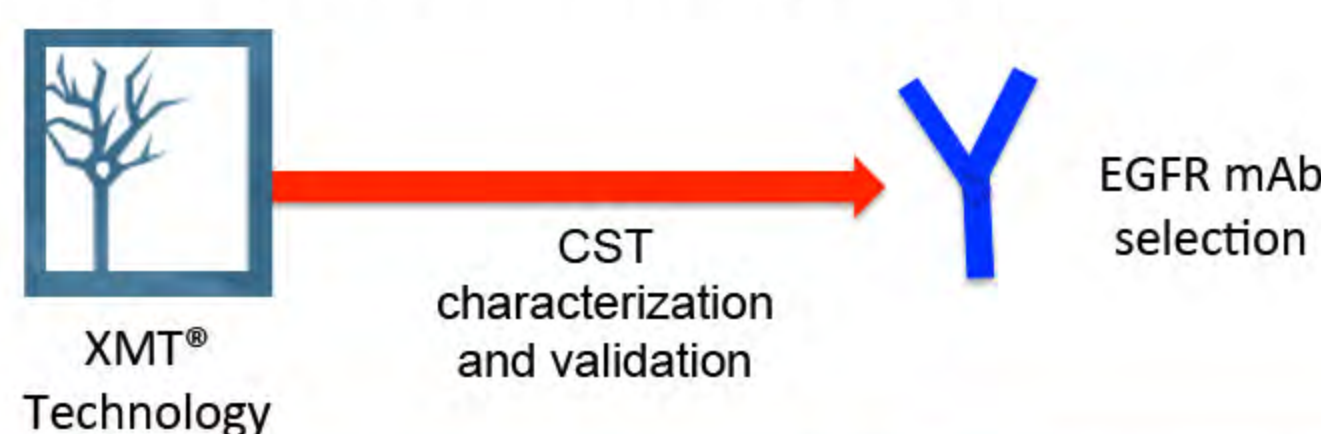


Figure 1C: Monoclonal Antibody Creation



Assessing Specificity

D1D4J was tested according to validation procedures at CST. D1D4J was validated and optimized for WB, IF, FC, and IP.

Figure 2A: D1D4J detects EGFR in human and mouse by WB and does not cross react with other ErbB/HER family members

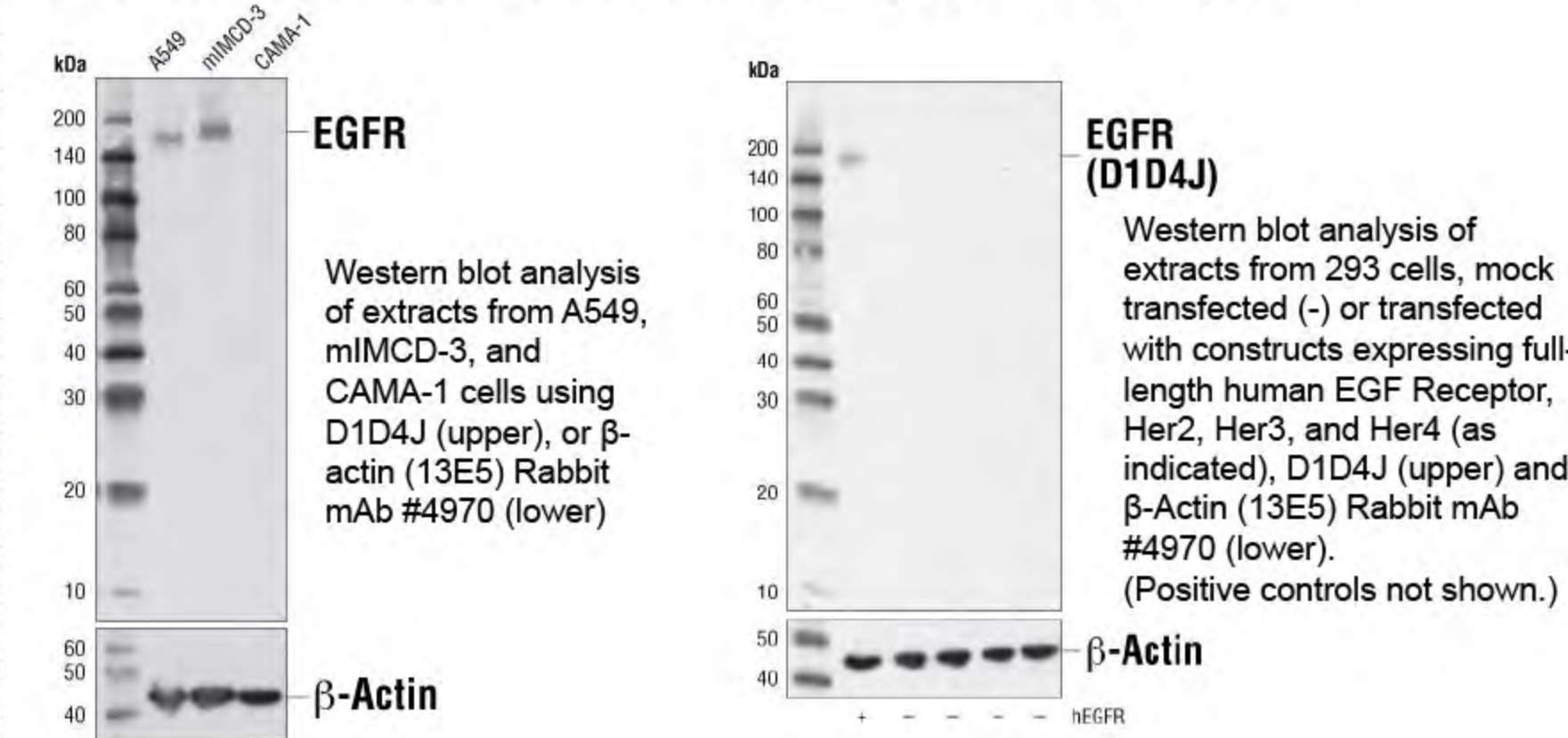
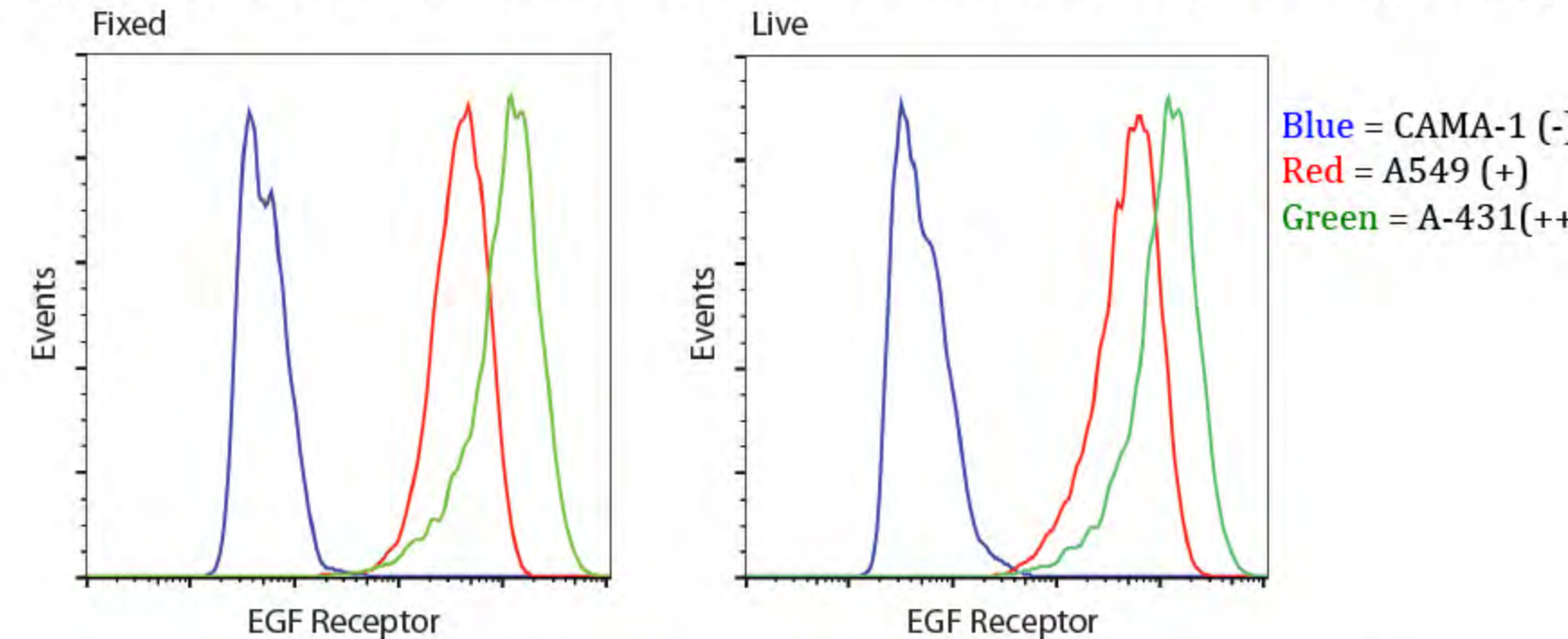
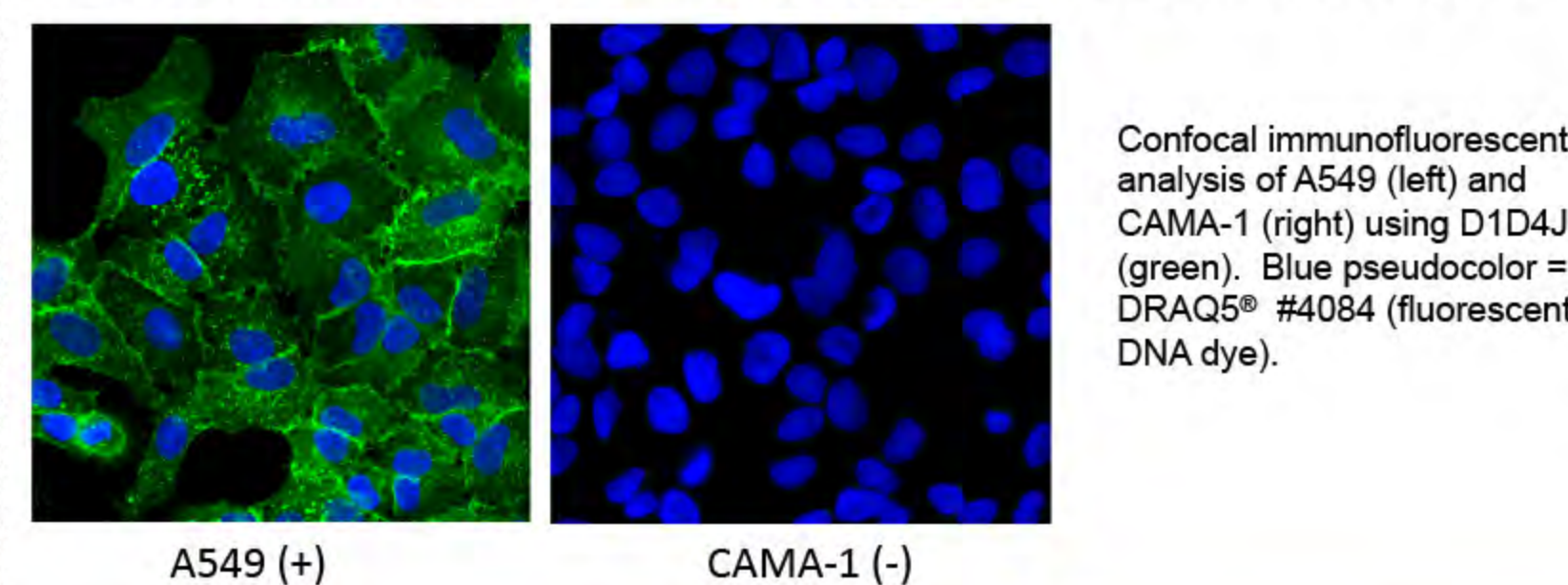


Figure 2B: D1D4J detects EGFR in fixed and live cells by Flow Cytometry

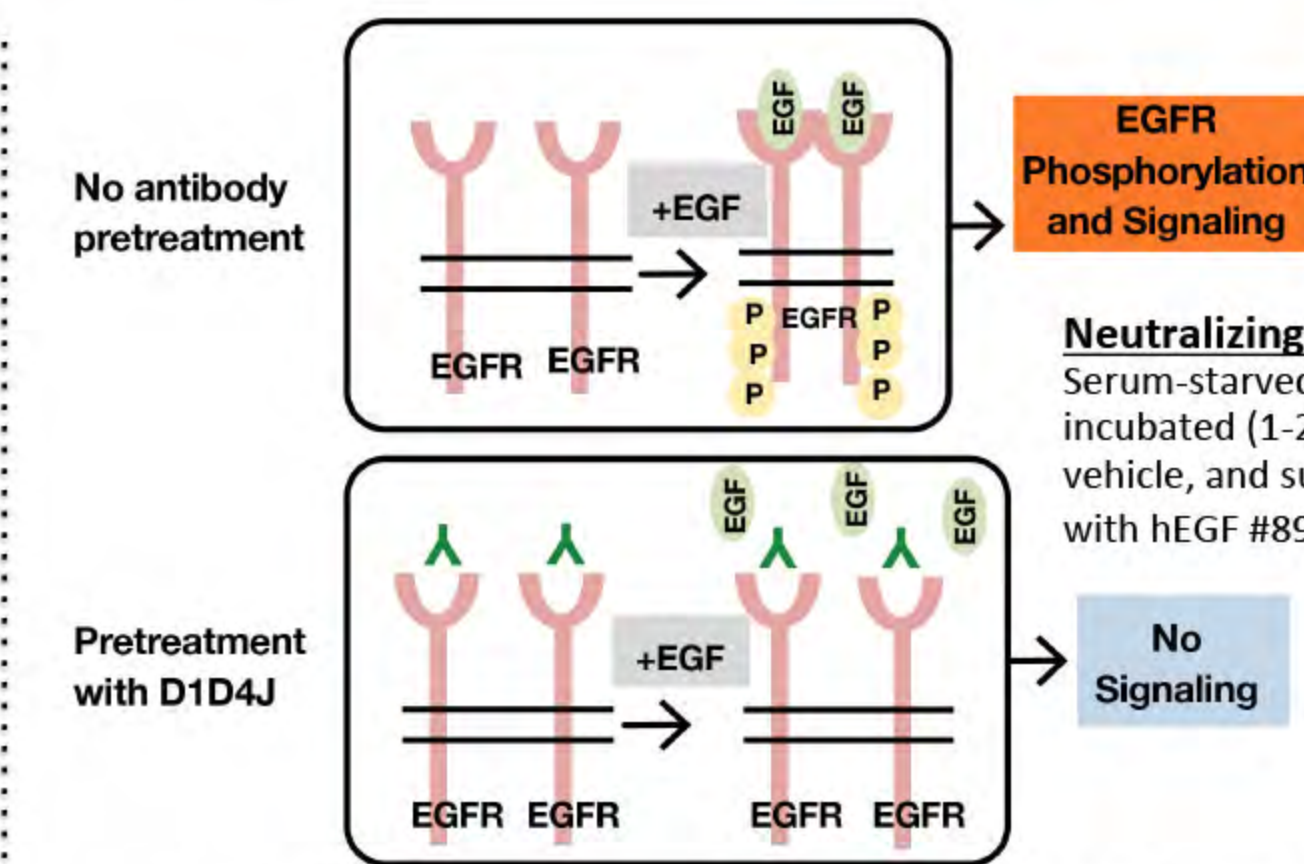


Flow cytometric analysis of fixed, unpermeabilized (left) or live (right) CAMA-1 cells (blue), A549 cells (red), and A-431 cells (green) using D1D4J. Anti-rabbit IgG (H+L), F(ab')₂ Fragment (Alexa Fluor[®] 488 Conjugate) #14412 was used as a secondary antibody.

Figure 2C: D1D4J detects EGFR by Immunofluorescent analysis

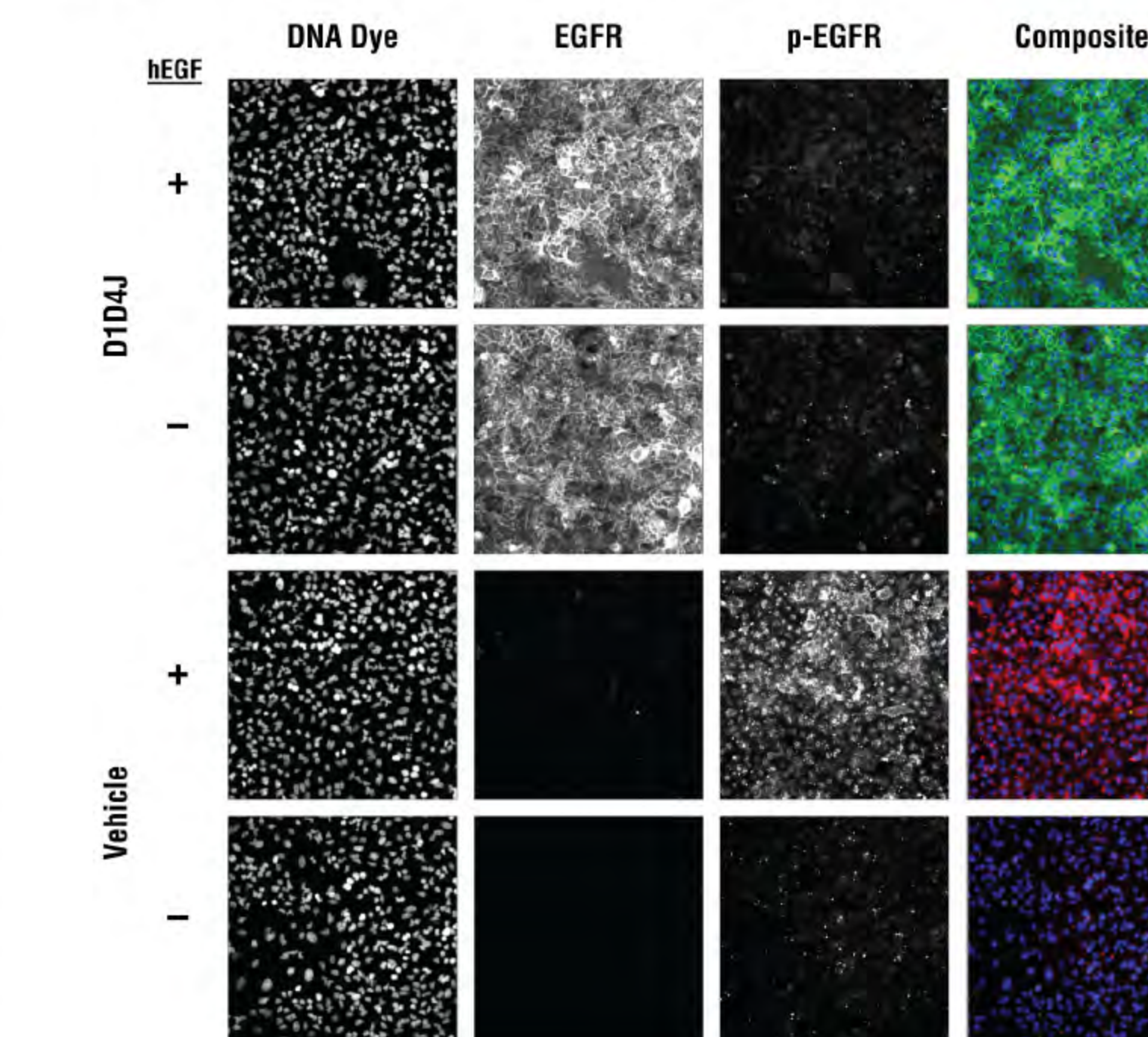


Assessing Neutralizing Activity

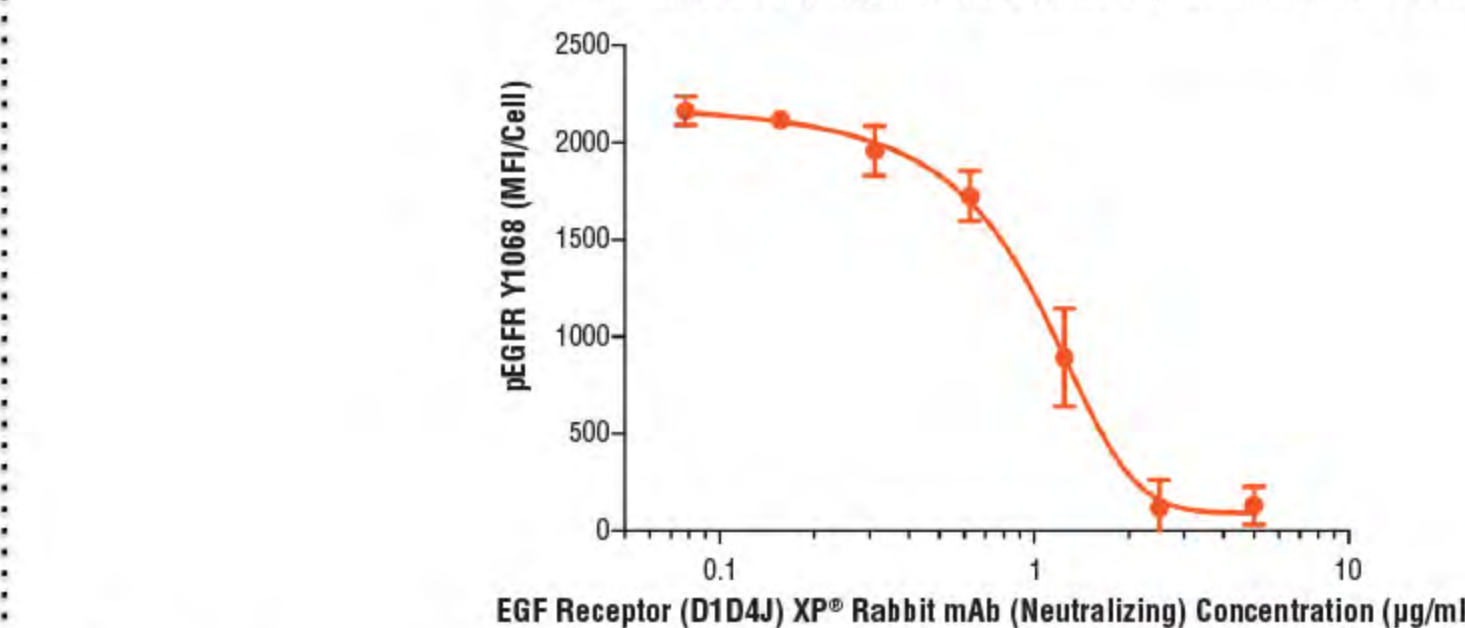


Neutralizing Treatment Method:
Serum-starved A549 cells were pre-incubated (1-2hr, 5ug/ml) with D1D4J, or vehicle, and subsequently stimulated with hEGF #8916 (15min, 100ng/ml; +).

Figure 3A: D1D4J neutralizing activity confirmed by IF analysis

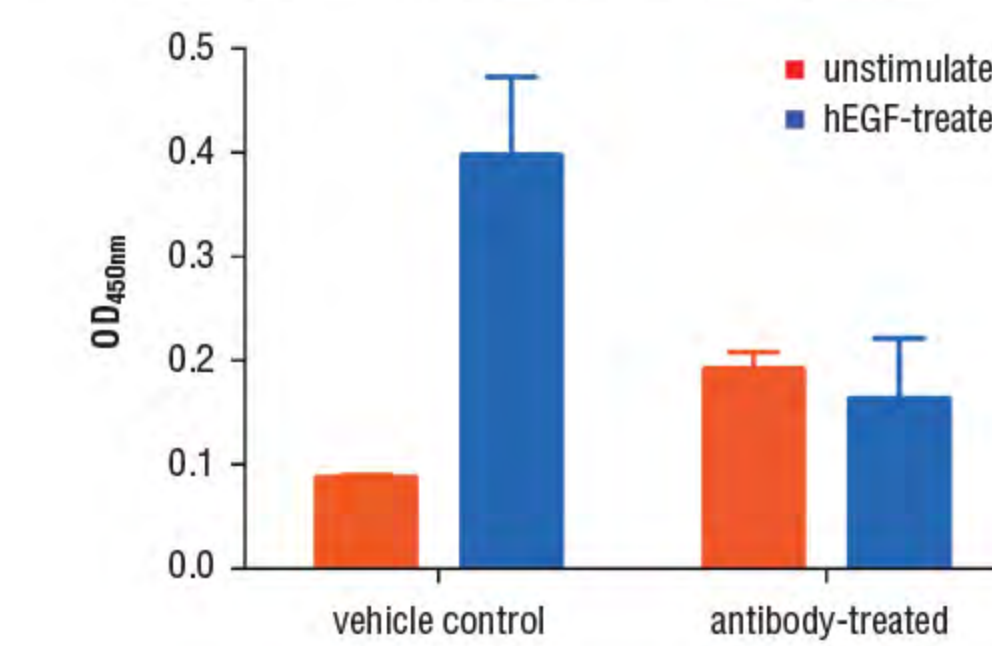


Inhibition of EGFR phosphorylation by antibody treatment in A549 cells



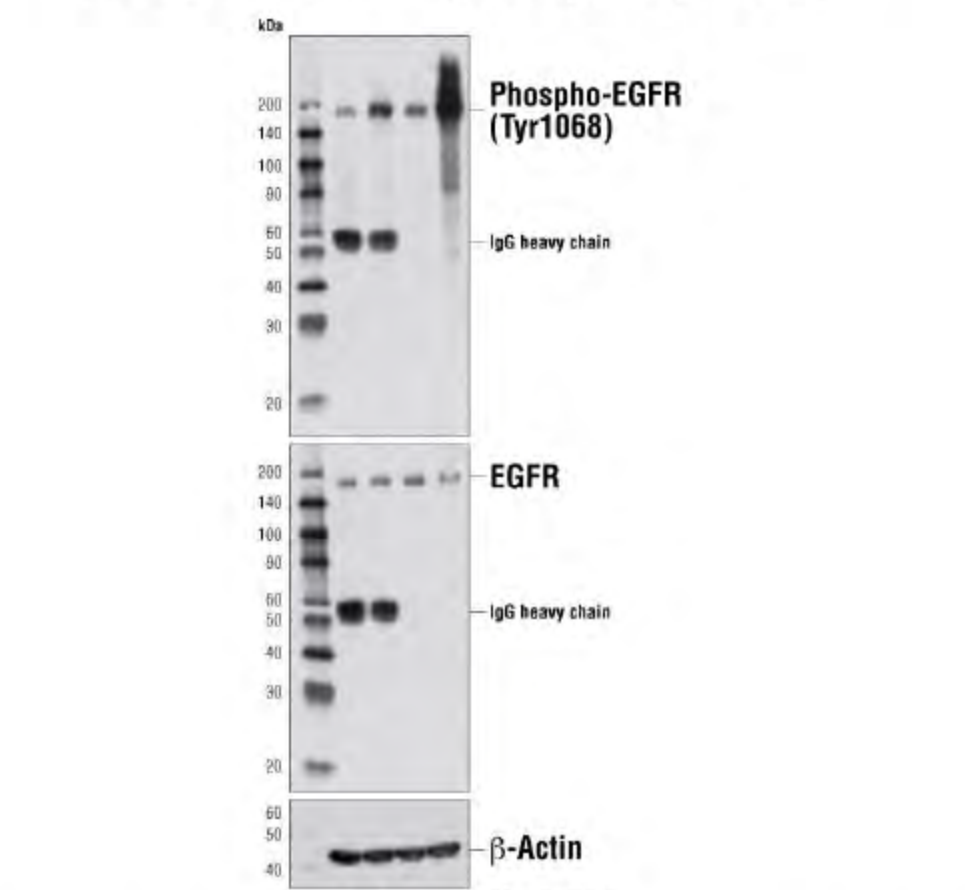
IF images (upper) of treated A459 cells pre-incubated with D1D4J (5ug/mL, 1 hr) or vehicle (PBS); and subsequently stimulated with hEGF #8916 0.....10 (15 min, 100 ng/ml) as shown. Cells were then fixed and stained for activated EGFR (p-EGFR Tyr1068 in red); D1D4J was also labeled post-fixation (in green). Blue pseudocolor = Hoechst 33342 (fluorescent DNA dye). Phospho-EGFR signal was quantified for the full titration range of D1D4J treatment using an ArrayScan[™] VTI HCS Reader (bottom).

Figure 3B: D1D4J neutralizing activity confirmed by PathScan[®] ELISA analysis



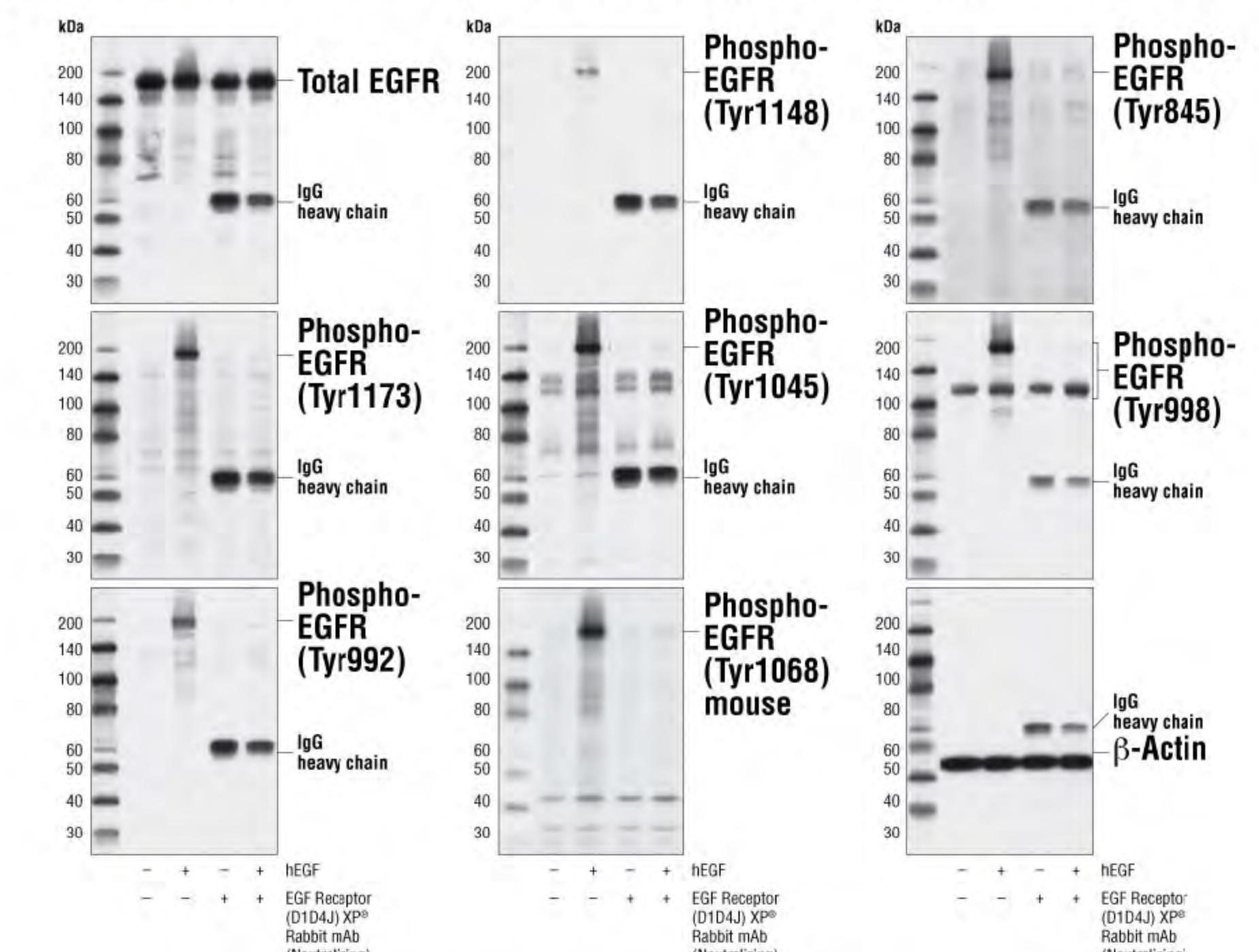
ELISA analysis of cell extracts from treated or control A549 cells. The extracts were subsequently left unstimulated (red) or stimulated with hEGF #8916 (15 min, 100 ng/ml; blue), using PathScan[®] Phospho-EGF Receptor (panTyr) Sandwich ELISA Kit #7911.

Figure 3C: D1D4J pre-treatment reduces EGFR phosphorylation



WB analysis of extracts from treated (+) or control (-) A549 cells using p-EGFR (Tyr1068) (D7A5) #3777 (upper), D1D4J (middle), or beta-actin (13E5) Rabbit mAb #4970 (lower).

Figure 3D: D1D4J neutralizes EGFR phosphorylation at numerous tyrosine residues



WB analysis of extracts from treated (+) or control (-) A549 cells using: Total EGFR (D38B1) #4267 (left, upper), p-EGFR (Tyr1173) #4407 (left, middle), p-EGFR (Tyr992) #2235 (left, lower), p-EGFR (Tyr1148) #4404 (center, upper), p-EGFR (Tyr1045) #2237 (center, middle), p-EGFR (Tyr1068) (1H12) Mouse mAb #2236 (center, lower), p-EGFR (Tyr845) (D63B4) #6963 (right, upper), p-EGFR (Tyr998) (C24A5) #2641 (right, middle), beta-actin (13E5) #4970 (right, lower).