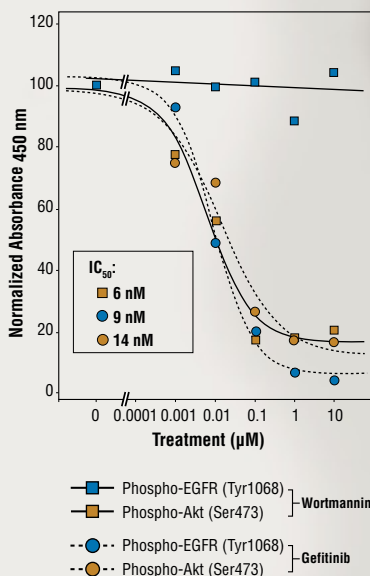


### Analysis of intracellular EGFR and Akt Signaling using PathScan® Sandwich ELISA Kits



# PathScan® Sandwich ELISA Products

from Cell Signaling Technology

## Unparalleled product quality, validation, and technical support

Our line of PathScan® Sandwich ELISA products provides researchers with a wide selection of assays to measure endogenous levels of critical regulatory proteins from cell lysates. In-house development, production, and validation ensure the highest possible product quality and technical support. Custom ELISA products are available upon request.

- **Over 170 Modification-state and Total Protein Sandwich ELISA Kits** available to cover a broad spectrum of signaling pathways.
- **Chemiluminescent Sandwich ELISA Kits** provide the broadest dynamic range and assay sensitivity while requiring half the sample size.
- **Antibody Pairs** provide an economical alternative to our complete PathScan® ELISA Kits.
- **Control Cell Extracts** provide the appropriate positive and negative controls, and also allow the standardization of signal obtained from different plates.

**Analysis of intracellular EGFR and Akt Signaling using PathScan® Sandwich ELISA Kits:** A-431 cells were exposed to varying concentrations of gefitinib (2 hr) or wortmannin (1 hr), then treated with 100 ng/ml EGF (5 min). With increasing concentrations of gefitinib, a significant decrease (~25 fold and ~6.25 fold, respectively) in phospho-EGFR and phospho-Akt1 signals were detected using the PathScan® Phospho-EGFR (Tyr1068) Sandwich ELISA Kit #7240 and the PathScan® Phospho-Akt1 (Ser473) Sandwich ELISA Kit #7160. With increasing concentrations of wortmannin, a significant decrease (~5 fold) in phospho-Akt1 signal was detected. Wortmannin had no effect on phospho-EGFR levels.

for quality products you can trust...

[www.cellsignal.com](http://www.cellsignal.com)

 **Cell Signaling**  
TECHNOLOGY®