

# MSD® EGFR Family Whole Cell Lysate Kit

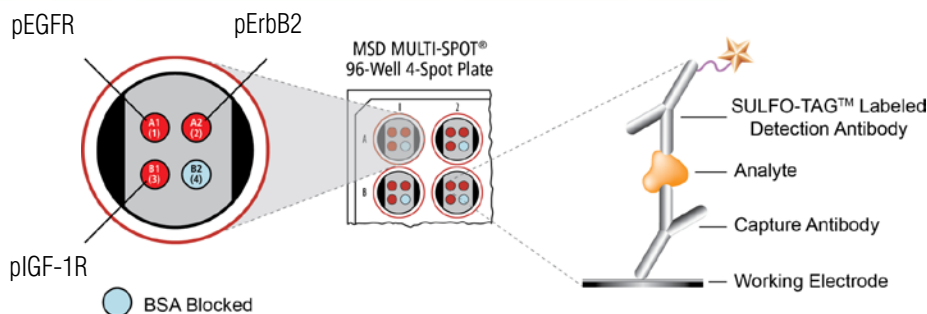
For quantitative determination of phospho-EGFR, phospho-ErbB2, and phospho-IGF-1R in human whole cell lysate samples



Alzheimer's Disease  
BioProcess  
Cardiac  
**Cell Signaling**  
Clinical Immunology  
Cytokines  
Hypoxia  
Immunogenicity  
Inflammation  
Metabolic  
Oncology  
Toxicology  
Vascular

## Catalog Numbers

| EGFR Family Whole Cell Lysate Kit |           |
|-----------------------------------|-----------|
| Kit size                          |           |
| 1 plate                           | K15106D-1 |
| 5 plates                          | K15106D-2 |
| 20 plates                         | K15106D-3 |

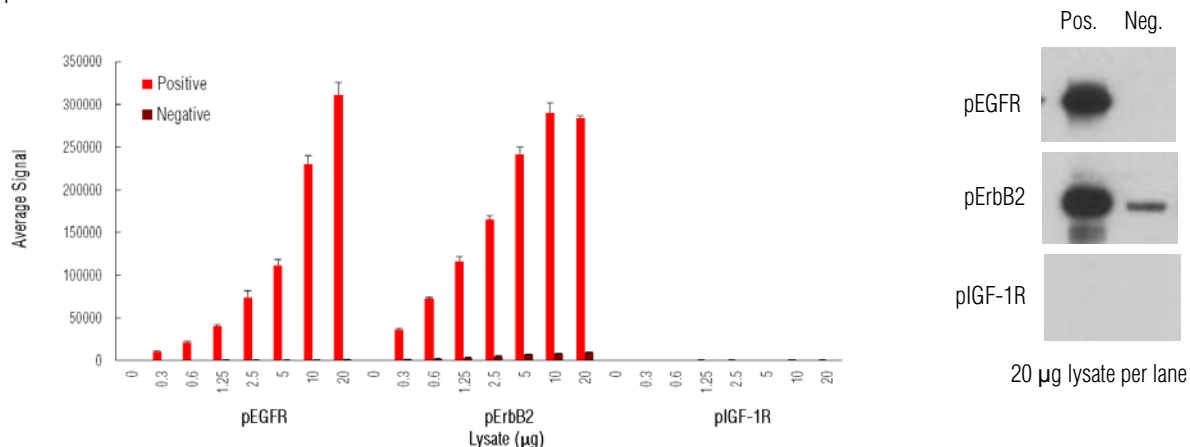


**Epidermal Growth Factor Receptor (EGFR), Human Epidermal Growth Factor Receptor 2 (Her2/Neu/ErbB2), and Insulin like Growth Factor I Receptor (IGF-1R)** are all receptor tyrosine kinases with extracellular ligand binding domains and intracellular tyrosine kinase domains. In normal cells these receptors play key roles in growth and development whereas IGF-1R also plays a critical role in metabolism and determination of overall nutrient state of the organism. All three of these receptors also signal through the PI3K-Akt signaling pathway and play a role in both normal as well as cancer cell development. EGFR and ErbB2 are oncoproteins and are involved in carcinogenesis, tumor growth/progression, and metastasis.<sup>1</sup> Many different cancers (such as breast cancer, prostate cancer, GI cancers, ovarian, neuroblastoma, and others) show increased expression of IGF-1R as well as IGF-1 and the level of expression tracks with tumor progression and cancer stage.<sup>2</sup> EGFR and ErbB2 are also involved in many different types of cancers, such as head, neck, colon, breast, pancreas, and lung cancers.<sup>3</sup> These important intracellular signaling receptors have been the subject of many basic research and drug development programs.

The MSD EGFR Family Assay is available on 96-well 4-Spot plates. This datasheet outlines the performance of the assay.

## Typical Data with SKOV3 Cells: Control Lysates for Phospho-EGFR and Phospho-ErbB2

Representative results for the EGFR Family are illustrated below. The signal and ratio values provided below are example data; individual results may vary depending upon the samples tested. Serum deprived SKOV3 cells were treated with Compound 56 and AG825 (1  $\mu$ M each, 2.5 hours) (negative) or pretreated with sodium vanadate (1 mM, 4 hours) and stimulated with EGF (100 ng/mL, 10 minutes) (positive). Whole cell lysates were added to MSD MULTI-SPOT® 4-Spot plates coated with anti-total EGFR, anti-total ErbB2, and anti-total IGF-1R antibodies on three of the four spatially distinct electrodes within a well. Phosphorylated EGFR, ErbB2, and IGF-1R were detected with an anti-phosphotyrosine antibody conjugated with MSD SULFO-TAG™ reagent. Western blot analyses of each lysate type were performed with phospho-EGFR, phospho-ErbB2, and phospho-IGF-1R antibodies and are shown below for comparison.



**Fig. 1:** Sample data generated with MULTI-SPOT EGFR Family. Increased signals for pEGFR and pErbB2 were observed with only the positive cell lysate. Signals for phospho-IGF-1R remained low throughout the titration of positive lysate. Signals for negative lysate were also low throughout the titration for all assays.

## Company Address

MESO SCALE DISCOVERY®  
A division of  
Meso Scale Diagnostics, LLC.  
9238 Gaither Road  
Gaithersburg, MD 20877 USA

[www.mesoscale.com](http://www.mesoscale.com)®

For Research Use Only.  
Not for use in diagnostic  
procedures.

# MSD Phosphoprotein Assays

## Lysate Titration

Data for positive and negative SKOV3 cell lysates using the MULTI-SPOT EGFR Family are presented below.

|         | Lysate<br>( $\mu$ g) | Positive       |        |      | Negative       |        |      | P/N |
|---------|----------------------|----------------|--------|------|----------------|--------|------|-----|
|         |                      | Average Signal | StdDev | %CV  | Average Signal | StdDev | %CV  |     |
| pEGFR   | 0                    | 81             | 26     | 32.1 | 74             | 13     | 17.2 |     |
|         | 0.31                 | 10782          | 779    | 7.2  | 141            | 27     | 19.4 | 76  |
|         | 0.63                 | 21653          | 981    | 4.5  | 195            | 9      | 4.6  | 111 |
|         | 1.3                  | 40238          | 2499   | 6.2  | 239            | 13     | 5.6  | 169 |
|         | 2.5                  | 74001          | 7543   | 10.2 | 295            | 16     | 5.4  | 251 |
|         | 5.0                  | 111613         | 6754   | 6.1  | 348            | 34     | 9.7  | 321 |
|         | 10                   | 230708         | 8906   | 3.9  | 706            | 63     | 8.9  | 327 |
|         | 20                   | 311189         | 14667  | 4.7  | 1166           | 183    | 15.7 | 267 |
| pErbB2  | 0                    | 139            | 29     | 21.2 | 123            | 17     | 13.9 |     |
|         | 0.31                 | 36591          | 1301   | 3.6  | 1152           | 52     | 4.5  | 32  |
|         | 0.63                 | 72742          | 1514   | 2.1  | 2117           | 46     | 2.1  | 34  |
|         | 1.3                  | 116381         | 5075   | 4.4  | 3448           | 130    | 3.8  | 34  |
|         | 2.5                  | 165277         | 5082   | 3.1  | 4818           | 413    | 8.6  | 34  |
|         | 5.0                  | 241421         | 8847   | 3.7  | 7184           | 231    | 3.2  | 34  |
|         | 10                   | 290486         | 11243  | 3.9  | 8151           | 255    | 3.1  | 36  |
|         | 20                   | 284703         | 2170   | 0.8  | 9648           | 110    | 1.1  | 30  |
| pIGF-1R | 0                    | 93             | 8      | 8.2  | 103            | 16     | 15.7 |     |
|         | 0.31                 | 503            | 60     | 11.9 | 131            | 21     | 15.8 | 3.8 |
|         | 0.63                 | 614            | 24     | 3.9  | 154            | 6      | 3.8  | 4.0 |
|         | 1.3                  | 784            | 18     | 2.3  | 167            | 5      | 3.0  | 4.7 |
|         | 2.5                  | 963            | 81     | 8.5  | 167            | 12     | 7.3  | 5.8 |
|         | 5.0                  | 592            | 36     | 6.1  | 133            | 40     | 29.8 | 4.4 |
|         | 10                   | 1102           | 29     | 2.6  | 181            | 4      | 2.2  | 6.1 |
|         | 20                   | 813            | 53     | 6.5  | 193            | 17     | 8.8  | 4.2 |

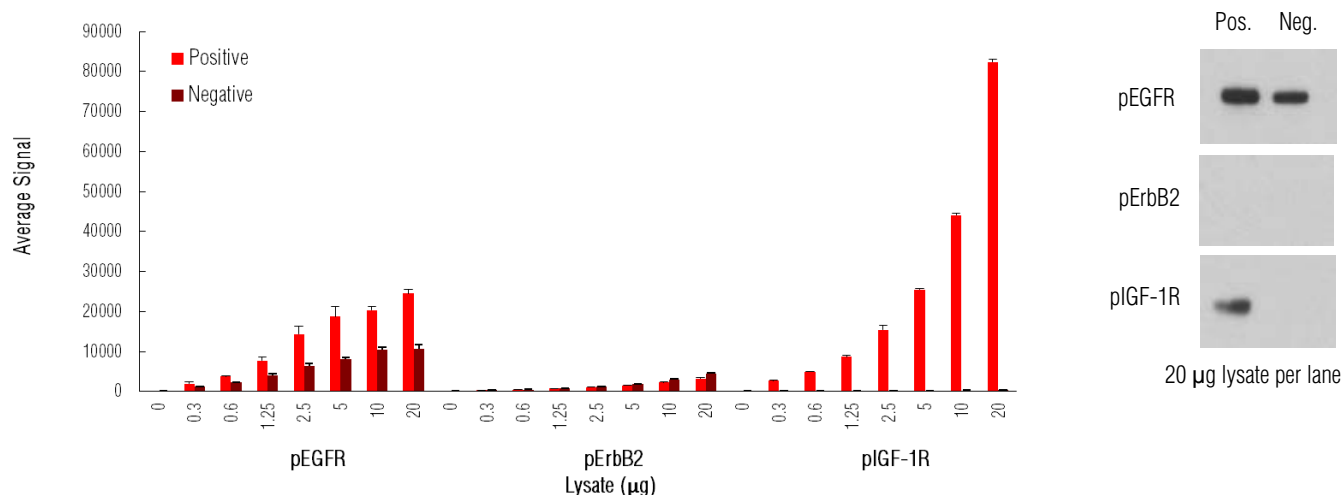
For Research Use Only. Not for use in diagnostic procedures.



# MSD Phosphoprotein Assays

## Typical Data with A431 Cells: Control Lysates for Phospho-IGF-1R

Representative results for the EGFR Family are illustrated below. The signal and ratio values provided below are example data; individual results may vary depending upon the samples tested. Serum deprived A431 cells (negative) were pretreated with sodium vanadate (1 mM, 4 hours) and stimulated with IGF-1 (100 nM, 10 minutes) (positive). Whole cell lysates were added to MSD MULTI-SPOT 4-Spot plates coated with anti-total EGFR, anti-total ErbB2, and anti-total IGF-1R antibodies on three of the four spatially distinct electrodes per well. Phosphorylated EGFR, ErbB2, and IGF-1R were detected with anti-phosphotyrosine antibody conjugated with MSD SULFO-TAG reagent. Western blot analyses of each lysate type were performed with phospho-EGFR, phospho-ErbB2, and phospho-IGF-1R antibodies and are shown below for comparison.



**Fig. 2:** Sample data generated with MULTI-SPOT EGFR Family. Increased signals for pEGFR and pIGF-1R were observed with the positive cell lysate. Signals for phospho-ErbB2 were relatively low.

### Lysate Titration

Data for positive and negative A431 cell lysates using the MULTI-SPOT EGFR Family are presented below.

|       | Lysate (µg) | Positive       |        |      | Negative       |        |      | P/N |
|-------|-------------|----------------|--------|------|----------------|--------|------|-----|
|       |             | Average Signal | StdDev | %CV  | Average Signal | StdDev | %CV  |     |
| pEGFR | 0           | 72             | 7      | 10.0 | 70             | 15     | 20.9 |     |
|       | 0.31        | 1933           | 406    | 21.0 | 1174           | 35     | 3.0  | 1.6 |
|       | 0.63        | 3720           | 129    | 3.5  | 2289           | 80     | 3.5  | 1.6 |
|       | 1.3         | 7733           | 907    | 11.7 | 3972           | 462    | 11.6 | 1.9 |
|       | 2.5         | 14358          | 2046   | 14.2 | 6346           | 672    | 10.6 | 2.3 |
|       | 5.0         | 18697          | 2466   | 13.2 | 8113           | 471    | 5.8  | 2.3 |
|       | 10          | 20289          | 1016   | 5.0  | 10537          | 627    | 5.9  | 1.9 |
|       | 20          | 24540          | 1068   | 4.4  | 10558          | 1230   | 11.6 | 2.3 |

For Research Use Only. Not for use in diagnostic procedures.



# MSD Phosphoprotein Assays

|         | Lysate<br>( $\mu$ g) | Positive       |        |      | Negative       |        |     | P/N |
|---------|----------------------|----------------|--------|------|----------------|--------|-----|-----|
|         |                      | Average Signal | StdDev | %CV  | Average Signal | StdDev | %CV |     |
| pErbB2  | 0                    | 110            | 13     | 12.1 | 111            | 1      | 1.0 |     |
|         | 0.31                 | 297            | 23     | 7.7  | 312            | 19     | 6.1 | 0.9 |
|         | 0.63                 | 480            | 15     | 3.1  | 483            | 29     | 5.9 | 1.0 |
|         | 1.3                  | 729            | 22     | 3.0  | 740            | 17     | 2.3 | 1.0 |
|         | 2.5                  | 1027           | 54     | 5.2  | 1177           | 29     | 2.4 | 0.9 |
|         | 5.0                  | 1500           | 95     | 6.4  | 1807           | 20     | 1.1 | 0.8 |
|         | 10                   | 2254           | 96     | 4.3  | 2970           | 122    | 4.1 | 0.8 |
|         | 20                   | 3228           | 226    | 7.0  | 4506           | 231    | 5.1 | 0.7 |
| pIGF-1R | 0                    | 109            | 7      | 6.3  | 93             | 5      | 5.4 |     |
|         | 0.31                 | 2682           | 178    | 6.6  | 152            | 12     | 7.7 | 18  |
|         | 0.63                 | 4975           | 82     | 1.6  | 165            | 4      | 2.6 | 30  |
|         | 1.3                  | 8623           | 373    | 4.3  | 196            | 10     | 4.8 | 44  |
|         | 2.5                  | 15432          | 1163   | 7.5  | 228            | 7      | 3.1 | 68  |
|         | 5.0                  | 25444          | 297    | 1.2  | 222            | 22     | 9.7 | 115 |
|         | 10                   | 44107          | 563    | 1.3  | 267            | 11     | 4.1 | 165 |
|         | 20                   | 82362          | 649    | 0.8  | 299            | 21     | 7.0 | 276 |

## MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample amounts of 25  $\mu$ g/well or less without compromising speed or performance
- **Large dynamic range:** Linear range of up to five logs enables the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions
- **Minimal background:** The stimulation mechanism (electricity) is decoupled from the signal (light)
- **Simple protocols:** Only labels near the electrode surface are detected, enabling no-wash assays
- **Flexibility:** Labels are stable, non-radioactive, and conveniently conjugated to biological molecules
- **High sensitivity and precision:** Multiple excitation cycles of each label enhance light levels and improve sensitivity

For a complete list of products, please visit our website at [www.mesoscale.com](http://www.mesoscale.com)

## References:

1. Gazdar AF. Activating and resistance mutations of EGFR in non-small-cell lung cancer: role in clinical response to EGFR tyrosine kinase inhibitors. *Oncogene*. 2009 Aug;28 Suppl 1:S24-31.
2. Li R, Pourpak A, Morris SW. Inhibition of the insulin-like growth factor-1 receptor (IGF-1R) tyrosine kinase as a novel cancer therapy approach. *J Med Chem*. 2009 Aug 27;52(16):4981-5004.
3. Wheeler DL, Dunn EF, Harari PM. Understanding resistance to EGFR inhibitors—impact on future treatment strategies. *Nat Rev Clin Oncol*. 2010 Sep;7(9):493-507. Epub 2010 Jun 15.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, SECTOR, SECTOR HTS, SECTOR PR, 4-SPOT (DESIGN) and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.  
© 2011 Meso Scale Diagnostics, LLC. All rights reserved.

For Research Use Only. Not for use in diagnostic procedures.

