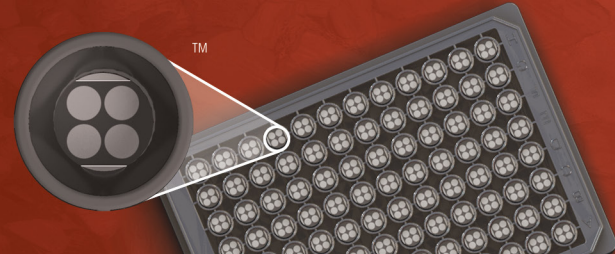


MSD® Phospho(Tyr1173)/Total EGFR Assay Whole Cell Lysate Kit

For quantitative determination 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

Phospho (Tyr1173)/Total EGFR Whole Cell Lysate Kit	
Kit size	
1 plate	K15104D-1
5 plates	K15104D-2
20 plates	K15104D-3

Phospho-EGFR Whole Cell Lysate Set	
200 µg	C11C1-1

Ordering information

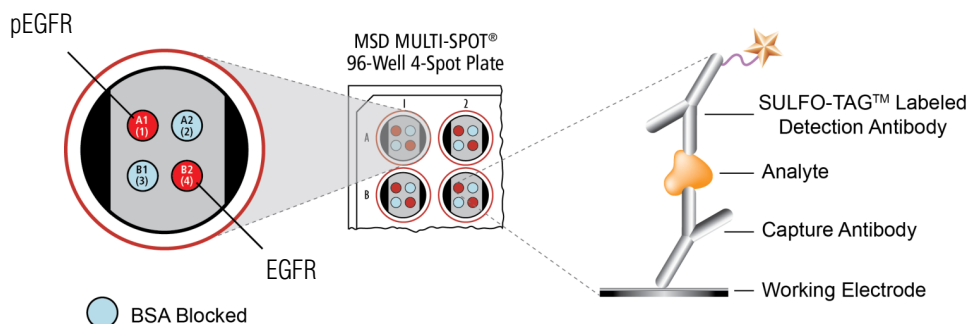
MSD Customer Service
Phone: 1-301-947-2085
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com

Company Address

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

www.mesoscale.com®

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



EGFR (Epidermal Growth Factor Receptor) is a 170 kDa transmembrane receptor tyrosine kinase, consisting of a ligand-binding extracellular domain, a single transmembrane domain, an intracellular protein-tyrosine kinase catalytic domain, and a tyrosine-containing cytoplasmic tail. EGFR (ErbB1/HER1) is one of a family of four ErbB/HER (1-4) receptor tyrosine kinases, each essential to embryonic survival. Upon binding its ligand EGF, the EGFR forms hetero- or homodimers. Dimerization results in the activation of its intrinsic tyrosine kinase activity and the phosphorylation of multiple tyrosines in the cytoplasmic domain, including Tyr992, Tyr1068, Tyr1086, Tyr1148, and Tyr1173. The phosphorylated tyrosines are binding sites for proteins containing SH2-domains. These binding events activate many intracellular signaling pathways including MAPK/ERK, PI-3K, PKC, and p38, controlling cell growth, survival, cell cycle arrest, and transformation. Due to its central role in many cellular physiological processes, EGFR overexpression and aberrant signaling is associated with many types of cancer, making EGFR an attractive target for chemotherapeutic drug development

The MSD Phospho(Tyr1173)/Total EGFR Assay is available on 96-well 4-Spot plates. This datasheet outlines the performance of the assay.

Typical Data with A431 cells treated with EGF or Compound 56

Representative results for the Phospho(Tyr1173)/Total EGFR Assay 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 were treated with compound 56 (5 nM; 3 hours) (negative), or with EGF (100 ng/mL; 10 minutes) (positive). Whole cell lysates were added to MSD MULTI-SPOT® 4-Spot plates coated with anti-phospho-EGFR (Tyr1173) and anti-total EGFR antibodies on spatially distinct electrodes within a well. Phosphorylated and total EGFR were detected with anti-total EGFR antibody conjugated with MSD SULFO-TAG™ reagent. Western blot analyses of each lysate type were performed with phospho-EGFR (Tyr1173) and total EGFR antibodies and are shown below for comparison.

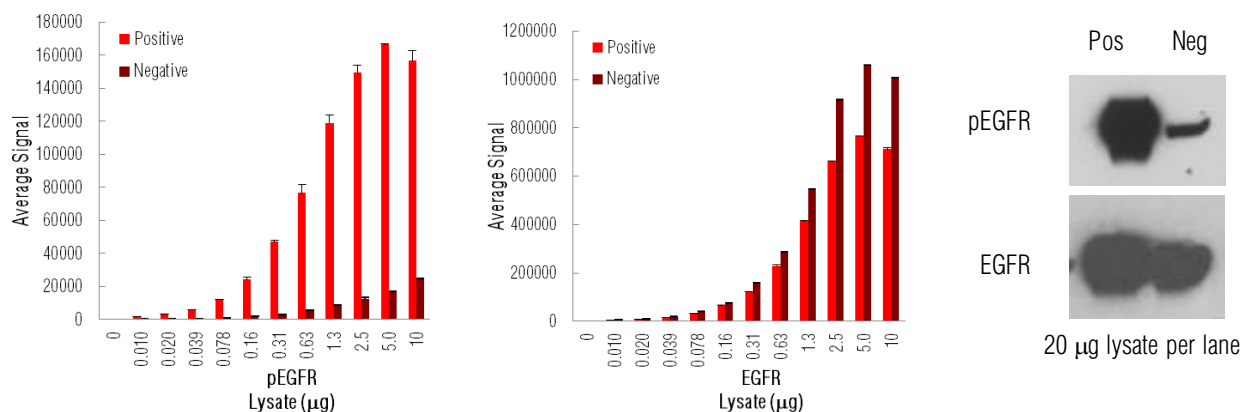


Fig. 1: Sample data generated with MULTI-SPOT Phospho(Tyr1173)/Total EGFR Assay. Increased signal for phosphorylated EGFR was observed with pEGFR positive cell lysate. Total EGFR signal increased throughout the titration of both pEGFR positive and negative cell lysates. The Phospho(Tyr1173)/Total EGFR Assay provides a quantitative measure of the data obtained with the traditional Western blot.

MSD Phosphoprotein Assays

Lysate Titration

Data for positive and negative A431 cell lysates using the MULTI-SPOT Phospho(Tyr1173)/Total EGFR Assay are presented below.

	Lysate (μ g)	Positive			Negative			P/N
		Average Signal	StdDev	%CV	Average Signal	StdDev	%CV	
pEGFR	0	83	37	45.4	83	37	45.4	
	0.010	1535	28	1.8	149	2	1.4	10
	0.020	3093	272	8.8	242	11	4.7	13
	0.039	5902	12	0.2	445	46	10.3	13
	0.078	11857	193	1.6	913	53	5.8	13
	0.16	24147	1647	6.8	1688	121	7.2	14
	0.31	47152	573	1.2	3174	29	0.9	15
	0.63	76705	4835	6.3	5215	573	11.0	15
	1.3	118913	4691	3.9	8261	431	5.2	14
	2.5	149480	4329	2.9	12501	725	5.8	12
	5.0	166478	311	0.2	16809	272	1.6	9.9
10	156688	6253	4.0	24320	398	1.6	6.4	
EGFR	0	184	52	28.1	184	52	28.1	
	0.010	3861	18	0.5	4958	325	6.5	0.8
	0.020	7637	180	2.4	9914	670	6.8	0.8
	0.039	15669	236	1.5	18656	1650	8.8	0.8
	0.078	30948	817	2.6	39656	276	0.7	0.8
	0.16	64429	4687	7.3	76377	2477	3.2	0.8
	0.31	121570	292	0.2	157668	5238	3.3	0.8
	0.63	228249	9044	4.0	285751	19146	6.7	0.8
	1.3	413168	675	0.2	544371	2655	0.5	0.8
	2.5	659744	7092	1.1	914856	23540	2.6	0.7
	5.0	766507	63037	8.2	1058852	3263	0.3	0.7
10	711199	1256	0.2	1006322	14843	1.5	0.7	

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



MSD Phosphoprotein Assays

Typical Data with COS7 cells treated with EGF

Representative results for the Phospho(Tyr1173)/Total EGFR are illustrated below. The signal and ratio values provided below are example data; individual results may vary depending upon the samples tested. Serum deprived COS-7 cells (negative) were treated with EGF (100 ng/mL, 10 minutes) (positive). Whole cell lysates were added to MSD MULTI-SPOT 4-Spot plates coated with anti-phospho-EGFR (Tyr1173) and anti-total EGFR antibodies on spatially distinct electrodes within a well. Phosphorylated and total EGFR were detected with anti-total EGFR antibody conjugated with MSD SULFO-TAG reagent. Western blot analyses of each lysate type were performed with phospho-EGFR (Tyr1173) and total EGFR antibodies and are shown below for comparison.

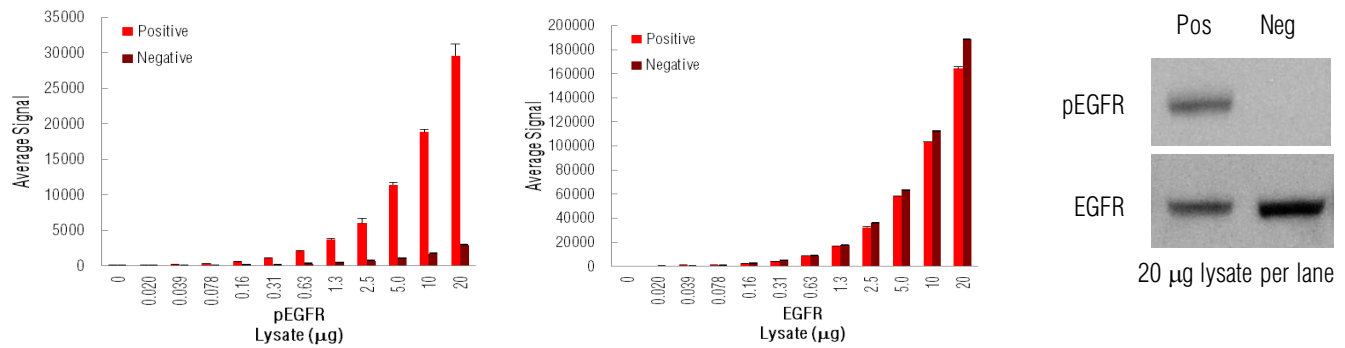


Fig. 2: Sample data generated with MULTI-ARRAY Phospho-EGFR (Tyr1173) Assay. Increased signal for phosphorylated EGFR was observed with pEGFR positive cell lysate. Total EGFR signal increased throughout the titration of both pEGFR positive and negative cell lysates. The Phospho(Tyr1173)/Total EGFR Assay provides a quantitative measure of the data obtained with the traditional Western blot.

Lysate Titration

Data for positive and negative COS7 cell lysates using the MULTI-SPOT Phospho(Tyr1173)/Total EGFR are presented below.

	Lysate (µg)	Positive			Negative			P/N
		Average Signal	StdDev	%CV	Average Signal	StdDev	%CV	
pEGFR	0	74	13	17.2	74	13	17.2	
	0.020	124	12	9.7	74	1	1.9	1.7
	0.039	226	23	10.3	81	8	9.7	2.8
	0.078	319	19	6.0	95	8	8.2	3.4
	0.16	606	6	0.9	161	6	4.0	3.8
	0.31	1083	44	4.0	214	5	2.3	5.1
	0.63	2068	119	5.7	311	32	10.2	6.7
	1.3	3742	95	2.6	424	8	2.0	8.8
	2.5	5986	717	12.0	705	40	5.6	8.5
	5.0	11415	305	2.7	1098	21	1.9	10
	10	18828	373	2.0	1712	34	2.0	11
	20	29591	1651	5.6	2935	38	1.3	10

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



MSD Phosphoprotein Assays

	Lysate (μ g)	Positive			Negative			P/N
		Average Signal	StdDev	%CV	Average Signal	StdDev	%CV	
EGFR	0	117	17	14.5	117	17	14.5	
	0.020	382	1	0.4	431	38	8.9	0.9
	0.039	1046	536	51.2	700	9	1.3	1.5
	0.078	1115	81	7.2	1268	59	4.6	0.9
	0.16	2199	111	5.0	2489	115	4.6	0.9
	0.31	4421	203	4.6	4866	172	3.5	0.9
	0.63	8729	93	1.1	9086	222	2.4	1.0
	1.3	16657	489	2.9	17907	643	3.6	0.9
	2.5	32427	1176	3.6	35960	1232	3.4	0.9
	5.0	58184	924	1.6	63150	2567	4.1	0.9
	10	103256	4692	4.5	112130	465	0.4	0.9
	20	164260	7383	4.5	188770	4751	2.5	0.9

MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample amounts of 25 μ 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

References:

1. Rogers SJ, Box C, Chambers P, Barbachano Y, Nutting CM, Rhys-Evans P, Workman P, Harrington KJ, Eccles SA. Determinants of response to epidermal growth factor receptor tyrosine kinase inhibition in squamous cell carcinoma of the head and neck. *J Pathol.* 2009 May;218(1):122-30.
2. Cao L, Yu Y, Darko I, Currier D, Mayeenuddin LH, Wan X, Khanna C, Helman LJ. Addiction to elevated insulin-like growth factor I receptor and initial modulation of the AKT pathway define the responsiveness of rhabdomyosarcoma to the targeting antibody. *Cancer Res.* 2008 Oct 1;68(19):8039-48.
3. Martin SE, Jones TL, Thomas CL, Lorenzi PL, Nguyen DA, Runfola T, Gunsior M, Weinstein JN, Goldsmith PK, Lader E, Huppi K, Caplen NJ. Multiplexing siRNAs to compress RNAi-based screen size in human cells. *Nucleic Acids Res.* 2007;35(8):e57. Epub 2007 Mar 28.
4. Gowan SM, Hardcastle A, Hallsworth AE, Valenti MR, Hunter LJ, de Haven Brandon AK, Garrett MD, Raynaud F, Workman P, Aherne W, Eccles SA. Application of meso scale technology for the measurement of phosphoproteins in human tumor xenografts. *Assay Drug Dev Technol.* 2007 Jun;5(3):391-401.

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.

