

Bio-Plex™ suspension array system

tech note 5301

Target Expression and Target Modulation Studies in Patient-Derived Tumor Xenografts Using a Bank of Protein Lysates

Martina Maurer, Andreas Wortmann, André Korrat, Heike Legrand, Julia Schüler, Sandra Kissel, Niko Bausch, Thomas Metz, and Heinz-Herbert Fiebig, Oncotest GmbH, Institute for Experimental Oncology, Am Flughafen 12-14, D-79108 Freiburg, Germany

Introduction

In this tech note, we present an efficient method that facilitates the preclinical analysis of antitumor compounds. By using specialized human tumor models and multiplex bead suspension assays, we are able to decide which tumor models to use for preclinical profiling of anticancer agents as well as identify biomarkers for specific drug response or resistance.

Methods

Protein Tumor Lysate Bank

By directly transplanting patient tumors subcutaneously into immune-compromised nude mice, more than 400 solid tumors, including all major human tumor types, have been established in serial passage (Fiebig et al. 1992, Fiebig and Burger 2001). At Oncotest, we have generated a protein tumor lysate bank consisting of 150 tumor models (Table 1).

Table 1. Xenografts selected for the protein tumor lysate bank (n = 150).

Bladder	4	Lymphoma	4	Pancreas	4
Colon	19	Mammary	14	Pleuroesothelioma	2
Gastric	4	Melanoma	12	Prostate	5
Head and neck	4	Non-small cell lung cancer (NSCLC)	33	Sarcomas	4
Kidney	6	Small cell lung	6	Uterus	5
Leukemias	5	Ovarian	8	Others	8
Liver	3				

Multiplex Bead Suspension Assays

Bead suspension assays are flexible bioassay systems that allow the parallel detection and quantitation of many targets in a single sample. For example, we can measure up to 28 cytokines and chemokines or 7 phosphorylated signal transduction proteins simultaneously. In a typical multiplex bead suspension assay, a target protein is captured from a crude cell lysate, tissue lysate, or serum with bead-bound antibodies. The total target protein amount and/or its phosphorylation status is quantitated with a secondary biotinylated antibody followed by an incubation step with streptavidin-phycoerythrin to complete a sandwich immunoassay. In this study, we used Bio-Rad's Bio-Plex multiplex suspension array system and Oncotest xenograft lysates (Figure 1).

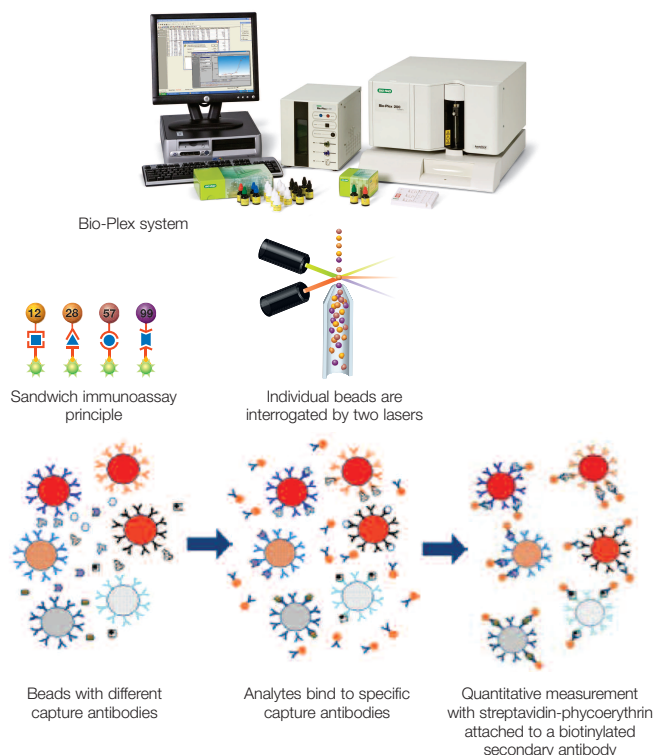


Fig. 1. Multiplex bead suspension assay test principle. The Bio-Plex multiplex system and Oncotest xenograft lysates (n = 150) were used to determine total target amount and phosphorylation status.

Results and Discussion

Expression Profiling

Multiplex bead suspension assays (Table 2) were used to determine the expression and phosphorylation state of a multitude of molecular markers involved in signal transduction, cell cycle, apoptosis, and so on. The expression profiles allowed the selection of optimal tumor models for the in vivo testing of novel anticancer agents (Figure 2). Furthermore, we used the complete molecular profiles to identify biomarkers predictive for drug response or resistance.

Target Monitoring

Another application of the multiplex bead suspension assays was the monitoring of target expression and activity (e.g., phospho-EGFR) during therapy experiments in tumor-bearing nude mice (Figure 3).

Table 2. Multiplex bead assays available at Oncotest as of February 2006. Highlighted assays are developed and manufactured at Bio-Rad.

Phosphorylated Signal Transduction Proteins	Phospho-STAT3 (Ser ⁷²⁷)	Apoptosis-Associated Proteins	Cytokines and Chemokines
Phospho-Akt/PKB (Ser ⁴⁷³)	Phospho-STAT5A/B (Tyr ^{694/699})	Active caspase 3	Eotaxin
Phospho-ATF-2 (Thr ⁷¹)	Phospho-Tau (Ser ¹⁹⁹)	Bcl-2	GM-CSF
Phospho-c-Jun (Ser ⁶³)	Phospho-Tau (Thr ¹⁸¹)	Cleaved PARP	G-CSF
Phospho-c-Kit	Phospho-TrkA (Tyr ⁴⁹⁰)	Single-stranded DNA	IL-1 α
Phospho-c-Met	Total Signal Transduction Proteins	Transcription Factors	IL-1 β
Phospho-CREB (Ser ¹³³)	Total Akt/PKB	AP-2	IL-2
Phospho-EGF receptor	Total ATF-2	CREB	IL-3
Phospho-ERK/MAP kinase 1/2 (Thr ¹⁸⁵ /Tyr ¹⁸⁷)	Total active β -catenin	EGR	IL-4
Phospho-GSK-3 α/β (Ser ²¹ /Ser ⁹)	Total c-Kit	HIF	IL-5
Phospho-HSP27 (Ser ⁷⁸)	Total c-Jun	NF1	IL-6
Phospho-I κ B- α (Ser ³²)	Total CREB	NF- κ B	IL-7
Phospho-IRS-1	Total EGF receptor	NFAT	IL-8
Phospho-JNK/SAPK1 (Thr ¹⁸³ /Tyr ¹⁸⁵)	Total ERK/MAP kinase 1/2	PPAR	IL-10
Phospho-Jun (Ser ⁷³)	Total ERK2	SRE	IL-12 (p40)
Phospho-Lck	Total HSP27	YY1	IL-12 (p70)
Phospho-NF- κ B (Ser ⁵³⁶)	Total I κ B- α	Matrix Metalloproteins	IL-13
Phospho-p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	Total IRS-1	MMP-1	IL-15
Phospho-p53 (Ser ¹⁵)	Total JNK/SAPK1	MMP-2	IL-17
Phospho-p70 S6 kinase (Thr ⁴²¹ /Ser ⁴²⁴)	Total Lck	MMP-3	Interferon- α 2
Phospho-p90RSK (Thr ³⁵⁹ /Ser ³⁶³)	Total p38 MAPK	MMP-7	Interferon- γ
Phospho-PDGFR α	Total p53	MMP-8	IP-10
Phospho-PRAS40 (Thr ²⁴⁵)	Total p70 S6 kinase	MMP-9	MCP-1
Phospho-Rb (Ser ²⁴⁹ /Thr ²⁵²)	Total p90RSK	MMP-12	MIP-1 α
Phospho-Rb (Thr ⁸²¹)	Total Rb	MMP-13	MIP-1 β
Phospho-RSK1/MAPKAP kinase 1a (Ser ³⁸⁰)	Total STAT1	Death Receptors	RANTES
Phospho-STAT1 (Tyr ⁷⁰¹)	Total STAT3	DR5	TNF- α
Phospho-STAT2	Total Tau	TNF-RI	TNF- β
Phospho-STAT3 (Tyr ⁷⁰⁵)		TNF-RII	

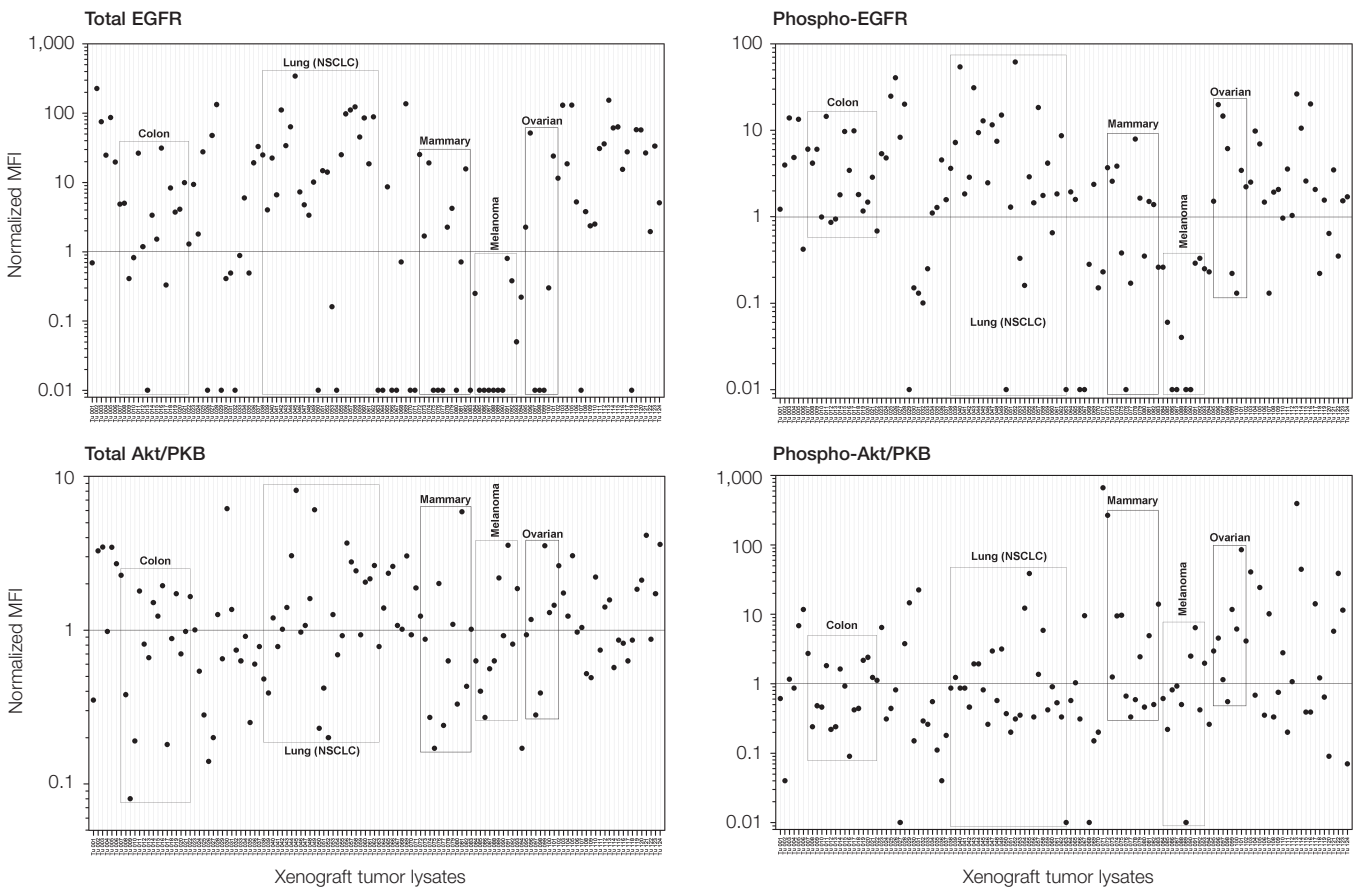


Fig. 2. Expression profiling of total proteins and phosphoproteins in 124 human tumor models. Upper panels, total and phospho-EGFR amounts; lower panels, Akt/PKB amounts. Each parameter was measured with the Bio-Plex system, and the black line shows the geometric mean value over all tumors. Tumors above the line exhibit enhanced protein expression or phosphorylation, respectively. These xenografts are suitable tumor models for target-directed in vivo studies.

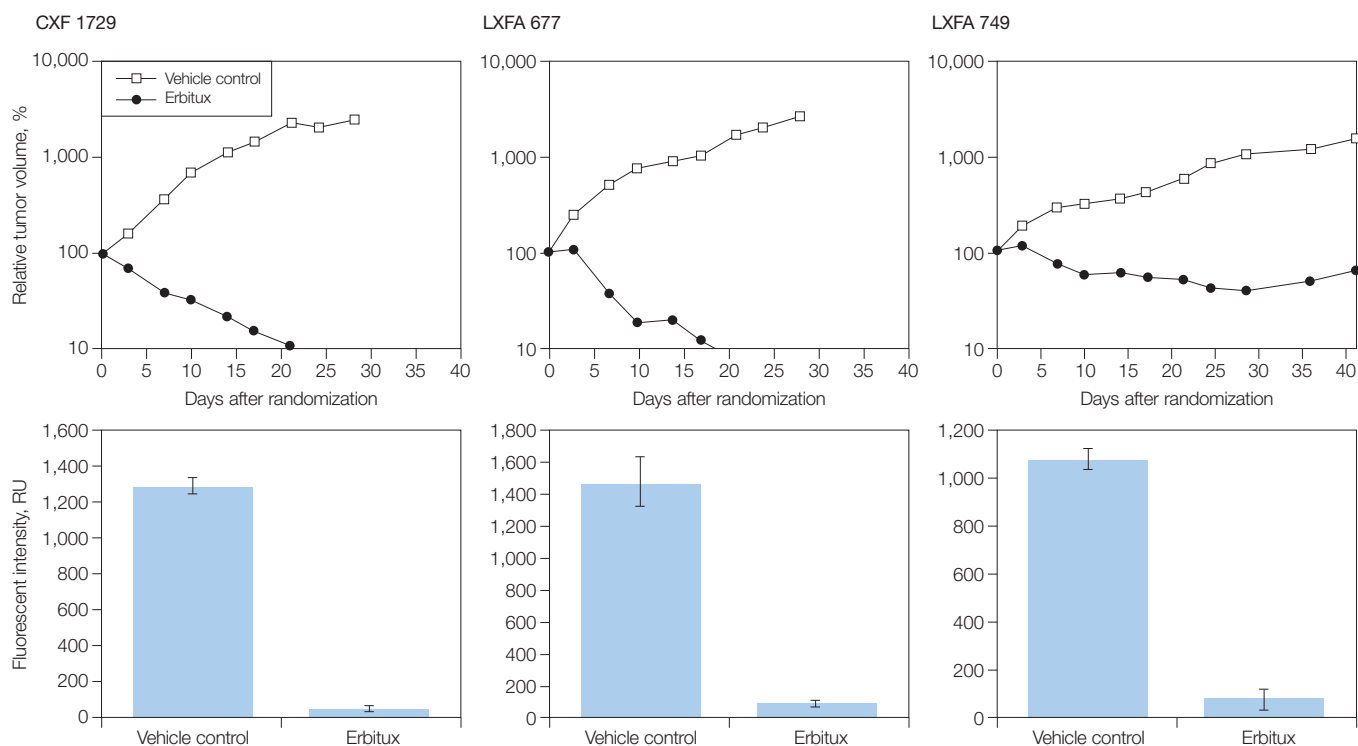


Fig. 3. Target monitoring for EGFR-inhibitory monoclonal antibody Erbitux in three sensitive tumor xenografts. Upper panels show tumor volume inhibition over time after intraperitoneal treatment with Erbitux on days 1, 8, and 15 at 30 mg/kg. Lower panels show the drop of EGFR phosphorylation 24 hr after administration of a single dose of Erbitux. CXF = colon carcinoma xenograft Freiburg; LXFA = lung adenocarcinoma xenograft Freiburg.

Conclusions

The detailed molecular characterization of Oncotest's tumor collection allows an optimal choice of tumor models for the preclinical profiling of promising new anticancer agents. This approach accelerates and improves the preclinical evaluation of novel target-directed compounds. In addition, the identification of biomarkers specific for drug response or resistance can be used as a diagnostic tool to determine the most promising individual therapy for cancer patients.

Acknowledgement

We thank Nicole Gollmitzer (Bio-Rad, Germany) for excellent technical advice.

References

- Fiebig H-H et al., Combined in vitro/in vivo test procedure with human tumor xenografts for new drug development, *Contrib Oncol* 42, 321-351 (1992)
- Fiebig H-H and Burger AM, Human tumor xenografts and explants, pp 113-137 in Teicher BA (ed), *Tumor Models in Cancer Research*, Humana Press, Totowa, NJ (2001)

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

Erbitux is a trademark of ImClone Systems, Inc.

Information in this tech note was current as of the date of writing (2005) and not necessarily the date this version (rev A, 2006) was published.



**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300/5013478 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 55 5200 05 20 **The Netherlands** 0318 540666 **New Zealand** 64 9415 2280
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 095 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 0861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189/2578 7241 **United Kingdom** 020 8328 2000