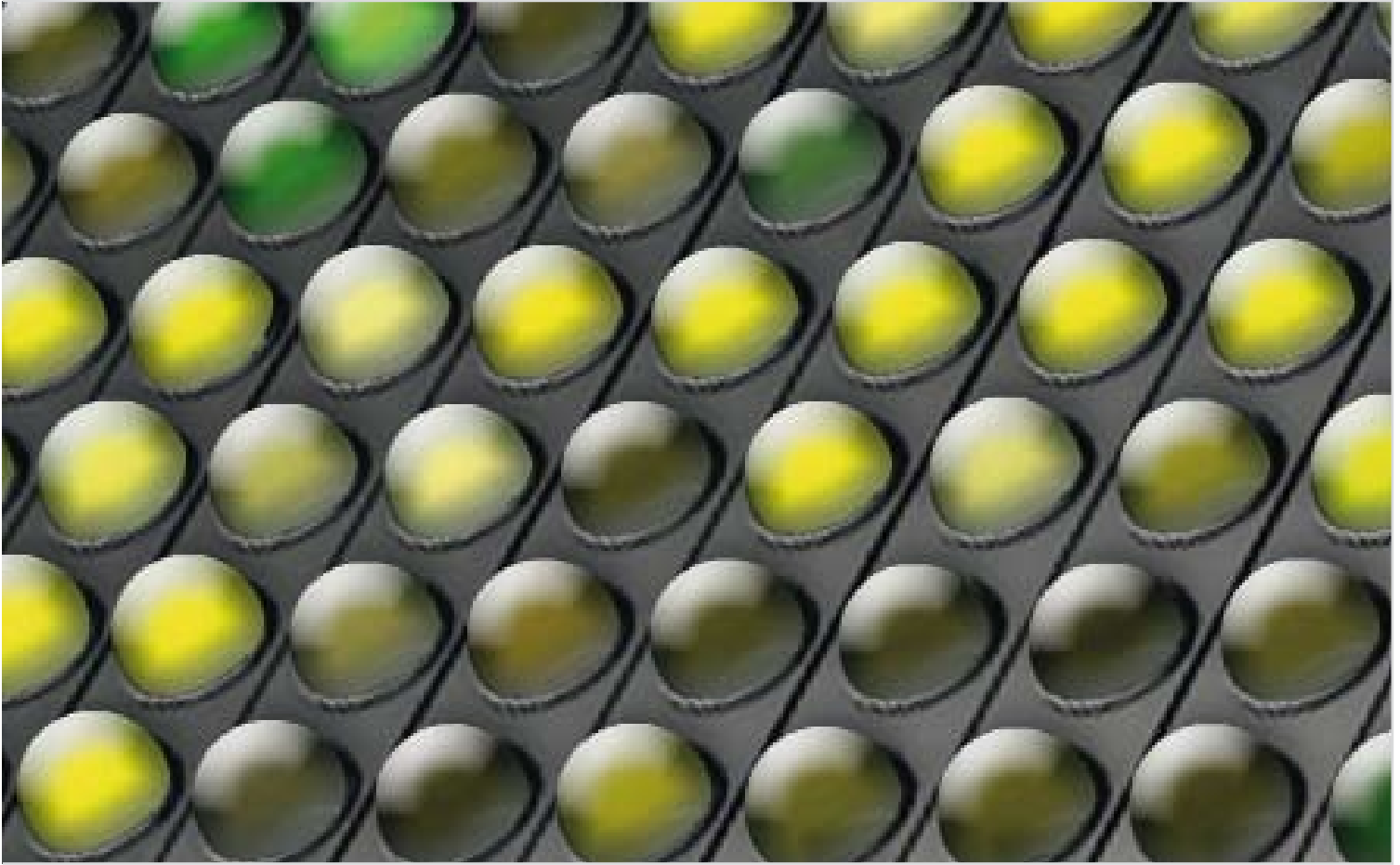


# ANTIBODY MICROARRAYS



EXTEND YOUR KNOWLEDGE ABOUT THE EXPRESSION OF  
SEVERAL PROTEINS IN ONE ASSAY

# PATHWAY & PHOSPHORYLATION ANTIBODY MICROARRAYS

Antibody MicroArray provides a high-throughput platform for efficient and accurate profiling of protein expressions. Our Pathway Focused Antibody MicroArrays and Phosphorylation Antibody MicroArrays are designed for researchers to study highly relevant proteins in their specific research fields. The antibodies are covalently immobilized on high quality glass surface coated with a proprietary 3-D polymer materials to ensure high binding efficiency and specificity.



Arrays utilize fluorescent detection and can be scanned on all MicroArray scanners that are compatible with 76 x 25 x 1 mm slides.

- > The antibodies are printed on standard-size microscope slides.
- > Each slide consists of an array of well-characterized antibodies with six replicates.
- > Multiple positive and negative controls are present.
- > Each antibody array kit comes with two identical slides.

## Applications

- > Qualitative Protein Expression Profiling
- > Measure Changes in protein expression level (normal vs. diseased; normal vs. treated)
- > Identify Candidate Biomarkers

## Specifications



- > Spot diameter: 260-280  $\mu\text{m}$
- > Slide dimensions: 76 x 25 x 1 mm
- > Spot-to-Spot distance 450-600  $\mu\text{m}$
- > Barcode label: 23 x 9 mm
- > Storage conditions: 4  $^{\circ}\text{C}$  (6 months)

## Key Features

- > Each array includes carefully selected antibodies highly relevant for a specific pathway.
- > High-throughput expression profiling using proteins from cell extracts, tissue lysates or treated samples.
- > 3D-Polymer coated glass substrates
- > High specificity
- > Low background
- > Sensitive fluorescent detection (Streptavidin – Cy<sup>®</sup>3 (Cy 5<sup>®</sup>/Alexa<sup>®</sup>532/Alexa<sup>®</sup>647)
- > Six replicates for increased data consistency
- > Suitable for detection on all fluorescent MicroArray scanners compatible with 76 x 25 x 1 mm slides

- > Optimised Antibody Array Assay Kit provides optimised reagents and buffers for protein extraction, labelling, conjugation and detection.
- > GenePix Array List (GAL) files are provided for convenient and accurate data analysis.

**Phospho-specific antibodies are available separately (50-150  $\mu\text{g}$ ). Please inquire for reference and price information.**

## COMPREHENSIVE ANTIBODY ARRAY

Description	Reference	# Abs	# Slides
Explorer™ Antibody Array	AR-ASB600	656	2
Phospho Explorer™ Antibody Array	AR-PEX100	1318	2
Tyrosine phosphorylation proarray AA	AR-PST228	228	2

## PATHWAY-SPECIFIC ARRAYS

Angiogenesis Antibody Array	AR-AGE1-25	127	2
Apoptosis Antibody Array	AR-APPO-69	73	2
Cancer Marker Antibody Array	AR-ACM0-83	86	2
Cell Cycle Antibody Array	AR-ACCO-58	60	2
Cytokine Related Antibody Array	AR-ACK0-75	77	2
Hematopoiesis Antibody Array	AR-AHT0-49	49	2
Hormone Antibody Array	AR-AHM0-30	30	2
Signal Transduction Antibody Array	AR-AST1-60	165	2
Stem Cell Antibody Array	AR-ASCO-48	58	2
Explorer Antibody Array	AR-ASB600	656	2

## PHOSPHO-SPECIFIC ARRAYS

AKT Pathway Phospho Antibody Array	AR-PAA1-37	137	2
AKT/PKB Signaling Phospho Antibody Array	AR-PAB2-16	216	2
Cancer/Apoptosis Phospho Antibody Array	AR-PAC1-55	155	2
AMPK Signaling Antibody Array	AR-PAM1-74	174	2
Apoptosis Phospho Antibody Array	AR-PAP2-47	247	2
Cell Cycle Phospho Antibody Array	AR-PCC0-76	76	2
Cell Cycle Control Phospho Antibody array	AR-PCC2-38	238	2
Cytoskeleton Phospho Antibody Array	AR-PCPO-95	95	2
Cytoskeleton II Phospho Antibody Array	AR-PCPI-41	141	2
CREB Antibody Array	AR-PCR1-74	174	2
Chromatin/Transcription Phospho Antibody Array	AR-PCT1-73	173	2
PDGF Antibody Array	AR-PDG1-95	195	2
EGF Pathway Antibody Array	AR-PEG2-14	214	2
Erk Signaling Antibody Array	AR-PEK2-08	208	2
ErbB/HER Signaling Antibody Array	AR-PER2-39	239	2
Phospho Explorer Antibody Array	AR-PEX1-00	1318	2
p53 Signaling Antibody Array	AR-PFT1-96	196	2
FGF Pathway Antibody Array	AR-PGF1-56	156	2
GPCR Signaling to MAPK/ERK Antibody Array	AR-PGP1-93	193	2
Insulin/Glucose Phospho Antibody Array	AR-PIG0-85	85	2
Insulin Receptor Phospho-Specific Antibody Array	AR-PIG2-19	219	2
IGF-IR Signaling Antibody Array	AR-PIR2-45	245	2
Jak/Stat Pathway Phospho Antibody Array	AR-PJS0-42	42	2
Jak/Stat II Phospho Antibody Array	AR-PJS2-02	202	2
MAPK Pathway Phospho Antibody Array	AR-PMK1-85	185	2
mTOR Phospho Antibody Array	AR-PMT1-38	138	2
NF-kB Pathway Phospho Antibody Array	AR-PNK0-63	63	2
NF-kB II Signaling Phospho Antibody Array	AR-PNK2-15	215	2
Nuclear/Membrane Receptor Phospho Antibody Array	AR-PNR0-52	52	2
Neuroscience Phospho Antibody Array	AR-PNS0-32	32	2
Tyrosine Phosphorylation ProArray Antibody Array	AR-PST2-28	228	2
T-Cell Receptor Signaling Antibody Array	AR-PTC1-88	188	2
TGF-beta Signaling Antibody Array	AR-PTG1-76	176	2
Tyrosine Kinase Adaptor Phospho Antibody Array	AR-PTK0-98	98	2
VEGF Pathway Antibody Array	AR-PVE1-85	185	2

## ANTIBODY ARRAY ASSAY KITS

- > The Antibody Array Assay Kits are designed for easy and reliable processing of antibody arrays.
- > Each kit comes with sufficient material and reagents to perform protein extract, labelling, conjugation and detection for two slides or twenty slides.
- > The buffers and reagents have been optimised to maximize protein extraction yield, labelling efficiency, and to produce high signal-to-noise ratio.

### ANTIBODY ARRAY ASSAY KITS

Antibody Array Assay Kit, 2 Reactions	AR-KAS0-02
Antibody Array Assay Kit, 20 Reactions	AR-KAS0-20

### Key features

- > Complete and easy-to-use reagent kit
- > Provide major reagents required to process antibody Arrays
- > Optimised to work with Thematic, Phospho-specific and Explorer antibody MicroArrays

## COMPLETE ANTIBODY ARRAY ASSAY SERVICE

We perform complete antibody array assays with your samples. You can select from a wide selection of antibody microarrays to examine normal and treated samples for changes in protein expression and phosphorylation events. This convenient assay service offers quick turnaround, accurate and reliable results, saving you time and resources!

- > Data acquisition. (Data provided in Excel format)
- > Full report including protocol, data in Excel format and images
- > Turnaround time: usually 8-12 days (maybe longer if there are more than 6 samples)

### The protein expression profiling service includes



- > One antibody Array slide
- > Protein Extraction and labelling
- > Assay condition optimisation
- > Conjugation of Biotin Labelled Protein to Antibody Array
- > Detection using Cy<sup>3</sup>-Streptavidin
- > Scan the arrays on Axon GenePix scanners

### Starting material

For each sample you wish to assay, please provide cell or tissue lysates with 400 µg or more total protein at a minimum of 2 µg/ µl. Acceptable lysis buffers include Antibody Array Protein Extraction Buffer (AR-EXB-050), RIPA buffer, or other buffers suggested by commonly used lysis protocols. The use of Tris buffer and detergents should be avoided or limited. No more than 50 mM of Tris, 0.20 % SDS, or 1 % non-ionic detergents should be present in the sample. Commonly used protease and/or phosphatase inhibitors may be added to the lysis buffer to preserve the proteins. The proteins in the sample should remain in their native structure (not denatured).

In addition to cell or tissue lysates, other suitable samples are:

- > Frozen cell pellets with 5 million cells
- > Frozen tissues, 50-100 mg (usually 75 mg)

**All samples must be frozen and shipped to us on dry ice.**

## EUROPEAN OFFICES

### BELGIUM

EUROGENTEC HEADQUARTERS  
TEL.: +32 4 372 74 00  
FAX: +32 4 372 75 00  
info@eurogentec.com

EUROGENTEC BIOLOGICS  
TEL.: +32 4 372 74 00  
FAX: +32 4 365 16 04  
info.biologics@eurogentec.com

### UNITED KINGDOM

EUROGENTEC LTD.  
TEL.: +44 1794 511 411  
FAX: +44 1794 522 417  
info.uk@eurogentec.com

### FRANCE

EUROGENTEC FRANCE S.A.S.U.  
TEL.: +33 2 41 73 33 73  
FAX: +33 2 41 73 10 26  
info@eurogentec.com

### GERMANY

EUROGENTEC DEUTSCHLAND GMBH  
TEL.: +49 221 258 94 55  
FAX: +49 221 258 94 54  
info@eurogentec.com

### THE NETHERLANDS

EUROGENTEC NEDERLAND B.V.  
TEL.: +31 43 352 06 98  
FAX: +31 43 354 19 65  
info@eurogentec.com

### SWITZERLAND

EUROGENTEC S.A.,  
SUCCURSALE DE GENÈVE

**FRANÇAIS:**  
TEL.: +32 4 372 74 00  
FAX: +32 4 372 75 00

**DEUTSCH:**  
TEL.: +49 221 258 94 55  
FAX: +49 221 258 94 54  
info@eurogentec.com

### CENTRAL EUROPE

REPRESENTED BY  
EUROGENTEC HEADQUARTERS  
TEL.: +32 4 372 74 00  
FAX: +32 4 372 75 00  
info@eurogentec.com

## ANTIBODY ARRAY PROFILING SERVICES

Description	Reference
Angiogenesis Antibody Array Service	AR-AGE125-SE
Apoptosis Antibody Array Service	AR-APP069-SE
Cancer Marker Antibody Array Service	AR-ACM083-SE
Cell Cycle Antibody Array Service	AR-ACC058-SE
Cytokine Related Antibody Array Service	AR-ACK075-SE
Hematopoiesis Antibody Array Service	AR-AHT049-SE
Hormone Antibody Array Service	AR-AHM030-SE
Signal Transduction Antibody Array Service	AR-AST160-SE
Stem Cell Antibody Array Service	AR-ASC048-SE
Explorer Antibody Array Service	AR-ASB600-SE
AKT Pathway Phospho Antibody Array Service	AR-PAA137-SE
AKT/PKB Signaling Phospho Antibody Array Service	AR-PAB216-SE
Cancer/Apoptosis Phospho Antibody Array Service	AR-PAC155-SE
AMPK Signaling Antibody Array Service	AR-PAM174-SE
Apoptosis Phospho Antibody Array Service	AR-PAP247-SE
Cell Cycle Phospho Antibody Array Service	AR-PCC076-SE
Cell Cycle Control Phospho Antibody array Service	AR-PCC238-SE
Cytoskeleton Phospho Antibody Array Service	AR-PCP095-SE
Cytoskeleton II Phospho Antibody Array Service	AR-PCP141-SE
CREB Antibody Array Service	AR-PCR174-SE
Chromatin/Transcription Phospho Antibody Array Service	AR-PCT173-SE
PDGF Antibody Array Service	AR-PDG195-SE
EGF Pathway Antibody Array Service	AR-PEG214-SE
Erk Signaling Antibody Array Service	AR-PEK208-SE
ErbB/HER Signaling Service	AR-PER239-SE
Phospho Explorer Antibody Array Service	AR-PEX100-SE
p53 Signaling Antibody Array Service	AR-PFT196-SE
FGF Pathway Antibody Array Service	AR-PGF156-SE
GPCR Signaling to MAPK/ERK Antibody Array Service	AR-PGP193-SE
Insulin/Glucose Phospho Antibody Array Service	AR-PIG085-SE
Insulin Receptor Phospho-Specific Antibody Array Service	AR-PIG219-SE
IGF-IR Signaling Antibody Array Service	AR-PIR245-SE
Jak/Stat Pathway Phospho Antibody Array Service	AR-PJS042-SE
Jak/Stat II Phospho Antibody Array Service	AR-PJS202-SE
MAPK Pathway Phospho Antibody Array Service	AR-PMK185-SE
mTOR Phospho Antibody Array Service	AR-PMT138-SE
NF-kB Pathway Phospho Antibody Array Service	AR-PNK063-SE
NF-kB II Signaling Phospho Antibody Array Service	AR-PNK215-SE
Nuclear/Membrane Receptor Phospho Antibody Array Service	AR-PNR052-SE
Neuroscience Phospho Antibody Array Service	AR-PNS032-SE
Tyrosine Phosphorylation ProArray Antibody Array Service	AR-PST228-SE
T-Cell Receptor Signaling Antibody Array Service	AR-PTC188-SE
TGF-beta Signaling Antibody Array Service	AR-PTG176-SE
Tyrosine Kinase Adaptor Phospho Antibody Array Service	AR-PTK098-SE
VEGF Pathway Antibody Array Service	AR-PVE185-SE

## Ordering information

**By fax:** +32 (0)4 264 07 88

**By e-mail:** order@eurogentec.com

### Eurogentec

LIEGE Science Park  
Rue du Bois Saint-Jean, 5  
4102 Seraing - Belgium  
TVA- (BE)-0427.348.346  
RPM Liège



Eurogentec is a distributor of Full Moon BioSystems Products and Services

