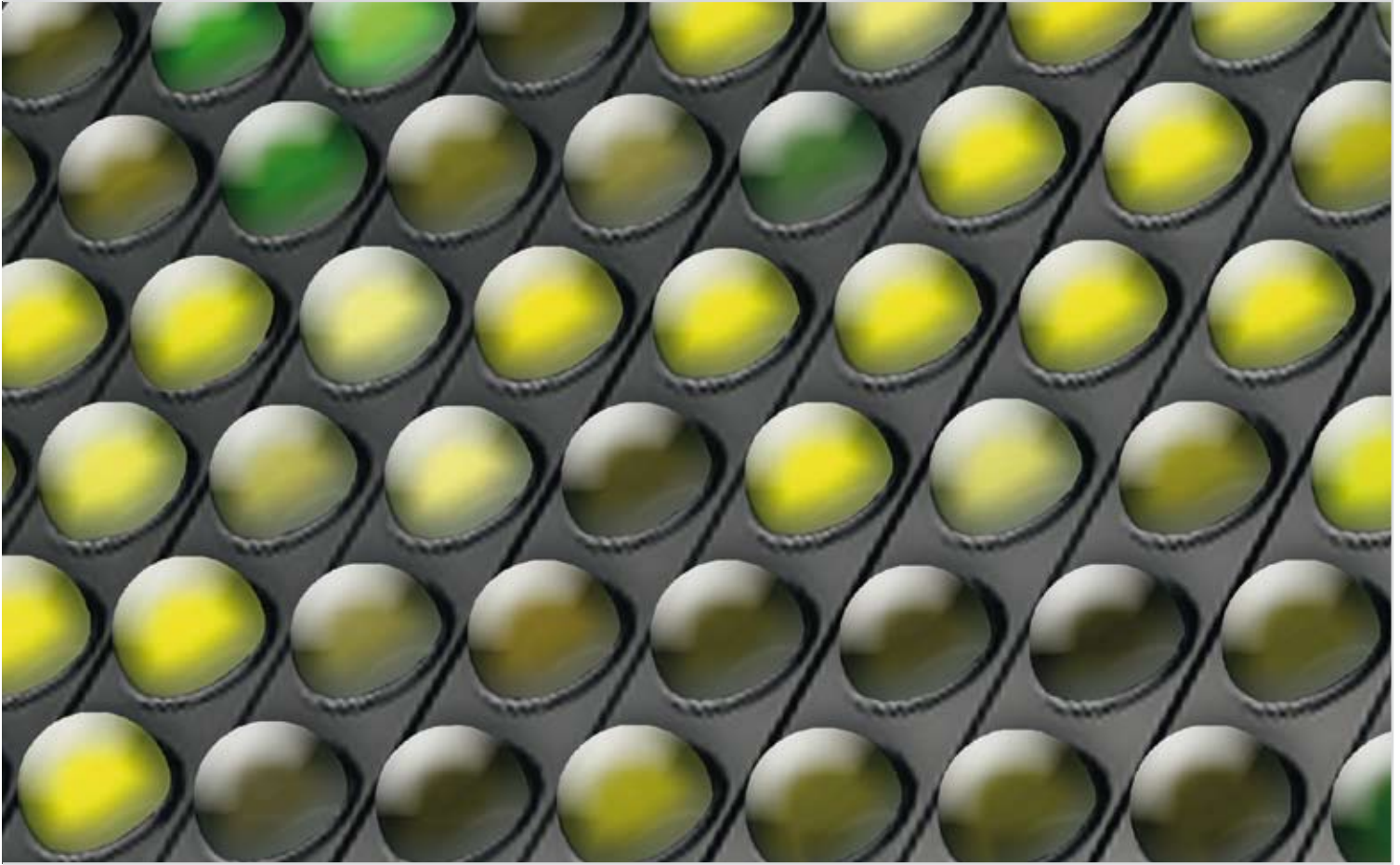


# ANTIBODY MICROARRAYS



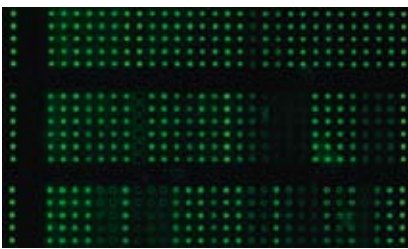
DON'T BE STUCK ANALYSING  
SINGLE PROTEIN EXPRESSION AT A TIME!



# 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.



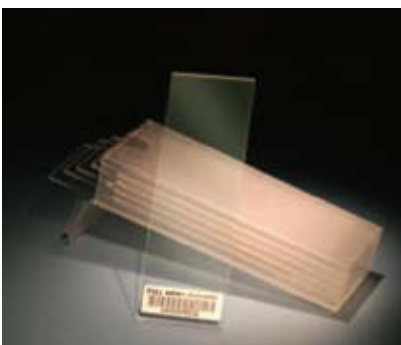
- > The antibodies are printed on standard-size microscope slides.
- > Each slide consists of an array of well-characterized antibodies with six replicates.
- > Multiple positive and negative controls are present.
- > Each antibody array kit comes with two identical slides.

## Applications :

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

## Specifications

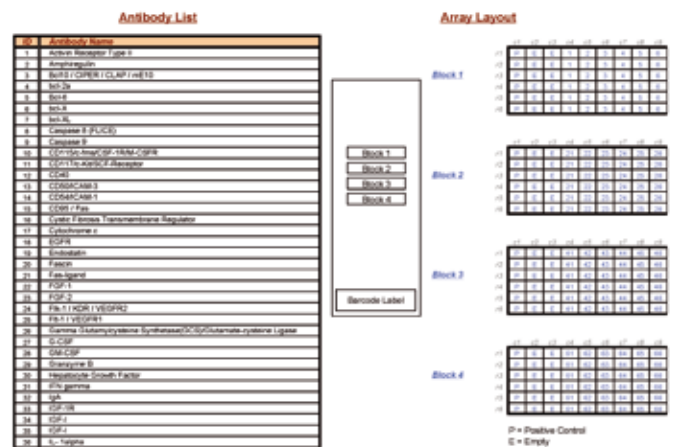
- > Spot diameter: 260-280 µm
- > Slide dimensions: 76 x 25 x 1 mm
- > Spot-to-Spot distance 450-600 µm
- > Barcode label: 23 x 9 mm
- > Storage conditions: 4 °C (6 months)



## 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®3 (Cy5®/Alexa®532/Alexa®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 µg). Please inquire for reference and price information.



Snapshot example of the layout of Antibody Arrays

The complete list of Antibodies printed on the focused Arrays is available on our website [www.eurogentec.com](http://www.eurogentec.com)

## 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.

### 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

### EXPLORER, THEMATIC & PHOSPHORYLATION ANTIBODY ARRAYS AND ASSAY KITS

Description	Reference	Species	Number of Ab	Number of slides	Price		
					EUR	CHF	GBP
Angiogenesis Antibody Array	AR-AGE1-25	Human	127	2	576.00	875.50	480.00
Apoptosis Antibody Array	AR-APP0-69	Human	71	2	428.80	651.80	357.30
Cancer Marker Antibody Array	AR-ACM0-83	Human	84	2	480.00	729.60	400.00
Cell Cycle Antibody Array	AR-ACC0-58	Human	58	2	390.40	593.40	325.30
Cytokine Related Antibody Array	AR-ACK0-75	Human	77	2	448.00	681.00	373.30
Hematopoiesis Antibody Array	AR-AHT0-49	Human	49	2	352.00	535.00	293.30
Hormone Antibody Array	AR-AHM0-30	Human	30	2	281.60	428.00	234.70
Signal Transduction Antibody Array	AR-AST1-60	Human	163	2	665.60	1011.70	554.70
Stem Cell Antibody Array	AR-ASC0-48	Human	48	2	384.00	583.70	320.00
AKT Pathway Phospho Antibody Array	AR-PAA1-37	Human, Mouse	137	2	595.20	904.70	496.00
Cancer/Apoptosis Phospho Antibody Array	AR-PAC1-55	Human, Mouse	155	2	633.60	963.10	528.00
Cell Cycle Phospho Antibody Array	AR-PCC0-76	Human, Mouse	76	2	441.60	671.20	368.00
Chromatin/Transcription Phospho Antibody Array	AR-PCT1-73	Human, Mouse	173	2	672.00	1021.40	560.00
Cytoskeletal Pathway Phospho Antibody Array	AR-PCP0-95	Human, Mouse	95	2	505.60	768.50	421.30
Insulin/Glucose Phospho Antibody Array	AR-PIG0-85	Human, Mouse	85	2	467.20	710.10	389.30
Jak/Stat Pathway Phospho Antibody Array	AR-PJS0-42	Human, Mouse	42	2	326.40	496.10	272.00
MAPK Pathway Phospho Antibody Array	AR-PMK1-85	Human, Mouse	185	2	704.00	1070.10	586.70
Neuroscience Phospho Antibody Array	AR-PNS0-32	Human, Mouse	32	2	288.00	437.80	240.00
NF-kappa B Pathway Phospho Antibody Array	AR-PNK0-63	Human, Mouse	63	2	403.20	612.90	336.00
Nuclear/Membrane Receptor Phospho Antibody Array	AR-PNR0-52	Human, Mouse	52	2	364.80	554.50	304.00
Tyrosine Kinase Adaptor Phospho Antibody Array	AR-PTK0-98	Human, Mouse	98	2	505.60	768.50	421.30
Explorer Antibody Array	AR-ASB600	Human	656	2	1018.00	1547.40	848.30
Antibody Array Assay Kit for Protein Extraction and Labelling (2 reactions)	AR-KAS0-02	-	-	-	382.70	581.70	318.90
Antibody Array Assay Kit for Protein Extraction and Labelling (20 reactions)	AR-KAS0-20	-	-	-	3187.20	4844.50	2656.00

## 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!

### 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

- > 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)

### 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.**

### ANTIBODY ARRAY PROFILING SERVICES

Description	Reference	EUR	CHF	GBP
Angiogenesis Antibody Array Profiling Service	AR-AGE125-SE	1951.70	2966.60	1626.40
Apoptosis Antibody Array Profiling Service	AR-APP069-SE	1797.90	2732.80	1498.20
Cancer Marker Antibody Array Profiling Service	AR-ACM083-SE	1797.90	2732.80	1498.20
Cell Cycle Antibody Array Profiling Service	AR-ACC058-SE	1797.90	2732.80	1498.20
Cytokine Related Antibody Array Profiling Service	AR-ACK075-SE	1797.90	2732.80	1498.20
Hematopoiesis Antibody Array Profiling Service	AR-AHT049-SE	1644.00	2498.90	1370.00
Hormone Antibody Array Profiling Service	AR-AHM030-SE	1644.00	2498.90	1370.00
Signal Transduction Antibody Array Profiling Service	AR-AST160-SE	2105.60	3200.50	1754.60
Stem Cell Antibody Array Profiling Service	AR-ASC048-SE	1644.00	2498.90	1370.00
Explorer Antibody Array Profiling Service	AR-ASB600-SE	2413.30	3668.20	2011.10
AKT Pathway Phospho Antibody Array Profiling Service	AR-PAA137-SE	1951.70	2966.60	1626.40
Cancer/Apoptosis Phospho Antibody Array Profiling Service	AR-PAC155-SE	2105.60	3200.50	1754.60
Cell Cycle Phospho Antibody Array Profiling Service	AR-PCC076-SE	1797.90	2732.80	1498.20
Chromatin/Transcription Phospho Antibody Array Profiling Service	AR-PCT173-SE	2105.60	3200.50	1754.60
Cytoskeletal Pathway Phospho Antibody Array Profiling Service	AR-PCP095-SE	1797.90	2732.80	1498.20
Insulin/Glucose Phospho Antibody Array Profiling Service	AR-PIG085-SE	1797.90	2732.80	1498.20
Jak/Stat Pathway Phospho Antibody Array Profiling Service	AR-PJS042-SE	1644.00	2498.90	1370.00
MAPK Pathway Phospho Antibody Array Profiling Service	AR-PMK185-SE	2105.60	3200.50	1754.60
Neuroscience Phospho Antibody Array Profiling Service	AR-PNS032-SE	1644.00	2498.90	1370.00
NF-kappa B Pathway Phospho Antibody Array Profiling Service	AR-PNK063-SE	1797.90	2732.80	1498.20
Nuclear/Membrane Receptor Phospho Antibody Array Profiling Service	AR-PNR052-SE	1797.90	2732.80	1498.20
Tyrosine Kinase Adaptor Phospho Antibody Array Profiling Service	AR-PTK098-SE	1797.90	2732.80	1498.20



## Arrays available for assay service:

### Pathway Antibody Arrays

Angiogenesis Array  
 Apoptosis Array  
 Cancer Marker Array  
 Cell Cycle Array  
 Cytokine Array  
 Explorer Array  
 Hematopoiesis Array  
 Hormones Array  
 Signal Transduction Array  
 Stem Cell Array

### Phospho-specific Antibody Arrays

Phospho AKT Pathway Array  
 Phospho Cancer/Apoptosis Array  
 Phospho Cell Cycle Array  
 Phospho Chromatin/Transcription Array  
 Phospho Cytoskeletal Pathway Array  
 Phospho Insulin/Glucose Array  
 Phospho JAK/STAT Pathway Array  
 Phospho MAPK Pathway Array  
 Phospho Neuroscience Array  
 Phospho NF- $\kappa$ B Pathway Array  
 Phospho Nuclear/  
 Membrane Receptors Array  
 Phospho Tyrosine Kinase Adaptors Array



## Ordering information

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

**By e-mail:** [order@eurogentec.com](mailto:order@eurogentec.com)

**By mail:**

**Eurogentec**

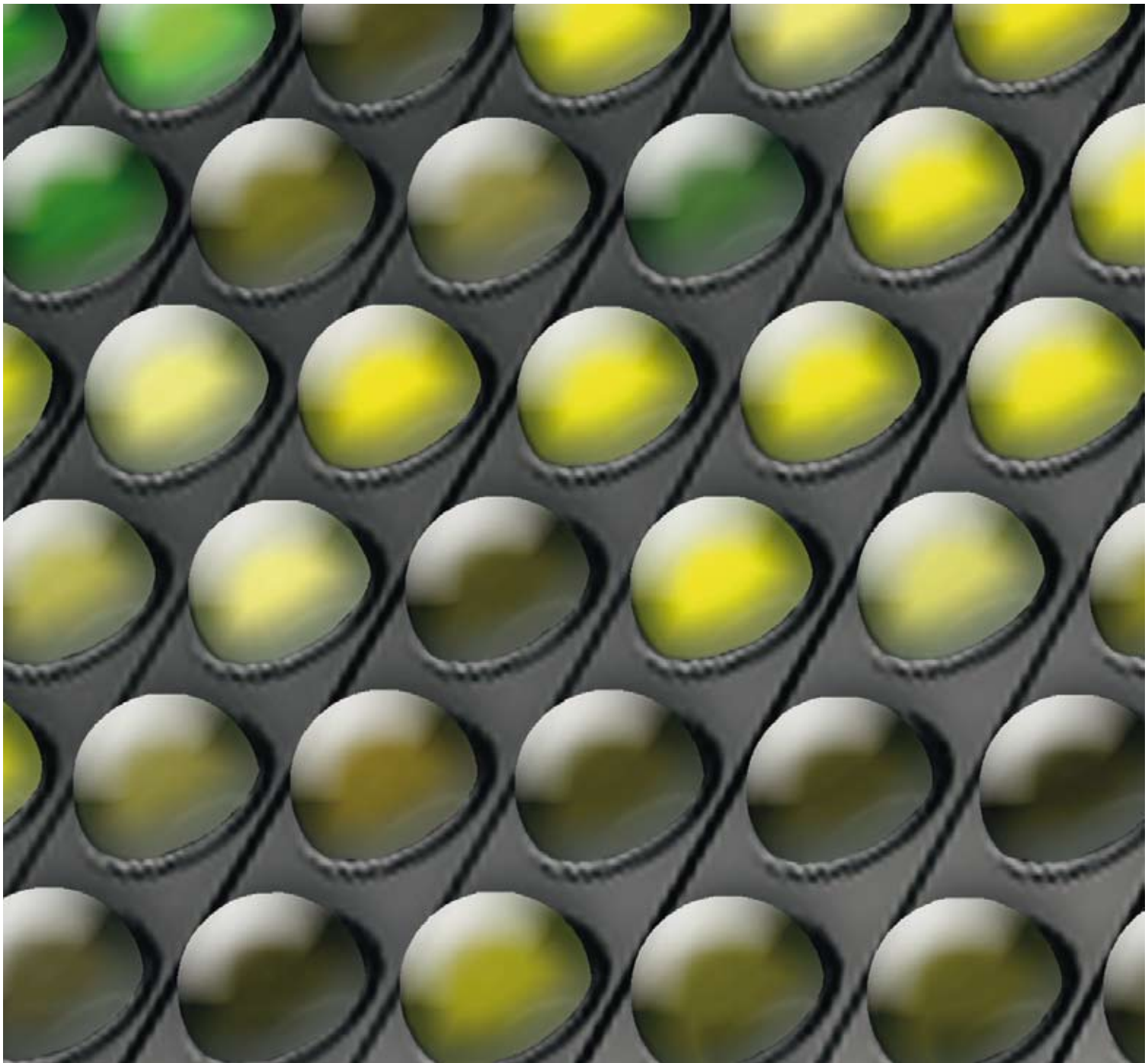
Liege Science Park  
 rue du Bois Saint-Jean, 5  
 4102 Seraing  
 Belgium  
 TVA- (BE)-0427.348.346  
 RPM Liège



**FULL MOON**  
*BioSystems*

Eurogentec is a distributor of Full Moon BioSystems Products and Services





GDA-FULLMOON-0409-V4

**EUROGENTEC S.A.**

Rue Bois St. Jean 5  
4102 Seraing, Belgium  
Tel: +32 (0)4 372 74 00  
Fax: +32 (0)4 372 75 00  
info@eurogentec.com  
www.eurogentec.com

 **Eurogentec**  
Experience true partnership