

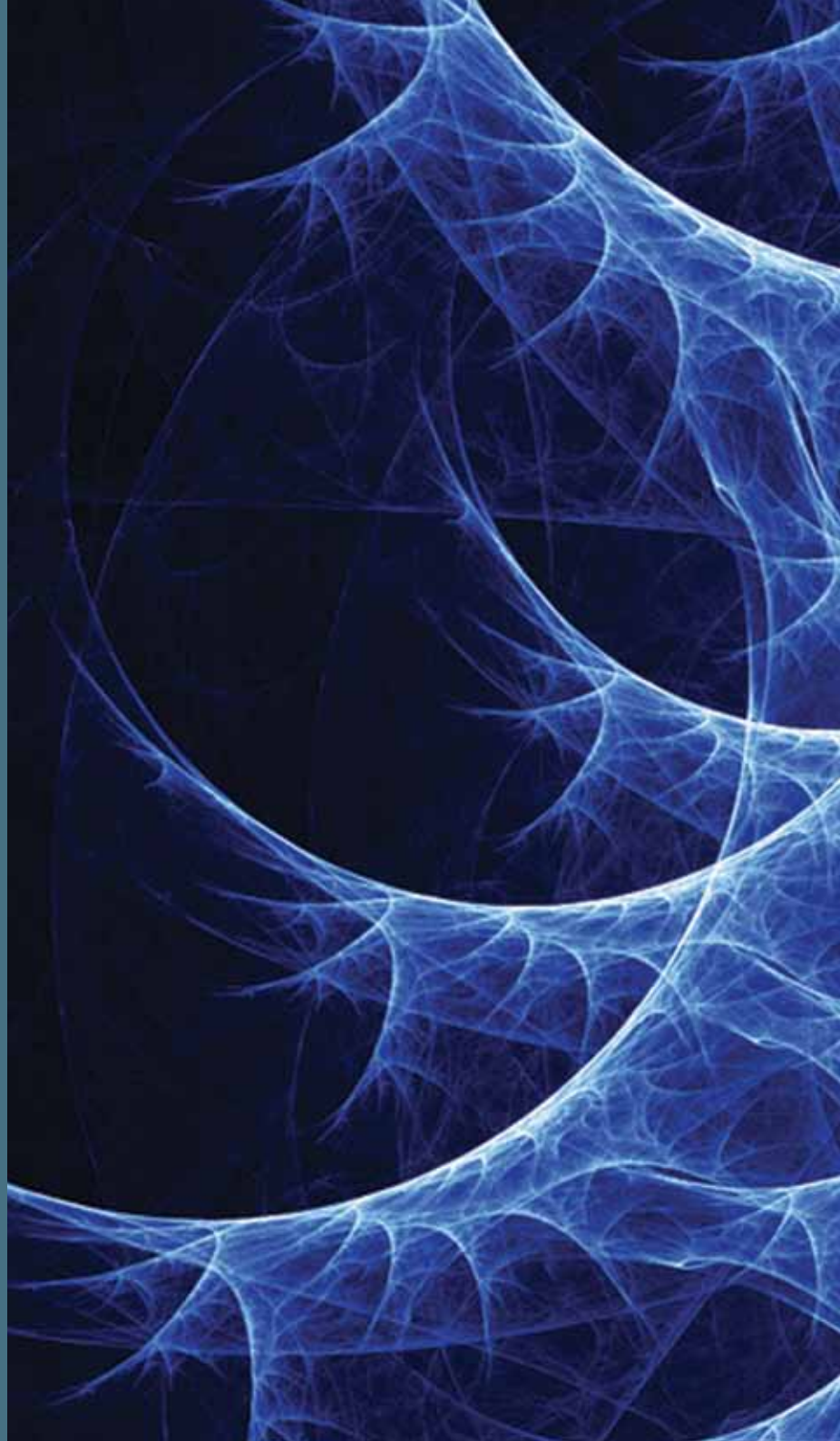
<b>1</b>	<b>General information</b>	<b>3</b>
<b>2</b>	<b>Genomic solutions</b>	
2.1	Oligonucleotides	20
2.2	Custom genes	112
2.3	Real-Time qPCR	116
2.4	Nucleic acid amplification	136
2.5	Gene cloning and expression in <i>E. coli</i>	150
2.6	Mammalian cell transfection reagents	168
2.7	Gene analysis	172
<b>3</b>	<b>Proteomic solutions</b>	
3.1	Custom polyclonal antibodies	180
3.2	Custom monoclonal antibodies	198
3.3	Fluorescent dyes	204
3.4	Catalogue antibodies	214
3.5	Catalogue peptides	236
3.6	Custom peptides	238
3.7	Sensolyte	246
3.8	Multiplex assay system	266
3.9	Custom protein expression service	272
3.10	Proteomic synthesis reagents	274
<b>4</b>	<b>Index</b>	<b>276</b>

# Table of contents

# 1

## General information

- 8 Ordering information
- 10 Company profile
- 12 Eurogentec offices
- 14 Registered trademarks & license statements
- 18 General conditions of sale



# 1 General information

<b>Table of contents</b>	<b>1</b>
<b>Ordering information</b>	<b>8</b>
<b>Company profile</b>	<b>10</b>
<b>Eurogentec offices</b>	<b>12</b>
<b>Registered trademarks &amp; license statements</b>	<b>14</b>
<b>General conditions of sale</b>	<b>18</b>

# 2 Genomic solutions

<b>2.1 Oligonucleotides</b>	<b>20</b>
<b>General Information</b>	<b>21</b>
<b>Life Science Oligonucleotides</b>	<b>24</b>
Oli&GO 1.2.3™	27
Custom Oligonucleotides	28
Backbone	28
Bases	28
Linkages	30

Modifications	31	2' O-Methyl RNA Molecular Beacons	55
Overview	31	Scorpions® Probes	56
Phosphate	32	LC Hybridisation probes	58
Biotins	32	Plexor™ Primers	58
Conjugates	33	NASBA®	58
Modifiers	35	MLPA®	60
Spacers	36	Multifinder	60
Non-Natural Bases	36	Real-Time qPCR	60
Degenerate Bases	39	RNAi oligonucleotides	62
Fluorophores	39	siRNA	62
Quenchers	43	Custom siRNA Duplexes	62
Purifications	46	Control siRNA Duplexes	63
SePOP desalting	46	miRNA	64
RP-Cartridge•Gold™	46	Clear-MiR™ miRNA Inhibitors	64
HPLC	47	Add-Mir™ miRNA Mimics	64
PAGE	47	Universal Primers	65
Dual HPLC	47	Unique Oligonucleotides	66
UltraPureGold™	48	Minimum Guaranteed Yields	67
Additional Services	48	Shipping	68
Design	48	Documentation	69
Quality Control (QC)	48	References	70
Format	49	<b>Clinical Lab Oligonucleotides (CLOs)</b>	
Packaging	50	<b>for European Lab Services</b>	<b>105</b>
Shipping	50	Increasing in-house molecular test	105
Real-time qPCR probes	51	CLOs benefits	105
Double-Dye Oligonucleotides	51	CLOs GMP manufacturing	105
Double-Dye Probes	51	<b>cGMP Oligonucleotides</b>	<b>106</b>
LNA® Double-Dye Probes	53	Experience true partnership with Eurogentec	106
Molecular Beacons	54		
Wavelength-Shifting Molecular Beacons	55		

Complete Custom Solutions: Benefits of a Eurogentec partnership	106
GMP & Regulatory Compliance	106
The Eurogentec Solution	107
To GMP or not to GMP	107
Manufacturing Processes & Options - Flexible Custom Solutions	108
Manufacturing Processes	108
Manufacturing Options	109
Classified clean rooms/ exceeding regulatory requirements	110
Risk Management & Quality Assurance	110
Partnering with Eurogentec: The Process	111

**2.2 Custom genes 112**

<b>Introduction</b>	<b>113</b>
<b>Case Study</b>	<b>113</b>
<b>Process Overview</b>	<b>114</b>

**2.3 Real-Time qPCR 116**

<b>Kits and Reagents</b>	<b>117</b>
Kits for every application	117
Kits in every format	117
Carry-over prevention	118
High performance enzyme	118
Custom Mixes	118
Quality Control	118
<b>qPCR kits &amp; platforms compatibility tables</b>	<b>119</b>
<b>qPCR Miscellaneous</b>	<b>123</b>

Uracil-N-glycosylase	123
Universal RT kit for qPCR	123
Fluorescein additive	123
ROX passive reference	123

**Controls 124**

Universal Exogenous qPCR Positive Control (IPC)	124
Housekeeping Internal Controls	125

**Plate and seals 126**

qPCR 96-well plates	127
qPCR 384-well plates	128
qPCR optical caps and seals	128

**Assay Dispensing Services 134**

**2.4 Nucleic Acid Amplification Reagents 136**

**DNA Polymerases 137**

PCR Enzyme Selection Guide	137
Diamond <i>Taq</i> <sup>®</sup> Family	138
Diamond <i>Taq</i> <sup>®</sup> Family benefits	138
Diamond <i>Taq</i> <sup>®</sup> Family specifications	138
Quality	139
Professional support for manufacturers of Diagnostic Assays	139
Diamond & Red Diamond <i>Taq</i> <sup>®</sup>	140
Key Benefits	140
Additional information	140
Performance test	140
Hot Diamond <i>Taq</i> <sup>®</sup>	141
Key Benefits	141
Additional information	141
Performance test	141

HGS Diamond <i>Taq</i> <sup>®</sup>	142
Key Benefits	142
Additional information	142
Performance test	142
DAP GoldStar <sup>®</sup> DNA Polymerase	143
Key Benefits	143
Applications	143
Additional information	143
SilverStar DNA Polymerase	144
Key Benefits	144
Applications	144
Additional information	144
Performance test	144
AccuStar <sup>™</sup> DNA Polymerase	145
Key Benefits	145
Additional information	145
Performance test	145
Booster Mix	145
Key Benefits	145
Activity	145
Authentik <sup>™</sup> DNA Polymerase	146
Key Benefits	146
Additional information	146
Tth DNA Polymerase	146
Additional information	146

**Ready-to-use PCR Mixes 147**

Key benefits	147
GoldStar <sup>®</sup> Mix / Red'y'GoldMix	147
HotGoldStar Mix / Red'y'StarMix	147

**Reverse Transcriptase 148**

Mu-MLV Reverse Transcriptase	148
------------------------------	-----

<b>Nucleotides</b>	<b>149</b>
dNTP Mix	149
dNTP Set	149

## 2.5 Gene cloning & expression in *E. coli* 150

<b>General Principle of the Staby™ Operating System</b>	<b>150</b>
Introduction	150
Staby™ Operating System application	150
Gene cloning	150
Protein expression	150

<b>Gene Cloning</b>	<b>152</b>
StabyCloning™Kit	152
Principle	152
Results	152
Vector map	153
Staby™ SubCloning Kits	154
Principle	154
Vector maps	154
GetStaby™ Kit	156
Description	156
Vector map	156
Benefits	156

<b>Protein Expression</b>	<b>157</b>
Cherry™Express Kit	157
Vector map	158
Cherry™Codon kit	159
Vector map	159
StabyExpress™ T7 kit	160
Vector Map	160
Staby™Codon T7 kit	162

Vector Map	162
<b>Auto-inducible medium</b>	<b>164</b>
Staby-Switch	164

<b>Competent bacteria</b>	<b>165</b>
For Staby™Technology	165
CYS21 Strain	165
SE1 strain	165
For General Cloning and Expression	166
DG1 Strain	166
BL21 strain	166

<b>Electroporation Cuvettes</b>	<b>167</b>
---------------------------------	------------

## 2.6 Mammalian cell transfection reagents 168

<b>DNA transfection: ICAFectin™ 441</b>	<b>169</b>
<b>siRNA transfection: ICAFectin™ 442</b>	<b>170</b>

## 2.7 Gene analysis 172

<b>Nucleic Acid Extraction Kits</b>	<b>173</b>
Kits for total RNA Extraction	173
Kits For Total RNA, DNA and Protein Extraction	173
Kits for DNA Extraction	173
<b>Mupid®-One electrophoresis system</b>	<b>174</b>
<b>Agarose, Molecular Biology Grade</b>	<b>175</b>
<b>Molecular Weight DNA Ladders</b>	<b>176</b>
SmartLadders	176
SmartLadder	176
SmartLadder SF	176
BioMarker®	176

Marker for Genotyping, RFLP, Tilling	177
MapMarker® 400	177
MapMarker® 1000	177

# 3 Proteomic solutions

## 3.1 Custom Polyclonal Antibodies 180

<b>General overview</b>	<b>181</b>
<b>Choosing your parameters</b>	<b>182</b>
Selecting your antigen	182
Antibody titer evolution	182
Antibody affinity evolution	182
Protein, Peptide or Genetic Immunisation	183
<b>Selecting your animal host</b>	<b>184</b>
<b>Selecting your immunisation protocol</b>	<b>185</b>
Speedy 28-day polyclonal	185
The classical 87-day protocol	187
<b>Ready to start an immunisation programme?</b>	<b>188</b>
Anti-peptide antibodies	188
Anti-protein antibodies	188
Construct your own programme	188
<b>Post Translational Modifications or Epigenomic Events detection</b>	<b>189</b>
Peptide immunisation protocols	190
<b>DNA immunisation</b>	<b>193</b>
<b>More for Your Polyclonal Programmes</b>	<b>195</b>
<b>Magnetic beads coupling services</b>	<b>196</b>

**3.2 Custom Monoclonal 198**

<b>Mouse Monoclonal Antibody Development</b>	<b>199</b>
Options	200
Additional quality control	200
<b>In vitro Monoclonals expression service</b>	<b>200</b>
<b>Mouse monoclonal antibody additional services</b>	<b>201</b>
Antibody labelling	201
Fragmentation	201
Assay development	201
Cell banking	201
<b>Magnetic beads coupling services</b>	<b>202</b>

**3.3 Fluorescent dyes & protein labelling kits 204**

<b>HiLyte Fluor™ Dyes</b>	<b>205</b>
Superior fluorescent labelling dyes spanning the full visible spectrum	205
HiLyte Fluor™ 488 Dye	206
HiLyte Fluor™ 555 Dye	207
HiLyte Fluor™ 594 Dye	208
HiLyte Fluor™ 647 Dye	209
HiLyte Fluor™ 680 Dye	210
HiLyte Fluor™ 750 Dye	210
<b>AnaTag™ protein labelling kits</b>	<b>211</b>
Easy labelling with AnaTag™ protein labelling kits	211
AnaTag™ HiLyte Fluor™ protein labelling kits	211
Biotins, Avidins and Streptavidins	212
Phycobiliproteins and Their Conjugates	212
Reactive fluorescent dyes	213

**3.4 Catalogue Antibodies 214**

<b>Catalogue Antibodies</b>	<b>215</b>
<b>Neuroscience antibodies</b>	<b>216</b>
Alzheimer's Disease Research	216
Anti-Tau Antibodies	216
Anti-β-Amyloid Antibodies and ELISA kits	218
BetaMark™ ELISA kits	219
Other Alzheimer's disease Related Antibodies	220
Multiple Sclerosis: Anti-MOG & Anti-PLP ELISA kits	221
Anti-MOG ELISA kit	221
Anti-PLP ELISA kit	221
Parkinson's Disease Research	222
Synuclein Antibodies	222
DJ-1 Antibodies	222
Parkin Antibodies	223
Prion Antibodies	223
Ubiquitous Antibodies	223
Apolipoprotein Antibodies	223
Ubiquitin Antibodies	223
Heat Shock Proteins Antibodies	224
Neuronal receptors antibodies	224
Neuronal markers antibodies	226
<b>Immunopathology products</b>	<b>229</b>
<b>Detection chemistries</b>	<b>230</b>
<b>Ancillaries</b>	<b>230</b>
<b>Epitope Tags and Fusion Protein Antibodies</b>	<b>231</b>
<b>Cellular components</b>	<b>232</b>
5-methylcytidine monoclonal antibody	232
Other components	232

**Labelled Secondary Antibodies 233**

Z-Fish™ Antibodies 234

**3.5 Catalogue peptides 236**

<b>Catalogue Peptides (Go™ Peptides)</b>	<b>237</b>
Amyloid Peptides	237
MOG Peptides	237
CEF Control Peptides	237
Histone H3 (1-21), Biotinylated	237
FRET protease substrates	237

**3.6 Custom peptides 238**

<b>Peptide Manufacturing Process</b>	<b>239</b>
<b>Custom Peptides modifications</b>	<b>241</b>
Conjugations-Labeling	241
Terminal Modifications	241
Structural Modifications	241
Amino acid Modifications	241
<b>Peptide purity levels &amp; applications</b>	<b>242</b>
Crude peptides / SePOP purified	242
Immunograde (> 70 %)	242
Biochemistry grade (> 85 %)	242
Very high purity (> 95 %)	242
<b>FRET peptides</b>	<b>243</b>
<b>ClearPoint™ peptides</b>	<b>244</b>
<b>Custom peptides librairies</b>	<b>245</b>

**3.7 Protein detection 246**

<b>Sensolyte® Assay kits</b>	<b>247</b>
Alzheimer disease	248
Oxidative stress	249
Epigenetics	250
Tissue Remodeling: MMPs	251
Apoptosis: Caspase	253
Bone-related proteolytic pathways & diseases	254
Vascular diseases	255
Viral diseases	256
ELISA & Gene expression assays	257
Cell proliferation, viability and cytotoxicity assays	259
Generic Protease Assay kits	260

<b>Duolink®</b>	<b>261</b>
Visualize and locate protein interactions	261

**3.8 Multiplex Assay Detection 266**

<b>Luminex®</b>	<b>267</b>
xMAP® Multiplexing assay development & testing services	267
Luminex® Epitope Mapping Service	269

<b>Full Moon Antibody Arrays</b>	<b>270</b>
Pathway & Phosphorylation Antibody MicroArrays	270
Complete Assay Service	271
Antibody Array Assay kit	271

**3.9 Custom Protein Expression 272****3.10 Peptide Synthesis Reagents 274**

<b>Resins</b>	<b>275</b>
<b>Reagents</b>	<b>275</b>
<b>Unusual Amino Acids</b>	<b>275</b>
<b>Amino Alcohols</b>	<b>275</b>
<b>Amino Acid Cartridges</b>	<b>275</b>
<b>Accessories</b>	<b>275</b>

**4 Index 276**

# Ordering information

## Products and services

From now on, you can order all your Oligonucleotides and Kits & Reagents via the Eurogentec Ordering System (EOS). You can continue to order all Products & Services via e-mail or fax as before.

### Oligonucleotides

Custom Oligonucleotides  
Real-Time qPCR Probes  
RNAi Oligonucleotides  
Universal Primers

### Kits & Reagents

Oli&GO 1.2.3™ Activations & Reloads  
Cloning & Expression Kits  
Transformation Products  
Transfection Reagents  
DNA Polymerases  
PCR Mixes  
Molecular Weight DNA Ladders  
Electrophoresis System  
Real-Time qPCR Kits  
Real-Time qPCR Plastics  
Catalogue Peptides  
Catalogue Antibodies

### Custom & Unique Services

Custom Genes  
Custom Peptides  
Custom Monoclonal Antibodies  
Custom Polyclonal Antibodies  
Custom Protein Expression



Unique Oligonucleotides  
Unique Peptides  
Unique Polyclonals



**Eurogentec**  
**Ordering System (EOS)**  
**www.eurogentec.com**



#### Continental Europe

✉ order@eurogentec.com  
☎ +32 4 264 07 88

#### United Kingdom

✉ order.uk@eurogentec.com  
☎ +44 1794 522 417

## Required information

When placing your order, please include the following information:

- ▣ client account number
- ▣ contract number / quotation number
- ▣ contact person
- ▣ delivery address
- ▣ invoice address
- ▣ telephone number
- ▣ product description and reference number
- ▣ quantity
- ▣ purchase order number

## Prices and sales terms

Prices are subject to change without prior notice. All prices are given excluding VAT and transport costs. At any time, please feel free to contact Eurogentec for the latest prices or the most recent price list.

Payment terms are 30 days net from the invoice date.

Eurogentec products conform, at the time of delivery and for 30 days thereafter, to the specifications specifically agreed by Eurogentec.

This warranty is limited to the replacement of the product.

No other warranty, express or implied, is provided by Eurogentec.

Eurogentec products are sold for research or laboratory use only and are not to be administered to humans or used for medical diagnostics.

## Bulk discounts and contracts

A number of products in this catalogue can be supplied in larger quantities than those listed. Please contact us for a price offer and delivery terms.

Special discounts are negotiable for annual volume contracts. Please contact our local sales representatives for the contract terms.

## Shipping

Upon receipt of an order before 14.00 h, the resales products in stock are shipped the same day. For custom products, production will start upon receipt of your order. Products are delivered by overnight parcel delivery.

Excepting special circumstances, products should reach our client within 24 hours of expedition.

No shipping fee is charged under certain conditions.

See our general conditions of sales

## Customer support

▣ If you need general information please contact us at our headquarters:

✦ Tel.: +32 4 372 74 00

✦ [info@eurogentec.com](mailto:info@eurogentec.com)

▣ If you need more precise or technical information please contact one of the following persons:

✦ Oligonucleotides

Oligo Centre

Tel.: 00 800 666 00 123

(European Toll Free number)

[oligo@eurogentec.com](mailto:oligo@eurogentec.com)

✦ Equipment and consumables

[customer.care@eurogentec.com](mailto:customer.care@eurogentec.com)

✦ Genomics and general technical information

Molecular Biology, Real-Time qPCR, PCR, Enzymes

[scientific.support@eurogentec.com](mailto:scientific.support@eurogentec.com)

✦ Proteomics & general technical information

Peptides, custom antibody

services (polyclonal & monoclonal, gene immunization), custom proteomics services:

[scientific.support@eurogentec.com](mailto:scientific.support@eurogentec.com)

✦ Custom Genes

[gene@eurogentec.com](mailto:gene@eurogentec.com)

# Company profile

Eurogentec was founded in 1985 as a spin-off from the University of Liège (Belgium). Eurogentec S.A. is a leading global supplier of innovative reagents, kits, specialty products and custom services to scientists in life science, biotechnology, diagnostic and pharmaceutical markets.

Experts in small and large-scale DNA, RNA and peptide synthesis, Eurogentec is an experienced Contract Manufacturing Organization (CMO), and is fully ISO 9001 and ISO 13485:2003 certified and authorised by Belgian ministry of health to manufacture GMP biologics.

It operates two full-service state-of-the-art GMP-compliant manufacturing facilities in Belgium that provide high value custom biologics and a wide range of oligonucleotide-based components for diagnostic and therapeutic applications.

Eurogentec's headquarters (administrative offices and production facilities) are located in the LIEGE Science Park. The company has subsidiaries in France, Germany, the UK, the Netherlands and Switzerland and production facilities in San Diego (USA), Fremont (USA) and in Toyama (Japan). The company is also present in Singapore thanks to a partnership with AIT Biotech Pte Ltd (previously Research Biolabs).

Eurogentec recently acquired AnaSpec, a privately held proteomics company based in Fremont, California. Founded in 1993, AnaSpec is a leading provider of integrated proteomics solutions for life science research and diagnostics with expertise in peptides synthesis, labelled peptides and antibodies, fluorescent dyes and enzyme activity assays.

In 2010 Kaneka and Eurogentec S.A. joined forces. This capital alliance with Kaneka will allow Eurogentec to reinforce and to accelerate its production capacity in the Biopharmaceuticals contract manufacturing unit as well as its capacity to develop and launch highly technical products that respond to the needs of the Life Science community, including the fast growing molecular Diagnostic market.

The EGT Group currently employs over 400 people, including around 40 PhDs.

## Products and Services

Eurogentec is composed of three inter-related departments:

- Life Science Research Products and Services
- Molecular Diagnostics Manufacturing Solutions
- GMP BioManufacturing

## Life Science Products and Services

**The Research department** offers to scientists innovative solutions in Genomics and Proteomics. Genomics activities include oligonucleotide synthesis with a large range of modifications, custom gene synthesis, Real-Time qPCR, microRNA, siRNA and proposes a wide range of kits and consumables. Proteomics gathers know-how in the design of peptides, either for antibody production or bioactive peptide purposes. Antibody production in a wide range of animal models as well as Peptide-, Antibody- and Protein arrays are also key areas of expertise.

## Molecular Diagnostics Manufacturing Solutions

**The Diagnostics department** provides exceptional cGMP quality products for DNA and RNA based Molecular Diagnostics Assay. Customized manufacturing processes take place in classified clean rooms and are fully compliant with European IVD directive 98/79 EC and U.S. FDA regulations for good Manufacturing Practice (FDA 21 CFR Part 820 QSR/cGMP).

The cGMP Oligonucleotides dedicated for the *in vitro* Diagnostic (IVD) are manufactured in classified clean rooms and use an ISO 13485-certified and GMP compliant quality Management System.

Although not required for compliance to either the cGMP guidelines or the ISO 13485 QMS, Eurogentec determined that the best approach for minimizing the environmental and oligonucleotide cross-contamination risks was to build a dedicated and self-contained cleanroom facility.

## GMP BioManufacturing Proteins, Plasmids, Vaccines

**“Eurogentec Biologics”** is a full-service Contract Manufacturing Organization (CMO) offering process development, scale-up and GMP manufacturing of proteins, plasmids, vaccines and conjugates for clinical trial and commercial use.

## Quality Management Systems

### ISO 9001 certification

Since 2000, Eurogentec S.A. has maintained a Quality Management System (QMS) that is compliant and certified for ISO 9001. This certification was first obtained for ISO 9002 and then renewed for ISO 9001:2000 and ISO 9001:2008.

The certification applies to following activities: “Development, production and sales of products and services for research and development in Life Sciences”.

In addition, Eurogentec North America, Inc., a wholly owned subsidiary of the EGT Group and serving the North, Central, and South American scientific community, has been certified in compliance with ISO 9001:2000 since September 2004. The registration was then renewed for ISO 9001:2008. Its scope applies to the: “Manufacture and sales of Oligonucleotides for research and development in Life Sciences”.

### ISO 13485:2003 certification and compliance to 21 CFR Part 820

As of 2004 we have observed an increasing demand for high quality large-scale cGMP oligos for the *In Vitro* Diagnostic industry. Therefore, the following global quality vision was formulated:

Eurogentec recognizes oligonucleotides as critical components of nucleic acid diagnostic medical devices.

Therefore our facility and QMS should be compliant to the ISO 13485:2003 Medical Device standard and FDA’s current Good Manufacturing Practices (cGMP) as outlined in the 21 CFR Part 820 quality system regulations.

In the last years our QC, QA and production staff has been very successful in fulfilling our global quality vision for the production of cGMP oligonucleotides.

In December 2008, as a result, Eurogentec (Liège production site) received ISO 13485:2003 certification for the “Production and sales of *in vitro* diagnostic oligonucleotides”.

At the moment, most of our cGMP oligonucleotides are still produced in our GMP facility in Belgium. We also established production facilities in North America and Japan, with complete harmonization in equipment, methods, procedures and QMS. This creates additional safety and a high quality standard for our diagnostic customers.

In 2009 our branch in Japan received ISO 13485:2003 certification for the “Production and sales of IVD oligonucleotides”. Our North American manufacturing site in San Diego received in 2010 ISO 13485:2003 certification for the “manufacture and sales of Oligonucleotides intended for *in Vitro* Diagnostic use”.

### Directive 2003/94/EC of the European Commission

Eurogentec is accredited by the Belgian Federal Agency for Medicines and Health Products for the manufacturing of clinical trial material. We manufacture biopharmaceuticals according to the European Commission’s Directive 2003/94/EC, EudraLex Volume 4 regulations as well as the FDA’s 21 CFR Part 210 and 211 cGMP manufacturing requirements for biopharmaceuticals.



# Eurogentec offices

## EUROPEAN OFFICES

### Belgium

#### Eurogentec Headquarters

LIEGE Science Park  
Rue du Bois Saint Jean 5  
4102 Seraing - Belgium  
Tel.: +32 4 372 74 00  
Fax: +32 4 372 75 00  
[info@eurogentec.com](mailto:info@eurogentec.com)

### Eurogentec Biologics

LIEGE Science Park  
Rue du Bois Saint Jean 14  
4102 Seraing - Belgium  
Tel.: +32 4 372 74 00  
Fax: +32 4 365 16 04  
[info.biologics@eurogentec.com](mailto:info.biologics@eurogentec.com)

### United Kingdom

#### Eurogentec Ltd.

Old Headmasters House  
Unit 1, Building 1  
Forest Business Centre  
Fawley Road, Fawley  
Southampton  
Hampshire SO45 1FJ - UK  
Tel.: +44 1794 511 411  
Fax: +44 1794 522 417  
[info@eurogentec.com](mailto:info@eurogentec.com)

### France

#### Eurogentec France SASU

Avenue Patton 152  
BP 80111  
49001 Angers CEDEX 01 - France  
Tel.: +33 2 41 73 33 73  
Fax: +33 2 41 73 10 26  
[info@eurogentec.com](mailto:info@eurogentec.com)

### Germany & Austria

#### Eurogentec Deutschland GmbH

Cäcilienstrasse 46  
50667 Köln - Germany  
Tel.: +49 221 258 94 55  
Fax: +49 221 258 94 54  
[info@eurogentec.com](mailto:info@eurogentec.com)

### The Netherlands

#### Eurogentec Nederland b.v.

Avenue Ceramique 223  
6221 KX Maastricht - The Netherlands  
Tel.: +31 43 352 06 98  
Fax: +31 43 354 19 65  
[info@eurogentec.com](mailto: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](mailto:info@eurogentec.com)

### Other European countries

#### Represented by Eurogentec Headquarters

LIEGE Science Park  
Rue du Bois Saint Jean 5  
4102 Seraing - Belgium  
Tel.: +32 4 372 74 00  
Fax: +32 4 372 75 00  
[info@eurogentec.com](mailto:info@eurogentec.com)

## US OFFICES

### North America

#### Anaspec, Inc. (Life Science Products)

34801 Campus Drive  
Fremont, CA 94555 - USA  
Tel.: +1 510 791 9560  
Toll-free: +1 800 452 5530  
Fax: +1 510 791 9572  
[info.anaspec@eurogentec.com](mailto:info.anaspec@eurogentec.com)

#### Eurogentec North America, Inc. (Diagnostics)

11111 Flintkote Avenue  
San Diego CA 92121-1222 - USA  
Tel.: +1 858 793 2661  
Toll-free: +1 877 387 6436  
Fax: +1 858 793 2666  
[ivd.na@eurogentec.com](mailto:ivd.na@eurogentec.com)  
[usa.eurogentec.com](http://usa.eurogentec.com)

## ASIAN OFFICES

### Japan

#### Nippon EGT Co., Ltd.

1-1-25, Arakawa  
Toyama 930-0982 - Japan  
Tel.: +81 76 411 02 77  
Fax: +81 76 452 03 99  
[info@n-egt.com](mailto:info@n-egt.com)

### Singapore

#### AIT Biotech Pte. Ltd.

Blk 73 Ayer Rajah Crescent  
#04-08/09 Ayer Rajah Industrial Estate  
Singapore 139952  
Tel.: +65 6778 9388 / 6778 6822  
Fax: +65 6778 5177

## WORLDWIDE DISTRIBUTION

### Australia

#### Integrated Sciences Pty. Ltd

Tel.: +61 2 9417 7866  
Fax: +61 2 9417 5066  
[tech@integratedsci.com.au](mailto:tech@integratedsci.com.au)  
[www.integratedsci.com.au](http://www.integratedsci.com.au)

### Denmark

#### Medinova Scientific A/S

Tel.: +45 3956 2000  
Fax: +45 3956 1942  
[medinova@medinova.dk](mailto:medinova@medinova.dk)  
[www.medinova.dk](http://www.medinova.dk)

### India

#### GeneX

Tel.: +91 44 2437 1711  
+91 44 6566 3944  
Fax: +91 44 2436 2638  
[info@genexindia.net](mailto:info@genexindia.net)  
[www.genexindia.net](http://www.genexindia.net)

### Israël

#### Synteza

Tel.: +972 2 586 7138  
Fax: +972 2 586 2605  
[info@synteza.com](mailto:info@synteza.com)  
[www.synteza.com](http://www.synteza.com)

### Italy

#### BioSense S.r.l.

Tel.: +39 02 612 59 11  
Fax: +39 02 612 59 33  
[contact@biosense.it](mailto:contact@biosense.it)  
[www.biosense.it](http://www.biosense.it)

**Diatech S.r.l.**  
Tel.: +39 0731 21 32 43  
Fax: +39 0731 21 32 39  
[info@diatech-meet.it](mailto:info@diatech-meet.it)  
[www.diatechpharmacogenetics.com](http://www.diatechpharmacogenetics.com)

### Japan

#### Nippon Gene CO., Ltd.

Tel.: +81 76 451 6548  
Fax: +81 76 451 6547  
[info@nippongene.com](mailto:info@nippongene.com)  
[www.nippongene.com](http://www.nippongene.com)

### Norway and Iceland

#### MedProbe AS

Tel.: +47 23 32 73 80  
Fax: +47 23 32 73 90  
[medprobe@medprobe.com](mailto:medprobe@medprobe.com)  
[www.medprobe.com](http://www.medprobe.com)

### Spain

#### Cultek, S.L.

Tel.: +34 91 729 03 33  
Fax: +34 91 358 17 61  
[cultek@cultek.com](mailto:cultek@cultek.com)  
[www.cultek.com](http://www.cultek.com)

### Sweden, Finland and Baltic countries

#### In vitro Sweden AB

Tel.: +46 8 30 60 10  
Fax+ 46 8 30 60 15  
[iv@invitro.se](mailto:iv@invitro.se)  
[www.invitro.se](http://www.invitro.se)

### Romania

#### Bio Zyme S.R.L.

Tel.: +40 264 523 281  
Fax.: +40 264 523 281  
[contact@biozyme.ro](mailto:contact@biozyme.ro)  
[www.biozyme.ro](http://www.biozyme.ro)

### Morocco

#### Univers Biotec

Tel.: +212 5 377 91 636  
Fax.: +212 5 377 91 636  
[infobiotech@yahoo.fr](mailto:infobiotech@yahoo.fr)  
[www.universbiotecmaroc.com](http://www.universbiotecmaroc.com)

# Trademarks labels

ABI Prism® 5700, 7000, 7500, 7700 and 7900 are registered trademarks of The Perkin-Elmer Corp.

Acrydite® is a registered trademark of Mosaic Technologies Corp.

Alexa Fluor® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

AnaTag™ is a trademark of Anaspec, Inc.

ArrayIt® is a registered trademark of TeleChem International, Inc.

Benzonase® is a registered trademark of Merck KGaA, Darmstadt, Germany.

BHQ-0®, BHQ-1®, BHQ-2®, BHQ-3® are registered trademarks of Biosearch Technologies, Inc.

BlackBerry® is a registered trademark of Berry & Associates, Inc.

Black Hole Quencher® is a registered trademark of Biosearch Technologies, Inc.

BODIPY® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

CAL Fluor® is a trademark of Biosearch Technologies, Inc.

Cascade Blue® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

CDP Star® is a registered trademark of Applera Corp.

Cherry™ is a trademark of Delphi Genetics S.A.

Cy® is a registered trademark of GE Healthcare Bio-Sciences Corp.

Cy5® Direct is a registered trademark of GE Healthcare Bio-Sciences Corp.

DHL™ is a trademark of Anaspec, Inc.

Diamond *Taq*® is a registered trademark of Eurogentec S.A.

DNA Engine Opticon® is a registered trademark of Bio-Rad Laboratories, Inc.

Dragonfly Orange™ is a trademark of Eurogentec S.A.

Eclipse® is a registered trademark of Epoch Biosciences, Inc.

Eppendorf® is a registered trademark of Eppendorf AG.

Excel™ is a trademark of Microsoft Corp.

Eurogentec® is a registered trademark of Eurogentec S.A.

Exiqon® is a registered trademark of Exiqon A/S

FLAG® is a registered trademark of Sigma-Aldrich Co

GeneAmp® is a registered trademark of Roche Molecular Systems, Inc.

GenePix® is a registered trademark of Astros Acquisition Sub II.

GenePix 4000® is a registered trademark of Astros Acquisition Sub II.

GetStaby™ is a trademark of Delphi Genetics S.A. Gold'n Oligo® is a registered trademark of Eurogentec S.A.

GO™ Peptide is a trademark of Anaspec, Inc.

GoldStar® is a registered trademark of Eurogentec S.A.

HiLyte™ is a trademark of Anaspec, Inc.

iCycler iQ® is a registered trademark of Bio-Rad Laboratories, Inc.

iCycler® is a registered trademark of Bio-Rad Laboratories, Inc.

I-MUPID® is a registered trademark of Advance Co., Ltd.

iQTM5 is a trademark of Bio-Rad Laboratories, Inc.

iQ-CHECK® is a registered trademark of Bio-Rad Laboratories, Inc.

LightCycler® is a registered trademark of Roche Diagnostics GmbH.

LNA® is a registered trademark of Exiqon A/S.

Lumavidin® is a registered trademark of Luminex Corp.

Luminex® is a registered trademark of Luminex Corp.

Luminex® 200™ is a trademark of Luminex Corp.

Luminex® XYP™ is a trademark of Luminex Corp.

Luminex® SD™ is a trademark of Luminex Corp.

Luminex® SeroMAP™ is a trademark of Luminex Corp.

MagPlex™ is a trademark of Luminex Corp.

MapMarker® is a registered trademark of BioVentures, Inc.

Marina Blue® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

MasterCycler® is a registered trademark of Eppendorf, Inc

MGB® is a registered trademark of Epoch Biosciences, Inc.

MGB Eclipse® is a registered trademark of Epoch Biosciences, Inc.

MicroPlex™ is a trademark of Luminex Corp.

Mupid® is a registered trademark of Advance Co., Ltd.

Mx3000® is a registered trademark of Stratagene

Mx3000P® is a registered trademark of Stratagene

Mx3005P® is a registered trademark of Stratagene

Mx4000® is a registered trademark of Stratagene

NASBA® is a registered trademark of bioMerieux

Oli&GO 1•2•3™ is a trademark of Eurogentec S.A.

OliGold® is a registered trademark of Eurogentec S.A.

Onco-Mir® is a registered trademark of Ambion Inc.

Opticon® is a registered trademark of MJ Research, Inc.

Oregon Green® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Oyster® is a registered trademark of Denovo Biolabels

Pacific Blue™ is a trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Plexor™ is a trademark of Promega Corporation

Pulsar™ is a trademark of Biosearch Technologies, Inc.

Quantica® is a registered trademark of Techne Ltd and Techne, Inc.

Quasar™ is a trademark of Biosearch Technologies, Inc.

Qiagen® and QiaQuick® are trademarks of Qiagen, Inc.

QXL™ is a trademark of Anaspec, Inc.

Rhodamine Red™ is a trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Rotor-Gene® is a registered trademark of Corbett Research Pty Ltd

ROX™ is a trademark of Applera Corp.

RP-Cartridge•Gold™ is a trademark of Eurogentec S.A.

Scorpions® is a registered trademark of DxS Ltd.

SELEX™ is a trademark of Gilead Sciences, Inc.

Smartcycler® is a registered trademark of Cepheid, Inc.

StabyCloning™ is a trademark of Delphi Genetics S.A.

StabyExpress™ is a trademark of Delphi Genetics S.A.

Staby™ Codon is a trademark of Delphi Genetics S.A.

StepOne® is a registered trademark of Applera Corp. or its subsidiaries

SuperScript® is a registered trademark of Invitrogen Corp.

SYBR® Green is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Taqman® is a registered trademark of Applied Biosystems, Inc.

Texas Red® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

TRIZOL® is a registered trademark of Molecular Research Center Inc.

UltraPureGold™ is a trademark of Eurogentec S.A.

VersArray® is a registered trademark of Bio-Rad Laboratories, Inc.

Vic® is a registered trademark of Applera Corp.

Word™ is a trademark of Microsoft Corp.

xPONENT® is a registered trademark of Luminex Corp.

xMAP® is a registered trademark of Luminex Corp.

Yakima Yellow® is a registered trademark of Epoch Biosciences, Inc.

# Notice to purchaser: limited license statements

## 1. Roche Enzyme Products Conveying Polymerase Patent Rights

- This Statement is to be included in all applicable product catalogs, product manuals, product inserts, and on product labels or packaging:

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,079,352, 5,789,224, 5,618,711, 6,127,155 and claims outside the US corresponding to US Patent No. 4,889,818. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim (such as the patented 5' Nuclease Process claims in US Patents Nos. 5,210,015 and 5,487,972), no right to perform any patented method, and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

- This "Short Statement" is provided solely for marketing materials such as advertisements and fliers, for use when space is limited:

Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights (such as 5' Nuclease Process patent rights) are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

## 2. Chemically-Modified Hot-Start Polymerase Reagents and Kits

- This Statement is to be included in all applicable product catalogs, product manuals, product inserts, and on product labels or packaging:

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,079,352, 5,789,224, 5,618,711, 6,127,155, 5,677,152, 5,773,258, and claims outside the US corresponding to US Patent No. 4,889,818. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim (such as the patented 5' Nuclease Process claims in US Patents Nos. 5,210,015 and 5,487,972) and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

- This "Short Statement" is provided solely for marketing materials such as advertisements and fliers, for use when space is limited:

Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights (such as 5' Nuclease Process patent rights) are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

## 3. Licensed dsDNA-Binding Dye Research Kits

- This Statement is to be included in all applicable product catalogs, product manuals, product inserts, and on product labels or packaging:

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,079,352, 5,789,224, 5,618,711, 6,127,155, 5,677,152, 5,773,258, 5,407,800, 5,322,770, 5,310,652, 5,994,056, 6,171,785, and claims outside the US corresponding to US Patent No. 4,889,818. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim (such as apparatus or system claims in US Patent No. 6,814,934) and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

- This "Short Statement" is provided solely for marketing materials such as advertisements and fliers, for use when space is limited:

Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

#### 4. Authorized 5' Nuclease Core Kit (i.e., without Licensed Probe)

- This Statement is to be included in all applicable product catalogs, product manuals, product inserts, and on product labels or packaging:

A license to perform the patented 5' Nuclease Process for research is obtained by the purchase of (i) both Authorized 5' Nuclease Core Kit and Licensed Probe, (ii) a Licensed 5' Nuclease Kit, or (iii) license rights from Applied Biosystems.

This product is an Authorized 5' Nuclease Core Kit. Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,079,352, 5,789,224, 5,618,711, 6,127,155, 5,677,152, 5,773,258, 5,407,800, 5,322,770, 5,310,652, 5,210,015, 5,487,972, and claims outside the US corresponding to US Patent No. 4,889,818. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. Separate purchase of a Licensed Probe would convey rights under the applicable claims of US Patents Nos. 5,538,848, 5,723,591, 5,876,930, 6,030,787, 6,258,569, 5,804,375 (claims 1-12 only), and 6,214,979, and corresponding claims outside the United States. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

- This "Short Statement" is provided solely for marketing materials such as advertisements and fliers, for use when space is limited:

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research when used with the separate purchase of Licensed Probe. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

#### 5. RNA interference

RNA interference is licensed under European Patents 1144623, 1214945 and foreign equivalents from Alnylam Pharmaceuticals, Inc., Cambridge, USA and is provided only for use in academic and commercial research (excluding the evaluation or characterization of this product as the potential basis for a siRNA-based drug) and not for any other commercial purposes. Information about licenses for commercial use (including discovery and development of siRNA-based drugs) is available from Alnylam Pharmaceuticals, Inc., 300 Third Street, Cambridge, MA 02142, USA.

#### 6. Cy<sup>5</sup> amidite, Cy<sup>3</sup> amidite, Cy<sup>3.5</sup> amidite, and Cy<sup>5.5</sup> amidite

The purchase of products with or in which Cy<sup>5</sup> amidite, Cy<sup>3</sup> amidite, Cy<sup>3.5</sup> amidite, and Cy<sup>5.5</sup> amidite may be used or combined, includes a limited non-exclusive sublicense under U.S. Patents 5,556,959 and 5,808,044 and foreign equivalent patents and other foreign and U.S. counterpart applications to use the amidites in such product to perform research. NO OTHER LICENSE IS GRANTED EXPRESSLY, IMPLIEDLY OR BY ESTOPPEL. Use of such product for commercial purposes is strictly prohibited without written permission from Amersham Biosciences Corp. For information concerning availability of additional licenses to practise the patented methodologies, please contact Amersham Biosciences Corp., Business Licensing Manager, Amersham Place, Little Chalfont, Bucks, HP79NA.

#### 7. Yakima Yellow<sup>®</sup> amidite and the Eclipse<sup>®</sup> Non-Fluorescent Quencher

The oligonucleotide probes and primers using the Yakima Yellow<sup>®</sup> amidite and the Eclipse<sup>®</sup> Non-Fluorescent Quencher are for research purposes only, and may not be used for commercial, clinical, diagnostic or any other use. These products or portions thereof are subject to proprietary rights of Epoch Biosciences, Inc. and are made and sold under license from Epoch Biosciences, Inc. under the patent rights WO 0142505 and other pending patent applications. There is no implied license for commercial use with respect to these products. A license must be obtained directly from Epoch Biosciences, Inc. with respect to any proposed commercial use of these products.

#### 8. Custom Molecular Beacon Probes

The use of Custom Molecular Beacon Probes employs the following patent rights licensed to Eurogentec S.A. from The Public Health Research Institute of the City of New York: (a) the claims of U.S. Patent Application Serial No. 08/439,819; (b) the claims of U.S. Patent Application Serial No. 08/990,176; and (c) all patent application claims and all patent claims, U.S. and foreign, that cover improvements in the foregoing claimed probes, kits and assays and whose subjects are inventions and discoveries that are made by the Public Health Research Institute of New York. The use of Custom Molecular Beacon Probes is limited to the field of use comprising the internal use by an end user of a product solely in assays of the end user (or in applications of the end user's customer, if the end user is performing contract research) in scientific research and development.

#### 9. LNA<sup>®</sup> oligonucleotides

LNA<sup>®</sup> oligonucleotides produced under license from Exiqon A/S for research use only. Not for resale or for therapeutic use or use in humans.

#### 10. UDG

Use of UDG employs U.S. Patents 5,035,996, 5,945,313, 5,683,896 and their foreign counterparts licensed to Eurogentec S.A. from Invitrogen Corporation.

**11. SYBR® Green I nucleic acid gel stain**

This product is provided under an agreement between Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corporation) and EUROGENTEC and the manufacture, use, sale or import of this product is subject to one or more of U.S. Patents and corresponding international equivalents, owned by Molecular Probes, Inc. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer, where such research does not include testing, analysis or screening services for any third party in return for compensation on a per test basis. The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, USA. Tel: (541) 465-8300, Fax: (541) 335-0354.

**12. Alexa Fluor®, BODIPY®, Cascade Blue®, Marina Blue®, Oregon Green®, Pacific Blue™, Rhodamine Red™ Texas Red® dyes**

This product is provided under an agreement between Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corporation), and Eurogentec, and the manufacture, use, sale or import of this product may be subject to one or more U.S. patents, pending applications, and corresponding non-U.S. equivalents, owned by Molecular Probes, Inc. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an

academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For information on purchasing a license to this product for any other use, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, USA, Tel: (541) 465-8300. Fax: (541) 335-0504.

**13. Scorpions® Probes**

The use of Scorpions® Probes for research purposes is covered by a license to Eurogentec S.A. from DxS Ltd. No rights to perform or offer diagnostic, commercial testing or other commercial services for money or money's worth are granted by the supply of this product. Should you wish to use this product for any other purpose not covered by this license, please contact DxS Limited at Manchester Incubator Building, 48 Grafton Street, Manchester, M13 9XX.

**14. Black Hole Quencher™**

Black Hole Quencher™, BHQ-0™, BHQ-1™, BHQ-2™, and BHQ-3™ are registered trademarks of Biosearch Technologies, Inc., Novato, California, U.S.A. ('BTI') Patents are currently pending for the BHQ technology and such BHQ technology is licensed by the manufacturer pursuant to an agreement with BTI and these products are sold exclusively for research and development use only. They may not be used for human or veterinary in vitro or clinical diagnostic purposes and they may not be re-sold, distributed or re-packaged. For information on licensing programs to permit use for human or veterinary in vitro or clinical diagnostic purposes, please contact Biosearch at [licensing@biosearchtech.com](mailto:licensing@biosearchtech.com).

**15. CAL Fluor™ and Quasar™**

CAL Fluor™ and Quasar™ dyes are trademarks of Biosearch Technologies, Inc., Novato, California, U.S.A. ('BTI'). Patents are currently pending for the CAL Fluor™ and Quasar™ dye technology and such CAL Fluor™ and Quasar™ dye technology is licensed pursuant to an agreement with BTI and these products are sold exclusively for research and development use only. They may not be used for clinical or diagnostic purposes and they may not be re-sold, distributed or re-packaged.

**16. Iso-base containing oligonucleotides**

Iso-base containing oligonucleotides are covered by US Patents 6,001,983, 6,037,120, and 6,140,496 and their foreign counterparts, and licensed to Eurogentec S.A. by EraGen Biosciences, Inc.

**17. C-5 Propyne and G-clamp modified oligonucleotides**

C-5 Propyne and G-clamp modified oligonucleotides are covered by patents or patents pending owned by Isis Pharmaceuticals, Inc. ('Isis'). Purchase of this product includes a limited license to use this product solely for internal research. This license specifically excludes (and you have no right to use this product for): (a) therapeutic or diagnostic applications (including products or services that incorporate this product), (b) any in vivo toxicity/safety study in support of an investigational new drug application (or foreign counterpart), (c) resale (including sale of any products or services that incorporate this product) or (d) gene functionalization activities (including products or services that incorporate data derived from gene functionalization activities) if such activities have commercial application, any and all of which require a separate license from Isis. Neither this product nor any product created through its use may be used in human clinical trials. In the event you have a separate agreement with Isis regarding this product which explicitly states that the foregoing is not applicable to you, your use of this product will be governed by the terms of such agreement. In no event does the limited license included with the purchase of this product expand or alter the scope of the license granted pursuant to such agreement.

**18. Testing services and contract assay development services**

Testing services and contract assay development services are covered by patents owned by Luminex Corporation. These services are sold exclusively for research use only.

**19. PNA**

All custom PNA oligomers sold by Eurogentec are manufactured by Panagene under an exclusive license and to be used for internal research only. And such custom PNA oligomers are not to be resold unless properly licensed to do so by separate license. All custom PNA oligomers are made and sold pursuant to license under one or more of US Patents Nos. US 5,773,571, 6,133,444, US 6,172,226, US 6,395,474, US 6,414,112, US 6,613,873, US 6,710,163 and US 6,713,602, or corresponding patent claims outside the US.

**20. Probes**

A license to perform the patented 5' Nuclease Process for research is obtained by the purchase of (i) both Licensed Probe and Authorized 5' Nuclease Core Kit, (ii) a Licensed 5' Nuclease Kit, or (iii) license rights from Applied Biosystems.

This product contains Licensed Probe. Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,538,848, 5,723,591, 5,876,930, 6,030,787, 6,258,569, and 5,804,375 (claims 1-12 only). The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. The right to use this product in the 5' Nuclease Process under the applicable claims of US Patents Nos. 5,210,015 and 5,487,972, and corresponding patent claims outside the United States, can be obtained through purchase of an Authorized 5' Nuclease Core Kit. Except under separate license rights available from Applied Biosystems, no right under any other patent claim, or to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, or to sublicense, repackaging with other products, or resell in any form, is conveyed

expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

# General conditions of sale

**1. Area of application**

- 1.1 These general conditions are applicable to all of our sales.
- 1.2 Every purchaser who gives us an order in any way whatsoever will be held to have accepted these general conditions prior to giving the order.
- 1.3 All other conditions provided for in an order form or in any other document emanating from purchaser will be regarded as not being applicable, with the exception of special conditions expressly accepted in writing by our company.

**2. Offer and order**

- 2.1 No order given by word of mouth will be acted upon until it is confirmed by a written order form.
- 2.2 Our agents and representatives have no representational powers. No offer, acceptance of order, or undertaking that these persons negotiate on our behalf will be binding on us until it has been accepted or confirmed in writing by a director or executive of our company duly authorized for this purpose.

**3. Price**

- 3.1 Our prices are quoted ex-VAT in Euros, in Pounds Sterling, or in Swiss franc. Any increase in the VAT rate or new tax imposed between the time of the order and the time of delivery will be borne by purchaser.
- 3.2 Unless otherwise provided in our special conditions, our prices are only for the supply of the products or services described in the special conditions to the exclusion of all other works and services.

**4. Payment**

- 4.1 Our invoices are payable at our principal office within thirty (30) calendar days from the date of invoices. Payments should mention the references indicated in our invoice.
- 4.2 Our agents and representatives have no authority to collect the amount of the invoice unless other expressly provided otherwise.
- 4.3 In the event of an invoice not being settled upon the expiry date, the entirety of our invoices to purchaser will become immediately payable.
- 4.4 Once the payment period has expired, the purchaser becomes automatically liable for the payment of interest for late payment of 12% per annum without formal notice of default being necessary.
- 4.5 All invoices that have not been settled by the due date will moreover be increased as of right and without formal notification of default by an automatic penalty of 10% of the outstanding sum by way of compensation and interests, with an absolute minimum charge of EUR 40,-.
- 4.6 Should the creditworthiness of purchaser deteriorate, we reserve the right, even after the partial completion of an order, to require the buyer to provide such guarantees that we deem fit with respect to the satisfactory performance of the undertakings that have been entered into. A refusal to provide such guarantees gives us the right to cancel all or part of the order.
- 4.7 All complaints regarding the invoiced amounts must reach us within fifteen (15) calendar days of the invoice date, in the absence of which the invoice will be held to be accepted irrevocably and in its entirety.

**5. Procedures for deliveries**

- 5.1 Except when provided otherwise in our special conditions, the delivery will be made Free Carrier at the address indicated by purchaser on his order form (FCA – LIEGE science park, B-4102 SERAING, BELGIUM) (Incoterms 2000). Nonetheless should purchaser so request, we can arrange the delivery, at the risk and expense of purchaser.

**6. Delivery times**

- 6.1 Except when express guarantee is given in our special conditions, the delivery time mentioned in our special conditions is not binding. We cannot be held liable unless the delay is significant and attributable to a grave error on our part.
- 6.2 Delays in delivery can in no case give rise to the cancellation of the order.

**7. Reservation of ownership**

- 7.1 The delivered products or services remain our property until the price has been paid in full, even in the event of the transformation or incorporation of the products into other goods.

**8. Approval**

- 8.1 Purchaser shall be deemed to have accepted the products or services fourteen (14) calendar days after the delivery except in the event of a precise and detailed report outlining the nonconformity of the services or the products sent to us by registered letter before the expiry of this period.
- 8.2 The approval will cover all apparent defects, that is to say all those defects that it is possible for the purchaser to discern at the time of the delivery and during the period of fourteen (14) calendar days following the delivery as a result of close and careful examination, in particular those defects relating to the specifications and the operation of the products or services sold.

**9. Limitation of liability**

- 9.1 Our sole liability will be, at our own option, the replacement of the non conforming products or services or the refund of the price paid for such non-conforming products or services.
- 9.2 After delivery, we cannot assume any other liability other than that provided for by Article 9.1.
- 9.3 In consequence we cannot be held liable for any damage or interest arising from accident sustained by persons, damage to goods distinct from the sold products, loss of earnings or any other loss arising directly or indirectly from defects in the products or services.

**10. Rescission of the sale**

- 10.1 We have the right to rescind any sale as of right and without formal notice of default, by notification to purchaser of our intention by registered letter, in the event of non-performance by purchaser of one or other of its contractual undertakings, in particular if purchaser is late in settling an invoice by more than thirty (30) calendar days, or if it should appear that purchaser will not perform or that there is reasonable likelihood that purchaser will not perform one or more of its main undertakings, even before purchaser is required to perform such obligation.
- 10.2 Upon rescission of the sale in application of the foregoing paragraph, purchaser will be liable for automatic damages and interest of 10% of the sale price.

**11. Applicable law**

- 11.1 Our sales shall be governed by and construed in accordance with the substantive laws of the country where our principal or registered office is situated.
- 11.2 All disputes directly or indirectly related to our contractual relations with the purchaser will be submitted exclusively to the Courts of the Judicial District of our principal office.