



Oligonucleotides

**About
Eurogentec**



↖
**DOWNLOAD THE
ONLINE BROCHURE**

10 000 m²
State-of-the-art facilities

1985
Since

**ISO 9001 &
ISO 13485**
Certified process

400+
Employees

3 Departments

Eurogentec's activities are spread over three business units to guarantee a smooth

continuum from research to higher demanding applications.



RESEARCH PRODUCTS



DIAGNOSTIC SERVICES



THERAPEUTIC SERVICES

6 Life sciences product lines

We produce various biomolecules for a large panel of applications.



OLIGONUCLEOTIDES



qPCR & PCR



ANTIBODIES



PEPTIDES



PROTEINS



GENES

Eurogentec is a **biotechnology** company providing life sciences products and personalized services to academic, medical and industrial scientists **worldwide since 1985**. We produce custom and catalog reagents, in different quality grades and scales.

We are also a trusted **CDMO** internationally recognized for the development and production of biopharmaceutical critical raw materials, vaccines and therapeutic molecules.

Based in **Belgium**, Eurogentec is a subsidiary of Kaneka Corporation.

08 — 15

Oligo grades

Research grade	11
Diagnostic services	12
Therapeutic services	14

16 — 47

Oligonucleotides overview

Custom Parameters	19
Backbones	20
Modifications	23
Scales	25
Purifications	26
Quality controls	28
Formats & Packaging	31
Specific Oligo Types	33
qPCR Probes	34
RNAi oligos	38
NGS	40
Aptamers	43
Catalog Reagents	45
Catalog products	46

48 — 74

Oligos ordering guidelines

Order settings	51
Order placement	57
Delivery	67
Handling instructions	71

Research grade	11
Diagnostic services	12
Therapeutic services	14

Oligo grades

We can provide research, diagnostic and therapeutic grade oligonucleotides based on your specific application and needs.



Since 1985, we have been providing high-quality reagents and custom-synthesized oligonucleotides to scientists around the world.

Research grade

Our oligos range from the routine research primers to highly modified sequences.

We have multiple synthesizers operating in parallel to ensure continuous production of any nucleic acid sequences. Our chemists are experts in complex oligo synthesis such as: siRNA, Double Dye Probes and long oligonucleotides.

Eurogentec oligonucleotides are produced with state-of-the-art equipment, using only the highest quality reagents from trusted suppliers.

Our team of experts is committed to delivering quality products in accordance with the customers' request.

FOCUS

225

From 2 to 225 bases

300+

Modifications

ISO 9001

Certified process

LARGE RANGE
of chemistries

CUSTOM
fill & finish

From lab to
LARGE SYNTHESIS SCALE



MORE INFO

Diagnostic services

FROM RESEARCH TO COMMERCIALIZATION

We manufacture diagnostic grade oligonucleotides with increased levels of quality, traceability and adapted batch records.

We offer a range of services designed to meet your specific needs, ensuring the perfect fit for your requirements.

Track™ Oligonucleotides are reasonably priced and include **documented traceability**, often required for demanding projects. They are delivered with a **quick turnaround time**, and their production process is certified to comply with ISO 9001.

Pre-diagnostic and **diagnostic** grade oligos are produced in a controlled environment within our ISO 7 and ISO 8 cleanrooms in Belgium. Their production process is certified to comply with ISO 9001 and ISO 13485 standards.

We are also a registered manufacturer of **Analyte Specific Reagent (ASR)** oligonucleotides for use in *in vitro* application.

Additional Services

IP-FRIENDLY

We offer IP-friendly solutions for Probes with our range of Access™ Dyes and Quenchers covering the full near blue to infrared spectrum.

INTEGRATED SOLUTION

Looking for an integrated solution for your diagnostic kit preparation? We have specified services for Molecular Diagnostic kit manufacturing or Lab Developed Test preparation.



ASK
A QUOTE

Oligo grade comparison & Development Process

Discovery	Feasibility	Design control	Validation	Commercialization
Research Oligos		Track™ Oligos		Diagnostic Oligos
		Pre-Diagnostic Oligos		

Oligonucleotide Grades	● Research	● Track™	● Pre-Diagnostic	● Diagnostic
Process				
Dedicated Project Management	Option	✓	✓	✓
Customized Fill & Finish	Option	Option	✓	✓
Quality Management				
ISO 9001 Certified Process	✓	✓	✓	✓
ISO 13485 Certified Process	-	-	✓	✓
Qualification/Validation [Equipment & Method]	-	Partial	✓	✓
Control				
Quantification	Single	Dual	Triple	Triple
Stringent QC Tests [Validated]	-	✓	✓	✓
Traceability	Partial	Documented	Documented	Full Documented
Batch Record [Archived for 5 years]	-	-	Partial	Full
Classified Cleanroom	-	-	✓	✓
Certificate of Analysis [CoA]	-	✓	✓	✓

Therapeutic services

FROM DEVELOPMENT TO CLINICAL PHASES

Our services cover every phase, from research and discovery, including oligo screening libraries up to larger quantities needed for preclinical and early clinical studies.

We provide **tailored and comprehensive solutions** for the development, manufacturing, and testing of oligonucleotides. Whether you need help with custom oligo design or large-scale production, our commitment to providing the right features, quantity, and quality is backed by our **extensive experience** as a CDMO for biotech and pharma companies.

Various chemistries and modifications

We specialize in creating and synthesizing various oligonucleotides, such as **antisense oligonucleotides (ASOs)**, **siRNA**, **aptamers**, and **CpG oligos**. Our extensive knowledge encompasses a wide array of chemistries (including 2'-O-Me RNA, 2'-O-MOE RNA) and modifications (such as fluorescent dyes, GalNAc, PEG, lipids), ensuring enhanced stability, binding affinity, specificity, and accurate site delivery of therapeutic oligonucleotides.

Equipped to provide quality

To achieve the required quality, we have a range of facilities up to **ISO7 and ISO8 cleanrooms** equipped with the latest technologies, to produce and purify your oligonucleotides. Moreover, for clinical oligonucleotides, we follow a validation approach that aligns with ICH guidelines, ensuring compliance with FDA, EMA, and other regulatory standards.

FOCUS

MORE THAN AN OLIGO SUPPLIER

Comprehensive project management

Our experienced project managers maintain clear communication, anticipate potential challenges, and implement strategic solutions to ensure smooth and efficient progress from start to finish.

Assistance through the testing process

We give you access to *in vivo* and *in vitro* preclinical studies prior to progress to clinical trials.

Administrative support

We assist you with the writing and review of CMC documents and provide comprehensive support for your IND file preparation, delivering documentation tailored to your specific needs.



ASK A QUOTE

Services Overview

	Research & Discovery	Preclinical development	Preclinical to clinical	Clinical trials
	Early Drug Discovery	<i>In vitro</i> testing	<i>In vivo</i> testing	Engineering batch Early Clinical Phases
Quality standards	● RUO		● GMP	
Examples of Applications	Early Drug Discovery steps incl. screening libraries to lead compound	Drug Discovery: <i>In vivo</i> & <i>in vitro</i> studies; other high demanding application	GLP tox/safety	Early Clinical Phases
Quantity	From micrograms to several grams	From hundred of milligrams to several grams	From hundred of milligrams to several grams	From several grams to hundred of grams
Agreements				
Confidentiality agreement	✓	✓	✓	✓
Master Service agreement	On request	On request	✓	✓
Quality agreement	-	-	✓	✓
Dedicated Project Management	✓	✓	✓	✓
Quality				
Quality systems	ISO9001	ISO9001	GMP	GMP
Incoming Raw material specification & testing	-	-	✓	✓
Optimization of analytical methods	-	On request	✓	✓
Equipments/methods qualification	-	-	✓	✓
Validation work	-	-	✓	✓
Classified clean rooms	-	-	✓	✓
Retained samples	-	-	✓	✓
Stability studies	-	On request	✓	✓
Documentation				
QA documentation	CoA	CoA, Summary batch record optional	Summary batch record or BMR	BMR
Change Control Notification	-	-	✓	✓

Custom Parameters	19
Backbones	20
Modifications	23
Scales	25
Purifications	26
Quality controls	28
Formats & Packaging	31
Specific Oligos Types	33
qPCR Probes	34
RNAi oligos	38
NGS	40
Aptamers	43
Catalog Reagents	45

Overview

We provide a large choice of oligonucleotides tailored to your specific application, ranging from unmodified primers to a diverse array of chemistries, modifications, specifications, and purification options.



Custom Parameters

A large choice of options including various backbones and over 300 modifications are available to help you design oligos with the appropriate properties.

We provide a wide range of modified backbones that grant different properties to your oligos.

Backbones

We offer classic DNA and RNA based oligos, but also modified backbones that grant different properties to your oligos.

LNA[®] backbones increase the affinity and specificity of the oligos. Indeed, their structure confers a very strong thermal stability towards complementary DNA and RNA template suitable for hybridization assays requiring high specificity and/or reproducibility.

Both **2'-O-Me RNA** backbone and **phosphorothioate linkages** increase the affinity and specificity for the target, while boosting the resistance to nucleases.

2'-O-MOE RNA backbones are used to increase resistance to nucleases.

We also offer **PNA backbones** well suited for **FISH studies**. This artificial DNA/RNA analogue confers a higher specificity and sensitivity to the oligo. PNA are also known to be resistant to enzymatic degradation and stable over a wide range of pH, temperature, and salt concentrations.

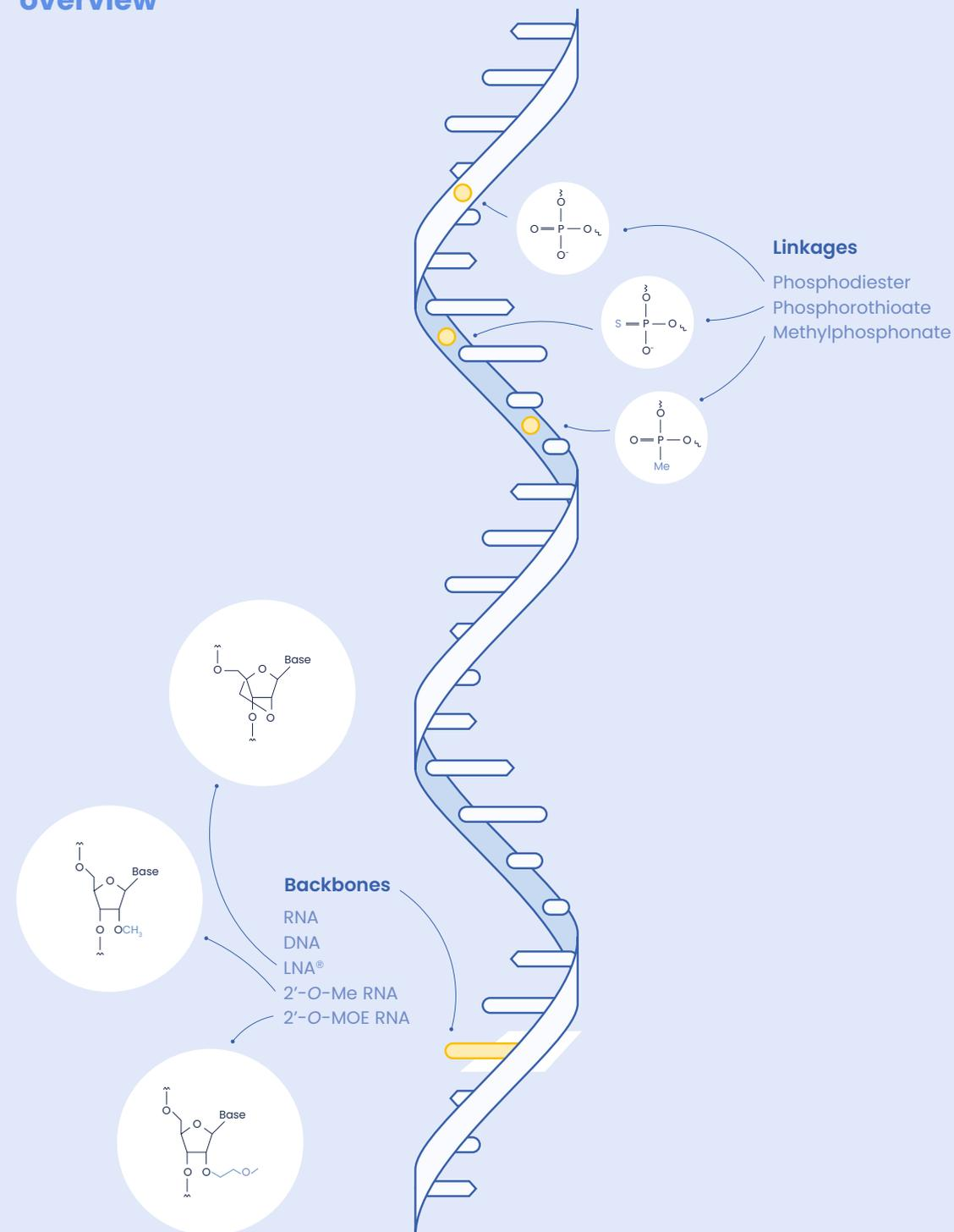
OUR PANEL OF LINKAGES INCLUDES

Phosphodiester bonds. They connect the 3' carbon atom and the 5' carbon of the sugar.

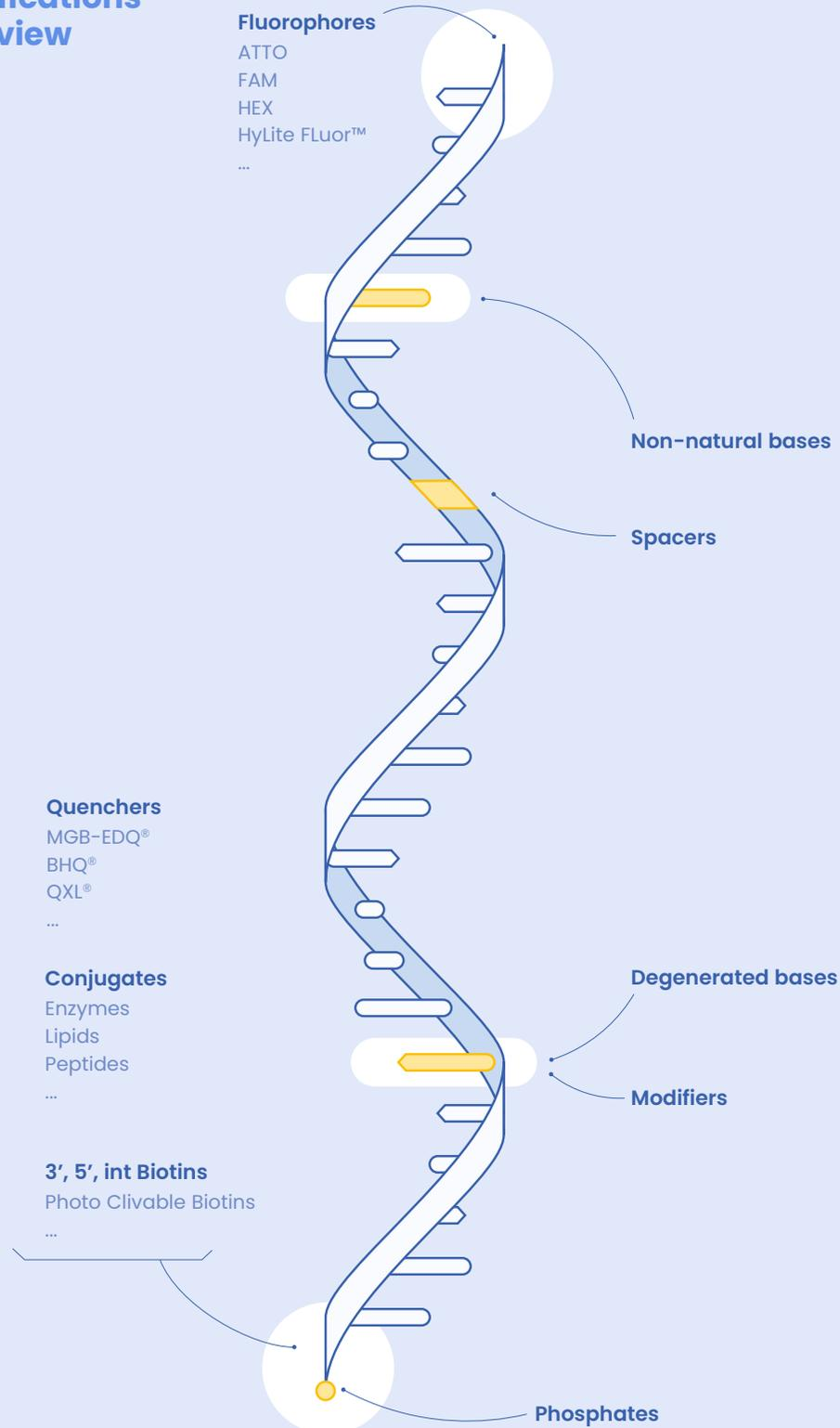
Phosphorothioate bonds possess an increased resistance against nucleases due to the substitution of non-bridging oxygen by sulfur.

Lipophilic bonds such as **methylphosphonate linkages** (only available for DNA bases) help the oligos cross the cell membrane. They are non-ionic nuclease resistant linkages. Methylphosphonate/RNA duplex are not recognized by RNaseH.

Backbones overview



Modifications Overview



Modifications

In addition to backbone chemistries, diverse additional modifications can also be incorporated to confer special characteristics to your oligos.

For instance, the ability to **cross cell membrane** can be obtained by adding **cholesterol** to the oligo, while increasing the **affinity and specificity** can be achieved with **C-5 methylated pyrimidine deoxy-nucleosides**, **2'fluoro RNA** or **methyl-dC**, among others.

Oligonucleotides can be modified by direct incorporation during the synthesis or by post-synthesis labeling.

Direct incorporation

Since oligonucleotide synthesis happens from 3' to 5', **3' direct incorporation** is not always possible. It depends on the availability of the solid support and compatibility with the chemistries used in the synthesis.

5' and internal modifications can be introduced using the phosphoramidites, but they need to support harsh cleavage-deprotection conditions.

Post-synthesis incorporation

Post-synthesis incorporations are used to introduce sensitive dyes or compounds that do not exist as phosphoramidites. It may influence the yield of the reaction. Indeed, a lower yield may result from poly-modifications and/or strong secondary structures.

All modifications are available in **research, diagnostic and therapeutic grades**.

300
Modifications

FOCUS

What are POCs?

POCs stand for **Peptide Oligonucleotide Conjugates** which serve many important roles as potential therapeutics.

Why using POCs?

POCs are frequently used for cell screening or *in vivo* studies thanks to the increase of the oligo membrane permeability. In therapeutics, one of the most interesting POCs applications is targeted drug delivery.

Our POCs can harbor specific modifications, either on the peptide or the oligonucleotide branch, to fit varied functionalities. Different conjugation strategies can be applied based on the sequences and customer requirements.



FULL LIST OF MODIFICATIONS



From small to large scales to fit the needs of your application.

Scales

We synthesize oligonucleotides from 10 nmol to 20 μmol scales according to the type of oligo, and even more on request.

When ordering an oligo, pay attention to the synthesis scale (*i.e.* the amount of raw material used to start the synthesis of oligonucleotides) and the yield of the final product recovered at the end of the manufacturing process.

The length, the sequence, the type/number of modifications and the purification, strongly influence the reaction yield. Based on that, Eurogentec has defined a minimum guaranteed yield in nmoles for all product categories.

[For more information consult p.52]

FROM
0.01 μMOL
TO
20 μMOL

FOCUS

Large scale synthesis

Does your project require larger quantities of oligonucleotides? Our large scale synthesis service is the answer for your production.

Scales range from micrograms to hundreds grams.

Custom service to guarantee the best production conditions.

Customized oligos with various modifications available.

Purity reached by HPLC method.

Format in bulk, plates or vials, dried or in solution.

Support with a dedicated project manager.

Production grades suitable for Research, Diagnostic, or Therapeutic use.

Choosing the appropriate purification method is key to tailoring each oligonucleotide for your application.

Purifications

The aim of any purification step is to remove the by-products resulting from the removal of the protecting groups and other synthesis by-products.

Whether you are engaged in cutting-edge research, diagnostic development, or therapeutic innovations, ensuring the right quality and purity of your oligonucleotides is essential for reliable and reproducible results.

Various parameters should influence your purification choice. Those parameters are the oligo length, the presence and the number of modifications, the intended applications and the expected yield.

[Consult the "Synthesis scale vs Guaranteed yield" Table p.52 to select the appropriate synthesis scale.]

IN VIVO LIKE OLIGONUCLEOTIDES

The solution for your project before entering preclinical testing.

The *in vivo* like process is recommended for the production of antisense oligonucleotides or siRNA testing at research level before entering pre-clinical testing. Therefore, they are an ideal solution for your animal studies.

In vivo like purification process includes:

- 1 Ion exchange HPLC
- 2 Desalting
- 3 0,2 µm filtration
- 4 Lyophilization

Thanks to this stringent purification process, endotoxins should remain undetectable using a chromogenic LAL endotoxin assay.

It will give more than 85 % purity for oligos shorter than 45 bases.

Please note that oligonucleotides manufactured with the *in vivo* like process are not endotoxins free nor sterile certified RUO oligos.



FOR MORE DETAILS

Purification methods



SEPOP DESALTING

Up to 70% purity for ≤79 bases oligos.

Differential precipitation to eliminate the largest part of contaminants (truncated material < 10 bases).

For common applications such as primers for standard PCR.



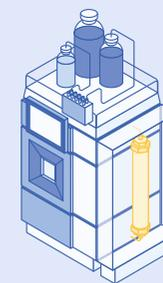
RP CARTRIDGE•GOLD

At least 70% purity for ≤45 bases oligos*.

Reverse Phase chromatography based on differences in hydrophobicity.

Best compromise for most applications such as PCR, mutagenesis or cell culture (absence of residues).

*Possibility to order oligos up to 59 bases, but 70% purity is not guaranteed.



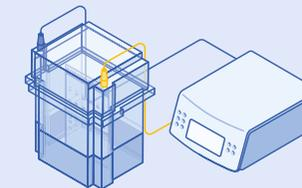
HPLC

At least 85% purity for ≤45 bases oligos*.

Reverse Phase (RP) or Ion Exchange (IEX) mode.

Default method for qPCR probes & antisense oligonucleotides.

* ≥ 80% for complex oligos shorter than 45 bases, ≥70% for longer oligos up to 59 bases



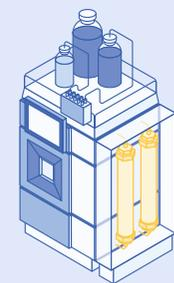
POLYACRYLAMIDE GEL ELECTROPHORESIS (PAGE)

At least 85% purity for ≤79 bases oligos*.

Separation based on the charge and mass of oligonucleotides, varying from only 1 base.

Recommended for long oligos starting from 60 bases.

* ≥ 75% for complex oligo shorter than 79 bases. Possibility to order oligos up to 139 bases, but 85% purity is not guaranteed.



DUAL HPLC

At least 85% purity for ≤45 bases oligos*.

Double purification: RP/RP-HPLC or RP/IEX-HPLC. The average purity will increase compared to simple HPLC.

Appropriate for antisense oligos or siRNAs, might be required for probes with post-synthesis coupling.

* ≥ 80% for complex oligo shorter than 45 bases, ≥70% for longer oligos up to 59 bases

Quality Controls

Quality Control (QC) is a crucial component of the oligonucleotide manufacturing process. All of our synthesizers are equipped with real-time trityl monitoring to ensure that each oligo meets our strict quality standards.

In addition to routine optical density (OD₂₆₀) measurement, which ensures you receive the correct quantity of oligos, we perform a range of automated quality control checks. These checks are tailored to the type and specifications of each oligo to ensure optimal performance and quality for your specific application.

Supplementary controls are available if required.

Default QCs

Oligo Type		MS	UHPLC
Custom Oligonucleotides	Unmodified	✓ ¹	
	Modified	✓ ²	
qPCR Probes	Double Dye Probes	✓	✓
	Molecular Beacons	✓	✓
	MGB Probes	✓	✓
RNAi Oligonucleotides	siRNA Duplexes	✓	
NGS Oligonucleotides		✓	✓
Complex Oligonucleotides		✓	✓ ³
Calibration Oligos		✓	✓
Universal Primers		✓	✓

TABLE LEGENDS & NOTES

MS Mass Spectrometry;
UHPLC Ultra High Performance Liquid Chromatography.

¹Randomized high throughput control.

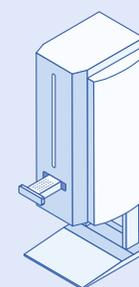
²Except for SePOP desalted oligonucleotides.

³Optional.

For technical reasons this general rules may be adapted to provide you with the most suitable accurate oligonucleotide.

QCs Overview

If the default QC is not sufficient for your application, you can request supplementary controls. We offer various options to perfectly meet your specific quality requirements.



MASS SPECTROMETRY

MALDI-TOF, ESI-TOF or LC-MS

The most precise information about the length, deprotection-product, and the presence of labels for modified oligonucleotides over a broad range of lengths.



CHROMATOGRAPHY

Ion-Exchange UHPLC (IEX-UHPLC) or Reverse Phase UHPLC (RP-UHPLC)

UHPLC will give you quantitative information about the purity level of oligonucleotides.



FLUORIMETRY ANALYSIS

Non destructive physical analysis method.

Qualitative information about fluorescent oligonucleotides.



Formats

According to your specifications, oligonucleotides may be provided in various formats.

Dried is the default format (except for SePoP unmodified oligonucleotides from 15 to 39 bases).

In solution: Select the type of reconstitution buffer (H₂O or TE (pH=8)), determine its volume (ranging from 50 to 1000 µl), and/or specify the final oligonucleotide concentration (ranging from 5 to 250 µM).

Annealed: Oligonucleotides are annealed by a short treatment at 94°C followed by a progressive cooling to room temperature. siRNA or cloning linkers are annealed by default.

Mixed: Similar amounts of different oligonucleotides can be combined in a single tube.

Packaging

We provide various packaging options based on your preferences.

2 mL tube: By default, each oligonucleotide is provided in individual 2 mL tube. Higher volume can be delivered on request (15 mL, 50 mL*).

96-well plates: Cluster tubes, well plates and deep well plates are available for oligos up to 59 bases.

Aliquoting: All the oligonucleotides in solution can be split on request in small aliquots of accurate volume (from 50 to 1000 µl).

We also offer customized packaging to meet the specific requirements of your project.

HIGH THROUGHPUT SCREENING & DISCOVERY

From Oligo screening for *in vitro* testing in cell models to lead optimization. We produce up to a thousand sequences in small scale and deliver them simultaneously.

DESALTED OR HPLC PURIFIED

IN VIVO LIKE PURIFICATION PROCESS AVAILABLE

DELIVERED IN PLATES OR TUBES

LYOPHILIZED OR IN SOLUTION

*Higher scales from 2.5 µmol may be delivered in 50 ml tubes by default.



Specific Oligo Types

We produce specialized oligonucleotides for specific applications, with preset parameters.

Various types of oligos such as custom probes, interfering RNAs, or oligos for NGS can be configured directly with our online configurators and even more upon request.

qPCR probes

We produce custom qPCR probes with a vast portfolio of fluorescent dyes and quenchers, compatible with every channel of any real-time thermocycler. Our fluorescent probes and quenchers are just as effective in digital PCR.

We have developed easy-to-use online configurators to order the custom probes you need. Watch the tutorial to get started or go directly to the appropriate configurator. [For more info consult p.59.]

Looking for a solution for commercial applications?

Our Access™ Dyes & Quenchers consist of IP-friendly efficient reporters & quenchers for dual labeled qPCR probes. Our selection of dyes for diagnostic tests covers the full visible spectrum, from blue to far-red-shifted wavelengths.

BHQ® QUENCHERS

Label your custom oligonucleotides and qPCR probes with the original BHQ® quenchers. They are license-free at Eurogentec for research, diagnostic and commercial use. BHQ® quenchers cover all qPCR channels and are compatible with most of the fluorescent dyes.

ATTO DYES

Improve your results with our extensive collection of IP-friendly ATTO dyes. Our selection covers the full visual spectrum, possesses strong absorption and high fluorescence quantum yield, and is well-suited for bioanalytical applications.

MOLECULAR BEACONS

Molecular Beacons are probes which contain a stem-loop structure, a fluorophore, and a quencher at their 5' and 3' ends, respectively.

We offer standard, wavelength-shifting and 2'-O-Me RNA Molecular Beacon probes.

Our custom Molecular Beacons are available with DABCYL or BHQ® non fluorescent quenchers, paired with compatible fluorochromes. Other dyes are available upon request.



CONFIGURE MB PROBES

NEED HELP TO DESIGN YOUR qPCR PROBES?

Our experts can help you design your probes and RNA oligonucleotides.

With our robust design software, we deliver optimized sequences, continuously updating it with the latest scientific advances and customers' feedback.



LEARN MORE

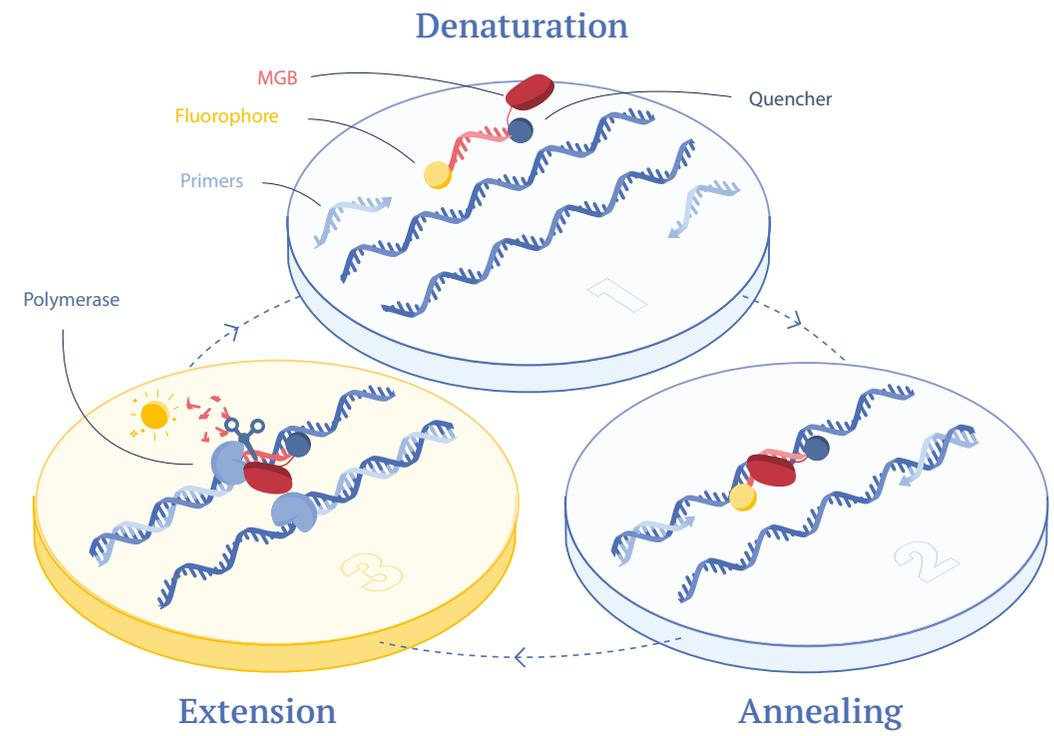
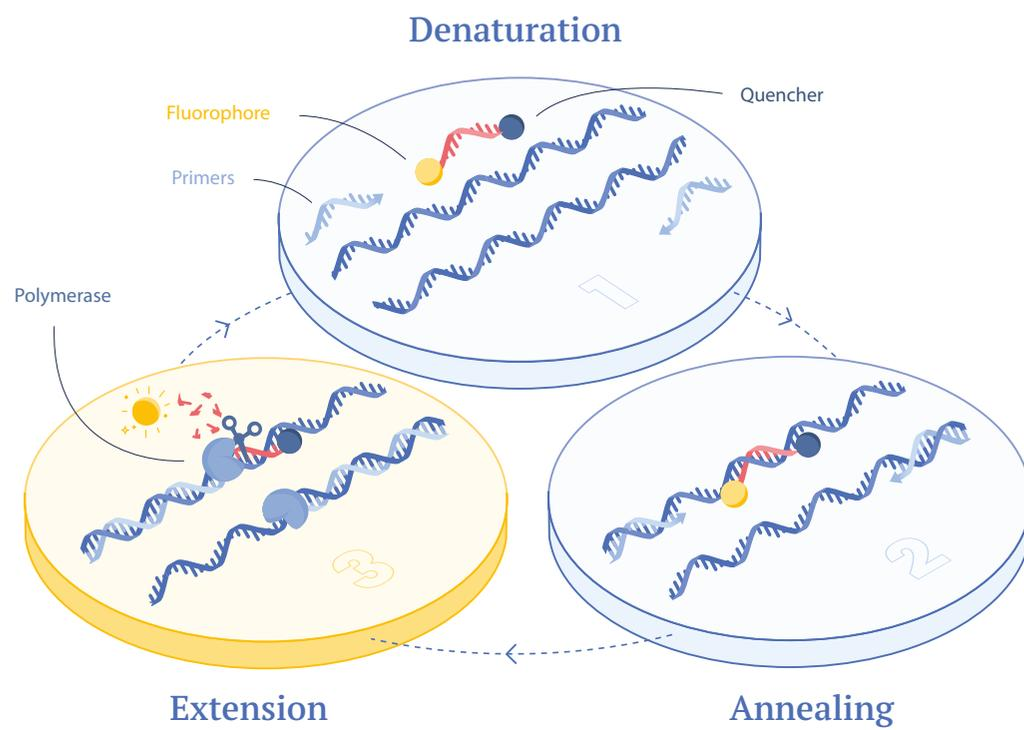
DON'T MISS OUT OUR OFFER!

For each probe configured online, receive two free primers!



← SCAN TO LEARN MORE





Double-Dye probes

Double-Dye probes are labeled with a fluorescent reporter dye and a quencher at their 5' and 3' ends, respectively.

The probes are available with classic DNA backbone, but also Locked Nucleic Acid (LNA[®]), MGB for higher thermal stability, specificity and reproducibility.

We have a large choice of fluorophore/quencher combinations. Various fluorescent dyes are available according to the selected quencher e.g. FAM/BHQ[®]-1, ROX/BHQ[®]-2, HEX/MGB-EDQ, and ATTO/BHQ[®] couples.

[View the full list on the online configurator].

ADVANTAGES OF (LNA[®]) TAQMAN[®] PROBES

- **THERMAL STABILITY**
Possibility to adjust T_m values of primers and probes in multiplexed assays.
- **SPECIFICITY**
- **REPRODUCIBILITY**

SPECIFICITIES

- Length**
from 8 to 45 bases
- Purification**
HPLC
- Quality Control**
MALDI-TOF MS + UHPLC
- Format**
Dried or in solution



CONFIGURE DOUBLE-DYES PROBES

MGB probes

MGB probes are dual-labeled probes incorporating a Minor Groove Binder at their 3' end.

MGB Eclipse[®] probes ordered at Eurogentec are no longer limited by any licence, and can be used for any applications all around the world!

With our portfolio of fluorescent dyes for MGB Eclipse[®] probes, our offer covers all qPCR channels. It will allow you to increase the specificity, sensitivity and efficiency of your assay.

SPECIFICITIES

- Length**
from 8 to 30 bases
- Delivered quantities**
6 - 20 - 50 nmol
- Purification**
RP-HPLC
- Quality Control**
MALDI-TOF MS + UHPLC
- Format**
In solution (H₂O or TE); dried or dried aliquoted in 10 nmol on request



CONFIGURE MGB PROBES

RNAi oligos

RNA Interference (RNAi) is a mechanism of gene silencing at the post transcriptional level, involving RNA molecules such as siRNAs and miRNAs.

siRNA

Synthetic siRNA duplexes are introduced into cells to cause RNA interference and inhibit the expression of a specific mRNA.

Custom siRNA synthesis

Typical siRNAs are 21-mer, with a core of 19 double-stranded RNA bases followed by 2 single-stranded DNA nucleotides in 3', usually Thymidine. Multiple options are available to customize your siRNAs.

Reach the level of purity needed by selecting the purification method among SePOP desalting, HPLC (RP or IEX), or our *in vivo like* purification process.

Negative and positive siRNA controls

No experiment is complete without proper controls. To monitor your experiment conditions, we provide you with siRNA control duplexes and kits including positive and negative siRNA controls. All siRNA control duplexes are IEX-HPLC purified, and 100 % MALDI-TOF Mass Spectrometry controlled. siRNA controls are annealed and shipped dried.

Positive controls: siRNAs targeting a range of endogenous and reporter genes. Each control contains one siRNA duplex.

Negative controls: siRNA with no homology with any known eukaryotic gene. The sequence is properly validated.

If needed, our experts will help you find the best design for your RNA oligonucleotides.



CONTACT OUR TECHNICAL SUPPORT

miRNA

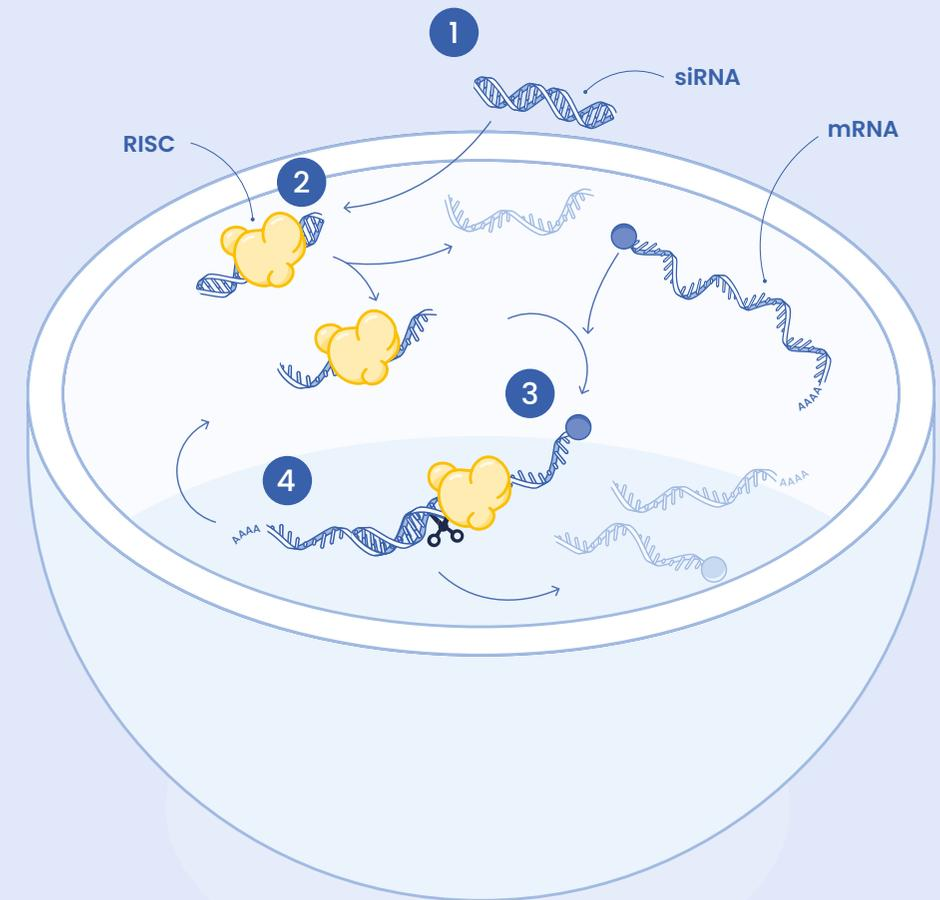
miRNAs (for microRNAs) are small non-coding RNAs forming short hairpins. They are involved in gene expression regulation and RNA silencing mostly by inducing the active degradation of the targeted mRNA or by preventing its translation.

We provide custom synthesized miRNA mimics and inhibitors available with different labels which can be linked to cholesterol to increase cellular uptake. Other modifications are available on request, like covalently-bond peptides, modified bases, dyes and more.



CONFIGURE YOUR siRNA DUPLEXES

siRNA-mediated RNA interference mechanism of action



siRNA-mediated RNA interference mechanism of action.

- 1 The siRNA duplex enters the cell.
- 2 The double stranded siRNA is incorporated into a multiprotein complex, forming the RISC (RNA-induced silencing complex). The guide strand is loaded into the RISC and binds complementary target mRNA. The siRNA passenger strand is discarded.
- 3 The target mRNA is cleaved.
- 4 The RISC starts a new cycle.

Eurogentec can support your sequencing projects by offering dedicated NGS grade oligonucleotides.

NGS oligos

A specific production process to ensure a low cross-contamination (<0.1%) for your Next-Generation Sequencing oligos.

Next-Generation Sequencing (NGS) is a high-throughput technology allowing the massive sequencing of nucleic acids following a DNA library preparation. The NGS technology requires that adapter sequences added to fragmented nucleic acids have a high level of purity (no by-products) and ultra-low cross-contamination.

Eurogentec can support your sequencing projects by offering dedicated NGS grade oligonucleotides. Specially produced to avoid cross contamination, they reduce barcode misalignment during multiplex next generation sequencing projects.

NGS SPECIFICITIES

Quality

Low cross contamination (<0,1%)

Length

from 20 to 85 bases

Quantity

10 nmol minimum delivered

5' Modifications

5' Phosphate / 5' Biotin-TEG

Purification

HPLC or Cartridge

Quality Control

100 % Quality Control checked by Maldi-TOF MS

Format

Dried in tubes

Bases Option

Phosphorothioate bond

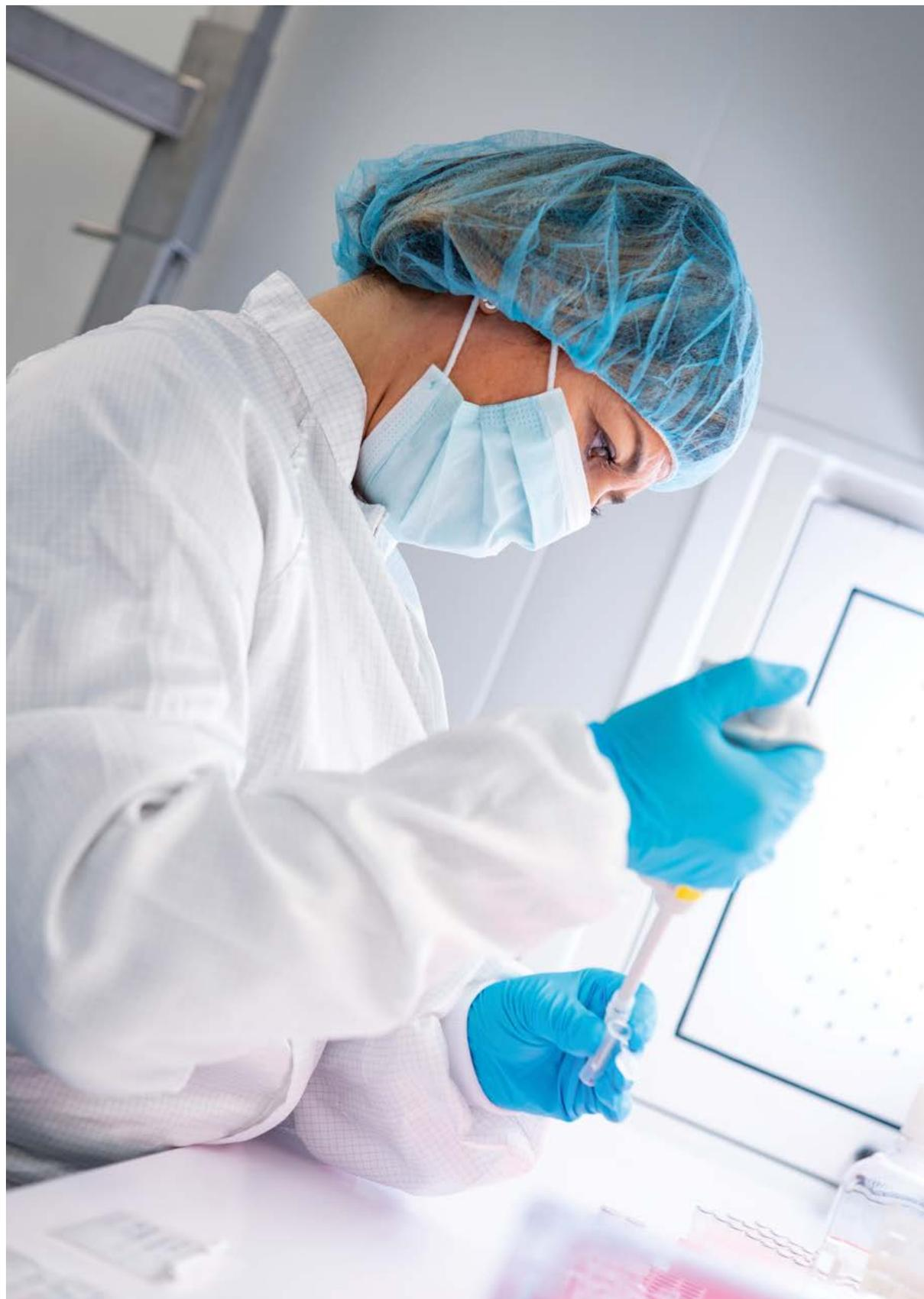
Wobble Bases

Available at no additional cost



CONFIGURE
NGS OLIGOS





We can produce subsequent large quantities of aptamers according to the setups defined during the development phase.

Aptamers

We collaborate with Novaptech, one of the world leaders in the development of oligonucleotide aptamers. They identify your ideal molecule, we synthesize it.

Aptamers are oligonucleotides that bind a specific target molecule, or molecule family, based on their tertiary structure. They are considered as **an alternative to antibodies**.

Novaptech identifies and optimizes aptamer candidates while Eurogentec is dedicated to synthesizing the final aptamer candidate, along with all oligonucleotides needed for your aptamer development process.

We provide high-value aptamer-related services which include the selection and production of custom optimized RNA and DNA aptamers for analytical, diagnostic, and therapeutic applications. Each aptamer can be provided in various quantities and delivered either unmodified or modified.

ANY TARGET – IN VARIOUS CONDITIONS

FAST & CONTROLLED PRODUCTION

GOOD ALTERNATIVE TO ANTIBODIES

Aptamers have shown similar, and even higher affinity and specificity compared to traditional monoclonal antibodies. They exhibit low immunogenicity and are chemically synthesized, requiring no animal or cell culture. Consequently, they overcome many limitations of antibodies and are well-suited for both diagnostic and therapeutic applications.



CONTACT OUR TECHNICAL SUPPORT



Catalog Reagents

We supply oligo synthesis reagents from Glen Research and ready-to-use nucleic acid sequences.

Easily order ready-to-use nucleic acid and oligo synthesis reagents from Glen Research.

Catalog oligos

In addition to our custom oligonucleotide synthesis service, we also provide ready to use nucleic acid sequences.

You can select **Universal primers** among our catalog. Those pre-designed primers are complementary to nucleotide sequences that occur very commonly in specific sets of DNA molecules and cloning vectors.

We also provide **Dye-labeled Calibration Oligos** used as a reference to calibrate real-time qPCR thermocyclers.

In addition, you will find in our catalog **PNA-FISH probes** for *in situ* hybridization and **negative and positive siRNA controls**.

No need to configurate those oligos. You can add your selected items directly to your cart.

GLEN RESEARCH SYNTHESIS REAGENTS

Eurogentec is **an exclusive distributor of Glen Research products** in most European countries. The Glen catalog includes many reagents and supports to synthesize DNA and RNA oligonucleotides.

- DNA & RNA nucleosides, analogs, and supports
- RNA synthesis
- Universal supports and synthesis supplies
- Modification and labeling
- Ancillary reagents
- Oligonucleotide purification supplies

Use the Fast Order option in your cart to order Glen Research product easily or contact us. [More infos about the Fast Order on p.62.]



LEARN MORE ON
CATALOG REAGENTS



DISCOVER OUR
GLEN RESEARCH
PRODUCT CATALOG



Order settings	51
Synthesis scale vs guaranteed yield	52
Compatibility cyclers channels vs fluorescent dyes	54
Determine the right synthesis scale for PCR	54
Order placement	57
Online configurators	59
Configure your oligos one by one	60
Order oligos by batch	61
Ask a quote for your complex oligos	62
Order a personal quote	63
Payment methods	64
Delivery	67
Shipping methods	68
Provided documentation	69
Handling instructions	71
Reconstitute your oligo	72
Storage	72
Quantify your oligo	73

Oligo ordering guidelines

We guide you through your ordering journey, from helping you select the best options for your oligos to ensuring their optimal use in your lab.



Order settings

We provide guidance to guarantee that your oligos meet the quantity and quality necessary for your application.

Compatibility cyclers channels vs fluorescent dyes

Table on the next page.

For complementary information, please refer to instrument manufacturer technical guide or contact us at scientific.support@eurogentec.com

Determine the right synthesis scale for PCR

*Please select in the minimum guaranteed yield table the synthesis scale corresponding to the desired minimum quantity.

FINAL OLIGO CONCENTRATION

50 nM 150 nM 300 nM 600 nM 900 nM

AVERAGE NUMBER OF REACTIONS
(TOTAL VOLUME 100µL)

MINIMUM QUANTITY
TO ORDER*

100	30	15	8	6	0.5 nmol
1000	300	165	80	55	5 nmol
5000	1650	830	415	275	25 nmol
10 000	3300	1660	830	555	50 nmol
100 000	33 300	16 660	8330	5555	500 nmol

LEGEND TABLE P.55

*perform a dye calibration for optimal results

Ch = Channel

YY = Yakima Yellow®

DFO = Dragonfly Orange™

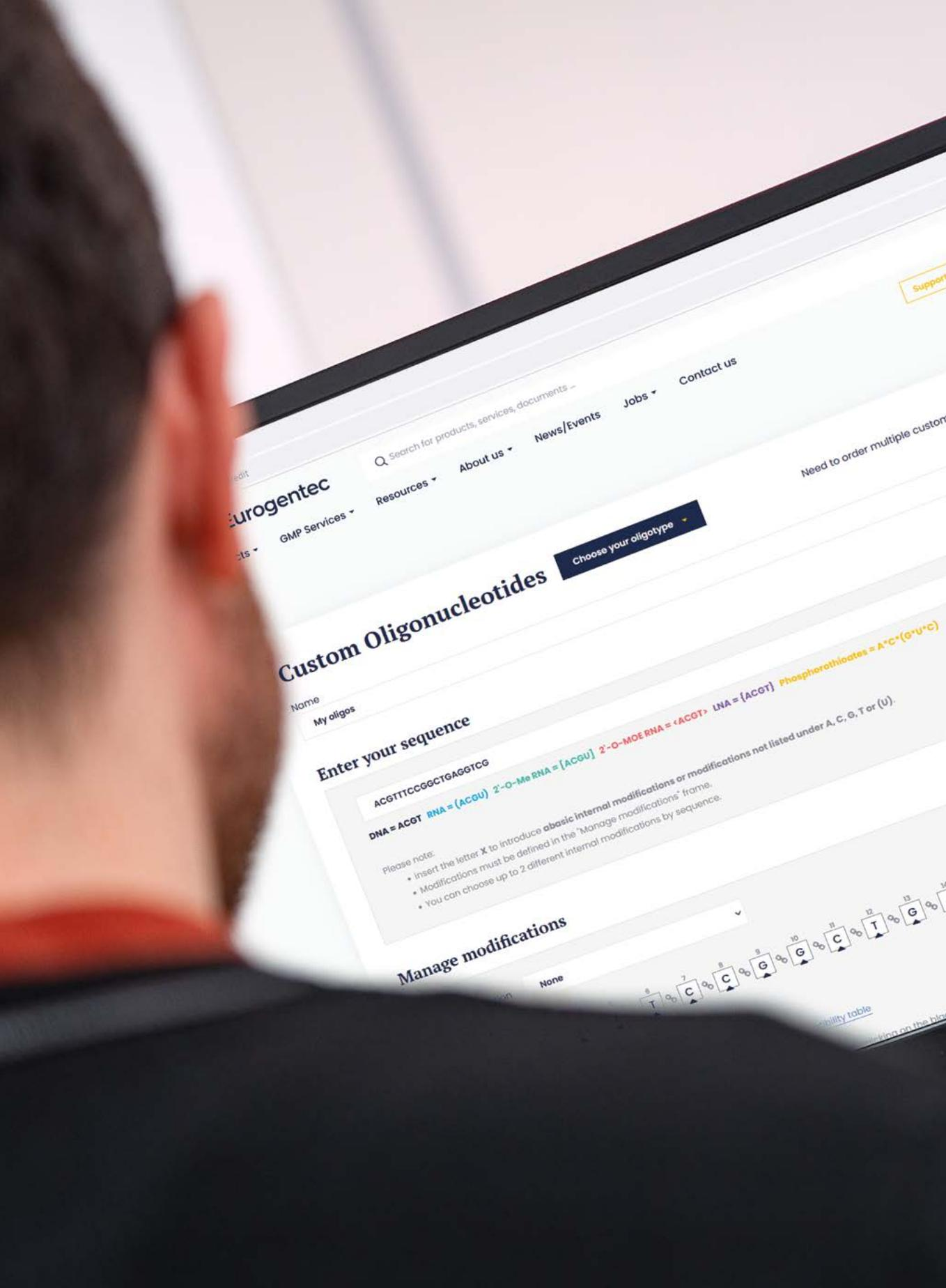
TR = Texas Red®

● In black = Recommended by Eurogentec

● In blue = Available at Eurogentec

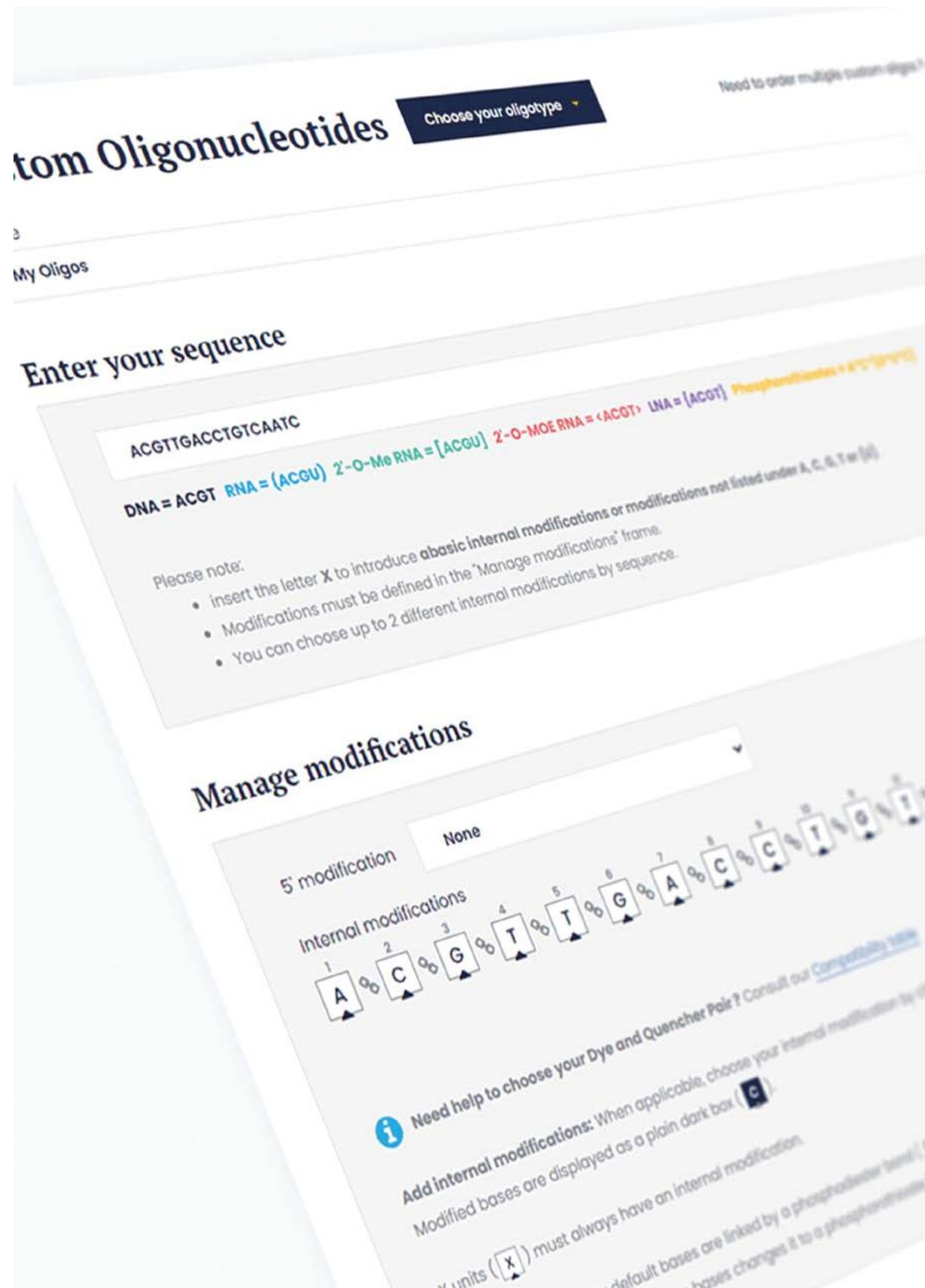
● In pale blue = Not available at Eurogentec

Thermocycler	Blue ch	Ch 1	Ch 2	Ch 3	Ch 4	Ch 5	Ch 6	Ch 7
Real-Time PCR Thermocyclers								
AriaMx & AriaDx (up to 6 filters)	ATTO 425	FAM	HEX	Gy ³	ROX	Gy ⁵		
qTOWER ³		FAM	YY/HEX/JOE/IC	DFO/TAMRA/NED	ROX/TR/Gy ^{3.5}	Gy ⁵	Gy ^{5.5}	
ABI Prism [®] SDS 7500		FAM	YY/HEX/JOE/IC/TET*	DFO/TAMRA/Gy ³ /NED	ROX/TR	Gy ⁵ /HL647		
ABI Prism [®] SDS 7900HT		FAM	YY/JOE/IC/TET*	DFO/TAMRA/NED	ROX			
QuantStudio [™] 1		FAM	YY/HEX/JOE/IC/TET	ROX/TR				
QuantStudio [™] 3		FAM	YY/HEX/JOE/IC/TET	DFO/TAMRA/Gy ³ /NED	ROX/TR			
QuantStudio [™] 5 (384-W) & 6 Flex/Pro		FAM	YY/HEX/JOE/IC/TET	DFO/TAMRA/Gy ³ /NED	ROX/TR	Gy ⁵ /HL647		
QuantStudio [™] 5 (96-W) & 7 Flex/Pro & 1.2k Flex		FAM	YY/HEX/JOE/IC/TET	DFO/TAMRA/Gy ³ /NED	ROX/TR	Gy ⁵ /HL647	Gy ^{5.5} /ATTO 700	
StepOnePlus		FAM	YY/HEX/JOE/IC/TET*	DFO/TAMRA/Gy ³ /NED	ROX			
ViiA7		FAM	YY/HEX/JOE/IC/TET	DFO/TAMRA/NED	ROX/TR*			
CFX Gained & CFX Duet		FAM	YY/HEX/JOE/IC/TET		ATTO 633/LIZ	ATTO 680/Alexa Fluor [®] 680		
CFX Opus & CFX96 Touch		FAM	YY/HEX/JOE/IC/TET	ROX/TR	Gy ⁵			
CFX384 Touch		FAM	YY/HEX/JOE/IC/TET	ROX/TR	Gy ⁵			
iCycler iQ [™]		FAM	YY/HEX/Gy ³ /TET	DFO/TAMRA/Gy ³ /NED	ROX/TR	Gy ⁵		
iQ [™] 5		FAM	YY/HEX/JOE/IC	DFO/TAMRA/Gy ³ /NED	ROX/TR	Gy ⁵		
QIAquant 96 & 384 5plex		FAM	YY/HEX/JOE/IC	ROX/TR	Gy ⁵	Gy ^{5.5}		
QIAquant 96 2plex		FAM	YY/HEX/JOE/IC					
RotorGene 6000 & Q		FAM	YY/HEX/JOE/IC/TET	TAMRA/ROX/Gy ^{3.5} /TR	Gy ⁵	ATTO 680		
LightCycler [®] 480 I	ATTO 425	FAM	YY/HEX/JOE/IC	ROX/TR/IC Red 610	ATTO 620/IC Red 640	Gy ⁵		
LightCycler [®] 480 II	ATTO 425	FAM	YY/HEX/JOE/IC	ROX/TR/IC Red 610	ATTO 620/IC Red 640	Gy ³ /Gy ^{5.5}		
LightCycler [®] 96		FAM	YY/HEX/IC	TR/IC Red 610	Gy ⁵			
Digital PCR Platforms								
Naica [®] Prism3		FAM	Gy ³ /HEX/IC	Gy ⁵				
Naica [®] Prism6		FAM	YY	ATTO 550	ROX	Gy ⁵	Gy ^{5.5} /ATTO 700	
Nico+		FAM	YY	ATTO 550	ROX	Gy ⁵	Gy ^{5.5}	DR-521-XL
QX One ddPCR		FAM	HEX/IC	Gy ⁵				
QX100 ddPCR		FAM	HEX/IC					
QX200 ddPCR		FAM	HEX/IC					
QuantStudio [™] Absolute Q		FAM	HEX/IC	Gy ⁵	Gy ^{5.5}	ROX	ATTO 590	
QIAcuity Four/Eight (5plex)		FAM	HEX/IC	TR/RBY	Gy ⁵ /LIN			
QIAcuity One (2plex)		FAM	HEX/IC	TAMRA	ROX	Gy ⁵		
Digital LightCycler	ATTO 425	FAM	HEX/IC	TR/CFR610/LC610	Gy ³ /A270/LC640	Gy ^{5.5}		



Order placement

We guide you step by step, simplifying your ordering journey to make it fast and effortless.



Order, save & manage your oligonucleotides easily with our online configurators.

Online configurators

Fast and easy for your research grade oligos.

Order, save, and manage your oligonucleotides easily with our online configurators for :

- Custom oligonucleotides
- Probes (Double-dye, MGB, Molecular Beacons)
- siRNAs
- NGS oligos

You can configure them one by one with our special tools, or order multiple oligos in a single batch thanks to our batch order.

NOMENCLATURE GUIDELINES TO WRITE YOUR SEQUENCE

CODE	BACKBONE
ACGTA	DNA
(ACGUA)	RNA
[ACGUA]	2'-O-Me RNA
<ACGTA>	2'-O-MOE
{ACGTA}	LNA®
A*C*(G*U)*A	Phosphorothioate links

IUB code

Mixed bases (also known as degenerate or wobble bases) follow the IUB codes :

IUB CODE	MIXED BASES
D	A/G/T
M	A/C
H	A/C/T
I	Inosine = Universal base
W	A/T
R	A/G
Y	C/T
V	A/C/G
S	C/G
K	G/T
N	A/G/C/T
B	C/G/T



WATCH TUTORIAL

Configure your oligos one by one



1 Connect to www.eurogentec.com



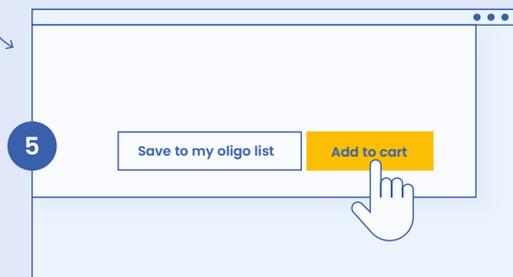
2 Go to the configurator page by clicking on the "Oligo configurators" in the "Resources" dropdown menu.



3 Select the oligonucleotide type via the dropdown menu available on the configurator page.



4 Fill in the configurator with your oligonucleotide specifications.



5 Add your oligonucleotide into the cart and finalize your order.



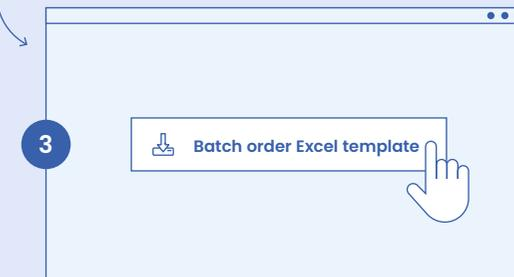
Order oligos by batch



1 Connect to www.eurogentec.com

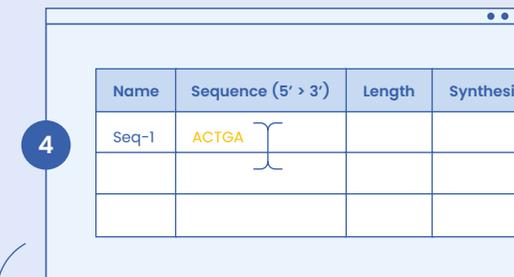


2 Go to the batch order section clicking on the "Oligo Batch order" link in the "Resources Menu".

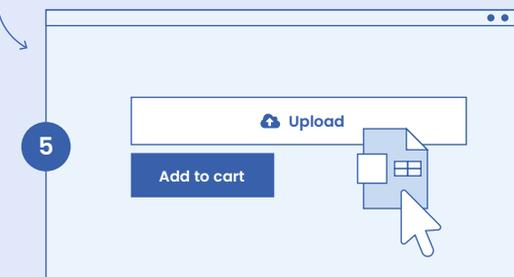


3 Download the batch order file.

NB : You need a spreadsheet software to edit the file.



4 Fill in the file with your oligonucleotides.



5 Upload the file and finalize your order.



We synthesize your highly modified oligonucleotides even with specifications out of our catalog offer.

Ask a quote for your complex oligos

Each request is analyzed by our experienced chemists prior to drawing up any quotation.

We manufacture complex oligonucleotides with specifications out of our catalog offer, such as atypical chemistries, rare modifications, small or large scales, personalized labels and packages...

Contact us through the form on our website and receive all the information regarding technical feasibility, pricing and turnaround time.

Our complex oligonucleotides service will bring you a tailor-made solution.

FAST ORDER

The Fast Order allows you to select products by entering either the product reference or name, adding them directly to your cart.

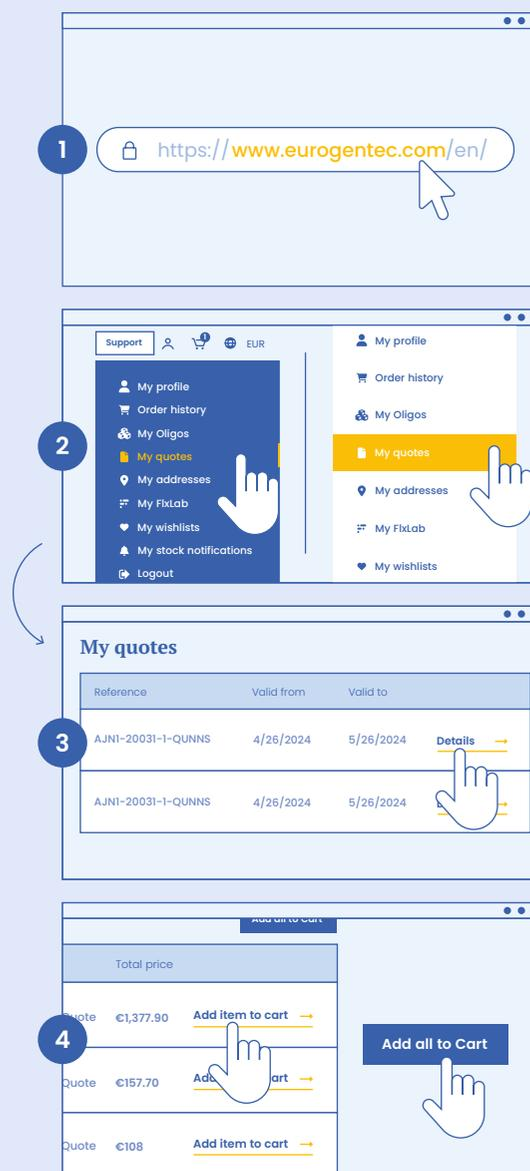
Here's how you can make the most of it :

- 1 Go to the Fast Order box in your cart.
- 2 Search for the desired product's reference or name in the Fast Order box.
- 3 Click on the product name, and it will be added to your cart.



LEARN MORE ABOUT FAST ORDER

Order a personal quote



1 Connect to www.eurogentec.com

2 You can easily order your quotes from your account, in "My quotes" section or *via* the dropdown menu from the account icon.

3 Select the desired quote. You will see the details of the available products.

4 You can add one item by one item or send all items directly to the cart.

NB : Please note that you can't add more than one quote to your cart per order and you have to go through this process to benefit the price from your quote. This is only available for catalog references.

MY OLIGOS TAB

In "My Oligos" section, you can create new oligonucleotides or select oligonucleotides in your saved sequences.

Each time you add an oligo to your cart, it will be added to your saved oligos in "My oligos tab" automatically.



ASK A QUOTE

Payment methods

Our standard methods of payment

We offer different payment methods to best suit your needs: Visa, MasterCard, Bank Transfers, and FlxLab. All these methods meet high-security standards to protect your payments.

Credit cards



Bank transfers (Payment on invoice)



FlxLab accounts



FLXLAB IS EUROGENTEC'S PREPAID SYSTEM

Place a defined amount on your FlxLab account, and use that amount over time. Please note that your account is credited only upon payment.

Only one invoice

A single invoice for multiple oligos orders over time.

Multiple accounts

One administrator can give restricted access to multiple users in a single FlxLab account.

Transactions history

Keep tracks of your budget and order history.



←
[LEARN MORE](#)

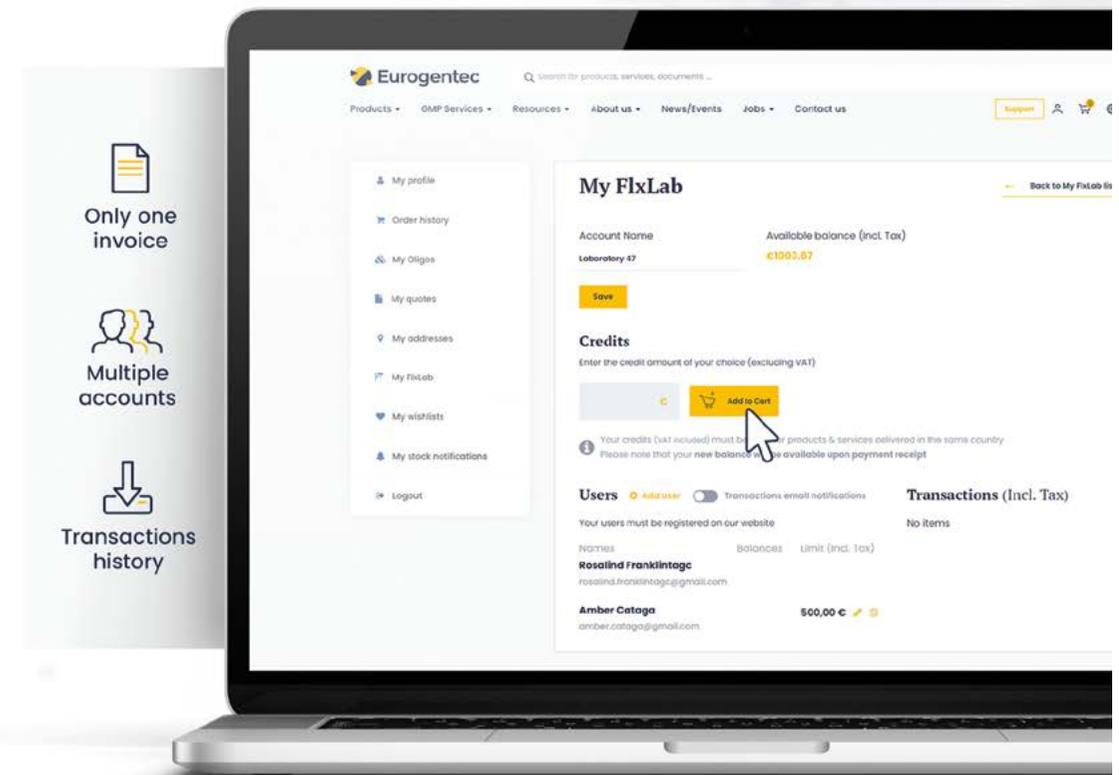
OUR FREE PAYMENT SERVICE



Manage your lab budget with less paperwork



←
[SCAN TO LEARN MORE](#)





Delivery

We focus on making the delivery process as simple and efficient as possible, ensuring that you receive everything you need on time.

Shipping methods

Express Delivery

Receive your oligonucleotides within 24 to 48 hours by international carrier.

Same day shipping Option

- Available on request (additional fees will be charged)
- For orders received before 10.00 AM (Central European Time)
- For custom oligonucleotides (max 24), 10/40 nmol scale, 5-30 DNA bases, unmodified, SePOP desalted or RP-Cartridge purified.

For large orders or unique oligonucleotides, please feel free to contact us at oligo@eurogentec.com to receive more details in terms of delivery schedules.

5'AP, BSA, HRP or SBP Conjugation: 3-5 Working Days Extra;

Additional Purification or Services: 2 Working Days Extra;

These values are indicative.

Delivery times (in working days)

Delivery time according to purification

Range	Product	Length (Bases)	SePOP	RP-Cartridge-Gold™	HPLC (RP or IEX)	PAGE	Dual HPLC
Custom Oligonucleotides	Non-Modified (DNA Only)	5-9	-	4-5	5	6	7
		10-39	2-3	4-5	5	6	7
		40-59	5	6	7	8	-
		60-79	-	6	-	8	-
		80-139	-	-	-	10	-
	Modified (including DNA, RNA, 2'-O-Me RNA, 2'-O-MOE RNA & LNA*)	10-39	5	7	7	8	9
Real-Time qPCR Probes	Double-Dye Probes	8-45	-	-	7	-	-
	Molecular Beacons	28-50	-	-	12-15	-	-
	MGB Taqman Probes	8-30	-	-	5-7	-	-
RNAi Oligonucleotides	siRNA Duplexes	21-27	5-7	-	9-10	10-11	-
NGS Oligonucleotides	-	20-85	-	4-6	5-7	-	-
Universal Primers	-	15-38	-	-	2-3	-	-
Unique Oligonucleotides	-	2-225	On Request				

Provided documentation

Each oligonucleotide is provided with a technical data sheet. Other documentations could be added depending on the oligonucleotide type. All the documents are sent as pdf files to your shipping email address.

	TDS	MS ¹	UHPLC ²
Custom Oligonucleotides	Unmodified	✓	
	Modified	✓	✓ ³
Real-Time qPCR Probes	✓	✓	✓
RNAi Oligonucleotides	siRNA Duplexes	✓	✓ ³
Universal Primers	✓	✓	
Unique Oligonucleotides	✓	✓	✓ ⁴
NGS Oligos	✓	✓	
Calibration Oligos	✓	✓	

TDS Technical Data Sheet; **MS** Mass Spectrometry; **UHPLC** Ultra High Performance Liquid Chromatography;

¹ Always provided up to 60 bases long Oligonucleotides.

² If applicable.

³ Except for SePOP desalted oligonucleotides.

⁴ Optional.

For technical reasons this general rule may be adapted to provide you with the most suitable and useful documentation.



Handling instructions

This resource offers all the information required to ensure the optimal use of your oligonucleotides.

Reconstitute your oligo

Refer to the dedicated technical data sheet for more information.

1. **Spin** the tube briefly to collect the pellet at the bottom of the tube.
2. **Add** the appropriate volume of recommended buffer.
3. Allow the tube to **stand** for a few minutes.
4. **Stir** the tube for 15 sec using a vortex shaker and **centrifuge** it briefly.

Oligo storage

Handling information

PRODUCTS	FORMAT	STORAGE ¹	STABILITY ²
Custom Oligonucleotides	Dried	RT	18 months
	TE Buffer (pH 8) ³ or dH2O	-20 °C	24 months
Real-Time qPCR Probes	Dried	RT	18 months
	TE Buffer (pH 8) ³ or dH2O	-20 °C	24 months
RNAi Oligonucleotides	Dried	RT	18 months
	RNase-free Buffer (pH 7.5)	-20 °C	24 months
Catalog Primers	Dried	RT	18 months
PNA FISH Probes / Custom PNA	Dried	RT	18 months

¹ Oligonucleotides and especially dye-labelled oligonucleotides should be protected from light for optimal stability. Tolerance -20°C ± 5° C. For custom PNA we suggest to prepare workable samples. Store these samples at 4°C for up to one month and the remaining at -20°C.

² Please note that depending on sequences and modifications, the stability of the oligos may vary substantially versus the given values, which should therefore be considered as indicative. Although it is generally recommended to avoid freeze-thaw cycles, we did not observe negative effects in qPCR tests after 10 freeze-thaw cycles.

³ Except for Cy® dye labelled oligonucleotides (pH 7)

Quantify your oligo

1. Concentration of your stock solution (c)

1. Prepare an aliquot of the resuspended oligonucleotides to 1ml final volume of distilled H₂O and spin briefly.
2. Measure the absorbance at 260 nm (A₂₆₀).
3. Calculate the concentration of your stock solution using this formula:

The formula is valid for absorbance A₂₆₀ ≤ 1.2

$$c(\mu\text{g/ml}) = A_{260} \times \text{dilution factor} \times \text{Weight per OD}_{260} \text{ of stock solution } (\mu\text{g/OD})$$

OD₂₆₀ measure

1 OD₂₆₀ (Optical Density) unit is defined as the amount of oligonucleotide which, when dissolved in a volume of 1.0 ml, results in an absorbance of 1.0 when measured at 260 nm in a 1 cm path-length quartz cuvette.

1 OD₂₆₀ unit corresponds to approximately 33 µg of single strand DNA. These relationships, however, can be inaccurate for short fragments of DNA, such as oligonucleotides. Base composition and even linear sequence will affect optical absorbance. Hence the precise value of the OD to mass relationship is unique for each oligonucleotide.

We carefully measure the OD value for your custom oligonucleotide by measuring the absorbance at 260 nm using UV spectrophotometer. This information is provided on the oligonucleotide Technical Data Sheet as the number of OD₂₆₀ units. The amount of oligo expressed in nanomoles and micrograms is derived from the OD measurement. →

TECHNICAL DATA SHEET (TDS)

You will find the precise data of your oligo on the Technical Data Sheet sent with each oligo.

2. Convert OD₂₆₀ into Nanomoles

This equation shows the relation between the oligo amount in nanomoles and the OD₂₆₀ value.

$$\text{Nanomoles} = (\text{OD}_{260} / \epsilon_{260}) \times 10^6$$

The molar extinction coefficient (ϵ_{260})

ϵ_{260} is the extinction coefficient at 260 nm.

$$\epsilon_{260} = 2 \times \left(\sum_1 \epsilon_{\text{Nearest Neighbour}}^{n-1} \right) - \sum_2 \epsilon_{\text{Individual}}^{n-1} + \sum_1 \epsilon_{\text{Modification}}^n$$

where $\epsilon_{\text{Nearest Neighbour}}$ is the nearest neighbour constant for a pair of bases, $\epsilon_{\text{Individual}}$ is the constant for an individual base, and n is the length of the oligonucleotide.

3. Convert Nanomoles into Micrograms

Use the following equation, using the nanomole value and the molecular weight of the oligonucleotide.

$$\text{Micrograms} = \text{Molecular Weight} \times \text{Nanomoles} \times 10^{-3}$$

The molecular weight (MW)

$$\text{Anhydrous MW (g/mol)} = \sum_{\text{Individual Base}} \text{MW} + \sum_{\text{Individual Mods}} \text{MW} - 63.98 + 2.016$$

For DNA bases: MW dA = 313.21; MW dC = 289.18; MW dG = 329.21; MW dT = 304.20; MW dU = 290.17; MW dI = 314.19

For RNA bases: MW DNA counterpart + 16

When determining the weight of Uracil (rU) start with dU and not dT

For LNA bases: MW DNA counterpart + 16 (+42 for dC)

For 2'-O-Methyl bases: MW DNA counterpart + 30.03

When determining the weight of mU start with dU and not dT

For phosphorothioated bases: MW DNA counterpart + 16.06

THE MOLAR EXTINCTION

COEF (ϵ_{260})

Example:

1 OD₂₆₀ unit of primer M13
Forward, 5'-GTA AAA CGA CGG
CCA GTG-3'
Molar extinction coefficient
(ϵ_{260}) = 182.800 L / (mole × cm)

Nanomoles = (1.0 / 182.800)
× 10⁶ = 5.47 nmoles

THE MOLECULAR WEIGHT

(MW)

Example:

1 OD₂₆₀ unit of primer M13
Forward,
5'-GTA AAA CGA CGG CCA
GTG-3'

Molecular Weight = 5558.7

Micrograms = 5558.7 × 5.47
× 10⁻³ = 30.4 µg

Licenses & trademarks

Alexa Fluor® is a registered trademark of Life Technologies Corp.

BHQ® is a registered trademark of Biosearch Technologies, Inc.

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

BODIPY® is a registered trademark of Life Technologies Corp.

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

Cascade Blue® is a registered trademark of Life Technologies Corp.

Cy® and CyDye® are registered trademarks of Cytiva.

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

Eclipse™ Dark Quencher is a trademark of ElitechGroup, Inc.

Eclipse™ Quencher is a trademark of ElitechGroup, Inc.

Excel™ is a trademark of Microsoft Corp.

Eurogentec™ is a trademark of Kaneka Eurogentec S.A.

Gold'N Oligo® is a registered trademark of Kaneka Eurogentec S.A.

Hilyte™ is a trademark of Anaspec, Inc.

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

Marina Blue® is a registered trademark of Life Technologies Corp.

MGB™ is a trademark of ElitechGroup, Inc.

MGB Eclipse® is a registered trademark of ElitechGroup, Inc.

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

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

Oregon Green® is a registered trademark of Life Technologies Corp.

Pacific Blue™ is a trademark of Life Technologies.

Corp.Plexor™ is a trademark of Promega Corporation.

Pulsar™ is a trademark of Biosearch Technologies, Inc.

QXL® is a registered trademark of Anaspec, Inc.

Rhodamine Red™ is a trademark of Life Technologies Corp.

ROX™ is a trademark of Applera Corp.

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

SYBR® Green is a registered trademark of Life Technologies Corp.

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

Taqman® is a registered trademark of Roche Diagnostics GmbH.

Texas Red® is a registered trademark of Life Technologies Corp.

Vic® is a registered trademark of Applera Corp.

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

Contact us

Headquarters

Belgium

+32 (0) 4 372 74 00

5, rue Bois Saint-Jean
4102 Seraing - Belgium

Europe

France

+33 (0) 2 41 73 33 73

152, Avenue Patton CS 80111
49001 Angers - France

United Kingdom

00 800 6660 0123

1 The Square Fawley
Southampton Hampshire
SO45 1DD - UK

Germany

+49 (0) 221 258 94 55

c/o debelux-Handelskammer
Friedrichstrasse, 42-44 50676
Köln - Deutschland

The Netherlands

+31 (0) 43 352 06 98

Avenue Ceramique 223
6221 KX Maastricht - The
Nederland

Switzerland

+32 (0) 4 372 74 00

Boulevard des Philosophes, 3
1205 Genève - Switzerland

America

USA

+1 (800) 452 5530

34801, Campus Drive
Fremont, CA 94555
USA

More details about our oligonucleotide services

Therapeutic services



DOWNLOAD FLYER

Diagnostic kits manufacturing



DOWNLOAD FLYER

Access™ Dyes & Quenchers



DOWNLOAD FLYER

Preclinical studies



DOWNLOAD FLYER

Takyon® Ultra



DOWNLOAD FLYER

MGB probes



DOWNLOAD FLYER



©Eurogentec 2024 - All rights reserved



www.eurogentec.com

Kaneka Eurogentec SA

5, rue Bois Saint-Jean
4102 Seraing - Belgium

T +32 (0) 4 372 74 00
F +32 (0) 4 372 75 00

RPM Liège
TVA (BE) 0427.348.346

PART OF
kaneka