



Track Oligonucleotides



Eurogentec is proud to introduce the **Track Oligonucleotides** as your right-priced solution. With over 25 years experience in Oligonucleotide Manufacturing, Eurogentec ensures high quality products & services covering research to commercialization.

Track Oligonucleotides offer the required traceability, including identification of the equipment used, the personnel involved and other important manufacturing specifications during the entire production process. And, all this, at an affordable price!

With the track oligos, you have the traceability required for most of the diagnostic market applications: Lab Services (Health, Agro, Veterinary, Environment), Hospital Lab, Agro Lab, MDx Companies...

Lab Services
Hospital & Clinical Labs

Diagnostic Companies

Discovery

Routine Assays

Contamination
Sensitive Assays

Discovery

Feasibility

Prototyping

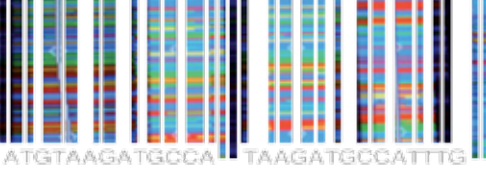
Validation

Commercialization

Oligonucleotide Grades	Research	Track	Pre-Diagnostic	Diagnostic
Process				
Dedicated Account Contact Person	Option	✓	✓	✓
Customized Fill & Finish	Option	Option	✓	✓
Quality Management				
ISO 9001 Certification	✓	✓	✓	✓
ISO 13485 Certification	-	-	✓	✓
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]	-	✓	✓	✓



- ⊕ **Custom Backbones:** DNA - RNA - 2' O-Me RNA - Phosphorothioate
- ⊕ **Modifications:** >250 Modifications Available incl. Proprietary Dyes (HiLyte™) and Quenchers (QXL™)
- ⊕ **Purification Chemistries:** RP/IE HPLC [Single or Dual]
- ⊕ **QC Validated Methods:** MS, HPLC and CGE
- ⊕ **Optional Services:** Additional QC - Customized Fill & Finish



Track Oligonucleotides



All products supplied are for Research and Development purposes only and may not be used for commercial, clinical, diagnostic, therapeutic, in humans or any other use for which an additional license is required. Customer will be responsible at its expenses for obtaining on behalf of Eurogentec all necessary consents and licenses with respect to the use of any third party intellectual property rights for process development and manufacturing of this project.

HilLyte™ dyes: Use of this product is covered by one or more of the following US patents owned by AnaSpec, Inc.: USP 7,465,810, USP 7,754,893, USP 7,820,783, USP 8,258,292, and divisionals, continuations, continuations-in-part, reissues, substitutes, and extensions thereof. The purchase of this product conveys to the buyer a limited, non-exclusive, nontransferable right (without the right to sell, repackage, or further sublicense) under these patent rights to use only the amount of product purchased for the purchaser's own internal research. No other license is granted to the buyer whether expressly, by implication, by estoppel or otherwise. In particular, the purchase of this product does not include nor carry any right or license to use, develop, or otherwise exploit this product commercially, and no rights are conveyed to the buyer to use the product or components of the product for any other purposes, including without limitation, provision of services to a third party, generation of commercial databases, or clinical diagnostics or therapeutics. This product is for research use only. For information on purchasing a license to this product for purposes other than research, contact licensing@eurogentec.com.

OXL® Quenchers: Use of this product is covered by one or more of the following US patents owned by AnaSpec, Inc.: USP 7,910,753, USP 8,093,411, and divisionals, continuations, continuations-in-part, reissues, substitutes, and extensions thereof. The purchase of this product conveys to the buyer a limited, non-exclusive, nontransferable right (without the right to sell, repackage, or further sublicense) under these patent rights to use only the amount of product purchased for the purchaser's own internal research. No other license is granted to the buyer whether expressly, by implication, by estoppel or otherwise. In particular, the purchase of this product does not include nor carry any right or license to use, develop, or otherwise exploit this product commercially, and no rights are conveyed to the buyer to use the product or components of the product for any other purposes, including without limitation, provision of services to a third party, generation of commercial databases, or clinical diagnostics or therapeutics. This product is for research use only. For information on purchasing a license to this product for purposes other than research, contact licensing@eurogentec.com.



More info



Please visit

www.eurogentec.com/track-oligos.html

Contact us



Please contact us at

track@eurogentec.com

Our specialists are available throughout the entire maturation of your project.