

Products & Services

For Life science Research

Eurogentec provides high-quality reagents for both proteomic and genomic applications. The company is known for its flexible and comprehensive valuable services related to the production and development of custom kits and reagents for research use. Eurogentec has a decades-long experience in the field. This ensures the guarantee of getting the products that best suit your needs.



Oligos



PCR/qPCR



Genes/DNA



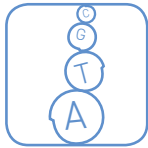
Antibodies



Peptides



Proteins



Oligos



- > Custom oligonucleotides
- > NGS oligonucleotides
- > RNAi oligonucleotides
- > Real-Time qPCR probes
- > Highly complex oligonucleotides
- > Catalogue oligonucleotides
- > Synthesis reagents



PCR/qPCR

Amplification technologies



- > 2X and 5X ready to use Takyon™ qPCR MMx
- > RT Stable lyophilised Takyon™ qPCR & qRT-PCR MMx
- > Takyon™ qPCR Core kits for adjustable setups
- > qPCR Controls
- > PCR Mixes and consumables
- > Plastics



SMART Gene Analysis



- > Custom genes
- > DNA/RNA Purification kits
- > DNA Extraction kits
- > Electrophoresis devices and consumables
- > Transformation products

SPECIFICATIONS

All types of oligonucleotides, from 2 to 225 bases

All chemistries: DNA, RNA, LNA®, 2'O-Me, PNA, 2'O-MOE...

More than 300 modifications

Synthesis scales from µg to grams

Various purification methods adapted to the application

Custom fill & finish

SPECIFICATIONS

From individual reagents and consumables to ready-to-use kits

Universal capabilities (end point or quantitative, SYBR®, Probes, Rox, HTP, One-Step converter...)

Suitable for multiplex

Clean detection of low copy targets

Maximum reproducibility of results

Efficient in small volume reaction

Custom formulation & OEM possibilities



Easy & user friendly web oligo configurators



1 Probe ordered
2 Free primers offered¹

¹ Valid for Double-Dye or Molecular Beacon Probe. Each primer must be 15-30 DNA bases, unmodified, purified by RP Cartridge and at the same synthesis scale as the probe (10, 40, 200 or 1000 nmol).



Select your qPCR kits online



First 150 RXNs sample for free
+ 50% on your first order²

² Sample for free offer is only valid on your first order. The samples are available for the MasterMix 2X and 5X



HIGHLIGHT

NGS Oligonucleotides: an optimised process ensuring a low cross contamination level <0,1%

Oli&GO™ prepaid system: Only one invoice for multiple orders

HIGHLIGHT

Takyon™ Dry: RT Stable qPCR/qRT-PCR MasterMixes for Probe Assays

Takyon™ 5X: 5X qPCR and qRT-PCR MasterMixes to maximise the template volume in the reaction mix

HIGHLIGHT

SMART Pure: for nucleic acid purification and extraction from various sample types

SMART Extract: for rapid (less than 10') column-less extraction from any type of tissue

Custom Genes: up to 50 Kbp sequence synthesis



Peptides



- > Custom peptides
- > Quantified peptides
- > Heavy isotope labelled peptides
- > FRET peptides
- > Cyclic peptides
- > Peptide-oligo conjugates
- > Catalogue peptides

SPECIFICATIONS

Length: from 2 up to 60 aa³

Quantity: 1 mg minimum to gram quantities (gross or net weight)

Purity: from crude up to >95%

³ depending on the peptide sequence



Check out the website to view the full list of catalogue peptides related to neurosciences, cancer, diabetes...

HIGHLIGHT

Quant-peptides: accurately quantified for high precision proteomics

Peptide-oligo conjugates: modified or not they constitute promising therapeutics



Antibodies



- > Custom monoclonal antibodies
 - Mouse hybridoma development
 - *In vitro* mAb production in bioreactors, roller bottles or CL1000
- > Custom polyclonal antibodies
 - Proprietary Speedy 28-Day programmes
 - Generation of anti-PTM antibodies
 - Genetic immunisation
- > Custom Ab fragments including VHH from Llama
- > Antibody purification, labelling and testing
- > Primary and secondary antibodies

SPECIFICATIONS

Stable and trusted antibody services since 1996

More than 1500 programmes every year

From ready-to-use catalogue antibodies to fully customised production programmes

From research antibodies to GMP production of antibody fragments

Highly skilled project managers dedicated to your project

HIGHLIGHT

PTM programmes: we can raise antibodies against phosphorylation, acetylation, ubiquitylation, SUMOylation....

Large project management: We are able to handle large demanding projects with remarkable flexibility to meet the customer's requirements and SOP's



Protein identification



- > Recombinant proteins manufacturing
- > Labelling service, kits and reagents
- > Protein analysis products (PAGE Gels, WB kits, devices and consumables)
- > iD Sensolyte® Assay kits for activity detection of specific proteases and other enzymes in biological samples



HIGHLIGHT

AnaTag™ protein labelling kits: get purified labelled proteins in only 4 steps

iD Blocking Kit: a 5-min blocking step for any membrane or solid surface of multi-well plates

www.eurogentec.com

Products & Services

Kaneka Eurogentec S.A.

Contact:

For more information, please contact
your Sales representative
or contact us at:
isales@eurogentec.com

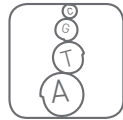
Toll free number: 00 800 666 00 123

© Eurogentec 2017

Follow us on



Life Science Research



Oligos



PCR/qPCR



Genes/DNA



Antibodies



Peptides



Proteins

DOWNLOAD

our White Papers on www.eurogentec.com



DISCOVER

the product range brochures on www.eurogentec.com

