

Oligonucleotides

# Aptamers

Ingeniously shaped to fit.



Eurogentec collaborates with Novaptech, Dr. Jean-Jacques Toulmé's French company, to provide high-value aptamer service.



## Custom RNA & DNA aptamer development

Novaptech identifies candidates by automated selection (SELEX) and generates optimized aptamers with strong binding affinity and high specificity for their targets.

Eurogentec synthesizes the final aptamer candidate and all the sequences required for the development process of your aptamers at Novaptech.

## Aptamer features

### From research to GMP therapeutic grade

Each aptamer sequence is produced with the appropriate grade (Research or GMP) according to the final applications (R&D, Diagnostics or Pre-clinical assays....).

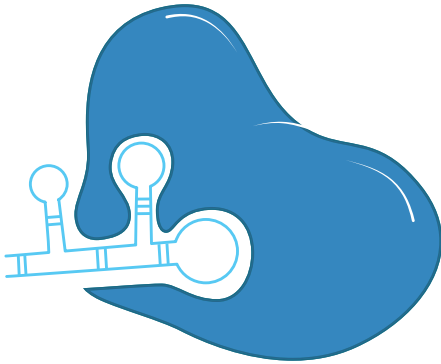
### From small to large scale – Any modification

For any upcoming production request Eurogentec will synthesize the oligonucleotide aptamer according to setups defined during the development process. Eurogentec will be able to deliver larger amounts of oligonucleotide aptamers with any chemical modification according to the customer's needs.

More info & contact:

[www.eurogentec.com/oligonucleotide-aptamers.html](http://www.eurogentec.com/oligonucleotide-aptamers.html)





## What is an aptamer?

Oligonucleotide aptamers are single stranded RNA or DNA sequences folded in a specific 3D conformation. Their tridimensional structure ensures complementarity with their target.

The perfect fit between both entities gives to the nucleic acid-based aptamers a high affinity and specificity for a wide range of molecules such as proteins/peptides, nucleic acids, small molecules, toxins, cells and viruses.

### GOOD ALTERNATIVE TO ANTIBODIES

- Similar and **even higher affinity** and specificity than traditional monoclonal antibodies.
- **Low** immunogenicity
- Chemical process requiring **no animal** or cell.
- **Small size** (10 times smaller than antibodies).

### FAST AND CONTROLLED PRODUCTION

- **Scalable** *in vitro* production.
- **Research and GMP** grades.
- **Batch to batch** reproducibility.

### ANY TARGET / VARIOUS CONDITIONS

- Effective even **against non-immunogenic** or toxic targets.
- **High stability** at room temperature and resistant to thermal denaturation.
- **Easy chemical modification** for improved efficiency, even in biological media.

Contact us

[www.eurogentec.com/oligonucleotide-aptamers.html](http://www.eurogentec.com/oligonucleotide-aptamers.html)