

INGENIOUSLY SHAPED TO FIT

Do you need Oligonucleotide Aptamers?



Eurogentec and Novaptech combine their expertise

We provide high-value services for the development and production of your nucleic acid-based aptamers.



Fast and controlled
production



Scalable
from µg to g



Large choice
of chemical modifications
for improved
performances



Various grades
for research, analytical,
diagnostic and
therapeutic applications